

# VWR® ULT Upright Freezers

## INSTRUCTION MANUAL

| North American Catalogue Number | European Catalogue Number | Model Number | Size (cu ft) | Size (liters) | Box | Voltage       | Temperature |
|---------------------------------|---------------------------|--------------|--------------|---------------|-----|---------------|-------------|
| <b>-86°C FREEZERS</b>           |                           |              |              |               |     |               |             |
| 10160-724                       | --                        | VWR24086A    | 13           | 368           | 240 | 120V/60Hz     | -86°C       |
| 10160-726                       | --                        | VWR24086D    | 13           | 368           | 240 | 208-230V/60Hz | -86°C       |
| --                              | 471-1140                  | VWR24086V    | 13           | 368           | 240 | 230/50Hz      | -86°C       |
| 10160-728                       | --                        | VWR32086A    | 17           | 490           | 320 | 120V/60Hz     | -86°C       |
| 10160-730                       | --                        | VWR32086D    | 17           | 490           | 320 | 208-230V/60Hz | -86°C       |
| --                              | 471-1141                  | VWR32086V    | 17           | 490           | 320 | 230/50Hz      | -86°C       |
| 10160-732                       | --                        | VWR40086A    | 23           | 651           | 400 | 120V/60Hz     | -86°C       |
| 10160-734                       | --                        | VWR40086D    | 23           | 651           | 400 | 208-230V/60Hz | -86°C       |
| --                              | 471-1142                  | VWR40086V    | 23           | 651           | 400 | 230V/50Hz     | -86°C       |
| 10160-736                       | --                        | VWR60086D    | 28           | 793           | 600 | 208-230V/60Hz | -86°C       |
| --                              | 471-1143                  | VWR60086V    | 28           | 793           | 600 | 230/50Hz      | -86°C       |
| <b>-40°C FREEZERS</b>           |                           |              |              |               |     |               |             |
| 10160-748                       | --                        | VWR24040A    | 13           | 368           | 240 | 120V/60Hz     | -40°C       |
| 10160-750                       | --                        | VWR24040D    | 13           | 368           | 240 | 208-230V/60Hz | -40°C       |
| --                              | 471-1135                  | VWR24040V    | 13           | 368           | 240 | 230V/50Hz     | -40°C       |
| 10160-752                       | --                        | VWR32040A    | 17           | 490           | 320 | 120V/60Hz     | -40°C       |
| 10160-754                       | --                        | VWR32040D    | 17           | 490           | 320 | 208-220V/60Hz | -40°C       |
| --                              | 471-1136                  | VWR32040V    | 17           | 490           | 320 | 230V/50Hz     | -40°C       |

**Important installer and user information:**

A redundant temperature sensing device has been included in this ULT freezer. This device is a type “T” thermocouple. For convenient access, the thermocouple (Figure 2-3) terminates in an interconnect jack (Figure 2-5) behind the base front cover. (May be located differently in chests. See Section 2.) It is strongly recommended that this thermocouple be attached to a redundant 24 hour 7 day monitoring system with alarm capabilities. Connecting the sensor to a monitoring and alarm system separate from the freezer provides the utmost in product safety, should the integral system fail.

| <b>Legal Address of Manufacturer</b>                | <b>Legal Address of Manufacturer</b>              |
|---|---|
| United States                                       | Europe  |
| VWR International, LLC                              | VWR International bvba                            |
| 100 Matsonford Rd                                   | Researchpark Haasrode 2020                        |
| Radnor, PA 19087                                    | Geldenaaksebaan 464                               |
| 800-932-5000  | B-3001 Leuven                                     |
| <a href="http://www.vwr.com">http://www.vwr.com</a> | + 32 16 385011                                    |
|   | <a href="http://be.vwr.com">http://be.vwr.com</a> |

Country of origin: United States

**Packing List**

| <b>Description</b>     | <b>Quantity</b> |
|------------------------|-----------------|
| Key                    | 2               |
| Neoprene Cap           | 2               |
| 1/4-20 x 5-1/2" Bolt   | 2               |
| Retaining Clip         | 1               |
| Remote Alarm Connector | 1               |

**MANUAL NUMBER 7085602**

|   |                   |          |  |     |
|---|-------------------|----------|--|-----|
| 8 | 31230             | 10/23/17 | Updated drawing 5602, 5702 bulkhead drawing  | bpg |
| 7 | 41728/41715       | 10/04/17 | Added door operation info/Removed condenser info, update acc chart, added schematics | bpg |
| 6 | 41370             | 8/14/17  | Updated BUS board information  | bpg |
| 5 | 41348/41159/40543 | 5/12/17  | Corrected intended use temp range, voltage rating, F-gas statement                   | bpg |
| 4 | 40797             | 4/7/16   | Updated Low Battery alarm to 8 hours, chart on pg 5-1                                | ccs |
| 3 | 40544             | 12/9/15  | BUS information (preventive maintenance and BUS control panel and operation)         | ccs |



**Important** Read this instruction manual. Failure to read, understand and follow the instructions in this manual may result in damage to the unit, injury to operating personnel, and poor equipment performance. ▲

**Caution** All internal adjustments and maintenance must be performed by qualified service personnel. ▲

Material in this manual is for information purposes only. The contents and the product it describes are subject to change without notice. VWR International makes no representations or warranties with respect to this manual. In no event shall VWR be held liable for any damages, direct or incidental, arising out of or related to the use of this manual.

**Intended Use:** The VWR freezers (see cover page for specific models) described in this manual are high performance units for professional use. These products are intended for use as cold storage in research use and as a general purpose laboratory freezer, storing samples or inventory at operating temperatures between -50°C and -86°C, or -10°C and -40°C, depending on model.

It is not considered a medical device and has therefore not been registered with a medical device regulatory body (e.g. FDA); that is, it has not evaluated for the storage of samples for diagnostic use or for samples to be re-introduced to the body.

**This unit is not intended for use in classified hazardous locations, not to be used for the storage of flammable inventory.**

©2015 VWR® International. All rights reserved.



Important operating and/or maintenance instructions. Read the accompanying text carefully.



Potential electrical hazards. Only qualified persons should perform procedures associated with this symbol.



Equipment being maintained or serviced must be turned off and locked off to prevent possible injury.



Extreme temperature hazards, hot or cold. Use special handling equipment or wear special, protective clothing.



This equipment is marked with the crossed out wheeled bin symbol to indicate that this equipment must not be disposed of with unsorted waste.

Instead it's your responsibility to correctly dispose of your equipment at lifecycle -end by handling it over to an authorized facility for separate collection and recycling. It's also your responsibility to decontaminate the equipment in case of biological, chemical and/or radiological contamination, so as to protect from health hazards the persons involved in the disposal and recycling of the equipment.

For more information about where you can drop off your waste of equipment, please contact your local dealer from whom you originally purchased this equipment.

By doing so, you will help to conserve natural and environmental resources and you will ensure that your equipment is recycled in a manner that protects human health.

- ✓ Always use the proper protective equipment (clothing, gloves, goggles, etc.)
- ✓ Always dissipate extreme cold or heat and wear protective clothing.
- ✓ Always follow good hygiene practices.
- ✓ Each individual is responsible for his or her own safety.

## **Do You Need Information or Assistance on VWR Products?**

The VWR Sales Group can provide information on pricing and give you quotations. We can take your order and provide delivery information on major equipment items or make arrangements to have your local sales representative contact you. Our products are listed on the internet and we can be contacted through our Internet home page.

Visit VWR's website at [vwr.com](http://vwr.com) for:

- Complete technical service contact information
- Access to VWR's Online Catalogue, and information about accessories and related products.
- Additional product information and special offers.
- Contact us: For information or technical assistance, contact your local VWR representative, or visit [vwr.com](http://vwr.com).

VWR can supply technical information about proper setup, operation or troubleshooting of your equipment. We can fill your needs for replacement parts or provide you with on-site service. We can also provide you with a quotation on our Extended Maintenance Program for your products.

Whatever VWR products you need or use, we will be happy to discuss your applications. If you are experiencing technical problems, working together, we will help you locate the problem and, chances are, correct it yourself...over the telephone without a service call.

When more extensive service is necessary, we will assist you with trained technicians or a qualified service organization for on-the-spot repair. If your service need is covered by the VWR Scientific Products warranty, we will arrange for the unit to be repaired at our expense and to your satisfaction.

# Table of Contents

|                  |   |            |
|------------------|---|------------|
| <b>Section 1</b> | <b>Product Specifications</b> .....             | <b>1-1</b> |
| <b>Section 2</b> | <b>Overview</b> .....                           | <b>2-1</b> |
|                  | Control Panel Keys, Displays & Indicators ..... | 2-3        |
|                  | Keypad Operation .....                          | 2-4        |
|                  | Install the Freezer .....                       | 2-5        |
|                  | Choose the Location .....                       | 2-5        |
|                  | Displays .....                                  | 2-5        |
|                  | Install the Shelves .....                       | 2-6        |
|                  | RS-232 Communications .....                     | 2-6        |
|                  | Install the Wall Bumpers .....                  | 2-6        |
|                  | Remote Alarm Contacts and Analog Output .....   | 2-7        |
|                  | Attach the Power Cord .....                     | 2-8        |
|                  | Connect Unit to Electrical Power .....          | 2-8        |
|                  | Freezer Start-Up .....                          | 2-9        |
|                  | Set the Operating Temperature .....             | 2-10       |
|                  | Set the High Temperature Alarm .....            | 2-10       |
|                  | Set the Low Temperature Alarm .....             | 2-11       |
|                  | Access Code .....                               | 2-11       |
|                  | Run Mode .....                                  | 2-11       |
| <b>Section 3</b> | <b>Calibrate</b> .....                          | <b>3-1</b> |
|                  | Calibrate the Control Probe .....               | 3-1        |
|                  | Calibrate the Optional Sample Probe .....       | 3-2        |
|                  | Temperature Stabilization Periods .....         | 3-2        |
| <b>Section 4</b> | <b>Configuration</b> .....                      | <b>4-1</b> |
|                  | High Alarm Test .....                           | 4-1        |
|                  | Low Alarm Test .....                            | 4-1        |
|                  | System Battery Test .....                       | 4-2        |
|                  | BUS Battery Test .....                          | 4-2        |
|                  | Display Temperature .....                       | 4-2        |
|                  | Clear High Stage Alarm .....                    | 4-3        |
|                  | Set Access Code .....                           | 4-3        |
|                  | RS485 Address .....                             | 4-4        |
|                  | Back-Up System Type .....                       | 4-4        |
|                  | Cold Excursion .....                            | 4-4        |
|                  | Warm Excursion .....                            | 4-4        |
|                  | Reset Excursion .....                           | 4-4        |

|                  |  |            |
|------------------|--|------------|
| <b>Section 5</b> | <b>Alarms</b> .....                                | <b>5-1</b> |
|                  | Lost Communication Alarm .....                     | 5-2        |
|                  | Micro Board Failure Alarm .....                    | 5-2        |
|                  | Multiple Alarms .....                              | 5-2        |
|                  | High Stage System Failure .....                    | 5-2        |
|                  | Error Messages .....                               | 5-3        |
| <b>Section 6</b> | <b>Maintenance</b> .....                           | <b>6-1</b> |
|                  | Clean Air Filter .....                             | 6-1        |
|                  | Clean the Condenser .....                          | 6-1        |
|                  | Defrost the Chamber .....                          | 6-3        |
|                  | Clean the Door Gasket .....                        | 6-4        |
|                  | Vacuum Relief Port .....                           | 6-4        |
|                  | Vacuum Relief Port Maintenance .....               | 6-5        |
|                  | Replace the Battery .....                          | 6-6        |
| <b>Section 7</b> | <b>Factory Installed Options</b> .....             | <b>7-1</b> |
|                  | Back-Up System (BUS) .....                         | 7-1        |
|                  | Install Vent Stack, Solenoid & Injection Asm. .... | 7-2        |
|                  | Install the Temperature Probe .....                | 7-3        |
|                  | Connect the Probe/Solenoid Harness .....           | 7-4        |
|                  | BUS Operation and Maintenance .....                | 7-5        |
|                  | Set Optional BUS Setpoing .....                    | 7-7        |
|                  | Test BUS Operation .....                           | 7-8        |
|                  | Cleat the Vent Stack .....                         | 7-8        |
|                  | Disconnect the Fitting Assembly .....              | 7-8        |
|                  | Chart Recorder .....                               | 7-9        |
|                  | Change the Recorder Temperature Range .....        | 7-9        |
|                  | Calibrate the Recorder .....                       | 7-10       |
| <b>Section 8</b> | <b>Warranty Information</b> .....                  | <b>8-1</b> |
| <b>Appendix</b>  | <b>Handling Liquid Nitrogen</b> .....              | <b>A-1</b> |
|                  | Introduction .....                                 | A-2        |
|                  | Handling Liquid Carbon Dioxide .....               | A-4        |
|                  | First Aid .....                                    | A-4        |

# Section 1 Product Specifications

|  |  |   |   |
|--|--|---|---|
| <b>Temperature Range</b>                     | -50°C (-58°F) to -86°C (-123°F) or -10°C (14°F) to -40°C (-40°F) in an 18°C to 28°C (64.4°F to 82.4°F) ambient |   |   |
| <b>Capacity</b>                              | 13.0 cu. ft. (368.1 liters)  |   |   |
|  | 17.3 cu. ft. (489.9 liters)  |   |   |
|  | 23.0 cu. ft. (651.3 liters)  |   |   |
|  | 28.0 cu. ft. (792.8 liters)  |   |   |
| <b>Refrigeration</b>                         | 2545 BTUH (Two compressors for -86C units), (One compressor for -40C units)                                    |   |   |
| <b>Insulation</b>                            | CFC-free, foamed-in-place urethane: minimum 5.0" (12.7cm) cabinet; 4.5" (11.4 cm) door                         |   |   |
| <b>Electrical -<br/>nominal voltage ±10%</b> | 208-230VAC, 60 Hz, 12 FLA  | 230VAC, 50 Hz, 12 FLA                                   | 120VAC, 60 Hz, 16 FLA                                   |
| <b>Breaker Requirements</b>                  | 15 Amp, Dedicated Circuit,<br>15 Amp Time Delay Breaker  | 15 Amp, Dedicated Circuit,<br>15 Amp Time Delay Breaker | 20 Amp, Dedicated Circuit,<br>20 Amp Time Delay Breaker |

## Certifications

Declaration of Conformity is available on request as a separate document. Contact VWR for information.

## Safety Specifications

Indoor Use Only

Altitude - Up to 2,000 meters

Temperature - 15°C to 32°C

Humidity - Maximum RH 80% for temperatures up to 31°C, decreasing linearly to 50% RH at 40°C

Mains Supply Fluctuations - Mains supply voltage fluctuations not to exceed ±10% of the nominal voltage

Installation Category II <sup>1</sup>

Pollution Degree 2 <sup>2</sup>

Class of Equipment



## Fluorinated Greenhouse Gases

Compliant with REGULATION (EU) No 517/2014 OF THE EUROPEAN PARLIAMENT AND OF THE COUNCIL on fluorinated greenhouse gases. This product contains fluorinated greenhouse gases in a hermetically sealed system. This product contains foam blown with fluorinated greenhouse gas, R-245fa. If a leak in the sealed system is detected, the operator shall repair without undue delay.

For ULT models rated 230V/50Hz, intended for import into the EU, the following sizes are designed for use with the following amounts of fluorinated greenhouse gases:

For -86°C rated units

| Size (ft <sup>3</sup> ) | 1st Stage   |             |       | 2nd Stage   |             |        |
|-------------------------|-------------|-------------|-------|-------------|-------------|--------|
|                         | Refrigerant | Amount (kg) | GWP   | Refrigerant | Amount (kg) | GWP    |
| <b>13</b>               | R-404a      | 0.723       | 3 922 | R-508b      | 0.306       | 13 396 |
| <b>17</b>               | R-404a      | 0.751       | 3 922 | R-508b      | 0.388       | 13 396 |
| <b>23</b>               | R-404a      | 0.751       | 3 922 | R-508b      | 0.425       | 13 396 |
| <b>28</b>               | R-404a      | 0.652       | 3 922 | R-508b      | 0.397       | 13 396 |

For -40°C rated units

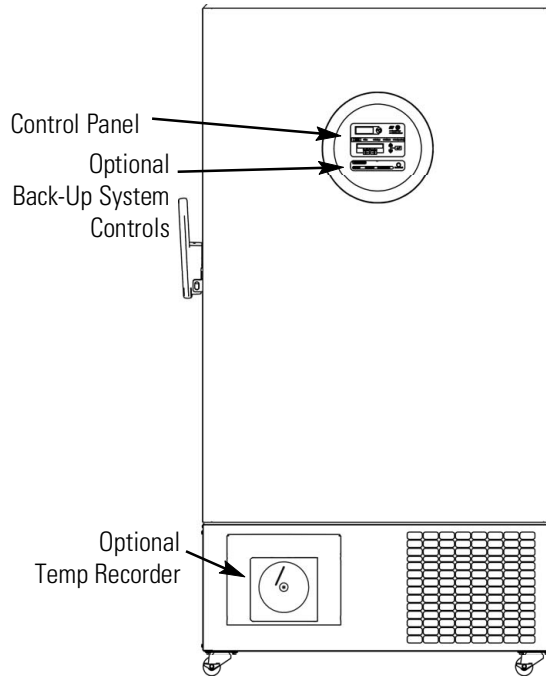
| Size (ft <sup>3</sup> ) | 1st Stage   |             |       |
|-------------------------|-------------|-------------|-------|
|                         | Refrigerant | Amount (kg) | GWP   |
| <b>13</b>               | R-404a      | 0.595       | 3 922 |
| <b>17</b>               | R-404a      | 0.624       | 3 922 |
| <b>23</b>               | R-404a      | 0.624       | 3 922 |

1 Installation category (overvoltage category) defines the level of transient overvoltage which the instrument is designed to withstand safely. It depends on the nature of the electricity supply and its overvoltage protection means. For example, in CAT II which is the category used for instruments in installations supplied from a supply comparable to public mains such as hospital and research laboratories and most industrial laboratories, the expected transient overvoltage is 2500V for a 230V supply and 1500V for a 120V supply.

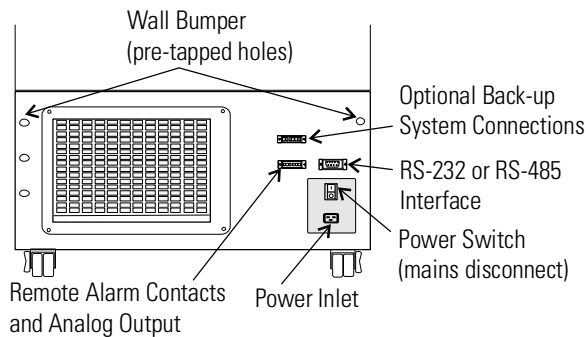
2 Pollution degree describes the amount of conductive pollution present in the operating environment. Pollution degree 2 assumes that normally only non-conductive pollution such as dust occurs with the exception of occasional conductivity caused by condensation.



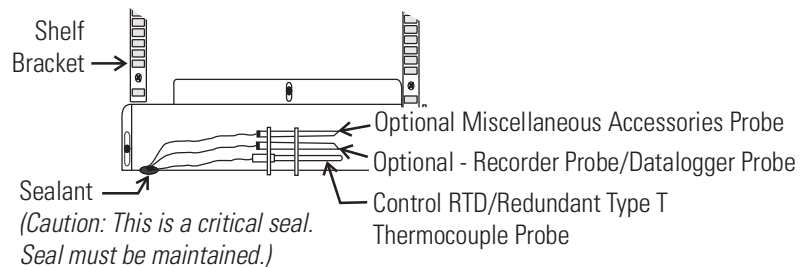
## Section 2 Overview



**Figure 2-1. Front View**



**Figure 2-2. Rear View**



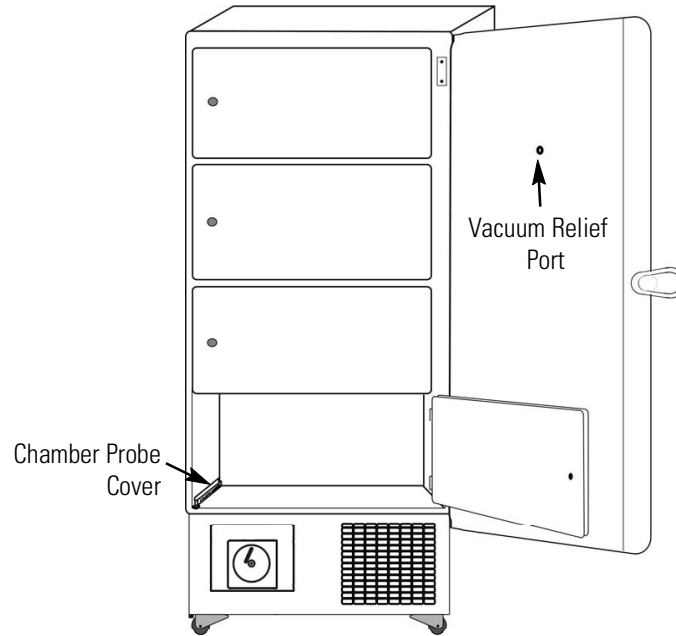
**Figure 2-3. Chamber Probes**

**Figure 2-1**

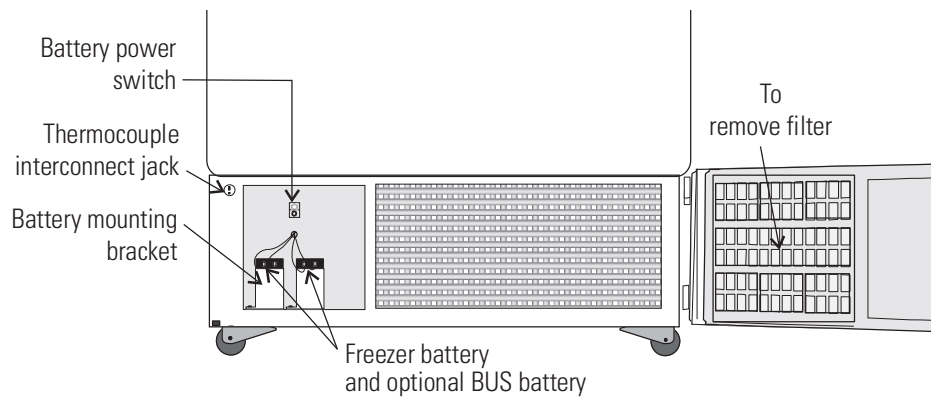
- Control Panel - keypad, displays and indicators
- BUS (Optional Back Up System) panel
- Optional temperature recorder (7 day, one pen)

**Figure 2-2**

- Remote alarm contacts and selectable analog output connection (0-1V, 4-20mA (default), 0-5V)
- Power inlet for power cord connection.
- Optional BUS connections for probe and solenoid
- RS-232 (default) or RS-485 interface
- Power switch (mains disconnect)



**Figure 2-4.** Vacuum Relief and Probe Cover Location



**Figure 2-5.** Battery(s) Location and Switch

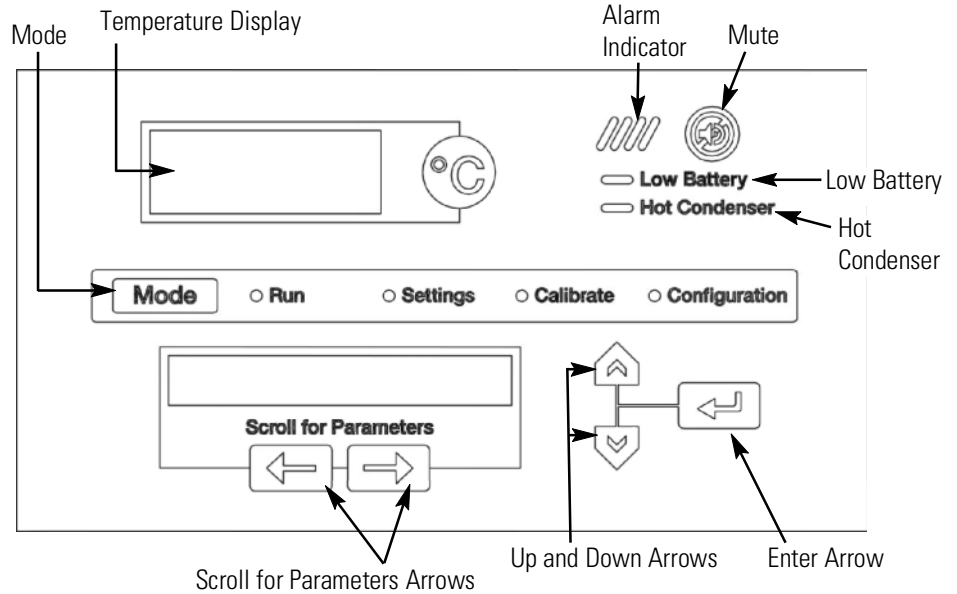
**Figure 2-4**

- Vacuum relief port - pressure equalization port
- Probe cover houses control, optional recorder, redundant alarm probes.

**Figure 2-5**

- Battery mounting bracket(s)
- Battery power switch (freezer and BUS)
- Freezer battery
- Optional BUS battery
- Freezer filter location

# Control Panel Keys, Displays & Indicators



**Figure 2-6.** Control Panel

- Mode Select Switch - Used to select Run, Settings, Calibrate and System Configuration Modes.  
Mode Select Indicators -  
Run: Run Menu  
Settings: Set Points Menu  
Calibrate: Calibrate Menu  
Configuration: Configuration Menu
- Temperature Display - Displays temperature in degrees Celsius.
- Alarm Indicator - Light pulses on/off during an alarm condition of the cabinet.
- Mute - Silences the audible alarm.
- Low Battery - indicates a low battery condition of the freezer battery.
- Hot Condenser - indicates a hot condenser condition.
- Message Center - displays system status and alarms.
- Scroll for Parameters Arrows - moves the operator through the choices of the selected mode.
- Up and Down Arrows - Increases or decreases values, toggles between choices.
- Enter Arrow - Stores the value into memory.

## Keypad Operation

This upright freezer has four basic modes which allow freezer setup: Run, Settings, Calibrate and Configuration.

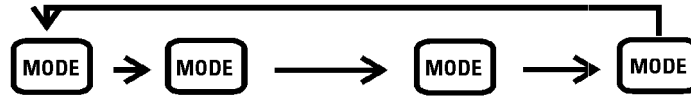
**Run** is the default mode for the freezer during normal operation.

**Settings** is used to enter system set points for freezer operation.

**Calibrate** is used to calibrate various system parameters.

**Configuration** allows for custom setup of various options.

The chart below shows the selections under each of the modes.



| Run                       | Settings                             | Calibrate                | Configuration             |
|---------------------------|--------------------------------------|--------------------------|---------------------------|
| Default Mode<br>System Ok | Control Set Point                    | Control Probe            | High Alarm Test           |
| Line Voltage              | High Alarm Set Point                 | Optional Sample<br>Probe | Low Alarm Test            |
| Compensated Voltage       | Low Alarm Set Point                  |                          | System Battery Test       |
| HSHX Temperature          | Optional Back Up<br>System Set Point |                          | BUS Battery Test          |
|                           |                                      |                          | Display Temperature       |
|                           |                                      |                          | Clear High Stage<br>Alarm |
|                           |                                      |                          | Set Access Code           |
|                           |                                      |                          | RS485 Address             |
|                           |                                      |                          | BUS type CO2 or LN2       |
|                           |                                      |                          | Cold Excursion            |
|                           |                                      |                          | Warm Excursion            |
|                           |                                      |                          | Reset Excursion           |

**Scroll for Parameters Arrows:** Steps the operator through the parameters of SETTINGS, CALIBRATE and CONFIGURATION Modes. The right arrow goes to the next parameter, the left arrow returns to the previous parameter.

**Up Arrow:** Increases or toggles the parameter value that has been selected in the SETTINGS, CALIBRATE, and CONFIGURATION Modes.

**Enter:** Must press Enter key to save to memory any changed values.

**Down Arrow:** Decreases or toggles the parameter values that have been selected in the SETTINGS, CALIBRATE and CONFIGURATION Modes.

**Mute Key:** Press to silence the audible alarm. See Section 4 for alarm ringback times.

## Displays

**Message Center:** Displays the system status (Mode) at all times. Displays SYSTEM OK during normal operation, or alarm messages if the system detects an alarm condition. See the Alarms section.

## Install the Freezer



**Caution** If tipped more than 45°, allow the unit to set upright for 24 hours before start up. ▲

To remove the freezer from the pallet, use the ½" wrench to remove all the bolts securing the shipping bracket to the pallet.

Remove the shipping bracket. Remove the ramp boards from the pallet and place the slotted end over the ramp brackets on the pallet. The support blocks on the ramps will be facing down. Before moving the freezer, make sure the casters are unlocked and moving freely. Align the caster with the ramp boards. Use adequate personnel to roll the freezer off the pallet.

The freezer can be easily pushed to the desired approved location, described below. If necessary, the doors and lower front panel may be opened to move the unit through tight openings. When the freezer is in position, set the front caster brakes.



**Caution** If the factory installed option water-cooled condenser is present, do not turn the freezer on without water connected and flowing. **Damage to the refrigeration system could occur within 5 minutes if water is not connected and flowing on unit start-up.** Refer to Section 6. ▲

## Choose the Location

Locate the freezer on a firm, level surface in an area with an ambient temperature between 18°C and 28°C. Provide ample room to reach the mains disconnect switch (power switch) located on the rear of the freezer.



**Caution** The freezer must not be moved with the product load inside. ▲

**Caution** For proper ventilation and airflow, a minimum clearance of 5" at the rear and top, and a clearance of 8" on the side of the freezer is required. Allow adequate space in front of the freezer for door opening. ▲

## Door Operation

Upright freezer models are equipped with an advanced assembly specifically designed for ultra-low temperature freezers.

Features include:

- One-hand operation
- A front-accessible lock
- Hasps for a standard padlock to provide additional security. Length of the shackle must be between ¾ inch (1.9 cm) and 1½ inch (3.8 cm).
- Durable construction for reliable operation and safe product storage.

## Install the Wall Bumpers

The parts bag, located inside the cabinet, contains the following parts.

| Quantity | Description        | Purpose       |
|----------|--------------------|---------------|
| 2        | 1/4-20x5-1/2" Bolt | Wall Bumper   |
| 2        | Neoprene Cap       | Cap Protector |

Install the bolts into the pre-tapped holes on the back of the compressor section. Install a neoprene cap on each bolt. Refer to Figure 1-2 for the location of the pre-tapped holes.

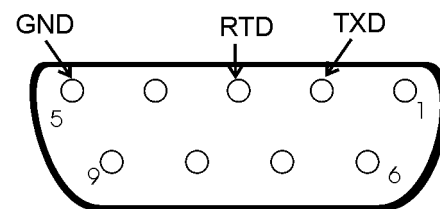
## Install the Shelves

Install the shelf clips into the shelf pilasters (front and back) at the desired shelf level. Install the shelves in the cabinet onto the clips.

**Note** On units having the optional 5 inner door option, refer to the instructions accompanying the inner door kit.

## RS-232 Communications

This upright freezer has a data communications interface. The factory default setting is RS-232.



**Figure 2-7.** Wiring Identification

The wiring identification for the interface is shown in Figure 2-7. One nine pin, sub "D" style connector is located on the back of the freezer. See Figure 1-2 for the location of the connector on the freezer.



## RS-232 Communications (continued)

The freezer transmits temperature information every 60 minutes. A standard DB9 serial extension cable can be used to connect the freezer to a serial device. Some serial devices may require a null modem adapter.

Data format:

Baud . . . . . 1200  
 Data bits . . . . . 8 (7 bit ASCII with leading zero)  
 Start bits . . . . . 1  
 Stop bits . . . . . 2  
 Parity . . . . . none

The data transfer sequence is transmitted in the following format. X refers to numerical temperature data.

(NUL) (-) XXX (SP) C (SP) (Error Message) (SP) (LF) (CR) (EOT) (SP)

In the event of a CNTRLFAIL, Er07, or the control probe is out of range error, the numerical temperature data (XXX) in the transmission would be replaced by T\_ERR.

If no alarm condition exists, spaces will be sent. A total of 20 characters will be sent.

|                           |                       |
|---------------------------|-----------------------|
| SP - Space                | LF - Line feed        |
| CR - Carriage return      | EOT - End of text (4) |
| NUL - Null character (00) |                       |

If an alarm condition does exist, “Error Message” in the protocol will be replaced by the following:

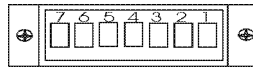
UNDERTEMP (temperature above the low alarm setpoint)  
 OVERTEMP (temperature below the high alarm setpoint)  
 PWRFAIL (AC power failure)  
 CNTRLFAIL (Control probe failure)  
 Er07 (micro failure)  
 HSHX FAIL (Heat exchanger failure)  
 HOT COND (Hot condenser)

## Remote Alarm Contacts and Analog Output

This upright freezer has remote alarm contacts and analog output. See Figure 2-2 for the location of the remote alarm contacts. The remote alarm connector is located in the parts bag provided with the manual. It must be installed if connecting the freezer to an alarm system. After installing the wiring from the alarm system to the connector, install the connector to the freezer microboard and secure with the two screws provided.

## Remote Alarm Contacts and Analog Output (cont.)

The remote alarm provides a NO (normally open) output, a NC (normally closed) output and COM (common). The contacts will trip on a power outage, high temperature alarm or low temperature alarm. They will also trip on high stage, control probe and microboard failures. Figure 2-8 shows the remote contacts in alarm state.



| REMOTE CONTACTS/ANALOG OUTPUT |                 |
|-------------------------------|-----------------|
| PIN# 1                        | Analog Output + |
| PIN# 2                        | Analog Output - |
| PIN# 3                        | Not Connected   |
| PIN# 4                        | Not Connected   |
| PIN# 5                        | Normally Closed |
| PIN# 6                        | Common          |
| PIN# 7                        | Normally Open   |

CONTACT RATING: 1A @ 30V  
CONTACTS IN ALARM STATE

**IMPORTANT USER INFORMATION**

CAUTION! Stored product should be protected by a redundant 24 hour/day monitoring system with alarm capability. An interconnect jack and thermocouple are installed for centralized monitoring, should on-board system fail.

**Figure 2-8.** Remote Alarm Contact Pins

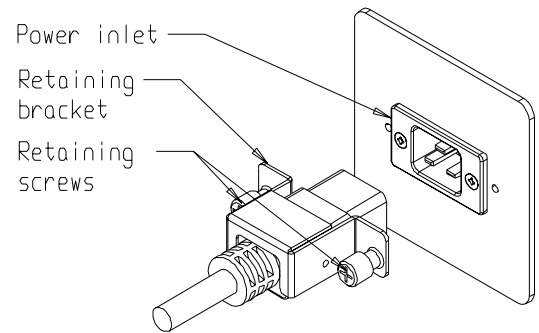
The analog output function allows the freezer to output signals representing the temperature of the freezer cabinet. The factory default setting is 4-20 mA. Refer to Figure 2-9 for output specifications.

|             | 4-20 mA       | 0-1V          | 0-5V          |
|-------------|---------------|---------------|---------------|
| Temperature | -100 to +50°C | -100 to +50°C | -100 to +50°C |

**Figure 2-9.** Specifications

## Attach the Power Cord

Insert the power cord into the power inlet module. Place the retaining bracket (P/N 195763) over the connector. Tighten retaining screws to secure.



**Figure 2-10.** Power Cord Connection

## Connect Unit to Electrical Power



**Caution** See the serial tag on the side of the unit for electrical specifications or refer to the electrical schematics in this manual. ▲

**Caution** If the factory installed option water-cooled condenser is present, do not turn the freezer on without water connected and flowing. **Damage to the refrigeration system could occur within 5 minutes if water is not connected and flowing on unit start-up.** Refer to Section 6. ▲

## Connect Unit to Electrical Power (continued)

The freezer should be operated on a dedicated grounded service. Check the voltage rating on the serial tag of the unit and compare it with the outlet voltage. Then, with the power switch turned off, plug the line cord into the wall outlet.

First turn on the freezer power switch. Then open the lower front door by grasping the bottom left corner. Locate the battery switch (Figure 1-5) and turn it to Standby mode (⏻). During initial freezer start-up, the system battery may require charging and the Low Battery message may appear in message center.



**Caution** Ensure the battery switch is turned to Standby mode (⏻). The rechargeable batteries require 36 hours to charge at initial start-up. A “Low Battery” alarm may occur until the batteries are fully charged. Should a power failure occur during the initial start-up period, the electronics will have limited operation. ▲

## Freezer Start-Up

With the freezer properly installed and connected to power, system setpoints can be entered. The following setpoints can be entered in Settings mode: Control temperature, high temperature alarm setpoint, low temperature alarm setpoint, and (optional) BUS setpoint. Default settings are shown in the table below.

**Table 2-2.** Default Settings

|                               |       |                               |       |
|-------------------------------|-------|-------------------------------|-------|
| <b>Control Set Point</b>      | -80°C | <b>Control Set Point</b>      | -40°C |
| <b>High Temperature Alarm</b> | -70°C | <b>High Temperature Alarm</b> | -30°C |
| <b>Low Temperature Alarm</b>  | -90°C | <b>Low temperature alarm</b>  | -50°C |
| <b>Optional BUS Set Point</b> | -60°C | <b>Optional BUS Set Point</b> | -30C  |



**Caution** If the setpoint is changed and the low temperature and high temperature alarms are set 10° from the set point, the alarm setpoints will be adjusted automatically to maintain a distance of at least 10° from setpoint. ▲



**Caution** If the factory installed option water-cooled condenser is present, do not turn the freezer on without water connected and flowing. **Damage to the refrigeration system could occur within 5 minutes if water is not connected and flowing on unit start-up.** Refer to Section 6. ▲

## Set the Operating Temperature

These upright freezers have an operating temperature range of  $-50^{\circ}\text{C}$  to  $-86^{\circ}\text{C}$ , or  $-10^{\circ}\text{C}$  to  $-40^{\circ}\text{C}$ , depending on model and ambient temperature. The freezer is shipped from the factory with a temperature set point of  $-80^{\circ}\text{C}$ , or  $-40^{\circ}\text{C}$ , depending on model. To change the operating temperature setpoint:

1. Press the Mode key until the Settings indicator lights.
2. Press right arrow until “SET PT = -XX” is displayed in message center.
3. Press the up/down arrow key until the desired temperature set point is displayed.
4. Press Enter to save the set point.
5. Press the Mode key until the Run indicator lights for Run mode or press the right/left arrow keys to go to next/previous parameter.

If no control keys are pressed, the freezer will automatically return to RUN mode after 5 minutes.

*For  $-40^{\circ}\text{C}$  units: At  $-20^{\circ}\text{C}$  operation, probe may need to be calibrated to ensure cabinet temperatures are within required range. Refer to Calibration section for procedure.*

## Set the High Temperature Alarm

The high temperature alarm will activate an audible/visual warning when the freezer chamber temperature has reached or exceeded the high temperature alarm set point. To set the high temperature alarm set point:

1. Press the Mode key until the Set indicator lights.
2. Press right arrow until “HI ALM = -XX” is displayed in message center.
3. Press the up or down arrow key until the desired high temperature alarm set point is displayed.
4. Press Enter to save the setting.
5. Press the Mode key until the Run indicator lights or press the right or left arrow to go to the next or previous parameter.

If no control keys are pressed, the freezer will automatically return to RUN mode after 5 minutes.



**Caution** The high alarm set point must be set at least  $5^{\circ}\text{C}$  from the control set point. At initial start-up, the high temperature alarm is disabled until the cabinet reaches set point, or 12 hours elapse. ▲

## Set the Low Temperature Alarm

The low temperature alarm will activate an audible/visual warning when the freezer chamber temperature has reached or decreased below the low temperature alarm set point. To set the low temperature alarm set point:

1. Press the Mode key until the Settings indicator lights.
2. Press the right arrow until “LO ALM = -XX” is displayed in the message center.
3. Press the up or down arrow key until the desired low temperature alarm set point is displayed.
4. Press Enter to save the setting.
5. Press the Mode key until the Run indicator lights or press the right or left arrow to go to the next or previous parameter.

If no control keys are pressed, the freezer will automatically return to RUN mode after 5 minutes.

**Note** The low alarm set point must be set at least 5°C from the control set point.

## Access Code

An access code of 000 is required to access the Settings, Calibrate or Configuration modes. If the access code is not set at the default ‘000’, a code must be entered to leave RUN mode. See Section 4 for instructions on modifying the access code.

## Run Mode

Run is the default mode for the freezer. The run mode will display the cabinet temperature on the temperature display and ‘SYSTEM OK’ on the message center under normal operating conditions. In addition, this mode allows display of the following information:

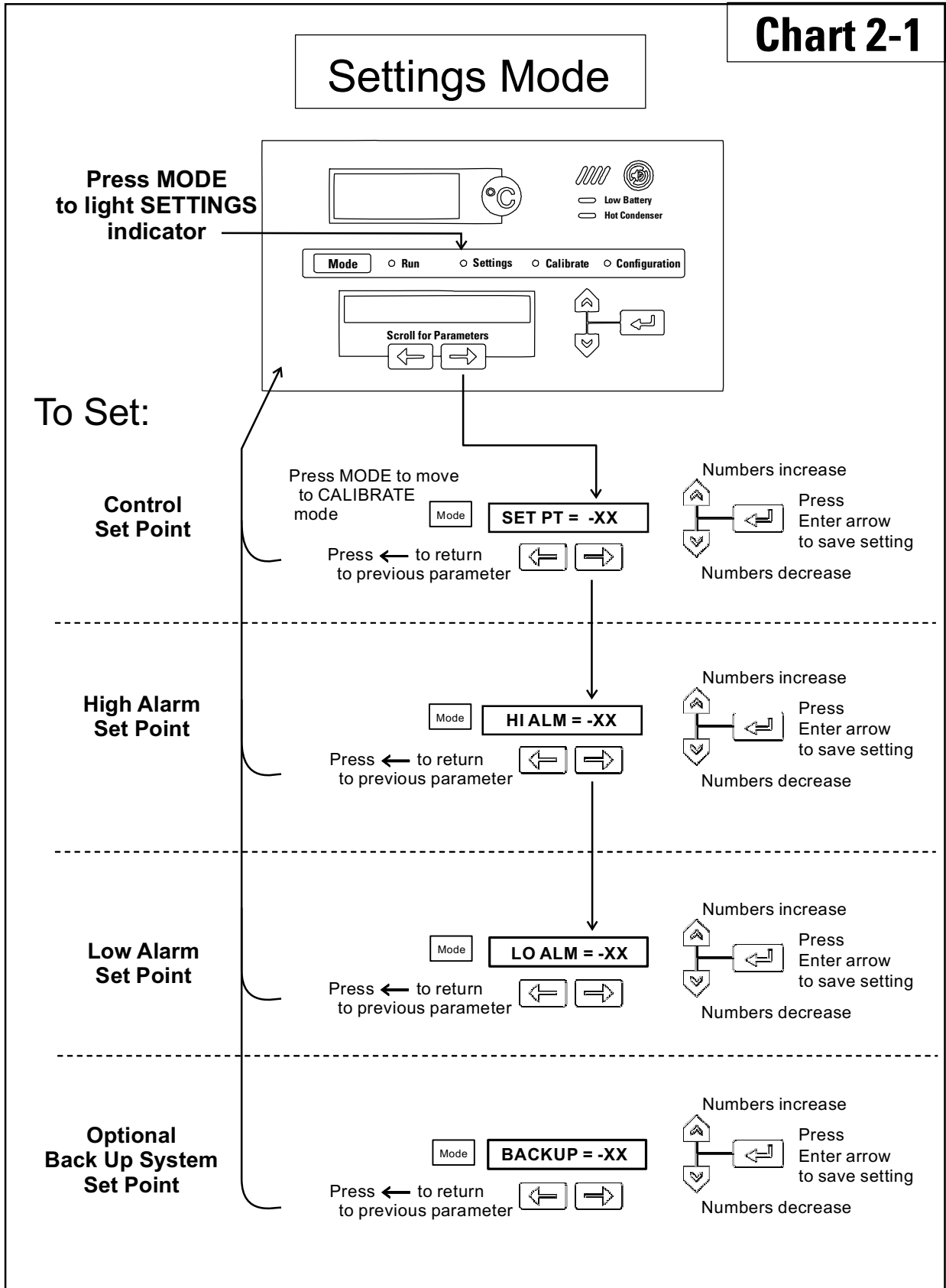
LINE VOLTAGE

COMPENSATED VOLTAGE

HSHX TEMPERATURE (heat exchanger temperature), - 86°C models

This information is scrolled individually by pressing the right arrow key. In each case, the message center returns to SYSTEM OK in 10 seconds if no keys are pressed.

**Chart 2-1**



## Section 3 Calibrate

Once the freezer has stabilized, the control or sample probe may need to be calibrated. Calibration frequency is dependent on use, ambient conditions and accuracy required. A good laboratory practice would require at least an annual calibration check. On new installations, all parameters should be checked after the stabilization period.

On -40°C models only: For -20°C operation, calibration may be needed to ensure cabinet air temperatures are within a specified range.



**Caution** Before making any calibration or adjustments to the unit, it is imperative that all reference instruments be properly calibrated. ▲

### Calibrate the Control Probe

Plug a type T thermocouple reader into the receptacle located inside the lower door (Figure 2-5). Compare the control temperature set point to the temperature of the measuring device. See Chart 3-1 at the end of this section for more detail.

1. Press the Mode key until the Calibrate indicator lights.
2. Press right arrow until “CONT T = -XX.X” appears in message center.
3. Press up/down arrow to match the display to calibrated instrument.
4. Press Enter to store calibration.
5. Press the Mode key to return to Run or the right/left arrow to go to next/previous parameter.

On -40°C models only: **Probe Calibration for -20°C Operation;** Note that if the peak variation is within  $\pm 5^{\circ}\text{C}$  of set value, then no calibration is required.

1. Set temperature to -20C.
2. Calibrate probe to -4°C (-2.5 for 13 ft units) per above procedure.
3. Allow unit to stabilize to -20°C per temperature stabilization period below.
4. Check peak variation after unit has achieved steady state operation. Probe may need to be calibrated a second time to achieve the required range.

## Calibrate the Optional Sample Probe

For freezers with the optional sample probe, place the calibrated instrument in the center of the sample bottle. The bottle should contain an appropriate medium and the measuring instrument should be centered in the bottle.

1. Press the Mode key until the Calibrate indicator lights.
2. Press the right arrow until "SAMP T = -XX.X" appears in the message center.
3. Press up/down arrow to match display to calibrated instrument.
4. Press Enter to store calibration.
5. Press the Mode key to return to Run or the right/left arrow to go to next/previous parameter.

See Chart 3-1 for calibration process functions.

## Temperature Stabilization Periods

### Temperature Stabilization Periods

Startup - Allow 12 hours for the temperature in the cabinet to stabilize before proceeding.

Already Operating - Allow at least 2 hours after the display reaches set point for temperature to stabilize before proceeding.

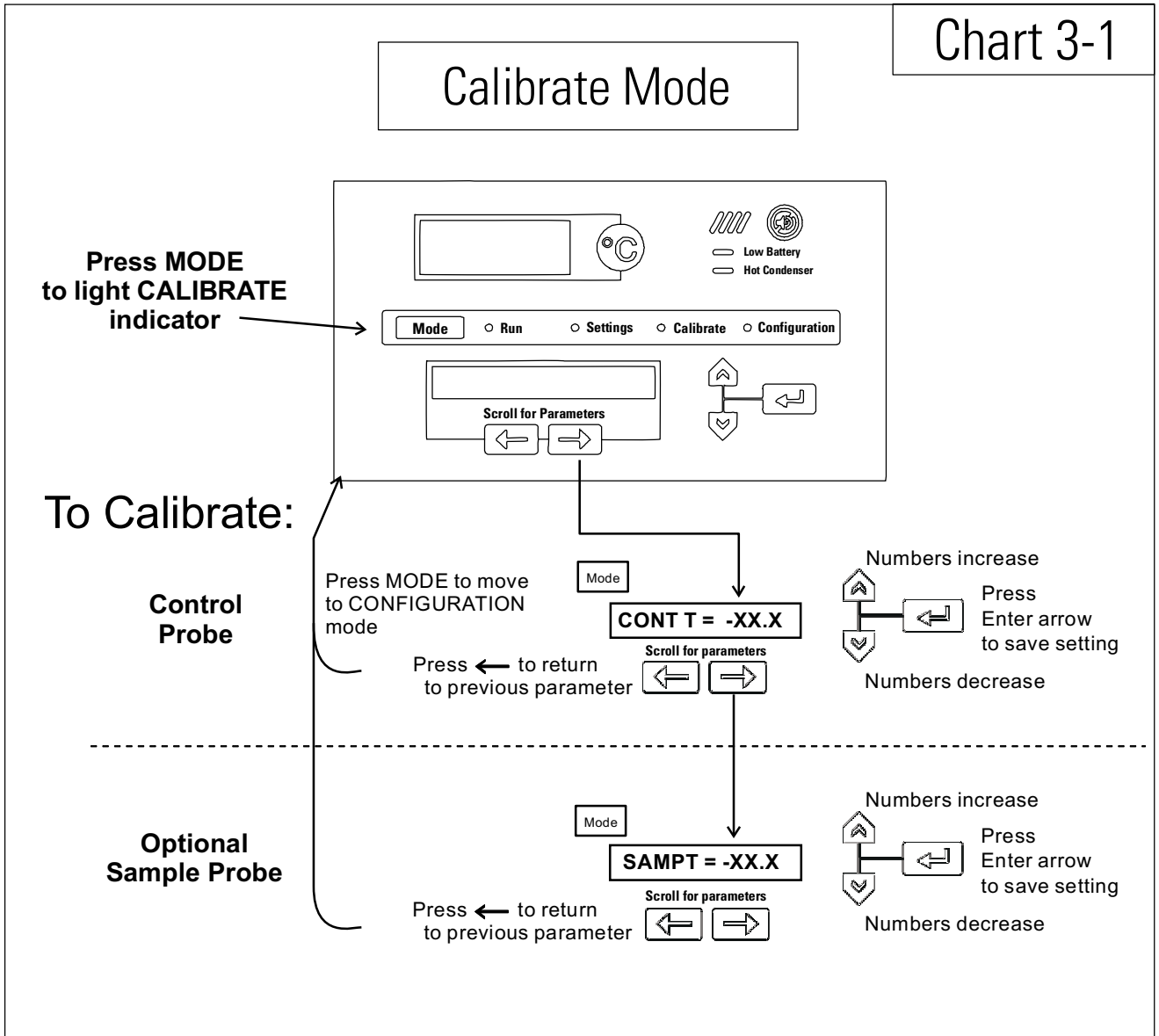


**Caution** During calibration, the temperature display is not available. ▲

If no keys are pressed for approximately five minutes while in calibration mode, the system will reset to Run mode.



Chart 3-1



## Section 4 Configuration

The Configuration Mode is used for testing and custom setup of the freezer. The configuration functions listed and described below may not be necessary in all applications, but are available if needed. See Chart 3-1 for more detail.

### High Alarm Test

The high alarm test is used to verify the high alarm will activate, should the freezer temperature equal or exceed the high alarm set point.

1. Press the Mode key until the Configuration indicator lights.
2. Press the right arrow until HI ALRM TEST is displayed in the message center.
3. Press Enter to initiate the test.

The temperature on the display will begin to increase until the high alarm set point has been reached. The audible alarm will sound and the alarm indicator will flash. Press the Mute key to silence the alarm.

### Low Alarm Test

The low alarm test is used to verify the low alarm will activate, should the freezer temperature equal or become less than the low alarm set point.

1. Press the Mode key until the Configuration indicator lights.
2. Press the right arrow until LO ALRM TEST is displayed in the message center.
3. Press Enter to initiate the test.

The temperature on the display will begin to decrease until the low alarm set point has been reached. The audible alarm will sound and the alarm indicator will flash. Press the Mute key to silence the alarm.

## System Battery Test

To test the charge of the freezer battery:

1. Press the Mode key until the Configuration indicator lights.
2. Press the right arrow until SYS BAT TEST is displayed in the message center.
3. Press Enter to initiate the test.

TESTING BATT displays during the testing period. Upon completion of the test, the message center displays BATT GOOD or BATT FAIL. When a test is failed, the audible alarm sounds, the alarm indicator and the Low Battery indicator light. Press the Mute key and the alarm indicator goes out. The Low Battery light stays on until a future battery test is performed and passed.

## BUS Battery Test

To test the charge of the BUS battery:

1. Press the Mode key until the Configuration indicator lights.
2. Press the right arrow until BUS BAT TEST is displayed in the message center.
3. Press Enter to initiate the test.

TESTING BATT displays during the testing period. Upon completion of the test, the message center displays BBAT GOOD or BBAT FAIL. When a test is failed, the audible alarm sounds, the alarm indicator and the Low Battery indicator lights. Press the Mute key. The audible alarm and alarm indicator go off. The Low Battery light stays on. If this test fails, it is recommended to replace the BUS battery.

## Display Temperature

This function, only available on freezers with the optional sample probe, allows the user to select which temperature is displayed in the temperature display window. The options are CONTROL or SAMPLE.

1. Press the Mode key until the Configuration indicator lights.
2. Press the right arrow until DISP CONTROL or DISP SAMPLE is displayed in the message center.
3. Press up/down arrow to toggle between the two display selections.
4. Press Enter to save.

If control probe is selected, the temperature display will be on continuously. If sample probe is selected, the temperature display will be preceded with a letter 'S'.

## **Clear High Stage Alarm**

Should a high stage alarm occurred, it may become necessary to clear the alarm condition after the condition has been corrected.

1. Press the Mode key until the Configuration indicator lights.
2. Press the right arrow until CLR HS ALARM is displayed in the message center.
3. Press Enter to clear the alarm.

## **Set Access Code**

To set the Access Code:

1. Press the Mode key until the Configuration indicator lights.
2. Press the right arrow until "SET ACC CODE" is displayed in the message center.
3. Press Enter.
4. The message center will display ACC CODE = 000. Press the up or down arrow key until the desired access code is displayed (000 - 999). Press the left or right arrow key to select digit 1, 2, 3.

**Note** The left and right arrow keys are used to move from the first through the third digits within the access code. ▲

5. Press Enter to save the setting
6. Press the Mode key until the Run indicator lights. A three digit Access Code can be entered to avoid unauthorized personnel from changing the set points, calibration, or configuration. A setting of 000 will bypass the access code. The factory setting is 000.

## RS485 Address

The freezer will need to have a unique identification address for data communications. This address is set through the Configuration mode.

1. Press the Mode key until the Configuration indicator lights.
2. Press the right arrow until RS485ADDR is displayed in the message center.
3. Press Enter. The message center will display 485 ADDR XX.
4. Press up or down arrow to select the appropriate address for the freezer (1 - 24).
5. Press Enter to save.

## Back-Up System Type

This function, which is only available on freezers with the optional BUS (back up system), allows the user to select which type of gas is injected into the freezer chamber. The options are CO2 and LN2.

1. Press the Mode key until the Configuration indicator lights.
2. Press the right arrow until BUS TYPE CO2 or BUS TYPE LN2 is displayed in the message center.
3. Press up/down arrow to toggle between the two display selections.
4. Press Enter to save.

## Cold Excursion

This function displays the coldest temperature recorded by the control probe.

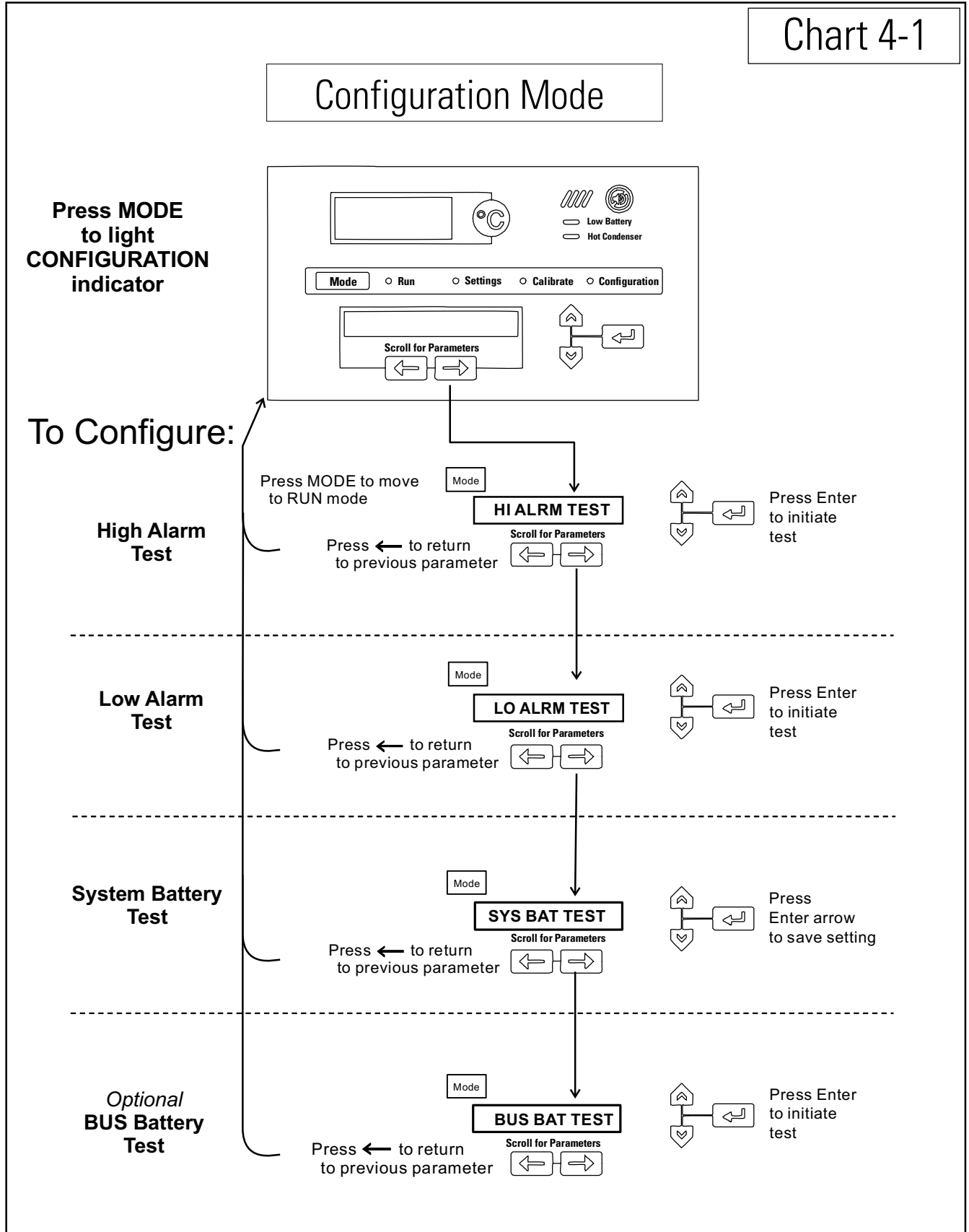
## Warm Excursion

This function displays the warmest temperature recorded by the control probe.

## Reset Excursion

This function resets the cold and warm excursions.

Chart 4-1



# Configuration Mode, Chart 4-1, Page 2 of 3

refer to previous page

To Configure:

Display Temperature

Mode

**DISP CONTROL**

Scroll for Parameters

Press ← to return to previous parameter

CONTROL probe  
Press Enter arrow to save the setting  
SAMPLE probe

Clear High Stage Alarm

Mode

**CLR HS ALARM**

Scroll for Parameters

Press ← to return to previous parameter

Press Enter arrow to clear the alarm condition

Access Code

Mode

**SET ACC CODE XXX**

Scroll for Parameters

Press Mode to exit.  
Press Enter to change.  
Press right or left arrow to move to next digit.

Numbers increase  
Press Enter to save the setting  
Numbers decrease

RS-485 Address  
*(if configured)*

Mode

**RS485 ADDR**

Scroll for Parameters

Press ← to return to previous parameter.  
Press Enter to change the address.

Scroll through the addresses 1 - 24  
Press Enter arrow to save the setting

Optional Back Up System Type

Mode

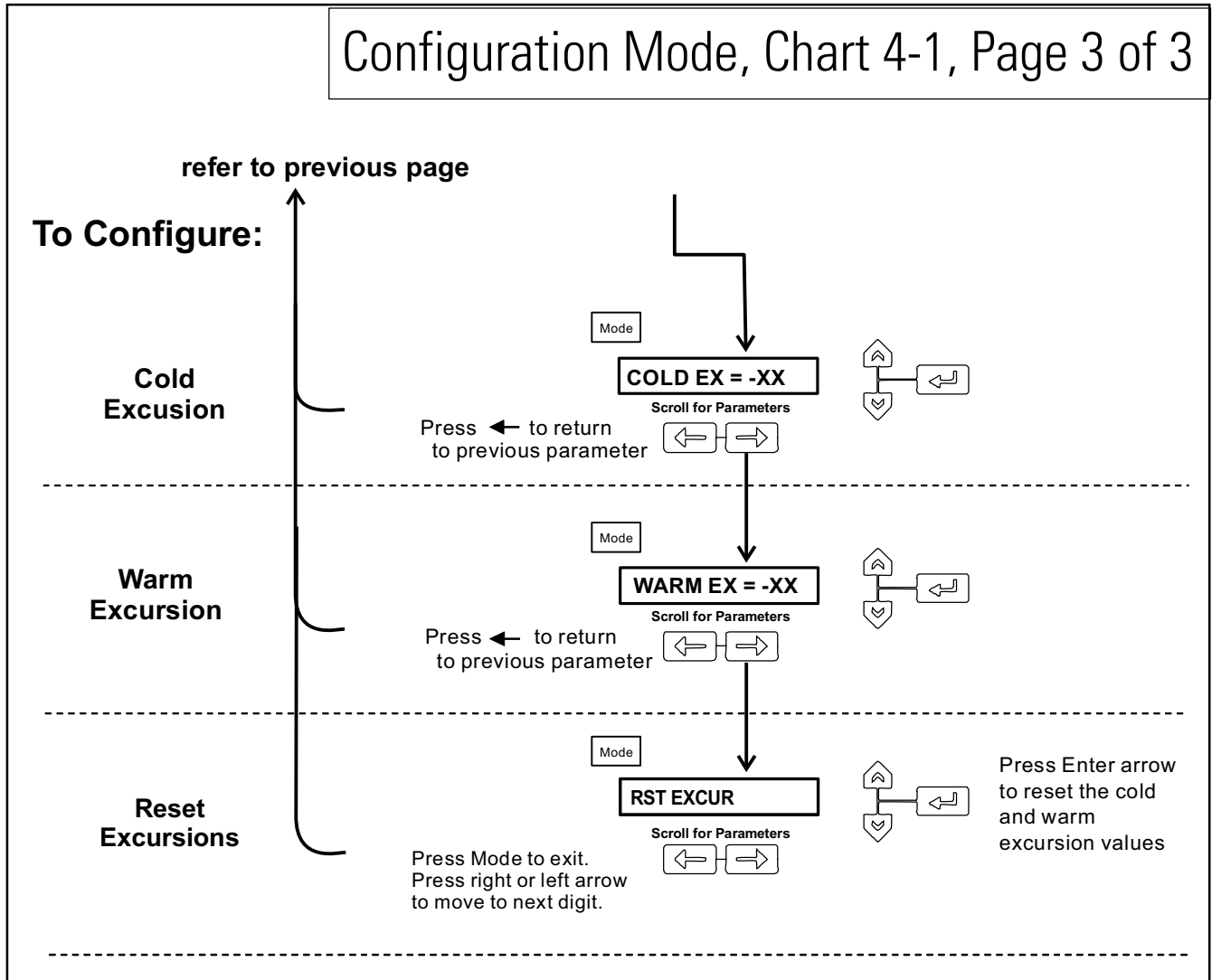
**BUS TYPE CO2**

Scroll for Parameters

Press ← to return to previous parameter

Type CO2  
Press Enter to save the setting  
Type LN2

# Configuration Mode, Chart 4-1, Page 3 of 3





## Section 5 Alarms

The upright freezer alarm system is shown in the table below. When an alarm is active, the message appears in the LED message center. Press the Mute key to silence the audible alarm for the ringback period. The visual alarm will continue until the freezer returns to a normal condition. The alarms are momentary alarms only. When an alarm condition occurs and then returns to normal, the freezer automatically clears the alarm condition and the message center.

| Description                                 | Message       | Delay  | Ringback | Relay |
|---|---------------|--------|----------|-------|
| No alarm condition exists                   | SYSTEM OK     | ----   | ----     | ----  |
| Power Failure                               | POWER FAIL    | 1 min. | 15 min.  | Yes   |
| High Temperature Alarm                      | TEMP IS HIGH  | 1 min. | 15 min.  | Yes   |
| Low Temperature Alarm                       | TEMP IS LOW   | 1 min. | 15 min.  | Yes   |
| Door Ajar                                   | DOOR IS OPEN  | 1 min. | 15 min.  | No    |
| Low Battery*                                | LOW BATTERY   | 1 min. | 8 hours  | No    |
| Low BUS Battery (optional)                  | LOW BUS BATT  | 1 min. | 15 min.  | No    |
| Control Probe Failure                       | PROBE 1 FAIL  | 1 min. | 15 min.  | Yes   |
| Heat Exchanger Probe Failure                | PROBE 2 FAIL  | 1 min. | 15 min.  | No    |
| Condenser Probe                             | PROBE 3 FAIL  | 1 min. | 15 min.  | No    |
| Sample Probe Failure (optional)             | PROBE 4 FAIL  | 1 min. | 15 min.  | No    |
| High Stage System Failure                   | HS SYST FAIL  | 1 min. | 15 min.  | Yes   |
| Condenser Hot Condition (-86°C models only) | HOT CONDENSER | 1 min. | none     | No    |
| Wrong Power                                 | WRONG POWER   | 0 min. | none     | Yes   |
| Micro Board Failure                         | MICRO FAIL    | 0 min. | 15 min.  | Yes   |

*All alarm delays and ringback times are  $\pm 30$  seconds.*

*\* The automatic battery test runs immediately on initial start-up, then every 8 hours thereafter.*

## **High Stage System Failure (-86°C models only)**

This condition is created when the high stage compressor and fans run for 30 minutes and are not capable of cooling the interstage heat exchanger to the proper temperature. Under this condition, the high stage compressor and fans will turn off after 30 minutes, and an audible and visual alarm will occur along with the "HS SYST FAIL" message in the LED message center.

## **Multiple Alarms**

When multiple alarm conditions occur, active messages are displayed in the message center one at a time, updating at 5 second intervals. Pressing Mute during multiple alarms causes all active alarms to be muted and to ring back in 15 minutes.

## **Micro Board Failure Alarm**

An internal communications failure has occurred with the micro board. During this alarm, the compressor(s) attempt to run continuously. However, with this type of failure, freezer operation becomes undependable.

## **Lost Communication Alarm**

Communication between the micro board and the display board has been lost. Under this condition, the visual alarm LED flashes along with dashes (----) in the temperature display. In addition, 'LOST COMM' flashes in the message center. Contact Technical Services.

## Error Messages

| Error | High End Message | Notes  |
|-------|------------------|--|
| Er00  | "INV. MODEL"     | <p><b>Name:</b> Improper model selected.</p> <p><b>Description:</b> Indicates that DIP SW3 has not selected a proper model or can't be accessed properly.</p> <p><b>Response:</b> Display shows "Er00" and will not start-up until a proper model is selected. Contact Technical Services.</p>   |
| ErA1  | " NO FREQUENCY"  | <p>This error condition will prevent peripherals (fans, compressors, etc.) from powering up with the incorrect voltage.</p> <p><b>Name:</b> Voltage/Frequency failure</p> <p><b>Description:</b> Indicates the measured RMS line voltage did not agree with the logic level sensed by the micros provided by the installed high voltage PCB; or the measured RMS voltage is not within a tolerable range (&lt;180VAC &lt; 270 for 230VAC unit / &lt;85 VAC &lt; 160 for 115VAC unit); or the frequency measured over 10 cycles was not within a tolerable range (55 Hz &lt; Freq &lt; 70 Hz for 60 Hz units / 40 Hz &lt; Freq &lt; 55 Hz for 50 Hz units)</p> <p><b>Response:</b> This condition is checked at power on reset and if it is active the unit will NOT power up. The unit will indefinitely display "Er_1" in the display and continue to monitor the frequency and voltage. Furthermore, the audible alarm will sound. Other startup error messages may be displayed prior to this message; however, the system will stop the startup sequence for this condition.</p> <p>ErA1 .. No pulses (zero crossings) detected to determine frequency (50 / 60 Hz)</p> <p>ErC1 .. Frequency detected is below 50 Hz</p> <p>ErD1 .. Frequency detected is above 60 Hz (Possible noise spikes on supply voltage)</p> <p>ErE1 .. Unit is 230V and the voltage detected is below the low limit (180VRMS)</p> <p>ErF1 .. Unit is 230V and the voltage detected is above the high limit (260VRMS)</p> <p>ErG1 .. Unit is 115V and the voltage detected is below the low limit (85VRMS)</p> <p>ErH1 .. Unit is 115V and the voltage detected is above the high limit (160VRMS)</p> |
| ErC1  | "FREQ <50Hz"     |  |
| ErD1  | "FREQ >60Hz"     |  |
| ErE1  | "VAC < 180V"     |  |
| ErF1  | "VAC > 260V"     |  |
| Erg1  | "VAC < 85V"      |  |
| ErH1  | "VAC > 160V"     |  |

**Section 5**

Alarms

| Error (cont.) | High End Message | Notes  |
|---------------|------------------|--|
| Er02          | "CNT PRB FLT"    | <p><b>Name:</b> Control (Cabinet) Sensor Failure</p> <p><b>Description:</b> This condition indicates that the control sensor has failed to produce a valid reading for <math>\geq 12</math> consecutive reads (~60 seconds).</p> <p><b>Response:</b> The unit will stage both compressors on (if necessary) and the unit will attempt to head to bottom out. If the sensor recovers, the system will begin to operate normally and respond to the temperature feedback. The remote alarm contacts will become active regardless of the key position for this mode of failure. 'Er02' will be added to the main display queue and the last valid cabinet temperature value will not be displayed</p>  |
| Er03          | "HSHX PRB FLT"   | <p><b>Name:</b> Heat Exchange Sensor Failure</p> <p><b>Description:</b> This condition indicates that the heat exchange sensor has failed to produce a valid reading for <math>\geq 12</math> consecutive reads (~60 seconds).</p> <p><b>Response:</b> The display will show "Er03" only when the button sequence to read the heat exchange sensor is depressed.</p>   |
| Er05          | N/A              | <p><b>Name:</b> Display Firmware Integrity Failure</p> <p><b>Description:</b> The display firmware has failed to pass its CRC CCITT checksum integrity test.</p> <p><b>Response:</b> The display performs this check at startup and the display board will fail to startup with out any error indication if it does not pass this at power on.</p>   |
| Er06          | N/A              | <p><b>Name:</b> Micro Firmware Integrity Failure</p> <p><b>Description:</b> The micro firmware has failed to pass its CRC CCITT checksum integrity test.</p> <p><b>Response:</b> This is checked at power on reset and the "Er06" will be displayed for ~10 seconds at startup if this condition exists.</p>   |
| Er07          | "MICRO FAIL"     | <p><b>Name:</b> Micro Fail - CS5521 SPI Failure / UISR Failure</p> <p><b>Description:</b> This condition indicates a micro board failure due to either the SPI bus is unable to communicate with the ADC device or a UISR event caused the microcontroller to be in an unstable state.</p> <p><b>Response:</b> The unit will try to recover from this fault three times by a hardware reset of the micro board. In the event that the system couldn't rectify the issue, the following sequence of events will occur:</p> <ol style="list-style-type: none"> <li>1. Remote alarm contacts will become active.</li> <li>2. Buzzer will annunciate audibly and will have a ringback of 15 minutes.</li> <li>3. "Seven segment" display will show "Er07".</li> <li>4. The system will have 10 minute staging between the high stage compressor and the low stage compressor activation.</li> <li>5. The system will go to bottom out temperatures.</li> </ol> |

| Error (cont.)                    | High End Message | Notes   |
|----------------------------------|------------------|---|
| Er09                             | N/A              | <p><b>Name:</b> Stuck Button</p> <p><b>Description:</b> This condition indicates that the display board has a stuck button.</p> <p><b>Response:</b> The Er09 will show on the display periodically.</p>   |
| Er11                             | "COND PRB FLT"   | <p><b>Name:</b> Condenser Probe Sensor Failure</p> <p><b>Description:</b> This condition indicates that the condenser probe sensor has failed to produce a valid reading for <math>\geq 12</math> consecutive reads (~60 seconds).</p> <p><b>Response:</b> The display shows "Er11".</p>          |
| N/A                              | "SMPL PRB FLT"   | <p><b>Name:</b> Sample Probe Sensor Failure</p> <p><b>Description:</b> This condition indicates that the sample probe sensor has failed to produce a valid reading for <math>\geq 12</math> consecutive reads (~60 seconds).</p> <p><b>Response:</b> The message center shows "SMPL PRB FLT".</p> |
| dErr                             | N/A              | <p><b>This is a general display error in which the value being displayed can not be represented within the characters provided.</b></p>   |
| (four dashes)<br>---- in display | N/A              | <p><b>Name:</b> Lost Communication</p> <p><b>Description:</b> Communication between the micro board and the display board has been lost. Under this condition, the visual alarm flashes along with dashes in the temperature display (----). Contact Technical Services.</p>                      |

## Section 6 Maintenance



**Warning** Avoid the excessive use of water around the control area due to the risk of electrical shock. Damage to the controls may also result. ▲

Wipe down the freezer exterior using soap and water and a general use laboratory disinfectant. Rinse thoroughly with clean water and dry with a soft cloth.



**Warning** If the unit has been in service, turn it off and disconnect the power cord connector before proceeding with any maintenance. ▲

### Clean Air Filter

Clean the air filter a minimum of four times per year.\*

1. Open the front lower door by grasping the bottom left corner.
2. Locate the grille on the door. See Figure 6-1. Grasp the middle of the grille material and gently pull out to remove.
3. Wash the filter material using water and a mild detergent.
4. Dry by pressing between two towels.
5. Install the filter back into the grille and attach the grille.

*\* The clean filter alarm occurs every three months as a reminder to clean the air filter. Depending upon environmental conditions, the filter may need to be cleaned or replaced more frequently. If the filter becomes torn or excessively dirty, a replacement can be purchased from VWR. Order part number 760203.*

### Clean the Condenser

Clean the condenser a minimum of once a year.\*

1. Open the front lower door by grasping the bottom left corner. See Figure 6-1.
2. Using a vacuum cleaner, exercising care to not damage the condenser fins, clean the condenser.

*\* Depending upon environmental conditions, the condenser may need to be cleaned more frequently.*

## Defrost the Chamber

1. Remove all product and place it in another freezer.
2. Turn the unit off and disconnect it from the power source.
3. Turn off the battery switch (O). See Figure 6-1.
4. Open all of the doors and place towels on the chamber floor.
5. Allow the frost to melt and become loose.
6. Remove the frost with a soft cloth.
7. After defrosting is complete, clean the interior with a non-chloride detergent. Rinse thoroughly with clean water and dry with a soft cloth.
8. Plug unit in and turn power switch on.
9. Turn the battery power switch to Standby mode (C)
10. Allow the freezer to operate empty overnight before reloading the product.

## Clean the Door Gasket

Clean the door gasket a minimum of once a month.\*

Using a soft cloth, remove any frost build-up from the gasket and door(s). The Clean Gasket alarm occurs every three months as a reminder to remove frost build-up from the gasket and door(s). Press the Silence key to disable the audible alarm.

*\*The door gasket may need to be cleaned more frequently if dirt or excessive frost build-up prevents the door from closing properly.*

The exterior door gasket provides an excellent seal that protects product, provides an energy efficient thermal barrier to keep cold air in and room temperature air out, and reduces frost build-up on the inner doors.

## Vacuum Relief Port

Because the door gasket seals so well, a vacuum can be created after a door opening. Warm air enters the cabinet, cools and contracts, creating a vacuum that pulls the door in tightly against the seal.

To equalize the pressure inside the cabinet after a door opening requires 1.5-3.0 cu.ft. of ambient air to be drawn into the cabinet. The amount of air required to equalize the pressure varies depending on the cabinet size, cabinet temperature, duration of door opening, inventory volume and the temperature/humidity of the ambient air. This unit is designed with a “vacuum relief port” that allows the pressure to be equalized.

## Vacuum Relief Port (cont)

The time required to draw 1.5-3.0 cu.ft. of air into the cabinet depends on two factors,

- a) size and number of paths available for air to enter cabinet, and
- b) pressure difference between internal cabinet and ambient room.

Cabinets with the vacuum relief port operating normally, (i.e. vacuum relief port is not iced over) will require a minimum of 30 seconds up to a maximum of 120 seconds for the cabinet to equalize. This is also a good indication that the exterior door is well sealed.

The vacuum relief port requires routine maintenance. It will ice over unless preventive measures are taken. If the vacuum relief port becomes iced over, the freezer will take several hours to equalize pressure.

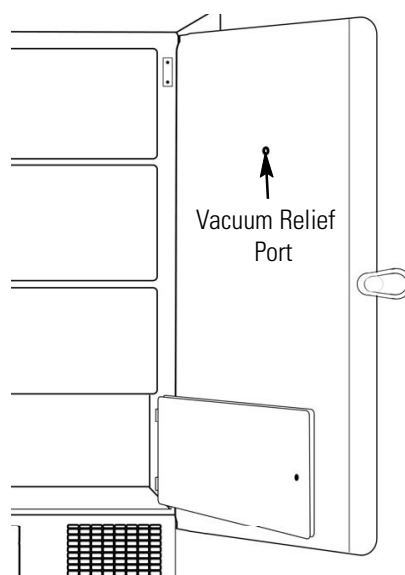


**Caution** Do not leave the freezer unattended while the door is unlatched. The vacuum could release at any time, resulting in the door opening and possible product loss. ▲

## Vacuum Relief Port Maintenance

Observe the inner side of port periodically for frost and ice build-up. Remove any frost with a soft dry cloth. If the tube should become clogged with ice, it must be cleaned. **Make sure during cleaning that the vacuum relief tube is completely free of ice to prevent rapid ice formation.**

Factors that can affect the vacuum relief port performance include: high ambient temperature, high humidity conditions and frequent door openings. Maintenance should be performed weekly or as needed.



**Caution** Failure to maintain the vacuum relief port may result in excessive ice build up inside the tube, clogging the port, and inability to open the door. The vacuum relief port may need to be cleaned more often with frequent door openings and high humidity environments. ▲

**Figure 6-2.** Port Location



## Check Battery

1. To gain access to the battery, open the lower door by grasping the bottom left corner. The battery is rectangular in shape, located on the front left corner of the compressor compartment and is secured in place by a mounting bracket.
2. Directly above the battery(s) is the battery power switch. Turn the battery power switch to the off position (O).
3. Cut the tie wrap securing the battery to the mounting bracket. Lift the battery out of the bracket.
4. Disconnect the red and black wires from the battery.
5. Use a voltmeter set to DC volts. Matching the wire colors, connect the meter to the battery.
6. If the voltage reads less than 10.8 volts, replace the battery. If above 10.8, re-install as previously.
7. Turn the battery power switch to Standby mode (⏻).
8. Close lower panel door.

## Replace the Battery

1. To gain access to the battery, open the lower door by grasping the bottom left corner. The battery is rectangular in shape, located on the front left corner of the compressor compartment and is secured in place by a mounting bracket.
2. Directly above the battery(s) is the battery power switch. Turn the battery power switch to the off position (O).
3. Disconnect the battery connections.
4. Remove the old battery and install the new battery.
6. Reconnect the battery (red to positive and black to negative).
7. Turn the battery power switch to Standby mode (⏻).
8. Close lower panel door.



**Warning** The % of charge can vary depending on the age, usage and condition of the battery. For a consistent and dependable charge, replace the battery every 2 years. Replacement batteries must be rechargeable and are available from VWR. Refer to the parts list for stock number and description of the replacement batteries. Dispose of the used batteries in a safe manner and in accordance with good environmental practices. ▲

## PREVENTIVE MAINTENANCE

### Tasks

Your equipment has been thoroughly tested and calibrated before shipment. Regular preventive maintenance is important to keep you and your food handling properly. The operator should perform routine cleaning and maintenance on a regular basis. For maximum performance and efficiency, recommendations for the unit is checked and calibrated periodically by a qualified service technician.

The following is a general list of preventive maintenance requirements. See the specific Section 6 of the operating manual for further details.

We have qualified service technicians, using NIST traceable instruments, available in many areas. For more information on Preventive Maintenance or Extended Warranties, please contact Technical Services.

Cleaning and calibration at shorter intervals are dependent upon use, environmental conditions and accuracy required.

### Tasks

- Pull an upright by starting at the bottom near the probe and pulling up on one side at a time. Allow freezer to recover to setpoint between cycles.
- Pull a chest by starting at the left side near the probe. Pulling with room temperature racks will result in a long pull-down time.
- Always make certain the vacuum relief port is free of frost and ice, making for timely pull-down. Use the reservoir as an opening.
- Fill unit with frozen product to be frozen, perform base frozen water tests, for example.

| Refer to Manual Section              | Action  | Monthly | Yearly | Every 2 Years |
|--------------------------------------|---|---------|--------|---------------|
| 4                                    | Verify alarm temperature control  | ✓       |        |               |
| 4                                    | Inspect door gasket for firm attachment   | ✓       |        |               |
| 4<br>4<br>4<br>4<br>4<br>4<br>4      | Check and seal drip cover, vacuum relief port, baskets, hoses and meter coils. Clean & log. | ✓       |        |               |
| 5                                    | Check air filter. Clean and replace as needed.  |         | ✓      |               |
| 4 & 5                                | Check alarm battery battery   | ✓       |        | Replace       |
| 4                                    | Check condenser fan motor for correct rotation & noise in VFD fan                           |         | ✓      |               |
| 4                                    | Verify alarm document calibration at the interval annual                                    |         | ✓      |               |
| 4                                    | Clean condenser compartment and area of condense  |         | ✓      |               |
| 4<br>5<br>5<br>5<br>5<br>5<br>5<br>5 | Verify Maintenance Tasks: Scales, Gas, Leak, Safety   | ✓       |        |               |

Qualified service technicians only.  
Disinfect properly according to local state and federal regulations.

To minimize ice buildup inside freezer:  
• Locate the freezer away from drafts or heat radiating vents.  
• Minimize the length of time doors are open.

• Reduce the number of door openings to a minimum.  
• Make sure doors are securely after opening.

## Section 7 Factory Installed Options

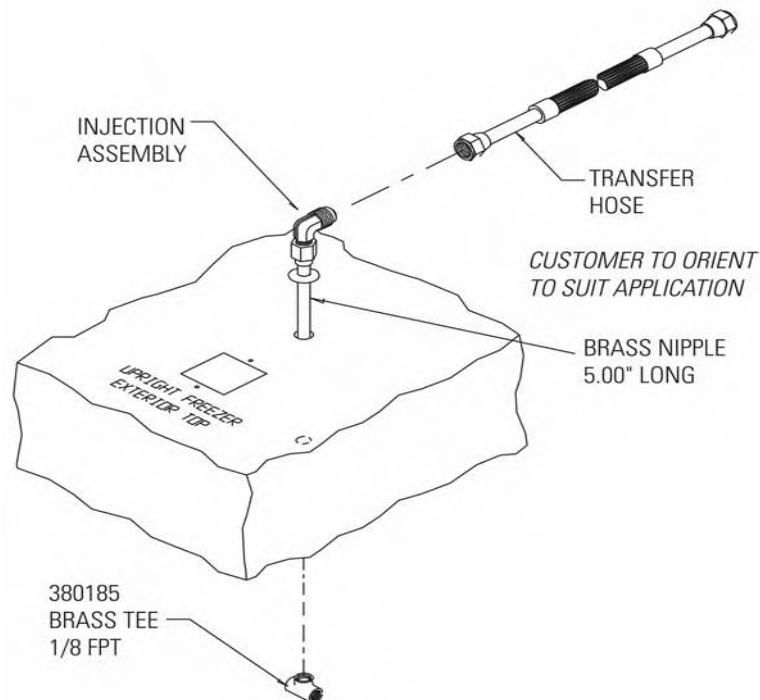
Options are available as factory or field installed, depending on the territory. Please contact your local VWR representative for details. Refer to the table in Appendix A-5 for the available field-installed options.

### Back-Up System (BUS)

**Warning** Before installation of BUS components, make sure the power to the freezer is disconnected, the battery switch is turned off (O) and the freezer has warmed to ambient temperature. ▲

The built-in BUS (back up system) will keep the freezer chamber temperature below the critical level in the event of a power or equipment failure. If power to the freezer fails, or temperature increases to the back up alarm set point, the BUS injects liquefied gas into the chamber to keep the chamber temperature within the specified range.

The BUS operates on an internal 12-volt, rechargeable battery which is kept charged during normal operation by the integral battery charger.



**Figure 7-1.** Injection

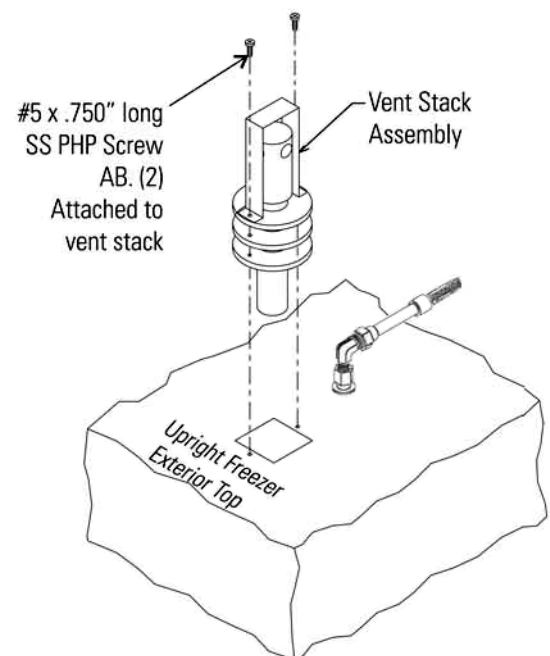
## Install Vent Stack, Solenoid & Injection Asm.

1. Install the injection assembly through the 1/2" pre-punched hole, directly behind the 2" vent stack hole in the center of the chamber ceiling.

**Note** Cover the open end of injection assembly with tape to keep insulation from entering the nipple. ▲

2. Slide 3/8" flatwasher over open end of nipple.
3. Insert the covered end of the injection assembly through the exterior hole.
4. Remove the tape covering from the end of the nipple and install the 1/8" NPT brass tee on the open end of the nipple. Place Permagum sealant between the brass tee and the interior top.
5. Remove the two Phillips head screws securing the metal bracket on the vent stack assembly.
6. Install the vent stack through the opening and secure it to the top of the freezer, using screws.
7. Go to the interior and seal around the end of the vent stack with Permagum.
8. Install the transfer hose connecting one end to the injection assembly, the other end to the solenoid valve. Install the solenoid valve to the supply source. The solenoid mounting bracket is not required and may be discarded.

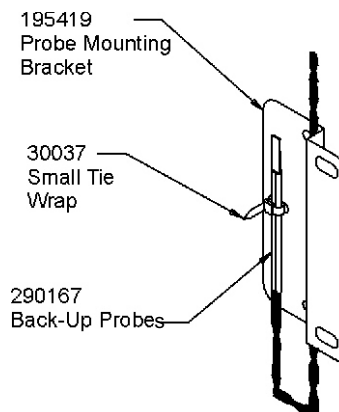
**Caution** When selecting a CO<sub>2</sub> supply cylinder, it must be equipped with a siphon tube. ▲



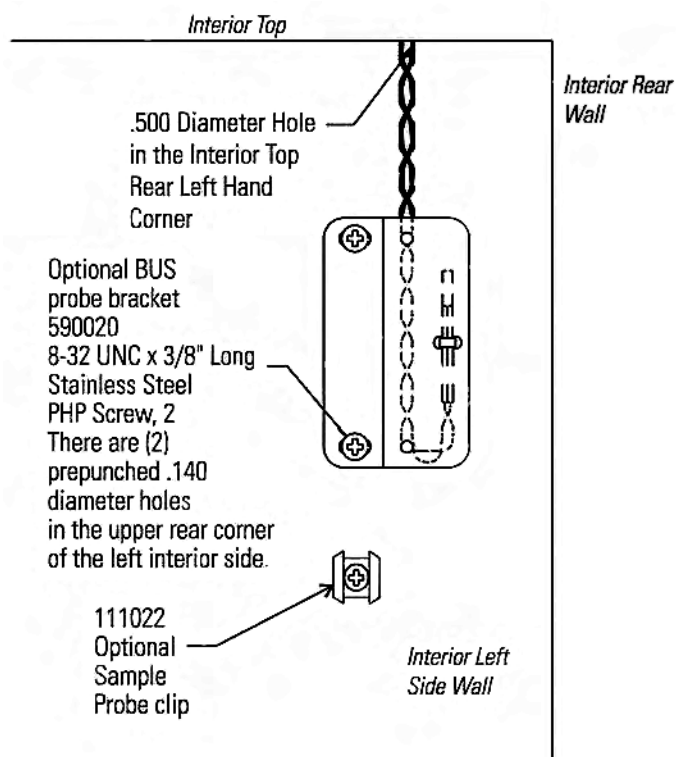
**Figure 7-2.** Vent Stack

## Install the Temperature Probe

1. Locate the 0.500" pre-punched hole in the upper left back corner of the chamber ceiling. Remove the tie wrap securing the coiled probe/solenoid harness. Uncoil the probe lead and run the probe tip (approximately 12") down through 0.500" porthole (Figure 7-4).
2. As in Figure 7-3, thread the small tie wrap through the openings in the front of the bracket. Secure the probe on the back of the bracket with the tie wrap.
3. Tap #8-32 the two pre-punched holes located on the interior left wall of the freezer. Mount the bracket. Figure 7-4 shows the Back-Up probe mounted on the interior left side wall of the freezer.

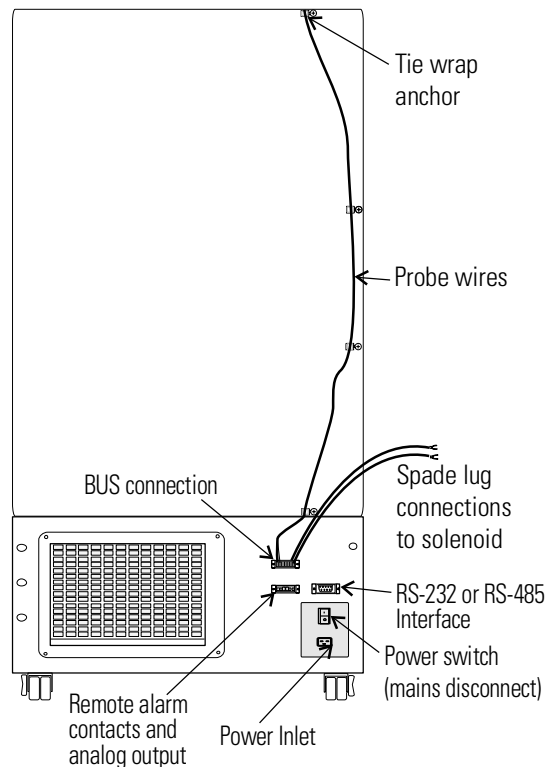


**Figure 7-3.** Probe



**Figure 7-4.** Probe Bracket

## Connect the Probe/Solenoid Harness



**Figure 7-5.** Connections

1. Remove the four screws on the freezer back panel and use them to mount the tie wrap anchors as shown in Figure 7-5. Secure the probe wire with tie wraps.
2. Plug the solenoid/probe connector into the BUS connection and secure with a screw on the right and left side. The connector is keyed.
3. Loosen the terminal screws on the solenoid. Slide the spade lug connectors under the screws and tighten to secure.
4. Connect power to the freezer. Turn the freezer On, with battery switch Off (O).
  - a. **Arming on Power On**

The arming function occurs only one time at initial power-on. It may occur in 3 ways; BUS timeout arming, both sensors fail and BUS temperature sensor arming.

At power-on, the BUS will perform a Hold-off period check. The BUS will not inject the refrigerant for a Hold-off period, calculated as follows:

## Connect the Probe/Solenoid Harness (cont.)

### Hold-off period:

- 12 hour is between 0 and +70C
- 12 hour is between 0 and (BUS Setpoint - Hysteresis)
- 0 hours is below (BUS Setpoint - Hysteresis)
- 0 hours is temperature Out Of Range or both probes fail.

b. The Low Battery indicator may also illuminate.

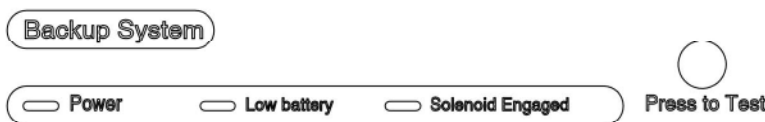
5. Turn the battery switch to Standby mode (⏻) to charge both batteries.

## BUS Operation and Maintenance

**Warning** When activated, this unit injects liquid nitrogen or carbon dioxide. Liquid nitrogen can cause serious freezing (frostbite) if it comes in contact with unprotected skin or eyes. Nitrogen and carbon dioxide gas suppresses oxygen levels and may cause suffocation if area is not well ventilated. Refer to Appendix A for the proper handling of liquid LN<sub>2</sub>. ▲

**Caution** Make sure the pressure relief valve on any LN<sub>2</sub> tank is adjusted to 30 PSI max blow-off. ▲

**Warning** Carbon dioxide gas suppresses oxygen levels and may cause suffocation if area is not well ventilated. Refer to “Handling Liquid CO<sub>2</sub> in Appendix B. ▲



**Figure 7-6.** Back-Up System (BUS) Control Panel

**Power** - indicates the unit has AC power.

1. Normal Operation

- Low Battery: Illuminates only after a battery test which occurs every 8 hours. The battery test is a loaded test and during the test, the battery voltage is monitored. When the battery voltage is less than a certain threshold, this light will illuminate and stay illuminated until the next battery test.
- Solenoid Engaged: Illuminates any time the controller is activating the solenoid. Illumination will only occur when the controller senses the solenoid is active and calling for injection.

## BUS Operation and Maintenance (cont.)

### 2. Fault Code

This same user interface will also permit fault codes to be easily viewed when the user presses and releases the Press to Test button in less than 2 seconds. When this occurs, both the Low Battery LED and Solenoid Engaged LED will blink twice ON/OFF quickly indicating that fault code display mode is active. For the next ~10 seconds, the fault codes will be displayed as shown in the table below. The display shows only the lower number fault code until that fault is resolved.

| Low Battery LED | Solenoid Engaged LED | Fault Code Number | Fault Code Definition   |
|-----------------|----------------------|-------------------|-------------------------|
| OFF             | OFF                  | 0                 | Normal Operation        |
| OFF             | Blinks 1 time        | 1                 | Primary Probe Failure   |
| OFF             | Blinks 2 times       | 2                 | Secondary Probe Failure |
| OFF             | Blinks 3 times       | 3                 | Solenoid Driver Failure |
| OFF             | Blinks 4 times       | 4                 | Low Tank                |

**Table 7-1.** BUS Fault Codes

Each blink will take one second, with half of the second interval being ON (illuminated) and the other half of the second interval being OFF. After the sequence concludes, both the Low Battery LED and the Solenoid Engaged LED will flash four times ON/OFF quickly to indicate a return to normal display.

### 3. BUS Injection Test

The user must press and hold the Press-to-Test button for  $\geq 2$  seconds before injection occurs. After the Press-to-Test button is pressed for  $>30$  seconds, the solenoid will turn off. This will avoid a button or button signal becoming stuck in an active condition.

During a Press-to-Test injection, if either RTD #1 or RTD #2 reads a temperature less than 20°C below BUS set point and it is determined to be a valid input, then injection will terminate immediately.

### 4. BUS Set point “Validation” Feedback

Whenever the BUS set point or working fluid changes, the BUS acknowledges the set point acceptance by causing the Low Battery LED and Solenoid Engaged LED to blink three times ON/OFF quickly. This does not imply that the BUS error-checked the values, other than that the value received was within the range of the BUS working correctly with prior programming.

**Note** Solenoid will not engage if door is open. ▲



## BUS Operation and Maintenance (cont.)

**Caution** The back-up system is designed to inject refrigerant (CO<sub>2</sub> or LN<sub>2</sub>). In the unlikely event of both probes failing, the back-up system will operate in PWM or Non-PWM mode based on jumper setting (Figure 7-2). ▲

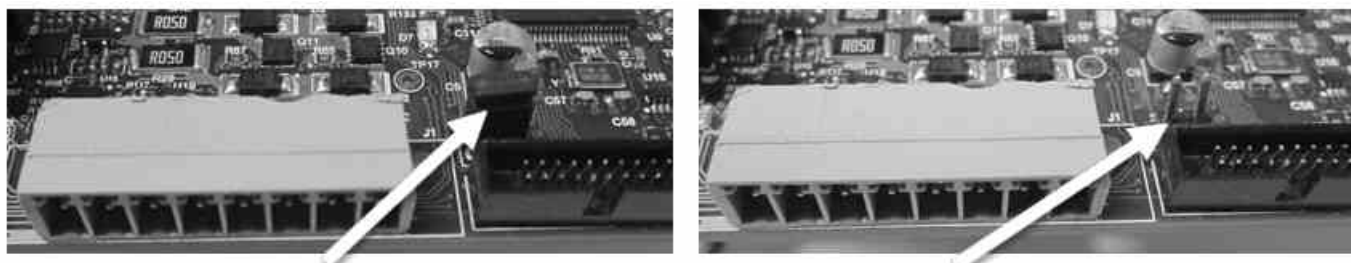
For Hysteresis & Dual probe fault mode settings (Figure 7-2), locate Jumper J17 near to solenoid/ sensor connector (J1), refer to table below for operation details:

| Jumper PIN        | Status                               | Hysteresis | Dual Probe Fault Mode |
|-------------------|--------------------------------------|------------|-----------------------|
| J17 PIN 1 & PIN 2 | Short Together/ Jumper Cap Installed | 5°C        | Non-PWM Mode          |
| J17 PIN 1 & PIN 2 | Open/No-Jumper Installed             | 9°C        | PWM Mode              |

**Table 7-2.** Probe Fault Modes

PWM mode: If both probes fail, the solenoid will activate periodically, allowing the flow of refrigerant (CO<sub>2</sub> or LN<sub>2</sub>).

Non-PWM mode: If both probes fail, the solenoid remains OFF and no refrigerant will flow



**Figure 7-7.** Jumper Cap for Hysteresis & Dual Probe Fault Mode Selection

### Set Optional BUS Setpoint

The optional back up system is designed to inject CO<sub>2</sub> or LN<sub>2</sub> into the freezer compartment if the temperature rises above back up system set point. To set the BUS set point:

1. Press the Mode key until the Settings indicator lights.
2. Press the right arrow until “BACKUP = -XX” is displayed in the message center.
3. Press the up or down arrow key until the desired BUS set point is displayed.
4. Press Enter to save the setting.
5. Press the Mode key until the Run indicator lights or press the right or left arrow to go to next or previous parameter.

If no control keys are pressed, the freezer automatically returns to RUN mode after 5 minutes.

**Set Optional BUS Setpoint (cont.)**

**Caution** The BUS setpoint cannot be set any colder than the high temperature alarm setpoint (see Section 2). If the back-up system is installed with CO<sub>2</sub>, then -65°C is the coldest BUS setpoint that can be used (if the cabinet setpoint is -75°C or colder).

Changing the operating temperature setpoint can affect the BUS setpoint. The BUS setpoint will self-adjust to maintain a temperature of at least 10°C above the operating temperature setpoint. ▲

**Test BUS Operation**

After the freezer has stabilized and both batteries are fully charged, the BUS can be tested to verify proper operation.

1. Disconnect the AC power to the freezer by turning the power switch off.
2. As the freezer warms up, verify the BUS injects at the desired temperature. Displayed temperature may vary by a few degrees from inject temperature due to the differences in probe locations.
3. Preventive Maintenance - Monthly maintenance action to check CO<sub>2</sub> or LN<sub>2</sub> back-up system operation, and battery voltage.  
  
- Check for fault codes on the BUS probe per Table 7-1. Also verify solenoid operation by performing a Press-To-Test.

**Clean the Vent Stack**

Routinely check the vent stack for frost or ice build-up. The type of frost that forms in the vent stack is generally very soft and may be easily removed with a bristle brush or soft cloth. If ice build-up has occurred, a complete defrost may occasionally be required. See Section 5 for freezer defrost instructions.

**Disconnect the Fitting Assembly and Transfer Hose**

To disconnect the freezer back-up from the gas supply:

1. Close the supply valve.
2. Depress the test button on the Back-Up System control box to remove the gas from the line.
3. Slowly disconnect the fitting assembly from the supply (in the event that any gas remains in the line).

## Chart Recorder

To install the chart paper in the recorder, follow the steps below.

1. Open the glass door of the recorder and press button #3 until the pen begins to move outward.
2. Unscrew the knob at the center of the chart and remove the paper.
3. Install the new chart paper, position the paper to the correct time line and replace the knob.
4. Remove the cap from the felt pen and press button #3.

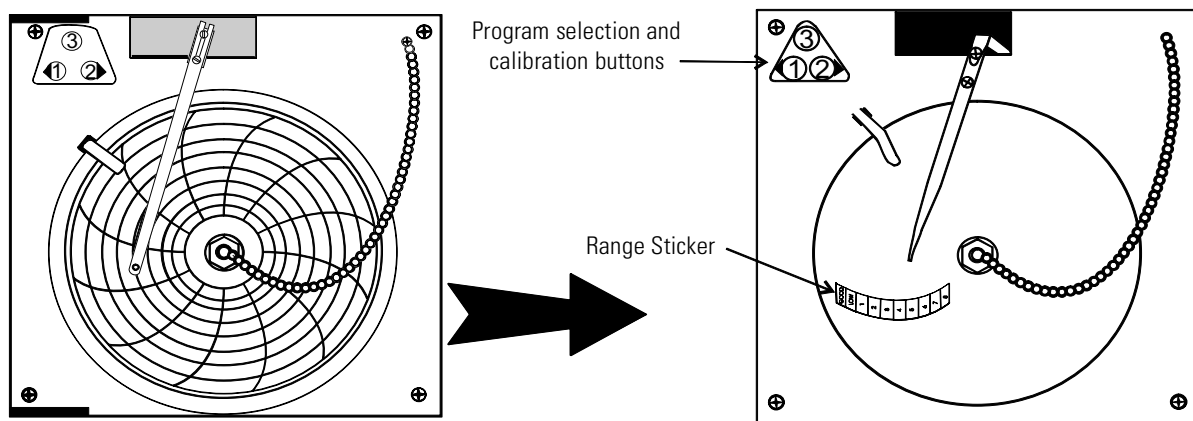


Figure 7-7. Recorder Details

## Change the Recorder Temperature Range

The chart recorder contains eight temperature ranges and is factory-programmed for the freezer. A list of the programs with temperature ranges follows.

1. Press and hold button #3 for one second, then let the pen move off the chart paper.
2. Press and hold for five seconds either button #1 or button #2.
3. Release the button and the green LED will begin to flash. Count the number of flashes to determine the present program setting.
4. To change the program setting, press the left or right arrows to increase or decrease the count.
5. When the desired program number is flashing, press button #3 to bring the pen arm back onto the chart. Recording will begin in the new program.

| Range | From | To    |
|-------|------|-------|
| 1     | -40  | 30°C  |
| 2     | 0    | 60°C  |
| 3     | -100 | 38°C  |
| 4     | -5   | 50°C  |
| 5     | 0    | 100°C |
| 6     | -100 | 200°C |
| 7     | -115 | 50°C  |
| 8     | -10  | 70°C  |

Table 7-3. Recorder Temp Range

## **Calibrate the Recorder**

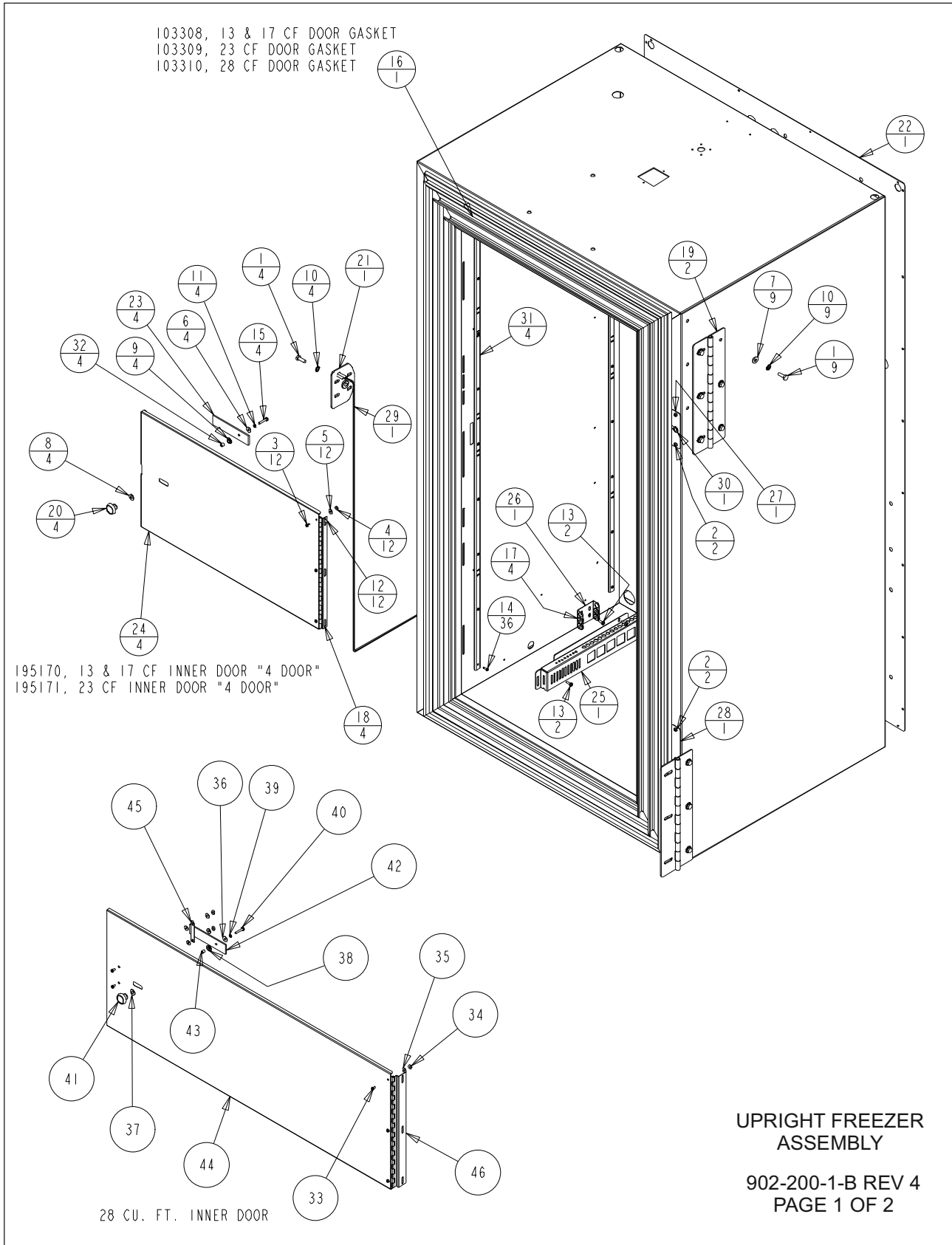
**Caution** The recorder must be in service for 24 hours before performing the following calibration procedure. ▲

1. Place an accurate thermometer in the chamber next to the recorder probe.
2. Temperature probes for the recorder are located in the left front corner of the freezer chamber (Figure 2-4).
3. After about three minutes, compare the thermometer reading with the chart recorder reading.
4. If an adjustment is necessary, press the #1 button to move the pen to the left or the #2 to move the pen to the right. The button must be held about five seconds before the pen begins to move. Release the button when the pen position matches the thermometer.

**Note** The felt-tip pen on the recorder requires periodic replacement. Usually the ink will appear to fade before replacement becomes necessary. Additional pen tips are available for purchase. ▲

**Section 6**

Factory Installed Options



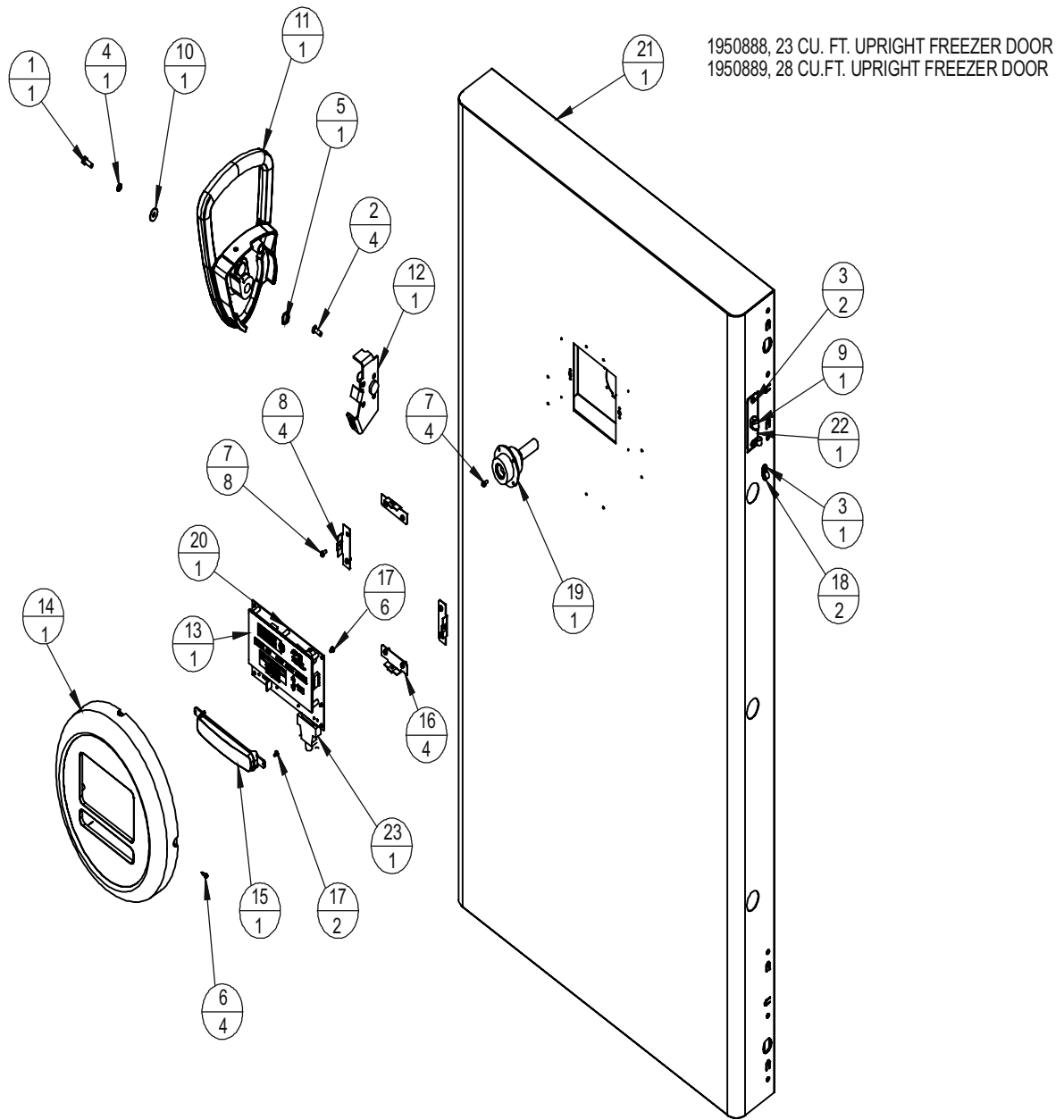
### BILL OF MATERIALS

| ITEM NO. | PART NO. | PART DESCRIPTION                      |
|----------|----------|---------------------------------------|
| 1        | 20003    | 1/4-20 X 3/4 SS HH CAP SCREW          |
| 2        | 22053    | #8-32 X 1/2 SS PHP SCREW              |
| 3        | 22115    | #6-32 X 1/4 SS PHP SCREW              |
| 4        | 23009    | #6-32 SS HEX NUT                      |
| 5        | 23020    | #6 SS FLAT WASHER                     |
| 6        | 23021    | #8 SS FLAT WASHER                     |
| 7        | 23023    | 1/4 SS FLAT WASHER                    |
| 8        | 23043    | NYLON FLAT WASHER                     |
| 9        | 23044    | 1/4" NYLON SHOULDER WASHER            |
| 10       | 23062    | 1/4 SS EXT TOOTH LOCKWASHER           |
| 11       | 23080    | #8 SS SPRING LOCKWASHER               |
| 12       | 24032    | #8-32 X 3/8 SS PHP SCREW F POINT      |
| 13       | 24042    | #8-32 X 1/2 SS PHP SCREW F POINT      |
| 14       | 25018    | 1/8 X 1/4 SS POP RIVET                |
| 15       | 59008    | #8-32 X 7/8 SS PHP SCREW              |
| 16       | 103308   | 13/17 CU.FT. SINGLE DOOR FRAME GASKET |
| 17       | 114020   | 5/8" X 1/2" ID GROMMET                |
| 18       | 116077   | FRONT PANEL HINGE                     |
| 19       | 116092   | EXTERIOR FREEZER DOOR HINGE           |
| 20       | 120400   | BLACK PLASTIC KNOB                    |
| 21       | 121099   | FRZ CAM LATCH STRIKE                  |
| 22       | 189921   | EXTERIOR BACK 13 & 17                 |
| 23       | 195169   | LATCH TAB                             |
| 24       | 195170   | 13/17 CU.FT. INNER DOOR               |
| 25       | 195866   | PROBE GUARD                           |
| 26       | 195867   | PROBE MOUNT                           |
| 27       | 195830   | UR DOOR WIREWAY COVER PLATE           |
| 28       | 195900   | MAGNETIC SWITCH                       |
| 29       | 30033    | RT. ANGLE STRAIN RELIEF               |
| 30       | 500177   | PILASTER STRIPS                       |
| 31       | 515083   | 1/4 DIA. X 1/4L SS SPACER             |
| 32       | 22115    | #6-32 X 1/4 SS PHP SCREW              |
| 33       | 23009    | #6-32 SS HEX NUT                      |
| 34       | 23020    | #6 SS FLAT WASHER                     |
| 35       | 23021    | #8 SS FLAT WASHER                     |
| 36       | 23043    | NYLON FLAT WASHER                     |
| 37       | 23044    | 1/4" NYLON SHOULDER WASHER            |
| 38       | 23080    | #8 SS SPRING LOCKWASHER               |
| 39       | 59008    | #8-32 X 7/8 SS PHP SCREW              |
| 40       | 120400   | BLACK PLASTIC KNOB                    |
| 41       | 195169   | LATCH TAB                             |
| 42       | 515083   | 1/4 DIA. X 1/4L SS SPACER             |
| 43       | 1950217  | 28 CU. FT. INNER DOOR                 |
| 44       | 1950218  | 28 CU.FT. INNER DOOR LATCH GUIDE      |
| 45       | 116090   | FRONT PANEL HINGE                     |
| 46       | 24030    | 8 X 1/2 ZP PHP SCREW TEKS             |

UPRIGHT FREEZER  
ASSEMBLY

902-201-1-B REV 4  
PAGE 2 OF 2

**Section 8**  
Parts



UPRIGHT FREEZER  
DOOR ASSEMBLY  
5602-201-1-B REV 9  
PAGE 1 OF 2

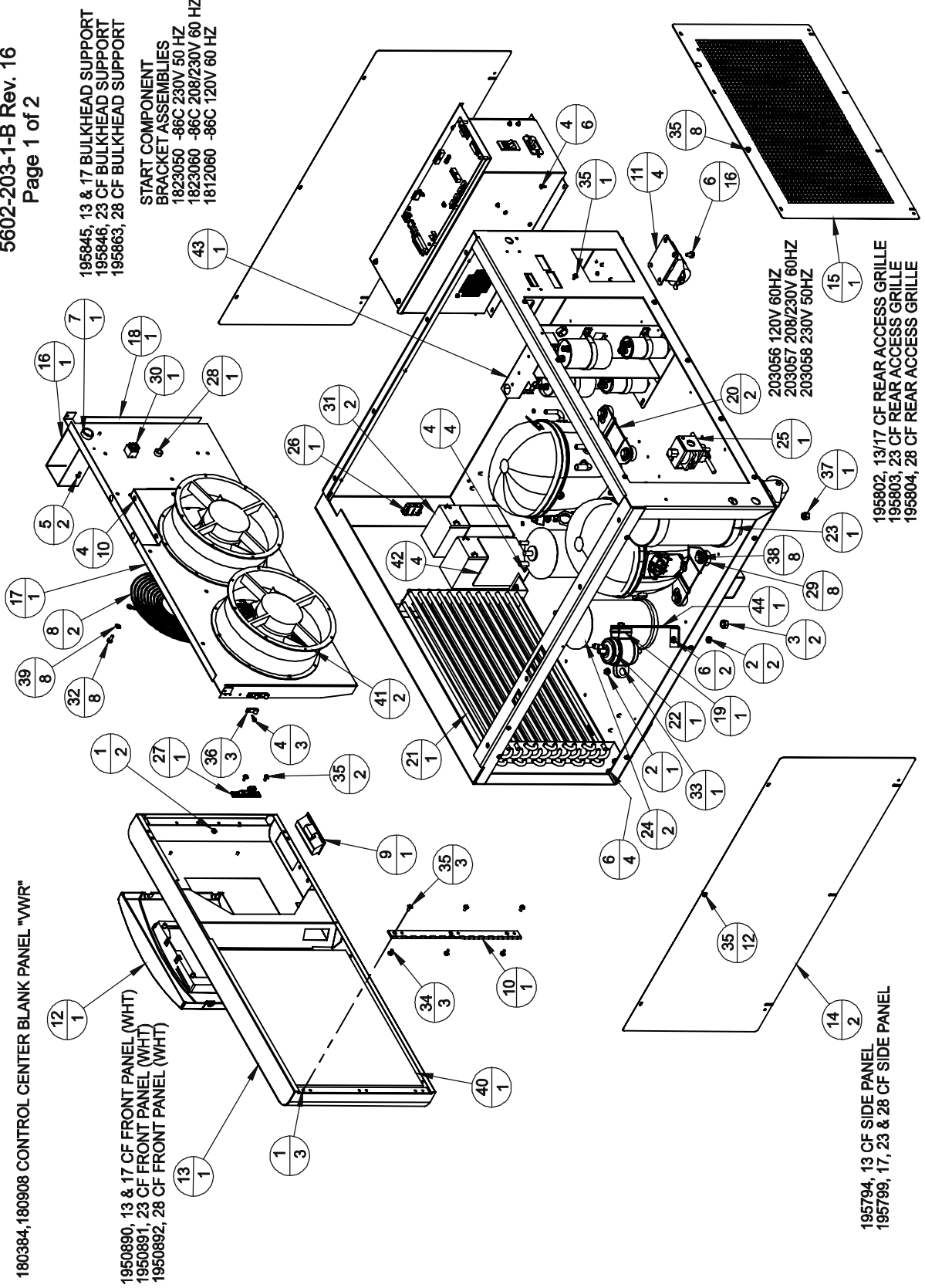
| BILL OF MATERIALS |          |                                      |
|-------------------|----------|--------------------------------------|
| ITEM NO.          | PART NO. | PART DESCRIPTION                     |
| 1                 | 20003    | 1/4-20 X 3/4 SS HH CAP SCREW         |
| 2                 | 20058    | #1/4-20 X 3/4 SS FHP UC SCREW        |
| 3                 | 22053    | #8-32 X 1/2 SS PHP SCREW             |
| 4                 | 23033    | 1/4 SS INTERNAL TOOTH LOCK WASHER    |
| 5                 | 23057    | 5/8 WAVE WASHER                      |
| 6                 | 24016    | #6 X 1/2" SS PHP SCREW AB POINT      |
| 7                 | 24032    | #8-32 X 3/8 SS PHP SCREW F POINT     |
| 8                 | 25040    | #6 U SPEED NUT STL. STL.             |
| 9                 | 30033    | RIGHT ANGLE STRAIN RELIEF            |
| 10                | 121092   | .856" OD FLAT WASHER                 |
| 11                | 121104   | FINISHED HANDLE VWR PMS 647 "BLU"    |
| 12                | 121102   | CAM LATCH MOUNT                      |
| 13                | 140474   | CONTROL PANEL                        |
| 14                | 180387   | VWR ROUND PLASTIC CONTROL BEZEL      |
| 15                | 180318   | VWR BACK-UP SYSTEM BLANK PANEL       |
| 16                | 1950895  | BRACKET MOUNTING                     |
| 17                | 590027   | #6-32 X 1/4 SS PHP EXT SEMS SCREW    |
| 18                | 600085   | 5/16 NYLON CABLE CLAMP               |
| 19                | 1950069  | HEATED VACUUM RELIEF PORT            |
| 20                | 191941   | FREEZER DISPLAY BOARD                |
| 21                | 1950887  | 13 & 17 CU. FT. UPRIGHT FREEZER DOOR |
| 22                | 195830   | UPRIGHT DOOR WIREWAY COVER PLATE     |
| 23                | 430336   | 15 FT, RS-232 CABLE 25 POS.          |

UPRIGHT FREEZER  
DOOR ASSEMBLY

5602-201-1-B REV 9  
PAGE 2 OF 2



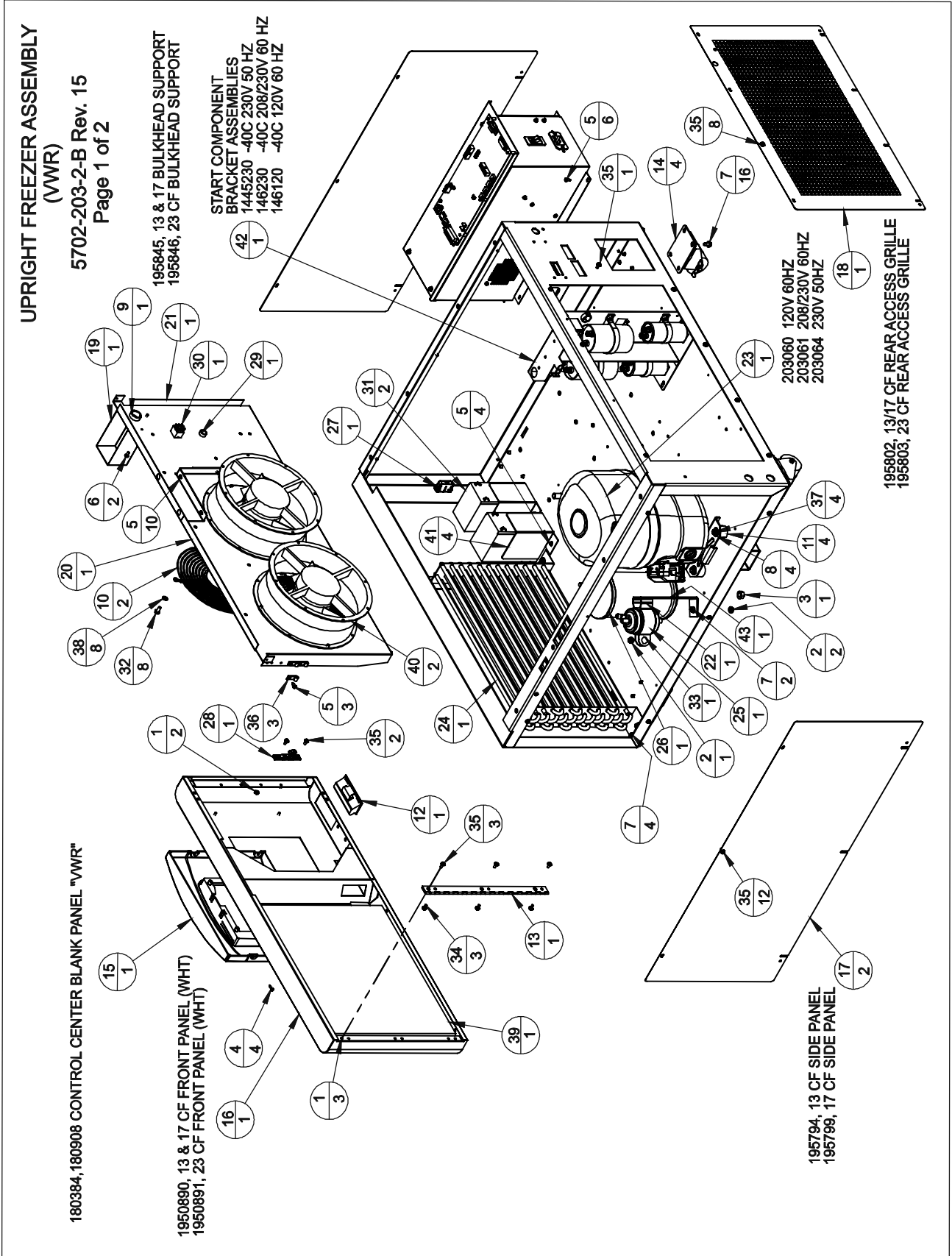
**UPRIGHT FREEZER ASSEMBLY  
(VWR)**  
5602-203-1-B Rev. 16  
Page 1 of 2



UPRIGHT FREEZER ASSEMBLY  
(VWR)  
5602-203-1-B Rev. 16  
Page 2 of 2

| BILL OF MATERIALS |  |
|-------------------|--|
| ITEM NO.          | PART DESCRIPTION                                 |
| 1                 | 23002 #8-32 ZP LKWASH HEX NUT                    |
| 2                 | 23011 1/4-20 ZP LKWASH HEX NUT                   |
| 3                 | 23013 3/8-16 ZP LKWASH HEX NUT                   |
| 4                 | 24030 #8 X 1/2" TEKS SCREW                       |
| 5                 | 24032 #8-32 X 3/8 SS PHP SCREW F POINT           |
| 6                 | 24038 1/4-20 X 1/2 SELF TAPPING SCREW            |
| 7                 | 30016 1" SNAP BUSHING                            |
| 8                 | 108020 10" WIRE FAN GUARD                        |
| 9                 | 115032 BLACK ABS PLASTIC PULL                    |
| 10                | 116115 FRONT PANEL HINGE                         |
| 11                | 120011 DUAL WHEEL CASTER                         |
| 12                | 180384 THERMO CONTROL CENTER BLANK PANEL         |
| 13                | 1950890 13&17 CU. FT. UR FRZ BASE FRONT PNL      |
| 14                | 195799 SIDE PANEL 17, 23 AND 28 CU. FT. UPRIGHT  |
| 15                | 195802 13/17 REAR ACCESS GRILLE                  |
| 16                | 195829 MULLION /DOOR SWITCH WIRE COVER           |
| 17                | 195844 UR FRZ FAN BULKHEAD                       |
| 18                | 195845 13&17 CU.FT. FAN BULKHEAD SUPPORT         |
| 19                | 200126 2" RIGID HANGER                           |
| 20                | 203057 UPRIGHT COMPRESSOR 208/230V 60HZ          |
| 21                | 204009 REFRIGERATION CONDENSER                   |
| 22                | 209020 LIQUID LINE FILTER DRYER WITH ACCESS PORT |
| 23                | 214018 10.000" H X 5.000" DIA. EXPANSION TANK    |
| 24                | 214023 OIL SEPARATOR                             |
| 25                | 220626 120V - 50/60 HZ SOLENOID VALVE            |
| 26                | 319388 LATCH CATCH, PART OF 319388 ASSEMBLY      |
| 27                | 319388 LATCH KEEPER, PART OF 319388 ASSEMBLY     |
| 28                | 330002 5/8" SNAP BUSHING                         |

| BILL OF MATERIALS |             |   |
|-------------------|-------------|---|
| ITEM NO.          | PART NO.    | PART DESCRIPTION                              |
| 29                | 350096      | RUBBER MOUNT FOOT EMBRACO                     |
| 30                | 360248      | MINI SNAP-IN POWER SWITCH                     |
| 31                | 400159      | SEALED LEAD ACID BATTERY - 12 VOLT - 7.2 Ah   |
| 32                | 510035      | #12-24 X 1/2 SS HH CAP SCREW                  |
| 33                | 550043      | 1/4-20 X 1" L ZP CARRIAGE BOLT                |
| 34                | 590020      | #8-32 X 3/8 SS PHP EXT SEMS SCREW             |
| 35                | 590029      | #8-32 X 3/8 SS PHP EXT SEMS SCREW W/PATCH     |
| 36                | 600080      | 1/4 ALUM CLAMP W/LINER                        |
| 37                | 610053      | 3/8-16 NYLON INSERT LOCK NUT                  |
| 38                | 680030      | 5/16-18 X 1.75 SS HH SHOULDER BOLT            |
| 39                | 730087      | #12 SS EXT TOOTH LOCKWASHER                   |
| 40                | 760203      | AIR FILTER                                    |
| 41                | 900113      | 10" TUBEAXIAL FAN, 115V                       |
| 42                | 1950074     | BATTERY MOUNTING BRACKET                      |
| 43                | 120230-01-2 | UPRIGHT FREEZER START COMPONENT BRACKET ASSY. |
| 44                | 1950850     | FILTER DRIER MOUNTING BRACKET                 |

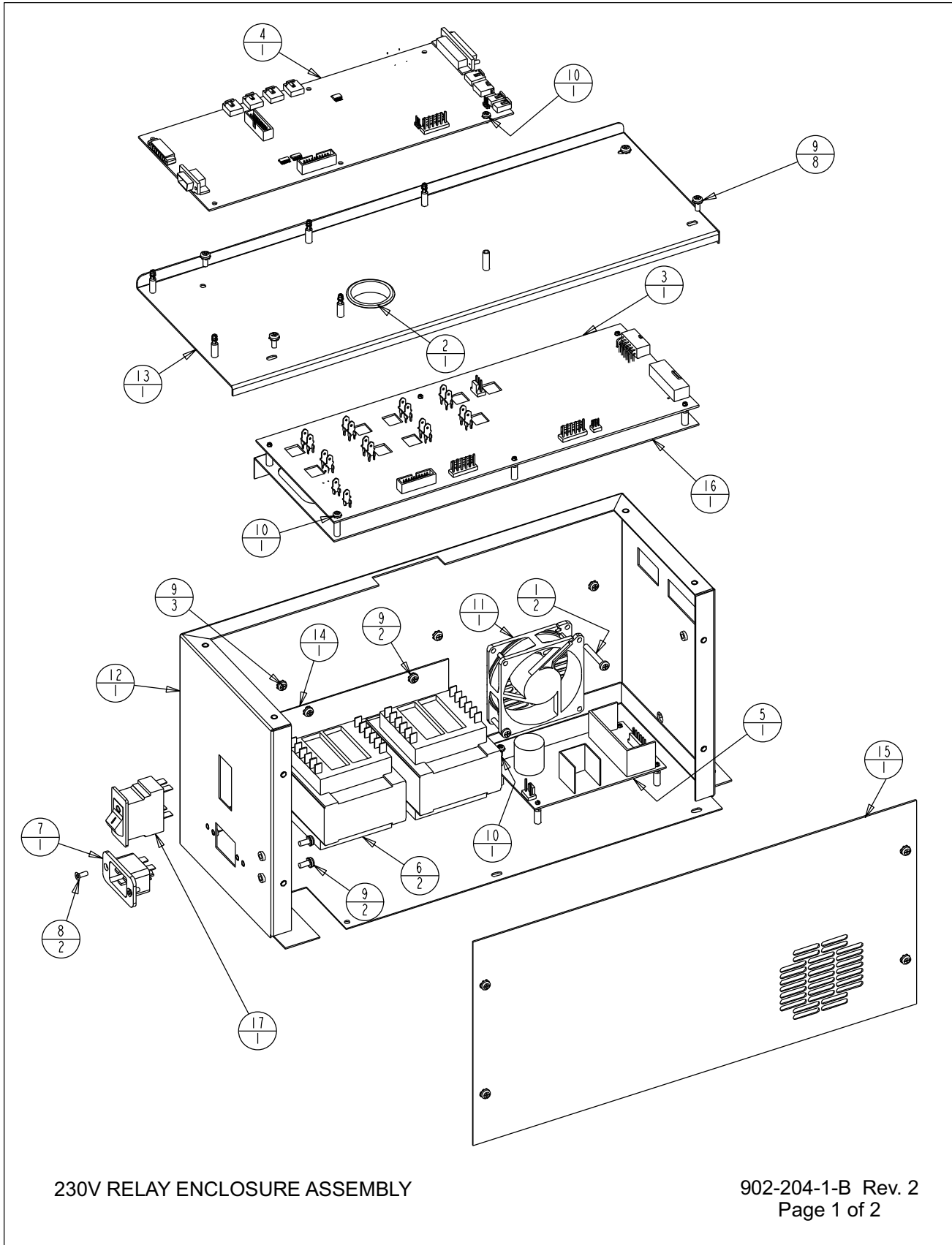


UPRIGHT FREEZER ASSEMBLY  
(VWR)  
5702-203-2-B Rev. 15  
Page 2 of 2

| BILL OF MATERIALS |  |
|-------------------|--|
| ITEM NO.          | PART NO. PART DESCRIPTION                        |
| 1                 | 23002 #8-32 ZP LKWASH HEX NUT                    |
| 2                 | 23011 1/4-20 ZP LKWASH HEX NUT                   |
| 3                 | 23013 3/8-16 ZP LKWASH HEX NUT                   |
| 4                 | 24016 #6 X 1/2" SS PHP SCREW AB POINT            |
| 5                 | 24030 #8 X 1/2" TEKS SCREW                       |
| 6                 | 24032 #8-32 X 3/8 SS PHP SCREW F POINT           |
| 7                 | 24038 1/4-20 X 1/2 SELF TAPPING SCREW            |
| 8                 | 24049 1/4 ZP FLAT WASHER                         |
| 9                 | 30016 1" SNAP BUSHING                            |
| 10                | 108020 10" WIRE FAN GUARD                        |
| 11                | 114034 COMPRESSOR MOUNTING SLEEVE                |
| 12                | 115032 BLACK ABS PLASTIC PULL                    |
| 13                | 116115 FRONT PANEL HINGE                         |
| 14                | 120011 DUAL WHEEL CASTER                         |
| 15                | 180384 THERMO CONTROL CENTER BLANK PANEL         |
| 16                | 1950890 13&17 CU. FT. UR FRZ BASE FRONT PNL      |
| 17                | 195799 SIDE PANEL 17, 23 AND 28 CU. FT. UPRIGHT  |
| 18                | 195802 13/17 REAR ACCESS GRILLE                  |
| 19                | 195829 MULLION /DOOR SWITCH WIRE COVER           |
| 20                | 195844 UR FRZ FAN BULKHEAD                       |
| 21                | 195845 13&17 CU.FT. FAN BULKHEAD SUPPORT         |
| 22                | 200126 2" RIGID HANGER                           |
| 23                | 203060 BRISTOL HIGH/LOW STAGE COMPRESSOR         |
| 24                | 204009 REFRIGERATION CONDENSER                   |
| 25                | 209020 LIQUID LINE FILTER DRYER WITH ACCESS PORT |
| 26                | 214023 OIL SEPARATOR                             |
| 27                | 319388 LATCH CATCH, PART OF 319388 ASSEMBLY      |
| 28                | 319388 LATCH KEEPER, PART OF 319388 ASSEMBLY     |

| BILL OF MATERIALS |   |
|-------------------|---|
| ITEM NO.          | PART NO. PART DESCRIPTION                                 |
| 29                | 330002 5/8" SNAP BUSHING                                  |
| 30                | 360248 MINI SNAP-IN POWER SWITCH                          |
| 31                | 400159 SEALED LEAD ACID BATTERY - 12 VOLT - 7.2 Ah        |
| 32                | 510035 #12-24 X 1/2 SS HH CAP SCREW                       |
| 33                | 550043 1/4-20 X 1" L ZP CARRIAGE BOLT                     |
| 34                | 590020 #8-32 X 3/8 SS PHP EXT SEMS SCREW                  |
| 35                | 590029 #8-32 X 3/8 SS PHP EXT SEMS SCREW W/PATCH          |
| 36                | 600080 1/4 ALUM CLAMP W/LINER                             |
| 37                | 680008 1/4-20 X 1-3/4 SELF TAPPING SCREW                  |
| 38                | 730087 #12 SS EXT TOOTH LOCKWASHER                        |
| 39                | 760203 AIR FILTER   |
| 40                | 900113 10" TUBEAXIAL FAN, 115V                            |
| 41                | 1950074 BATTERY MOUNTING BRACKET                          |
| 42                | 120230-01-2 UPRIGHT FREEZER START COMPONENT BRACKET ASSY. |
| 43                | 1950850 FILTER DRIER MOUNTING BRACKET                     |

**Section 8**  
Parts



230V RELAY ENCLOSURE ASSEMBLY

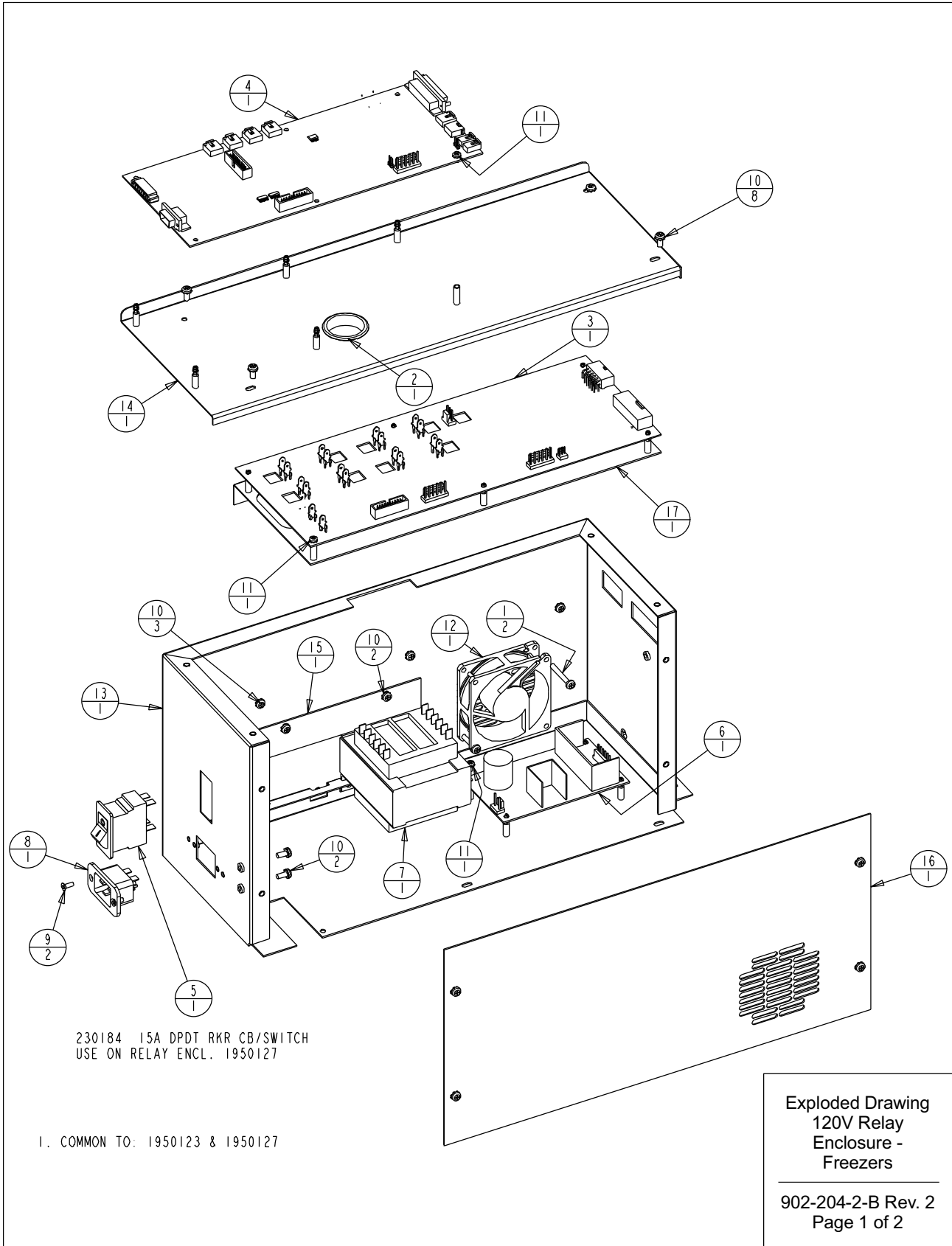
| BILL OF MATERIALS |             |                                       |
|-------------------|-------------|---------------------------------------|
| ITEM NO.          | PART NO.    | PART DESCRIPTION                      |
| 1                 | 22143       | #8-32 x 1-1/4 SS PHP SCREW            |
| 2                 | 30077       | 1-1/2" SNAP BUSHING                   |
| 3                 | 191923      | HIGH VOLTAGE BOARD 230V               |
| 4                 | 191989      | MICRO BOARD ( HIGH END)               |
| 5                 | 400165      | SWITCHER BOARD                        |
| 6                 | 420090      | 175V TRANSFORMER                      |
| 7                 | 460169      | POWER INLET, 16/20A                   |
| 8                 | 490009      | #6-32 X 3/8 SS FHP UC SCREW           |
| 9                 | 590020      | #8-32 X 3/8 SS PHP EXT SEMS SCREW     |
| 10                | 590027      | #6-32 X 1/4 SS PHP EXT SEMS SCREW     |
| 11                | 900134      | TUBEAXIAL FAN, 30 CFM, 12V            |
| 12                | 195631-16-1 | RELAY ENCLOSURE SPOTWELD SUB-ASSEMBLY |
| 13                | 195631-16-4 | RELAY ENCLOSURE COVER/191656 SUPPORT  |
| 14                | 195631-31-3 | TRANSFORMER HOLD DOWN                 |
| 15                | 195631-31-5 | RELAY ENCLOSURE COVER (MAIN)          |
| 16                | 195730-16-1 | 191658 SUPPORT BRACKET SUB-ASSEMBLY   |
| 17                | 230184      | 15A DPDT SWITCH/CIRCUIT BKR           |

I. COMMON TO: 1950124 & 1950128

230V RELAY ENCLOSURE ASSEMBLY

902-204-1-B Rev. 2  
Page 2 of 2

**Section 8**  
Parts



| BILL OF MATERIALS |             |                                       |
|-------------------|-------------|---------------------------------------|
| ITEM NO.          | PART NO.    | PART DESCRIPTION                      |
| 1                 | 22143       | #8-32 x 1-1/4 SS PHP SCREW            |
| 2                 | 30077       | 1-1/2" SNAP BUSHING                   |
| 3                 | 191932      | HIGH VOLTAGE BOARD 120V               |
| 4                 | 191989      | MICRO BOARD ( HIGH END)               |
| 5                 | 230184      | 15A DPDT SWITCH/CIRCUIT BKR           |
| 6                 | 400165      | SWITCHER BOARD                        |
| 7                 | 420065      | 175V TRANSFORMER                      |
| 8                 | 460169      | POWER INLET, 16/20A                   |
| 9                 | 490009      | #6-32 X 3/8 SS FHP UC SCREW           |
| 10                | 590020      | #8-32 X 3/8 SS PHP EXT SEMS SCREW     |
| 11                | 590027      | #6-32 X 1/4 SS PHP EXT SEMS SCREW     |
| 12                | 900134      | TUBEAXIAL FAN, 30 CFM, 12V            |
| 13                | 195631-16-1 | RELAY ENCLOSURE SPOTWELD SUB-ASSEMBLY |
| 14                | 195631-16-4 | RELAY ENCLOSURE COVER/191656 SUPPORT  |
| 15                | 195631-31-3 | TRANSFORMER HOLD DOWN                 |
| 16                | 195631-31-5 | RELAY ENCLOSURE COVER (MAIN)          |
| 17                | 195730-16-1 | 191658 SUPPORT BRACKET SUB-ASSEMBLY   |

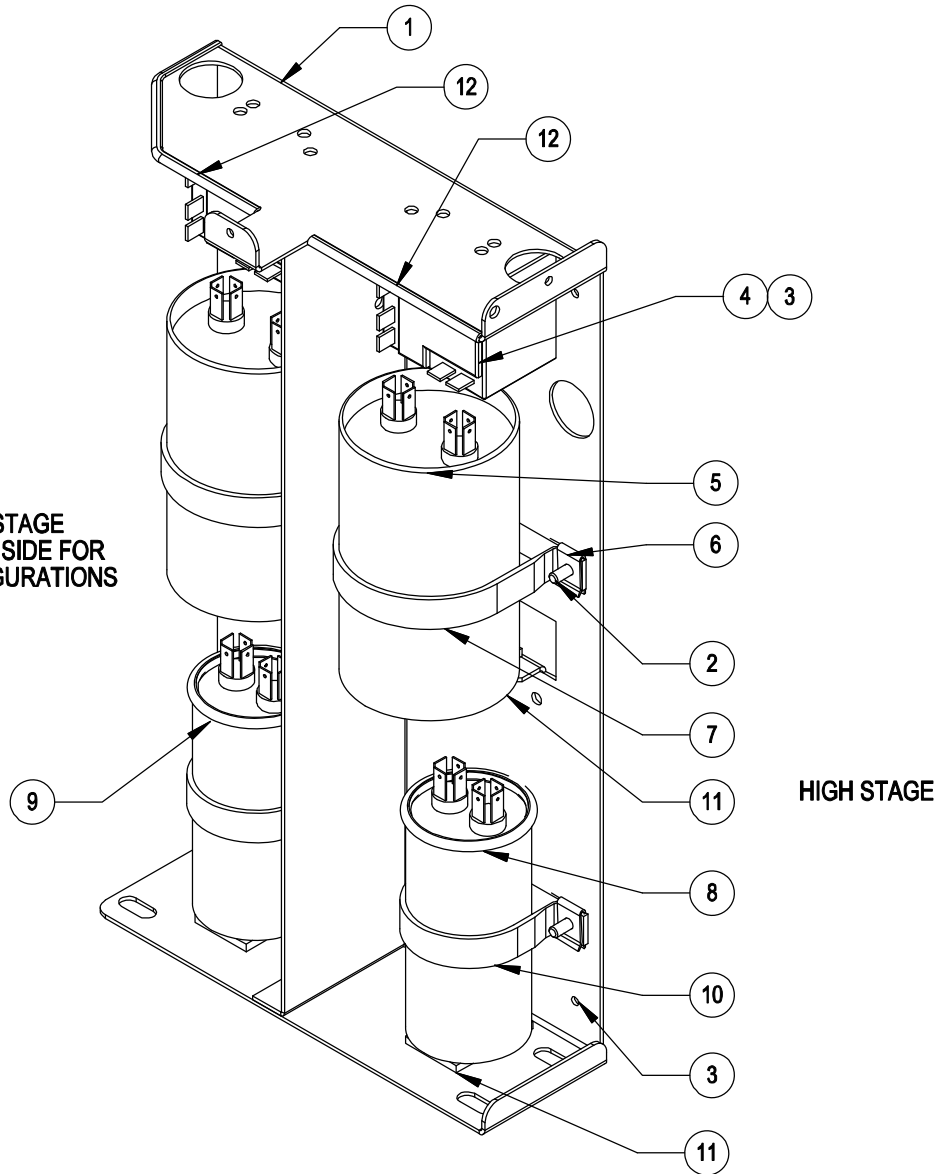
Exploded Drawing  
120V Relay  
Enclosure -  
Freezers

902-204-2-B Rev. 2  
Page 2 of 2



**Section 8**  
Parts

LOW STAGE  
USE THIS SIDE FOR  
-40 CONFIGURATIONS



HIGH STAGE

**NOTES:**

1. SEE PSS 8602-724-012 FOR ASSEMBLY WIRING AND OTHER DETAILS.
2. AFTER ATTACHING NOMEX COVER (250138) WITH SCREWS (590029), STICK ESD LABEL (220555) TO OUTSIDE OF COVER.
3. IDENTIFY EACH ASSEMBLY WITH LABELS ON THE BOTTOM AND BACK THAT INCLUDE THE FOLLOWING:  
 FIRST LINE - ASSEMBLY NUMBER AND REVISION  
 SECOND LINE - BAR CODE OF ASSEMBLY AND REVISION  
 THIRD LINE - UNIQUE SERIAL NUMBER INCLUDING DATE OF MANUFACTURE IN THE FOLLOWING FORMAT AND ORDER:  
 THREE LETTER MANUFACTURING SITE, TWO DIGIT YEAR MANUFACTURED, TWO DIGIT WEEK MANUFACTURED FOLLOWED  
 BY A UNIQUE FOUR DIGIT NUMBER.  
 FOURTH LINE - BAR CODE OF UNIQUE SERIAL NUMBER.  
 FIFTH LINE - ASSEMBLY DESCRIPTION IN THE FOLLOWING FORMAT: UPRIGHT, TEMPERATURE, VOLTAGE AND FREQUENCY.

Upright Freezer Start Component  
Enclosure Assembly

120230-01-2-D Rev 10  
Page 1 of 2

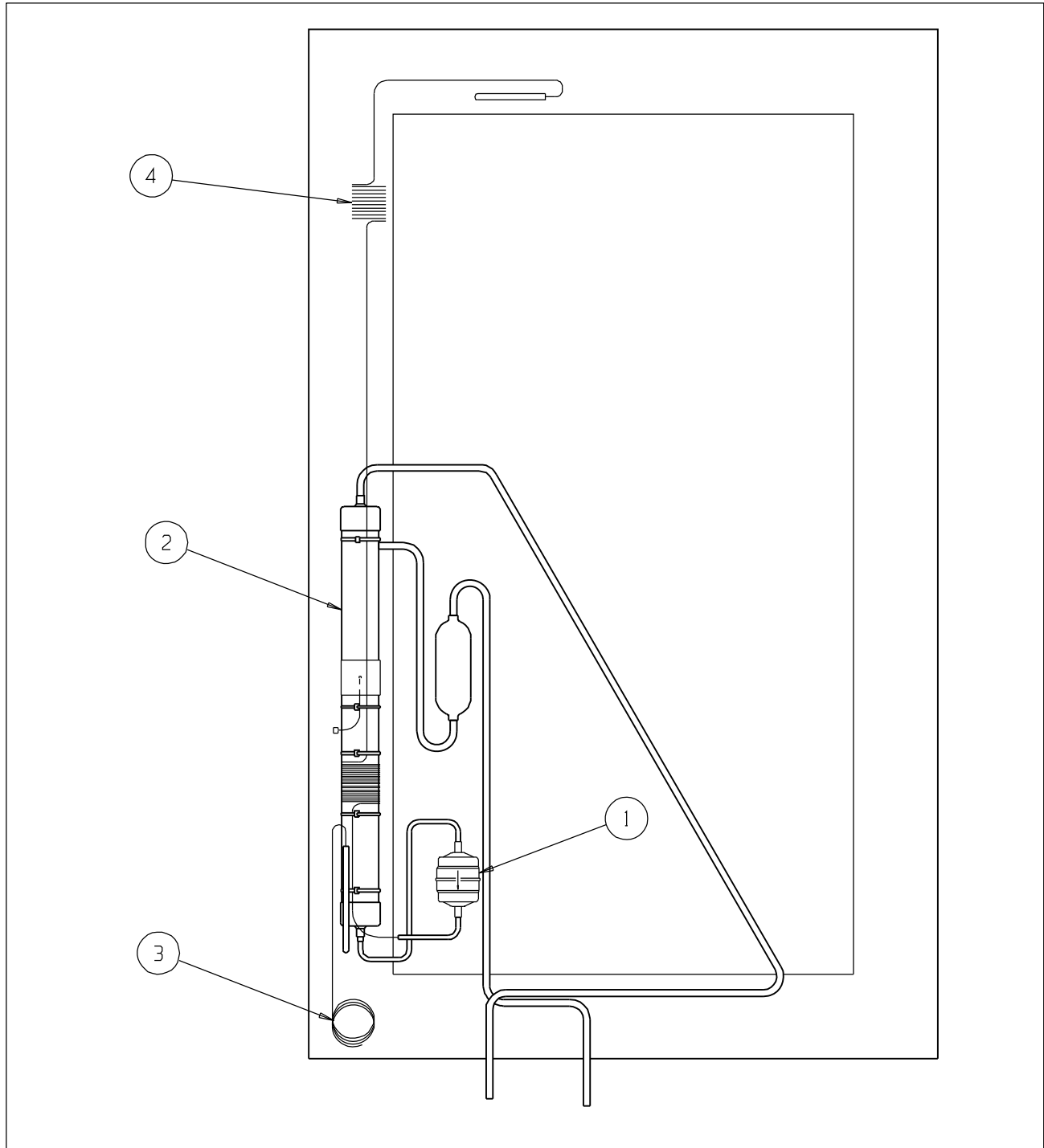
| UPRIGHT FREEZER START COMPONENT BRACKET ASSEMBLIES |                                 |                                      |                               |                            |                            |                              |                           |
|--|---------------------------------|--------------------------------------|-------------------------------|----------------------------|----------------------------|------------------------------|---------------------------|
| ITEM NO.   | DESCRIPTION                     | 1823050<br>-86C 230V 50 Hz           | 1823060<br>-86C 208/230V 60Hz | 1812060<br>-86C 120V 60 Hz | 1445230<br>-40C 230V 50 Hz | 146230<br>-40C 208/230V 60Hz | 146120<br>-40C 120V 60 Hz |
| 1  | COMPONENT BRACKET               | 120230                               | 120230                        | 120230                     | 120230                     | 120230                       | 120230                    |
| 2  | #8 X 1/2<br>PHP TYPE AB         | 24021 (4)                            | 24021 (4)                     | 24021 (4)                  | 24021 (2)                  | 24021(2)                     | 24021 (2)                 |
| 3  | 8-32 X 3/8 SCREW                | 590020 (4)                           | 590020 (4)                    | 590020 (4)                 | 590020 (2)                 | 590020 (2)                   | 590020 (2)                |
| 4  | START RELAY                     | 312286 (2)                           | 312186 (2)                    | 312086 (2)                 | 312287 (1)                 | 312187 (1)                   | 312087 (1)                |
| 5  | START CAPACITOR                 | 623050 (2)                           | 623060 (2)                    | 612060 (2)                 | 170236 (1)                 | 170154 (1)                   | 170154 (1)                |
| 6  | #8 SPEED NUT, S.S.              | 327675 (4)                           | 327675 (4)                    | 327675 (4)                 | 327675 (2)                 | 327675 (2)                   | 327675 (2)                |
| 7A   | 2.5" VINYL COATED<br>STRAP      | N/A                                  | N/A                           | 600106 (4)                 | N/A                        | N/A                          | 600106 (2)                |
| 7B   | 2" VINYL COATED STRAP           | 600108 (2)                           | 600108 (2)                    | N/A                        | 600108 (1)                 | 600108 (1)                   | N/A                       |
| 8  | RUN CAPACITOR                   | 223050                               | 223060                        | 212060                     | N/A                        | N/A                          | N/A                       |
| 9  | RUN CAPACITOR                   | 223050                               | 223050                        | 212060                     | 170097 (1)                 | 170226 (1)                   | 170224 (1)                |
| 10   | 1.75" VINYL COATED<br>STRAP     | 600107 (2)                           | 600107 (2)                    | SEE 7A                     | 600107 (1)                 | 600107 (1)                   | SEE 7A                    |
| 11   | NEOPRENE TAPE 1/8" X<br>1" X 2" | 28003 (4)                            | 28003 (4)                     | 28003 (4)                  | 28003 (2)                  | 28003 (2)                    | 28003 (2)                 |
| 12   | EDGE GUARD<br>(CUT TO LENGTH)   | 114012                               | 114012                        | 114012                     | 114012                     | 114012                       | 114012                    |
| NOT SHOWN *  | 13                              | ULT MAIN HARNESS<br>(REFERENCE ONLY) | 350092                        | 350092                     | 350092                     | 350093                       | 350093                    |
| NOT SHOWN  | 14                              | ENCLOSURE LOW STAGE<br>HARNESS       | 350108                        | 350108                     | 350108                     | 350108                       | 350108                    |
| NOT SHOWN  | 15                              | ENCLOSURE HIGH STAGE<br>HARNESS      | 350107                        | 350107                     | 350107                     | N/A                          | N/A                       |
| NOT SHOWN  | 16                              | NOMEX COVER                          | 250138                        | 250138                     | 250138                     | 250138                       | 250138                    |
| NOT SHOWN  | 17                              | ESD WARNING LABEL                    | 220555 (2)                    | 220555 (2)                 | 220555 (2)                 | 220555 (2)                   | 220555 (2)                |
| NOT SHOWN, SEE NOTE 3                              | 18                              | IDENTIFICATION LABEL                 | 823050 (2)                    | 823060 (2)                 | 812060 (2)                 | 423050 (2)                   | 423060 (2)                |

\* THE ULT MAIN HARNESS IS LISTED FOR INFORMATION ONLY. IT IS NOT PART OF THIS ASSEMBLY.

Upright Freezer Start Component  
Enclosure Assembly

120230-01-2-D Rev 10  
Page 2 of 2

**Section 8**  
Parts

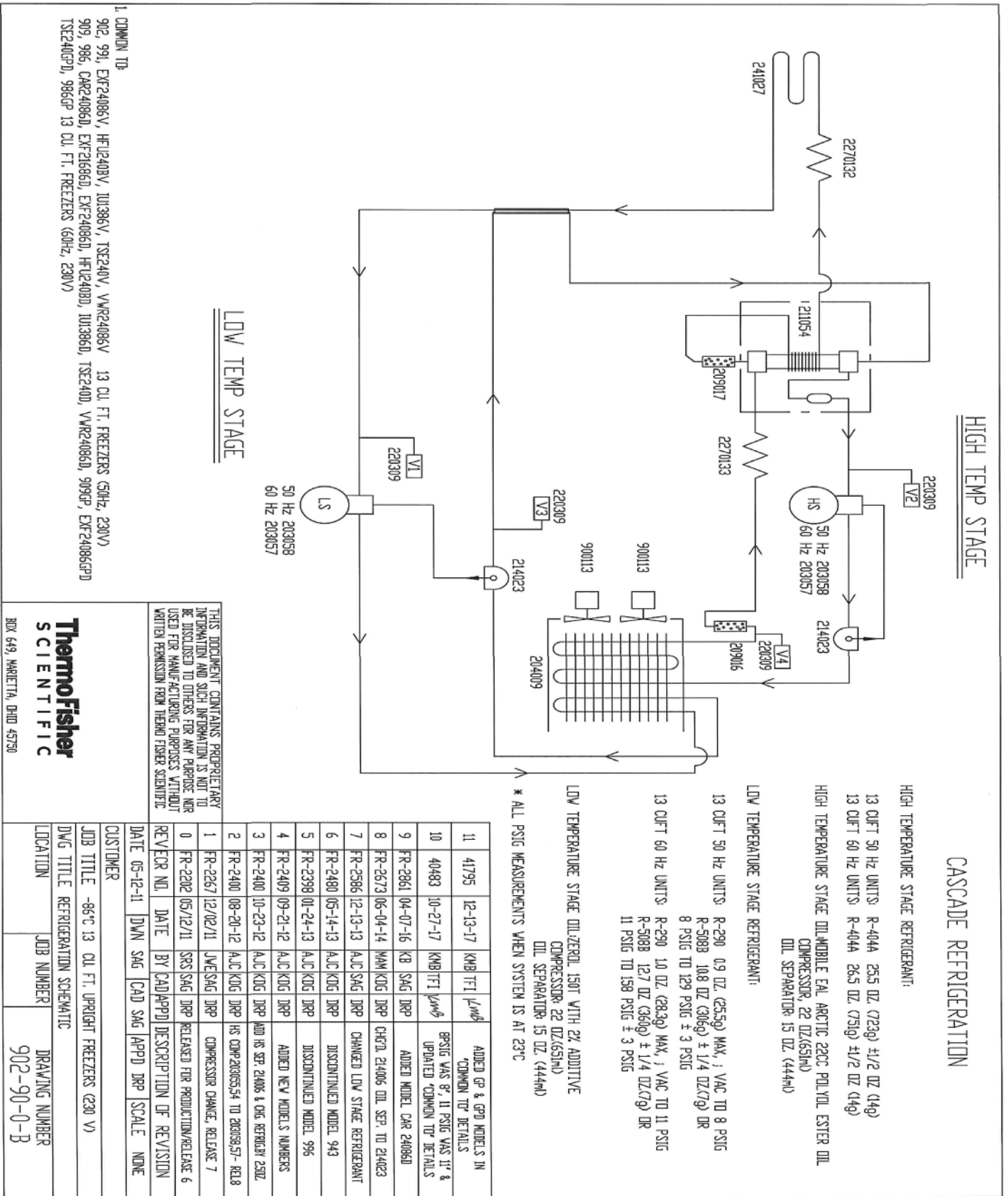


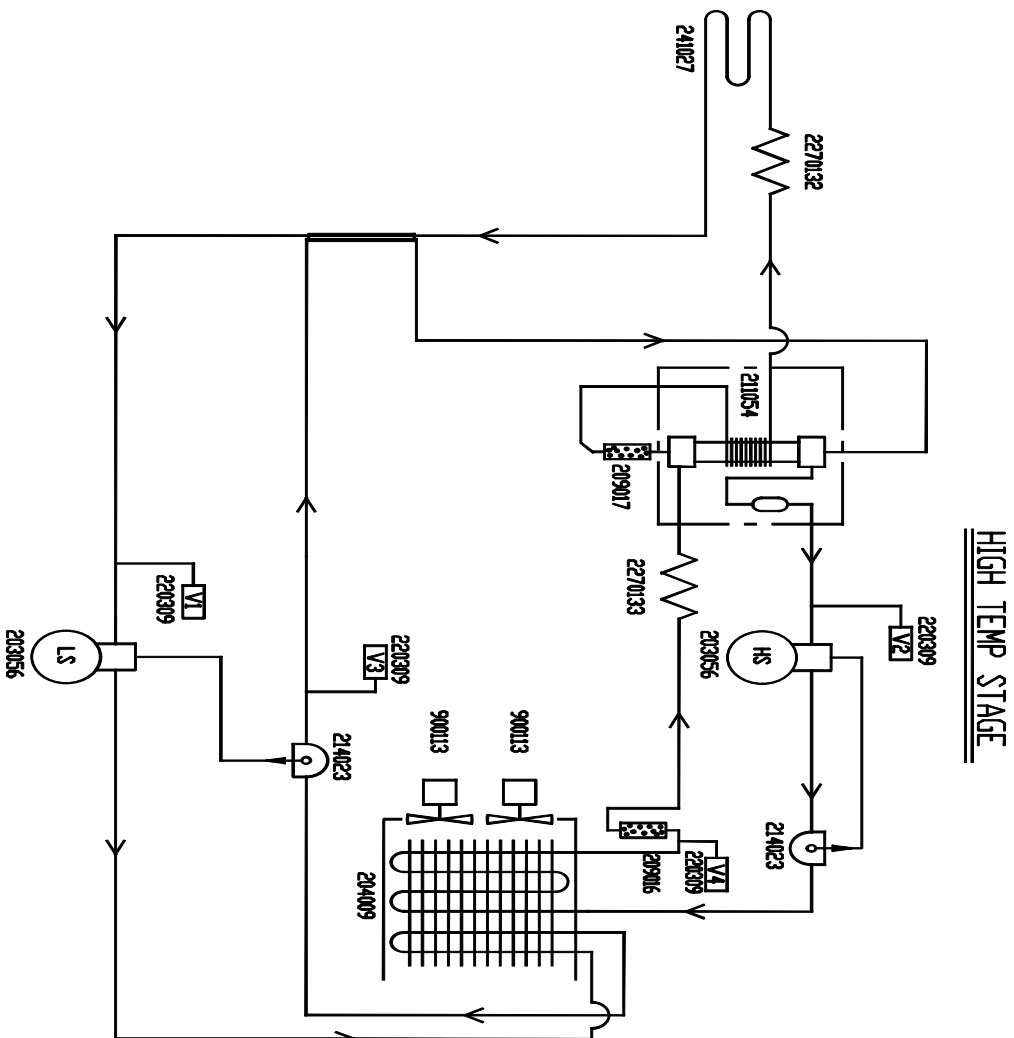
| BILL OF MATERIALS |          |                      |
|-------------------|----------|----------------------|
| ITEM NO.          | PART NO. | PART DESCRIPTION     |
| 1                 | 209017   | DRYER                |
| 2                 | 211050   | HEAT EXCHANGER       |
| 3                 | 227927   | HIGH STAGE CAP. TUBE |
| 4                 | 227928   | LOW STAGE CAP. TUBE  |

Exploded Drawing  
Heat Exchanger  
Upright Freezer

---

902-205-1-B Rev. 0  
Page 1 of 1





HIGH TEMP STAGE

LOW TEMP STAGE

CASCADE REFRIGERATION

HIGH TEMPERATURE STAGE REFRIGERANT:

13 CUFT UNITS R-404A 255 OZ (72g) 1/2 OZ (14g)

HIGH TEMPERATURE STAGE OIL/MODULE EAL ARCTIC 22CC POLYOL ESTER OIL  
COMPRESSOR 22 OZ (651ml)  
OIL SEPARATOR 15 OZ (444ml)

LOW TEMPERATURE STAGE REFRIGERANT:

13 CUFT UNITS R-290 08 OZ (227g) MAX, 1 VAC TO 8 PSIG  
R-5088 10.0 OZ (283.5g) ± 1/4 OZ (7g) DR  
8" TO 128 PSIG ± 3 PSIG

LOW TEMPERATURE STAGE OIL/ZERO L10T WITH 2% ADDITIVE  
COMPRESSOR 22 OZ (651ml)  
OIL SEPARATOR 15 OZ (444ml)

\* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C

THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC.

|   |            |          |      |      |                        |  |      |       |      |
|---|------------|----------|------|------|------------------------|--|------|-------|------|
| 7 | FR-2861    | 04-07-16 | KJ   | SAG  |                        | ADDED MODEL CAC24086A                  |      |       |      |
| 6 | FR-2673    | 06-04-14 | MM   | KDG  | DRP                    | CHG. 21406 OIL SEP. TO 214023          |      |       |      |
| 5 | FR-2586    | 12-13-13 | AJC  | SAG  | DRP                    | CHANGED LOW STAGE REFRIGERANT          |      |       |      |
| 4 | FR-2409    | 10-23-12 | AJC  | KDG  | DRP                    | ADDED NEW MODEL NUMBERS                |      |       |      |
| 3 | FR-2400    | 10-23-12 | AJC  | KDG  | DRP                    | ADD HS SEP. 21406 & OIL SEPARATOR 2102 |      |       |      |
| 2 | FR-2400    | 08-20-12 | AJC  | KDG  | DRP                    | HS COMP. 213053 TO 213056 - REL8       |      |       |      |
| 1 | FR-2267    | 12/02/11 | JVES | SAG  | DRP                    | COMPRESSOR CHANGE, RELEASE 7           |      |       |      |
| 0 | FR-2202    | 05-12-11 | SRS  | SAG  | DRP                    | RELEASED FOR PRODUCTION/RELEASE 6      |      |       |      |
|   | REVER. NO. | DATE     | BY   | APPD | DESCRPTION OF REVISION |  |      |       |      |
|   | DATE       | 05-12-11 | DVN  | SAG  | CAD                    | SAG                                    | APPD | SCALE | NONE |

CUSTOMER

JOB TITLE -86°C 13 CU. FT. UPRIGHT FREEZERS (220 V)

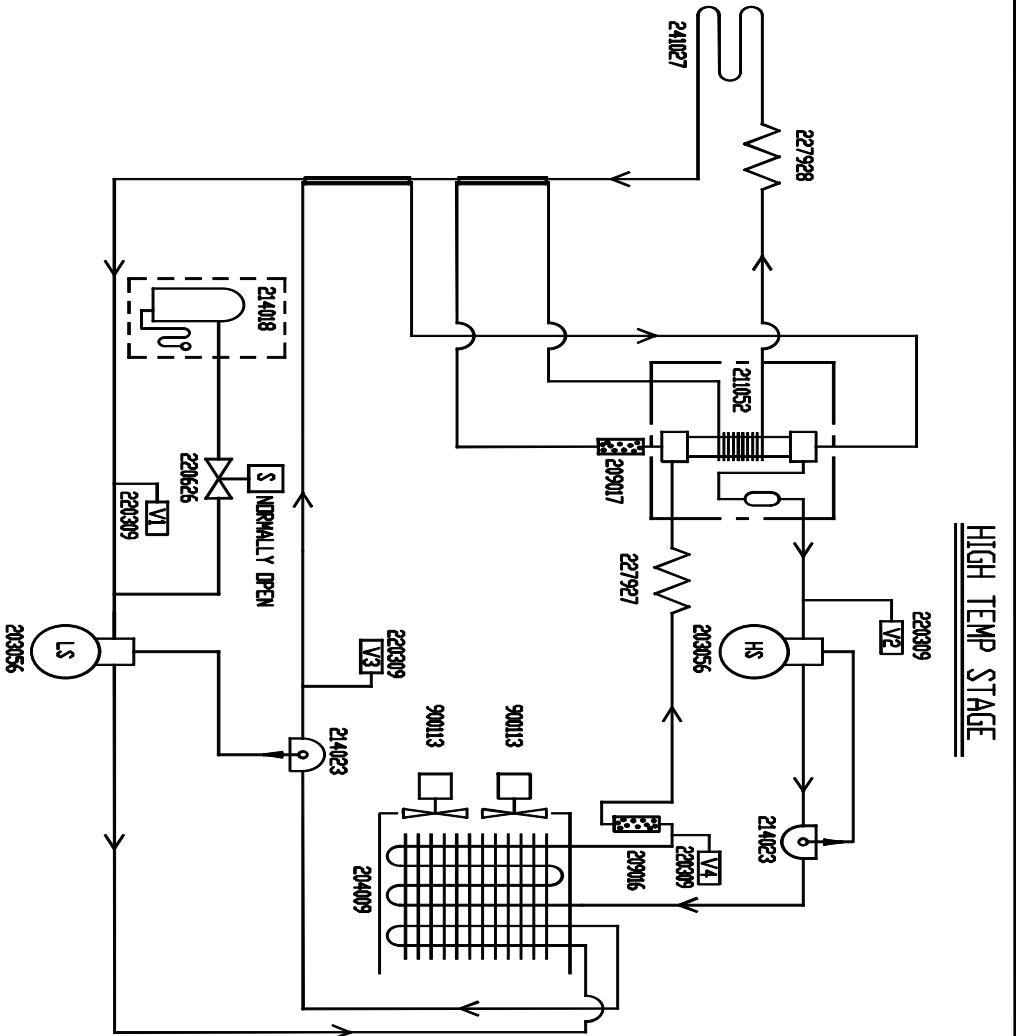
DWG TITLE REFRIGERATION SCHEMATIC

LOCATION

JOB NUMBER 903-90-0-B

**ThermoFisher**  
**SCIENTIFIC**  
BOX 649, WAUKESHA, WI 53186

1. COMMON TO: 903, 930, 992, 5703, 5603, 8922, 8933, 8955, 8967  
5903, 5992, 6422, 6433, 5930, CAC24086A 13 CU. FT. FREEZERS (220 V)



HIGH TEMP STAGE

LOW TEMP STAGE

**CASCADE REFRIGERATION**

HIGH TEMPERATURE STAGE REFRIGERANT:

17 CUFT UNITS R-404A 26.5 OZ (750g) 1/2 OZ (14g)

HIGH TEMPERATURE STAGE OIL/MOBILE EQ. ARCTIC 220C PL.VOL. ESTER OIL  
COMPRESSOR 22 OZ (650ml)  
OIL SEPARATOR 15 OZ (444ml)

LOW TEMPERATURE STAGE REFRIGERANT:

17 CUFT UNITS R-290 11 OZ (312g) MAX, 1 VAC TO 7 PSIG  
R-5088 14.4 OZ (408g) ± 1/4 OZ (7g) OR  
7 PSIG TO 128 PSIG ± 3 PSIG

LOW TEMPERATURE STAGE OIL/ZERO 150T WITH 2% ADDITIVE  
COMPRESSOR 22 OZ (650ml)  
OIL SEPARATOR 15 OZ (444ml)

\* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C

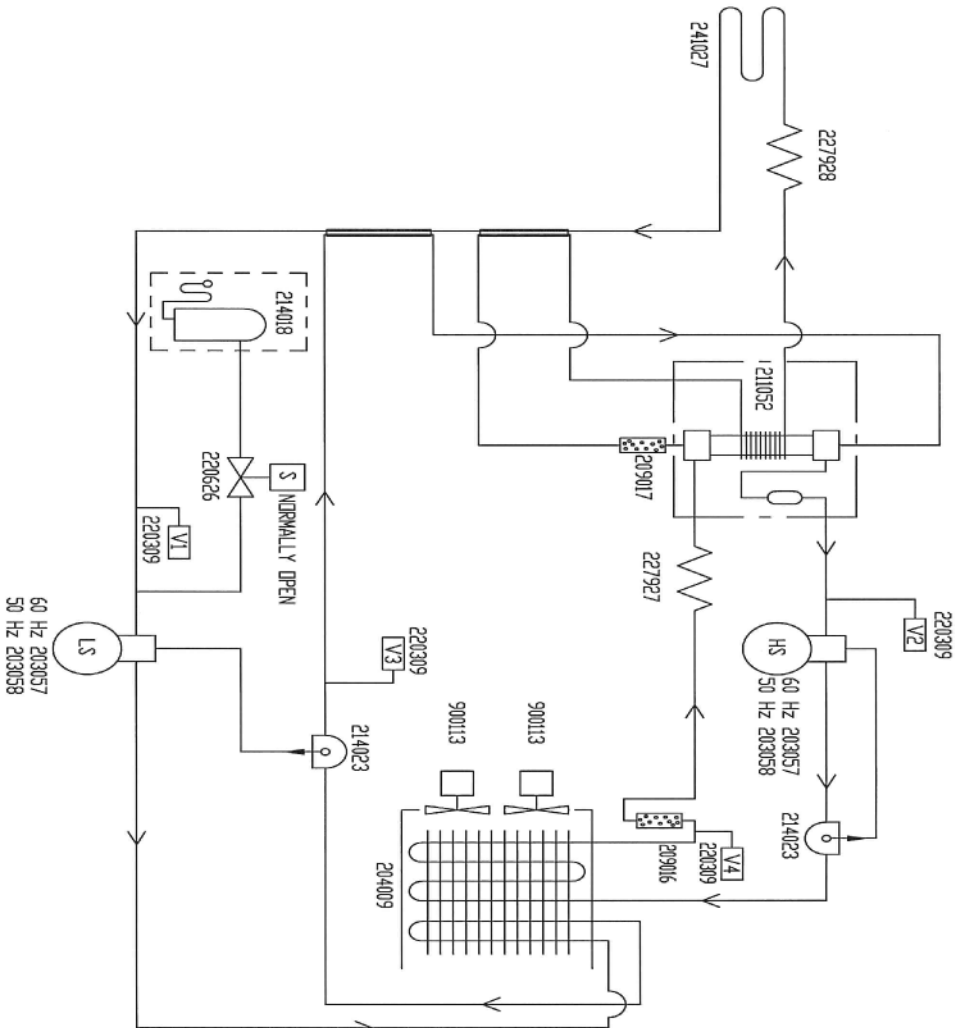
THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC

|  |          |          |     |     |      |                                     |     |       |      |
|--|----------|----------|-----|-----|------|-------------------------------------|-----|-------|------|
| 8  | FR-2861  | 04-07-16 | KB  | SAG |      | ADDED MODEL CAR22086A               |     |       |      |
| 7  | FR-2673  | 06-04-14 | MAN | KDG | DRP  | CHG. 21406 OIL SEP. TO 214023       |     |       |      |
| 6  | FR-2506  | 01-02-14 | AJC | SAG | DRP  | CHANGED LOW STAGE REFRIGERANT       |     |       |      |
| 5  | FR-2506  | 12-13-13 | AJC | SAG | DRP  | CHANGED LOW STAGE REFRIGERANT       |     |       |      |
| 4  | FR-2409  | 10-23-12 | AJC | KDG | DRP  | ADDED NEW MODEL'S NUMBERS           |     |       |      |
| 3  | FR-2400  | 10-23-12 | AJC | KDG | DRP  | ADD HS SEP. 21406 & OIL RETURN 2502 |     |       |      |
| 2  | FR-2400  | 08-20-12 | AJC | KDG | DRP  | HS COMP. 203056 TO 203056 - REL.8   |     |       |      |
| 1  | FR-2267  | 10-18-11 | LVE | SAG | DRP  | COMPRESSOR CHANGE/RELEASE 7         |     |       |      |
| 0  | FR-2202  | 05-12-11 | SRS | SAG | LMN  | RELEASED FOR PRODUCTION/RELEASE 6   |     |       |      |
| REV  | FR       | DATE     | BY  | CAD | APPD | DESCRIPTION OF REVISION             |     |       |      |
| DATE   | 05-12-11 | DVA      | SAG | CAD | SAG  | APPD                                | LMN | SCALE | NONE |
| CUSTOMER   |          |          |     |     |      |                                     |     |       |      |
| JOB TITLE -86°C 17 CU FT. UPRIGHT FREEZERS (220 V) |          |          |     |     |      |                                     |     |       |      |
| DVG TITLE REFRIGERATION SCHEMATIC                  |          |          |     |     |      |                                     |     |       |      |
| DRAWING NUMBER 904-90-0-B                          |          |          |     |     |      |                                     |     |       |      |
| LOCATION   |          |          |     |     |      |                                     |     |       |      |

1. COMMON TO: 904, 931, 993, 5704, 5604, 8923, 8934, 8956, 8968  
5904, 5993, 6423, 6434, 5931, CAR22086A 17 CU. FT. FREEZERS (220 V)

**ThermoFisher**  
**SCIENTIFIC**  
BOX 649, WAREHITA, OHIO 43750

HIGH TEMP STAGE



LOW TEMP STAGE

CASCADE REFRIGERATION

HIGH TEMPERATURE STAGE REFRIGERANT:

17 CUFT UNITS: R-404A 265 OZ. (75g) 1/2 OZ. (14g)

HIGH TEMPERATURE STAGE OIL: MOBILE EAL ARCTIC 220C POLYOL ESTER OIL  
COMPRESSOR: 22 OZ. (651ml)  
OIL SEPARATOR: 15 OZ. (444 ml)

LOW TEMPERATURE STAGE REFRIGERANT:

17 CUFT 50 HZ UNITS: R-290 11 OZ. (312g) MAX. 1 VAC TO 6 PSIG  
R-508B 13.7 OZ. (3884g) ± 1/4 OZ.(7g) DR  
6 PSIG TO 123 PSIG ± 3 PSIG

17 CUFT 60 HZ UNITS: R-290 11 OZ. (312g) MAX. 1 VAC TO 7 PSIG  
R-508B 14.5 OZ. (411g) ± 1/4 OZ.(7g) DR  
7 PSIG TO 128 PSIG ± 3 PSIG

LOW TEMPERATURE STAGE OIL: ZEROL 150T WITH 2% ADDITIVE  
COMPRESSOR: 22 OZ.(651ml)  
OIL SEPARATOR: 15 OZ. (444ml)

\* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C

THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC

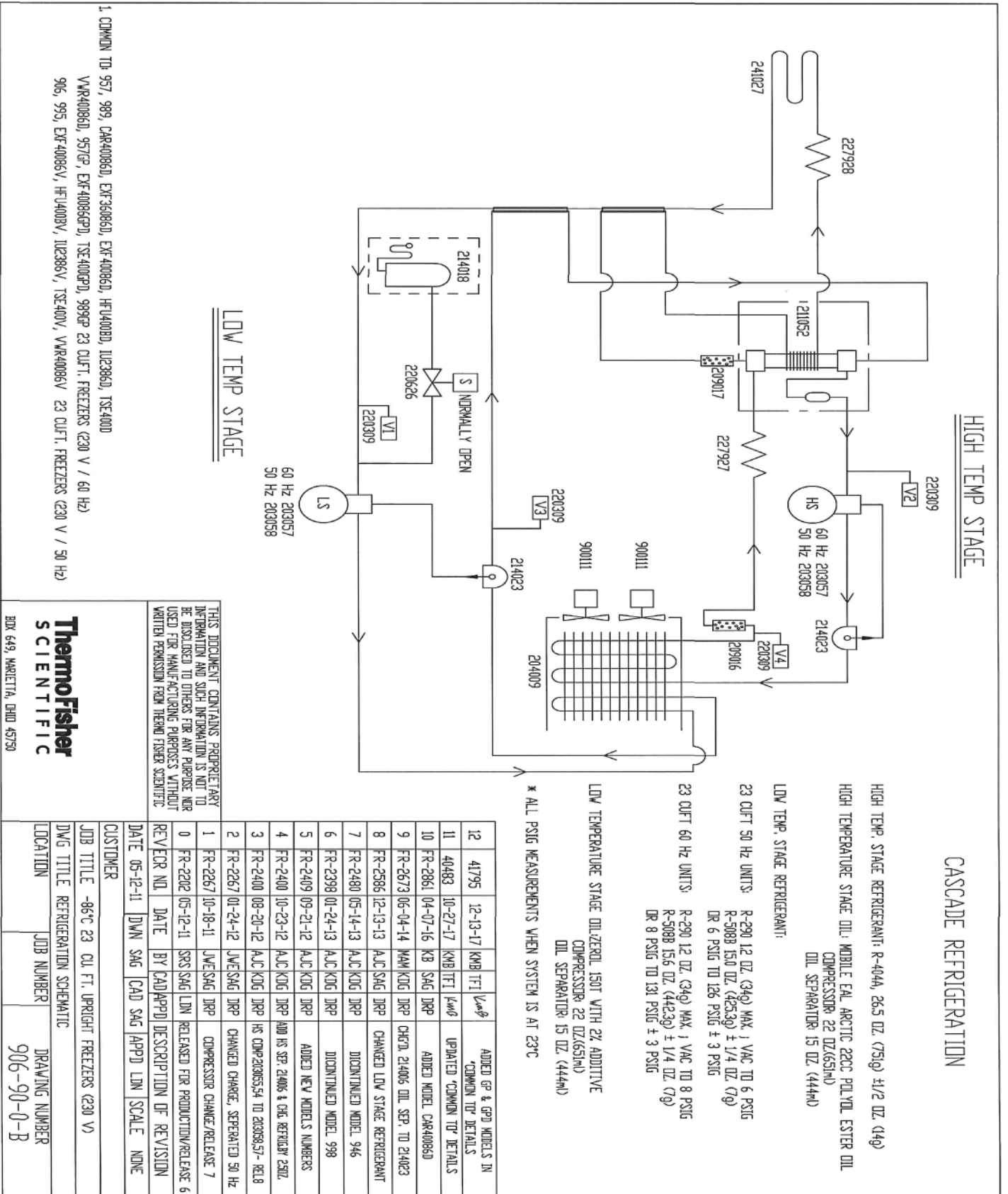
1. COMMON TO: 905, 994, EXF32086V, HFU3208V, IUL786V, TSE320V, VWR32086V 17 CU FT. FREEZERS (230 V / 50 HZ)

910, 988, CAR32086D, EXF28886D, EXF32086D, HFU3208D, IUL786D, TSE320D

VWR32086D, 910GP, EXF32086GPD, TSE320GPD, 988GP 17 CU FT. FREEZERS (230 V / 60 HZ)

**ThermoFisher**  
SCIENTIFIC  
BOX 649, MARLETTA, OHIO 45750

| DATE     | BY (APPD) | DESCRIPTION OF REVISION               |
|----------|-----------|---------------------------------------|
| 05-12-11 | DWN SAG   | CHD SAG (APPD) LHM SCALE NONE         |
| 05-12-11 | SRS SAG   | RELEASED FOR PRODUCTION/RELEASE 6     |
| 10-18-11 | JWE SAG   | COMPRESSOR CHANGE/RELEASE 7           |
| 10-23-12 | JWE SAG   | CHANGED CHANGE, SEPARATED 50 HZ       |
| 08-20-12 | AJC KDG   | HS COMP-203055-54 TO 203058-57- REL 8 |
| 10-23-12 | AJC KDG   | ADD HS SEP 24006 & CHG. REFRI. 250Z   |
| 10-23-12 | AJC KDG   | ADDED NEW MODEL NUMBERS               |
| 01-24-13 | AJC KDG   | DISCONTINUED MODEL 997                |
| 05-14-13 | AJC KDG   | DISCONTINUED MODEL 945                |
| 12-13-13 | AJC SAG   | CHANGED LOW STAGE REFRIGERANT         |
| 12-13-13 | AJC SAG   | CHANGED LOW STAGE REFRIGERANT         |
| 06-04-14 | MAN KDG   | CHGTL 214006 OIL SEP. TO 214023       |
| 04-07-16 | KB SAG    | ADDED MODEL CAR32086D                 |
| 10-27-17 | KMB TFI   | UPDATED COMMON TO* DETAILS            |
| 12-13-17 | KMB TFI   | COMMON TO* DETAILS                    |
| 4/7/95   |           | ADDED GP & GPD MODELS IN              |



1. COMMON TO: 957, 989, CAR40086D, EXF36086D, EXF40086D, HFTU4008D, ILC2386D, TSE400D  
 VWR40086D, 957GP, EXF40086GPD, TSE400GPD, 989GP 23 CU.FT. FREEZERS (230 V / 60 HZ)  
 906, 995, EXF40086V, HFTU4008V, ILC2386V, TSE400V, VWR40086V 23 CU.FT. FREEZERS (230 V / 50 HZ)

**LOW TEMP STAGE**

60 Hz 203057  
50 Hz 203058

**HIGH TEMP STAGE**

**CASCADE REFRIGERATION**

HIGH TEMP STAGE REFRIGERANT: R-404A, 26.5 OZ. (751g) 1/2 OZ. (14g)  
 HIGH TEMPERATURE STAGE OIL: MOBILE EAL ARCTIC 22CC POLYOL ESTER OIL  
 COMPRESSOR: 22 OZ.(651ml)  
 OIL SEPARATOR: 15 OZ. (444ml)

LOW TEMP STAGE REFRIGERANT:  
 R-290 12 OZ. (340g) MAX. 1 VAC TO 6 PSIG  
 R-5088 15.0 OZ. (425.3g) ± 1/4 OZ. (7g)  
 OR 6 PSIG TO 126 PSIG ± 3 PSIG

23 CU.FT. 50 HZ UNITS: R-290 12 OZ. (340g) MAX. 1 VAC TO 8 PSIG  
 R-5088 15.6 OZ. (442.3g) ± 1/4 OZ. (7g)  
 OR 8 PSIG TO 131 PSIG ± 3 PSIG

LOW TEMPERATURE STAGE OIL: ZEROL 150T WITH 2% ADDITIVE  
 COMPRESSOR: 22 OZ.(651ml)  
 OIL SEPARATOR: 15 OZ. (444ml)

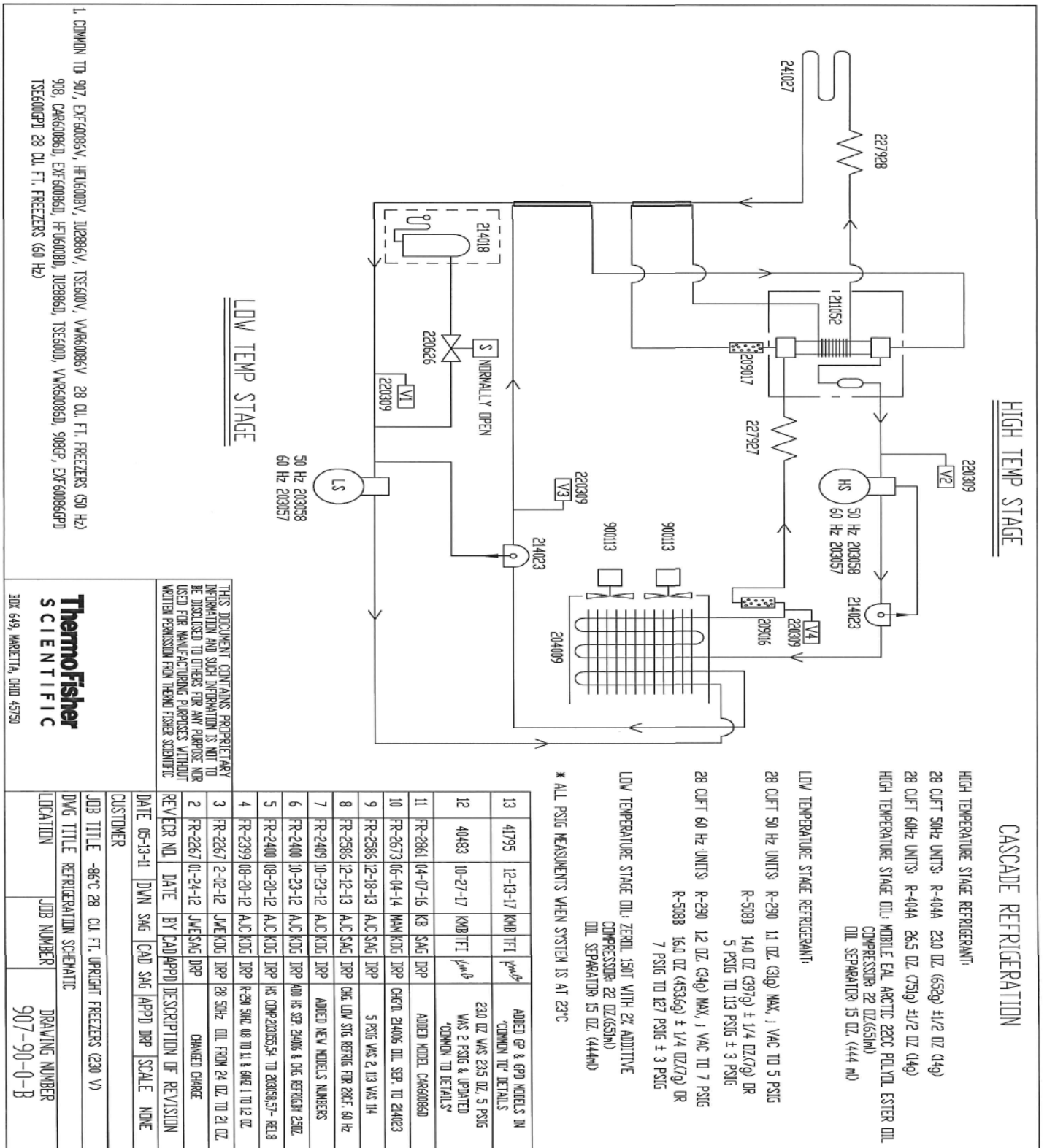
\* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C

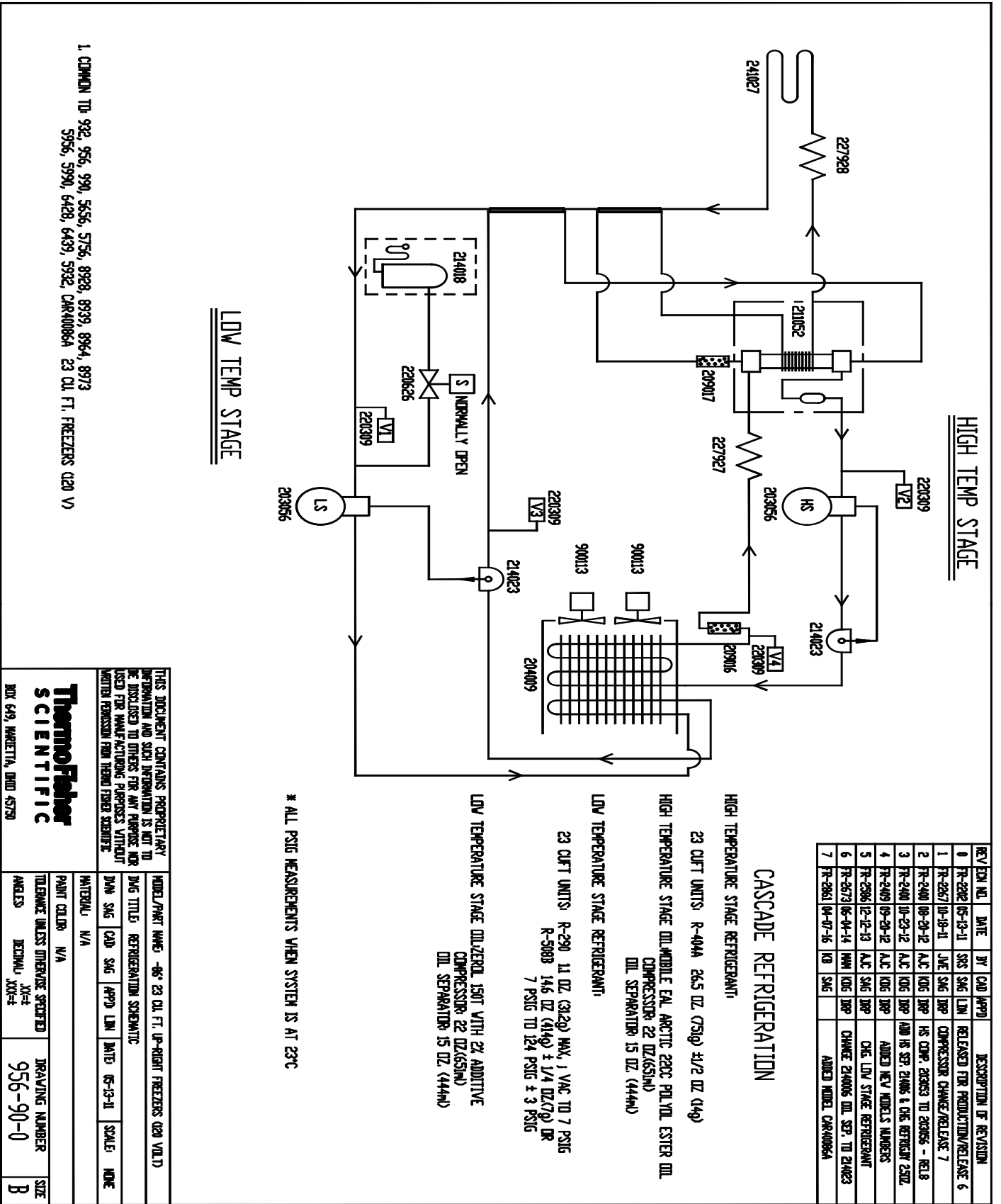
THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC



| NO. | DATE    | BY       | DESCRIPTION   |
|-----|---------|----------|---|
| 12  | 4/7/95  | 12-13-17 | KMB TFI Vamp ADDED GP & GPD MODELS IN COMMON TO DETAILS |
| 11  | 4/4/83  | 10-27-17 | KMB TFI Vamp UPDATED COMMON TO DETAILS                  |
| 10  | FR-2861 | 04-07-16 | KB SAG DRP ADDED MODEL CAR40086D                        |
| 9   | FR-2673 | 06-04-14 | MAM KDG DRP CHG'D 214006 OIL SEP. TO 214023             |
| 8   | FR-2585 | 12-13-13 | AJC SAG DRP CHANGED LOW STAGE REFRIGERANT               |
| 7   | FR-2480 | 05-14-13 | AJC KDG DRP DISCONTINUED MODEL 946                      |
| 6   | FR-2398 | 01-24-13 | AJC KDG DRP DISCONTINUED MODEL 998                      |
| 5   | FR-2409 | 09-21-12 | AJC KDG DRP ADDED NEW MODEL NUMBERS                     |
| 4   | FR-2400 | 10-23-12 | AJC KDG DRP ADD HS SEP. 21406 & CHG. REFRIG'R 230Z      |
| 3   | FR-2400 | 08-20-12 | AJC KDG DRP HS COMP:203055,54 TO 203058,57- REL:8       |
| 2   | FR-2267 | 01-24-12 | JWS SAG DRP CHANGED CHARGE, SEPERATED 50 HZ             |
| 1   | FR-2267 | 10-18-11 | JWS SAG DRP COMPRESSOR CHANGE/RELEASE 7                 |
| 0   | FR-2202 | 05-12-11 | SRS SAG LHM RELEASED FOR PRODUCTION/RELEASE 6           |





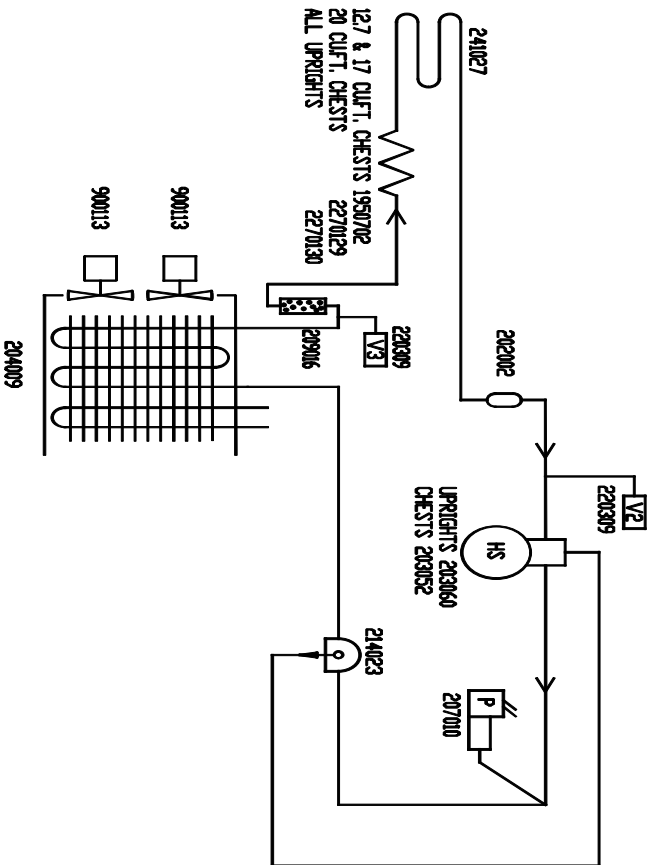


1. COMMON TO: 932, 956, 990, 5656, 5756, 8928, 8939, 8964, 8973  
 5956, 5990, 6428, 6439, 5932, CAR40086A 23 OIL FT. FREEZERS (20 V)

|  |  |
|--|--|
| <p>THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE AND USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM THERMO FISH SCIENTIFIC</p> |  |
| <p><b>ThermoFisher SCIENTIFIC</b><br/>                 BOX 649, WAUWATIGA, WI 53719</p>  | <p>MODEL/PART NAME: -66" 23 OIL FT. UP-RIGHT FREEZERS (20 VOL.)<br/>                 TAG TITLE: REFRIGERATION SCHEMATIC<br/>                 DWG SWS CAD SWS APPD LWN DATE: 05-13-11 SCALE: NONE<br/>                 MATERIAL: N/A<br/>                 PAINT COLOR: N/A<br/>                 TOLERANCE UNLESS OTHERWISE SPECIFIED: DECIMAL XX=± DRAWING NUMBER: 956-90-0 SIZE: B</p> |

| REV# | DATE            | BY  | CAUSE | DESCRIPTION OF REVISION              |
|------|-----------------|-----|-------|--------------------------------------|
| 0    | R-2302 05-13-11 | SBS | SWS   | RELEASED FOR PRODUCTION/RELEASE 6    |
| 1    | R-2367 09-08-11 | JVC | SWS   | COMPRESSOR CHANGE/RELEASE 7          |
| 2    | R-2400 08-29-12 | ALC | KDG   | HS COMP. 20265 TO 20266 - REL 9      |
| 3    | R-2400 08-29-12 | ALC | KDG   | ADD HS SEP. 21066 & OIL RETURN 2502. |
| 4    | R-2409 09-20-12 | ALC | KDG   | ADDED NEW MODEL'S NUMBERS            |
| 5    | R-2386 02-12-13 | ALC | SWS   | CHANGE 214006 OIL SEP. TO 214Q23     |
| 6    | R-2673 06-04-14 | MWH | KDG   | ADDED MODEL CAR40086A                |
| 7    | R-2861 04-07-16 | KD  | SWS   | ADDED MODEL CAR40086A                |

HIGH TEMP STAGE



REFRIGERATION

HIGH TEMPERATURE STAGE REFRIGERANT:

- 130 CUFT. UNITS R-404A 18.5 OZ. (525gr) 1/2 OZ (14gr)
- 170 CUFT. UNITS R-404A 22 OZ. (624gr) 1/2 OZ (14gr)
- 230 CUFT. UNITS R-404A 22 OZ. (624gr) 1/2 OZ (14gr)

HIGH TEMPERATURE STAGE OIL: ARCTIC 22CC POLYOL ESTER OIL  
COMPRESSOR, 40 OZ.(1183ml)

- 127 CUFT. UNITS R-404A 19 OZ. (538gr) 1/2 OZ (14gr)
- 170 CUFT. UNITS R-404A 21 OZ. (595gr) 1/2 OZ (14gr)
- 200 CUFT. UNITS R-404A 23 OZ. (652gr) 1/2 OZ (14gr)

HIGH TEMPERATURE STAGE OIL: ARCTIC 22CC POLYOL ESTER OIL  
COMPRESSOR, 40 OZ.(1183ml)  
OIL SEPARATOR, 15 OZ.(444ml)

THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC.

|    |         |          |         |     |                                   |
|----|---------|----------|---------|-----|-----------------------------------|
| 10 | FR-2761 | 01-14-15 | AJC SAG |     | CHG. 13 UPRIGHT 21 TO 18.5 CHANGE |
| 9  | FR-2761 | 01-07-15 | AJC SAG | DRP | CHANGED UPRIGHT CHARGES           |
| 8  | FR-2694 | 07-10-14 | AJC KOG | DRP | SPECIFIED OIL SEP. 15 OZ. OF OIL  |
| 7  | FR-2673 | 06-04-14 | MM/KOG  | DRP | CHG. 24006 OIL SEP. TO 214023     |
| 6  | FR-2499 | 07-30-13 | AJC SAG | DRP | OIL COMP. FROM 20365 TO 20360     |
| 5  | FR-2400 | 08-20-12 | AJC KOG | DRP | OIL COMP.20363 TO 20364-REL.8     |
| 4  | FR-2267 | 2/03/12  | JVE/KOG | DRP | UPRIGHT OIL FROM 24 OZ. TO 22 OZ. |
| 3  | FR-2267 | 01/16/12 | JVE/SAG | DRP | CHST OIL FROM 24 OZ. TO 40 OZ.    |
| 2  | FR-2267 | 11/07/11 | JVE/SAG | DRP | COMPRESSOR CHANGE/RELEASE 7       |
| 1  | FR-2202 | 08/02/11 | AJC SAG | DRP | ADDED CAP TUBE 2270120            |
| 0  | FR-2202 | 05/09/11 | SPS SAG | DRP | RELEASED FOR PRODUCTION/RELEASE 6 |

REVISOR NO. DATE BY (CAD/APPD) DESCRIPTION OF REVISION  
DATE (05/09/11) DWN SAG CAD SAG APPD DRP SCALE NONE

CUSTOMER  
JOB TITLE -40C ULTRA-LOW TEMPERATURE FREEZERS, 120V

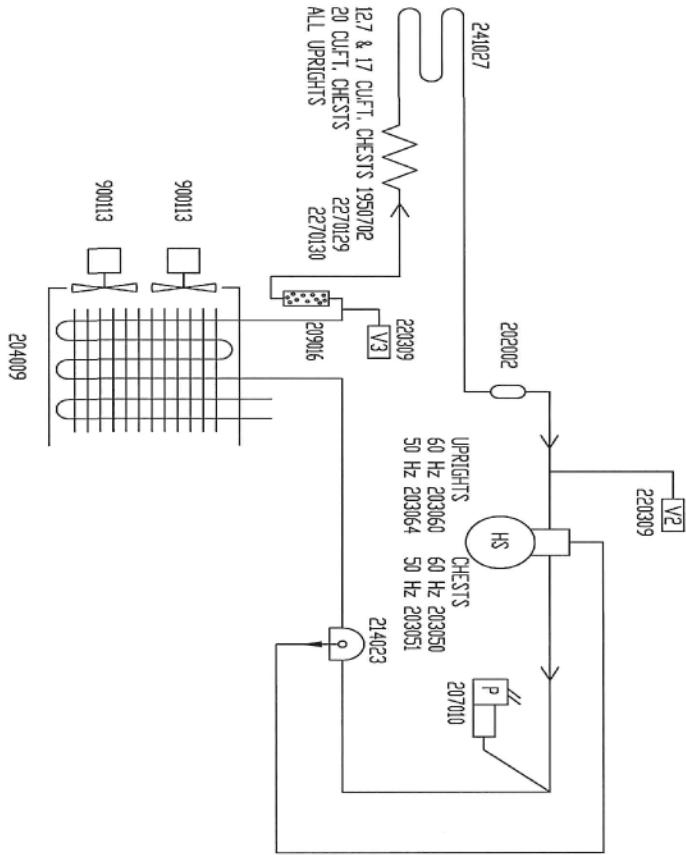
DWG TITLE REFRIGERATION SCHEMATIC

LOCATION JOB NUMBER DRAWING NUMBER  
FREEZERS 8800-90-0-B

1. COMMON TO: 5722, 8800, 8813, 8824, 8838 130 CUFT. UPRIGHT FREEZERS (120 V),  
5728, 8804, 8817, 8828, 8842 170 CUFT. UPRIGHT FREEZERS (120 V),  
8809, 8821, 8832, 8848 23 CUFT. UPRIGHT FREEZERS (120 V),  
5315, 8728, 8756, 8779 127 CUFT CHEST FREEZER (120V),  
5318, 8730, 8757, 8780 170 CUFT CHEST FREEZER (120V),  
5321, 8737, 8762, 8785 200 CUFT CHEST FREEZER (120V)

**ThermoFisher**  
**SCIENTIFIC**  
BOX 649, WAREHO, OHIO 45739

**HIGH TEMP STAGE**



1. COMMON TO
- 7240, DFC2440, FFC2440, TSM2440, VWC2440, Z240V, DFC2440V, FFC2440V, TSM2440V, VWC2440V, Z240V, DFC2440P, FFC2440P, TSM2440P, VWC2440P, Z240P, UPRIGHT FREEZERS (230 V)
  - 7230, DFC2340, FFC2340, TSM2340, VWC2340, Z230V, DFC2340V, FFC2340V, TSM2340V, VWC2340V, Z230V, DFC2340P, FFC2340P, TSM2340P, VWC2340P, Z230P, UPRIGHT FREEZERS (230 V)
  - 7400, DFC4040, FFC4040, TSM4040, VWC4040, Z400V, DFC4040V, FFC4040V, TSM4040V, VWC4040V, Z400V, UPRIGHT FREEZERS (230 V)
  - 7130, FFC1300, TSC1300, UL130-H-1, 713V, FFC1300V, TSC1300V, UL130-H-V, 7130P, TSC1300P, UL130-H-P, 127 CUFT CHEST FREEZER (230V)
  - 7170, FFC1700, TSC1700, UL170-H-1, 717V, FFC1700V, TSC1700V, UL170-H-V, 7170P, TSC1700P, UL170-H-P, 170 CUFT CHEST FREEZER (230V)
  - 7200, FFC2000, TSC2000, UL200-H-1, 720V, FFC2000V, TSC2000V, UL200-H-V, 7200P, TSC2000P, UL200-H-P, 200 CUFT CHEST FREEZER (230V)

**REFRIGERATION**

HIGH TEMPERATURE STAGE REFRIGERANT

HIGH TEMPERATURE STAGE REFRIGERANT

HIGH TEMPERATURE STAGE OIL/MOBILE CAL ARCTIC 220C PUL VOL. ESTER OIL

COMPRESSOR: 40 OZ(1183M)

OIL SEPARATOR: 15 OZ(444 M)

CHESTS

- 127 CUFT. UNITS: R-404A 19 OZ. (538gr) 1/2 OZ (14gr)
- 170 CUFT. UNITS: R-404A 21 OZ. (595gr) 1/2 OZ (14gr)
- 200 CUFT. UNITS: R-404A 23 OZ. (652gr) 1/2 OZ (14gr)

HIGH TEMPERATURE STAGE OIL/MOBILE CAL ARCTIC 220C PUL VOL. ESTER OIL

COMPRESSOR: 40 OZ(1183M)

OIL SEPARATOR: 15 OZ(444M)

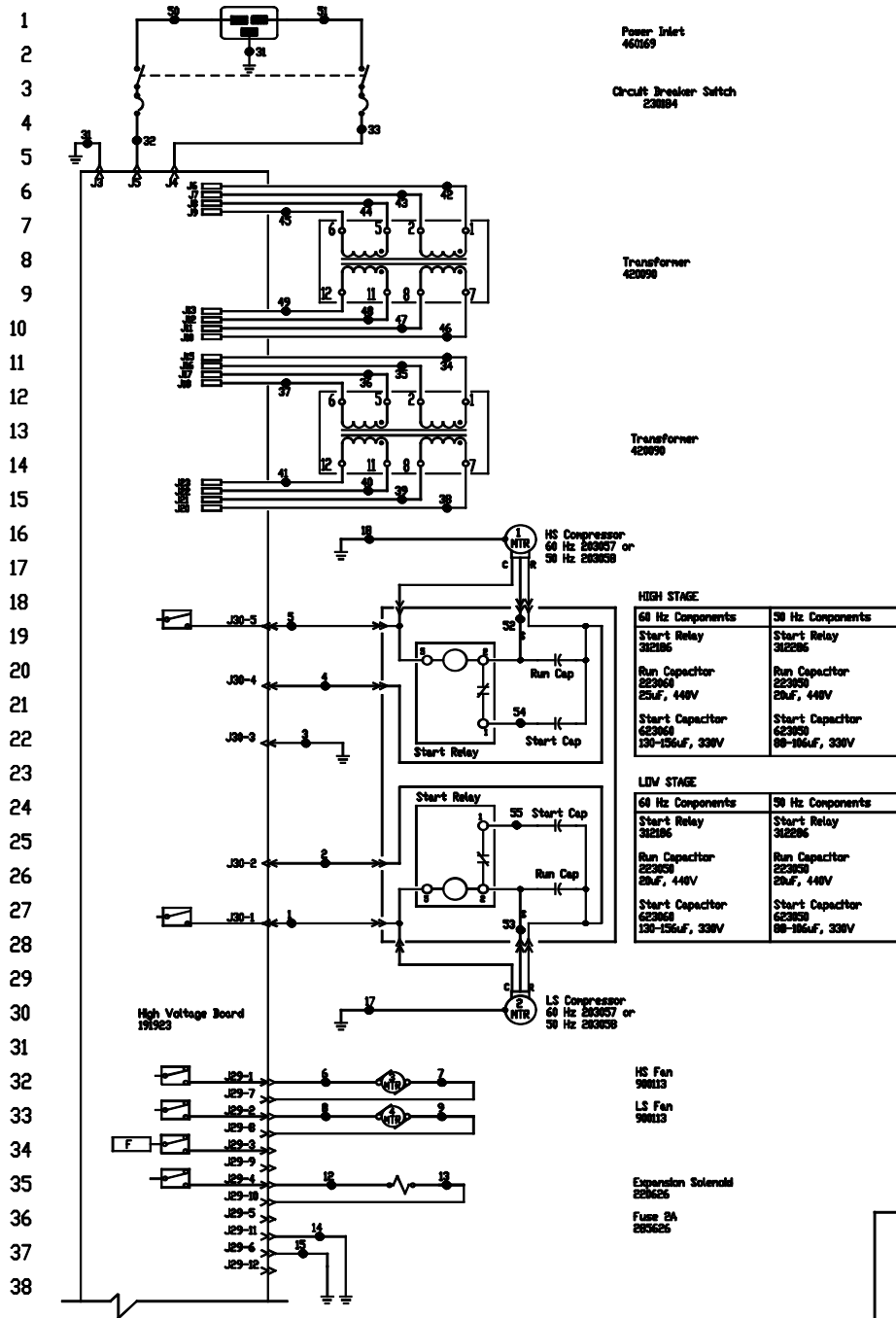
THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC

**ThermoFisher**  
**SCIENTIFIC**  
BOX 649, MALETTA, OHIO 45750

|    |            |   |                |          |  |
|----|------------|---|----------------|----------|--|
| 11 | 41795      | 12-13-17                                    | KWB TFI        | YMM      | ADDED P/N MODELS IN COMMON TO DETAILS                  |
| 10 | 40483      | 10-27-17                                    | KWB TFI        | YMM      | ADDED OIL SEPARATOR NOTE AND UPDATED COMMON TO DETAILS |
| 9  | FR-2761    | 01-07-15                                    | AJC SAG        | DRP      | CHANGED CHARGE ADDED 513 & 5726                        |
| 8  | FR-2694    | 07-10-14                                    | AJC KDG        | DRP      | SPECIFIED OIL SEP. 15 OZ. OF OIL                       |
| 7  | FR-2673    | 06-04-14                                    | MM/KDG         | DRP      | CHG'D. 214005 OIL SEP. TO 214023                       |
| 6  | FR-2499    | 07-30-13                                    | AJC SAG        | DRP      | CHANGED OIL COMPRESSORS & CHARGE                       |
| 5  | FR-2400    | 08-20-12                                    | AJC KDG        | DRP      | OR COMP.203064,55 TO 20307,50-R-6L,8                   |
| 4  | FR-2267    | 2/02/12                                     | JWE/KDG        | DRP      | UPRIGHT OIL FROM 24 OZ. TO 22 OZ.                      |
| 3  | FR-2267    | 01/16/12                                    | JWE/SAG        | DRP      | CHEST OIL FROM 24 OZ. TO 40 OZ.                        |
| 2  | FR-2267    | 11/07/11                                    | JWE/SAG        | DRP      | COMPRESSOR CHANGE/RELEASE 7                            |
| 1  | FR-2202    | 08/02/11                                    | AJC SAG        | DRP      | ADDED CAP TUBE 220020                                  |
| 0  | FR-2202    | 05/09/11                                    | SRS SAG        | LNK      | RELEASED FOR PRODUCTION/RELEASE 6                      |
|    | REVEER NLD | DATE  | BY             | CAD/APP/ | DESCRIPTION OF REVISION                                |
|    | DATE       | 05/09/11                                    | DWN SAG        | CAD SAG  | APPD LBN SCALE NONE                                    |
|    | CUSTOMER   |   |                |          |  |
|    | JOB TITLE  | -40° C ULTRA-LOW TEMPERATURE FREEZERS, 230V |                |          |  |
|    | DWG TITLE  | REFRIGERATION SCHEMATIC                     |                |          |  |
|    | LOCATION   | JOB NUMBER                                  | DRAWING NUMBER |          |  |
|    | FREEZERS   |   | 8801-90-0-B    |          |  |

**Section 10**  
Electrical Schematics

**POWER CONNECTION**  
230V/50HZ & 208/230V/60HZ, 16, 24, 32.0 FLA  
OPERATING RANGE 237-253



Power Inlet  
46069

Circuit Breaker Switch  
23894

Transformer  
42099

Transformer  
42099

1  
NTR

HS Compressor  
60 Hz 293957 or  
50 Hz 293958

HIGH STAGE

| 60 Hz Components                            | 50 Hz Components                          |
|---|---|
| Start Relay<br>32216                        | Start Relay<br>312206                     |
| Run Capacitor<br>22369<br>25uF, 440V        | Run Capacitor<br>22369<br>25uF, 440V      |
| Start Capacitor<br>62369<br>120-156uF, 330V | Start Capacitor<br>62369<br>60-96uF, 330V |

LOW STAGE

| 60 Hz Components                            | 50 Hz Components                          |
|---|---|
| Start Relay<br>32216                        | Start Relay<br>312206                     |
| Run Capacitor<br>22369<br>25uF, 440V        | Run Capacitor<br>22369<br>25uF, 440V      |
| Start Capacitor<br>62369<br>120-156uF, 330V | Start Capacitor<br>62369<br>60-96uF, 330V |

High Voltage Board  
19193

HS Fan  
96013

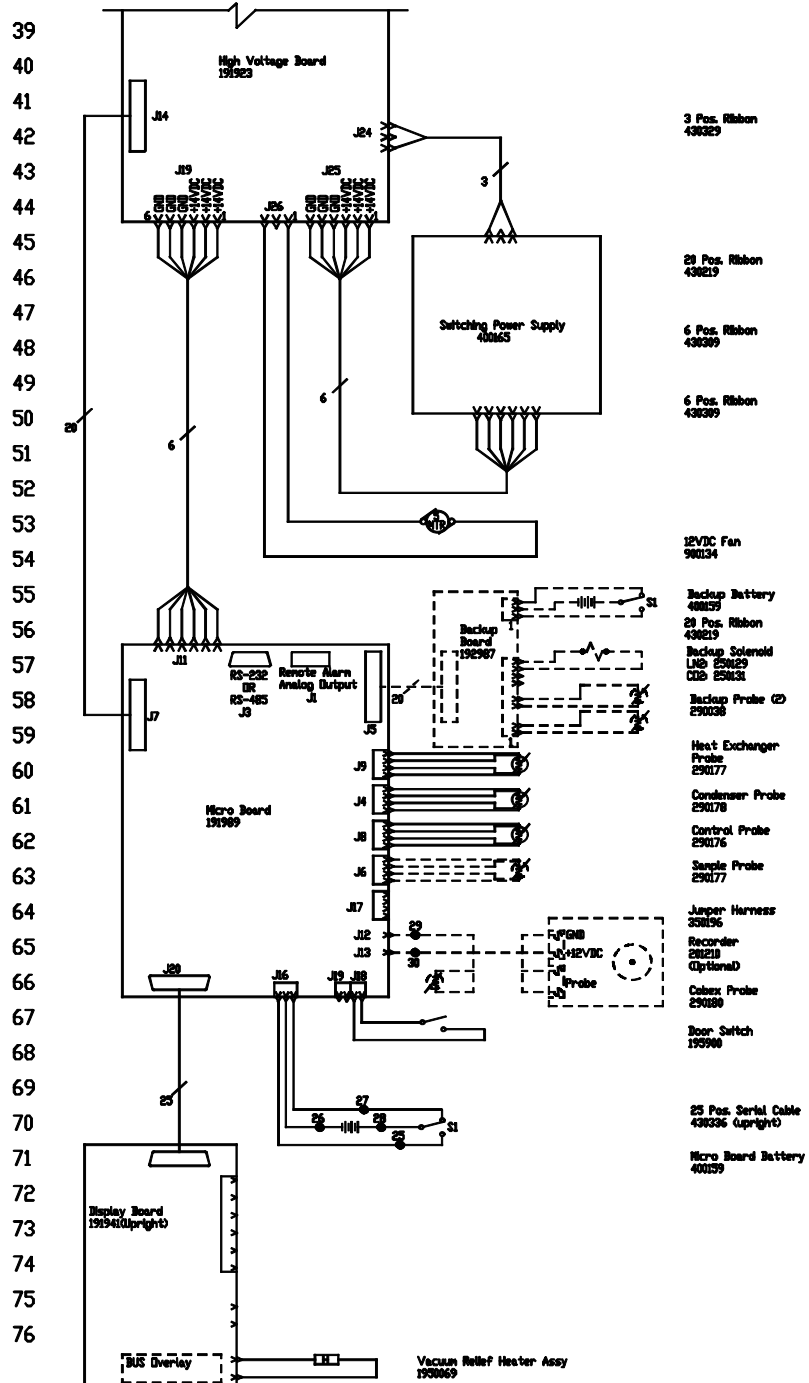
LS Fan  
96013

Expansion Solenoid  
23626

Fuse 2A  
24

Electrical Schematic  
Model:  
230V, 50 & 60Hz, -86°C  
ULT Upright Freezers

5602-70-2-0 REV 3  
Page 1 of 3



Electrical Schematic  
Model:  
230V, 50 & 60Hz, -86°C  
ULT Upright Freezers

77  
78  
79  
80  
81  
82  
83  
84  
85  
86  
87  
88  
89  
90  
91  
92  
93  
94  
95  
96  
97  
98  
99  
100  
101  
102  
103  
104  
105  
106  
107

| WIRE # | COLOR   | GAUGE | WIRE # | COLOR   | GAUGE |
|--------|---------|-------|--------|---------|-------|
| 1      | BLK     | 14    | 29     | BLK     | 22    |
| 2      | RED     | 14    | 30     | RED     | 22    |
| 3      | GRN/YEL | 14    | 31     | GRN/YEL | 14    |
| 4      | RED     | 14    | 32     | BLK     | 14    |
| 5      | BLK     | 14    | 33     | BLU     | 14    |
| 6      | BLK     | 18    | 34     | BLK     | 14    |
| 7      | BLK     | 18    | 35     | BLU     | 14    |
| 8      | BLK     | 18    | 36     | BLK     | 14    |
| 9      | BLK     | 18    | 37     | BLU     | 14    |
| 10     | -       | -     | 38     | BLK     | 14    |
| 11     | -       | -     | 39     | BLU     | 14    |
| 12     | BLK     | 18    | 40     | BLK     | 14    |
| 13     | BLK     | 18    | 41     | BLU     | 14    |
| 14     | GRN/YEL | 18    | 42     | BLK     | 14    |
| 15     | GRN/YEL | 18    | 43     | BLU     | 14    |
| 16     | -       | -     | 44     | BLK     | 14    |
| 17     | GRN/YEL | 14    | 45     | BLU     | 14    |
| 18     | GRN/YEL | 14    | 46     | BLK     | 14    |
| 19     | -       | -     | 47     | BLU     | 14    |
| 20     | -       | -     | 48     | BLK     | 14    |
| 21     | -       | -     | 49     | BLU     | 14    |
| 22     | -       | -     | 50     | BLK     | 14    |
| 23     | -       | -     | 51     | BLU     | 14    |
| 24     | -       | -     | 52     | YEL/WHT | 14    |
| 25     | WHT     | 18    | 53     | YEL     | 14    |
| 26     | BLK     | 18    | 54     | BLU/WHT | 14    |
| 27     | BLK     | 18    | 55     | BLU     | 14    |
| 28     | RED     | 18    |        |         |       |

UPRIGHT MODELS  
 13 CUFT. (230V) - VVR240863, VVR24086V  
 17 CUFT. (230V) - VVR220863, VVR22086V  
 23 CUFT. (230V) - VVR40863, VVR40863S3, VVR4086V  
 28 CUFT. (230V) - VVR60863, VVR6086V

RS-232 SPECIFICATION

BAUD = 1200  
 PARITY = N  
 BITS = 8  
 STOP BITS = 2

DS-232 POINT  
 PIN 2 TXD  
 PIN 3 RXD  
 PIN 5 GND

REMOTE CONTACTS/ANALOG OUTPUT  
 PIN 1 Analog Output 4  
 PIN 2 Analog Output -  
 PIN 3 Not Connected  
 PIN 4 Not Connected  
 PIN 5 Normally Closed  
 PIN 6 Common  
 PIN 7 Normally Open

CONTACT RATING 1A @ 30V  
 CONTACTS IN ALARM STATE

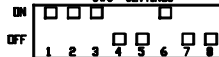
SV1 Settings for Analog Output



SV2 Settings for Communication Output



SV3 SETTINGS



- Expansion solenoid on 17, 23 & 28 cuft models only.
- Door switches shown in open position.
- Battery switch shown in the OFF position.
- Circuit breaker switch shown in OFF position.
- Options and accessories shown in dashed lines.

Electrical Schematic  
 Model:  
 230V, 50 & 60Hz, -86°C  
 ULT Upright Freezers

| REV | REV NO. | DATE     | BY  | CHK | APP'D | DESCRIPTION OF REVISION                |
|-----|---------|----------|-----|-----|-------|--|
| 3   | 42795   | 12-10-07 | KOB | WFL |       | UPDATED "UPRIGHT MODELS" DETAILS       |
| 2   | FR-0085 | 02-14-07 | KD  | KOB | KOB   | ISSUED "INCOMP" NUMBER TO FIELD        |
| 1   | FR-0250 | 02-07-06 | KOB | SAC | IMP   | DESIGN WAS CORRECT, DESIGN WAS CORRECT |
| 0   | FR-0250 | 02-28-04 | HW  | KOB | IMP   | RELEASED FOR PRODUCTION / RELEASE IS   |

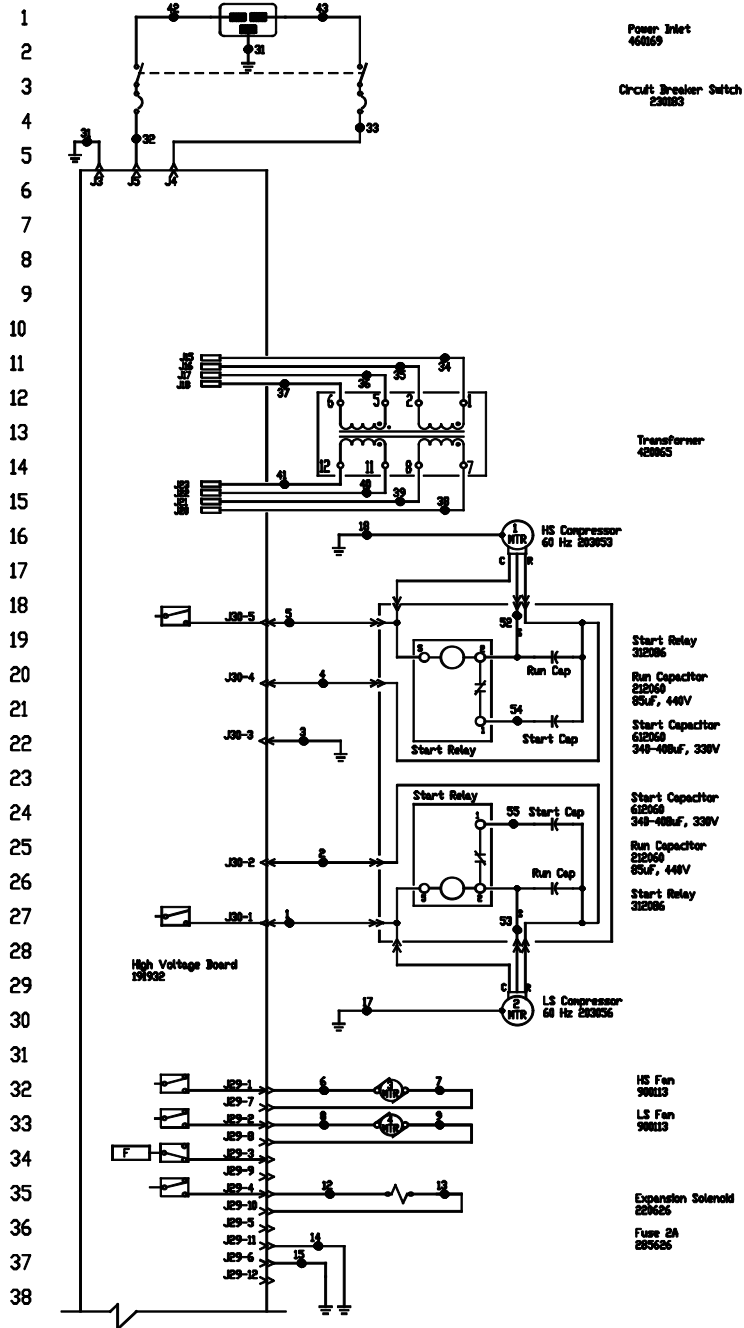


THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC

MODEL/PART NAME ULT -86C UPRIGHT FREEZERS, 230V, 50 & 60Hz  
 DWG TITLE UNIT SCHEMATIC  
 DRAWN BY CAD KOB APP'D DATED 12-10-04 SCALE NONE  
 MATERIAL:  
 PAINT COLOR:  
 TOLERANCE UNLESS OTHERWISE SPECIFIED  
 DIMENSIONS UNLESS OTHERWISE SPECIFIED  
 ANGLES DECIMALS 320-6 320-6  
 DRAWING NUMBER 5602-70-1  
 SIZE D

5602-70-1-0 REV 3  
 Page 3 of 3

POWER CONNECTION  
120V, 1Ø, 2W, 60Hz, 15A FLA  
OPERATING RANGE 100-130V

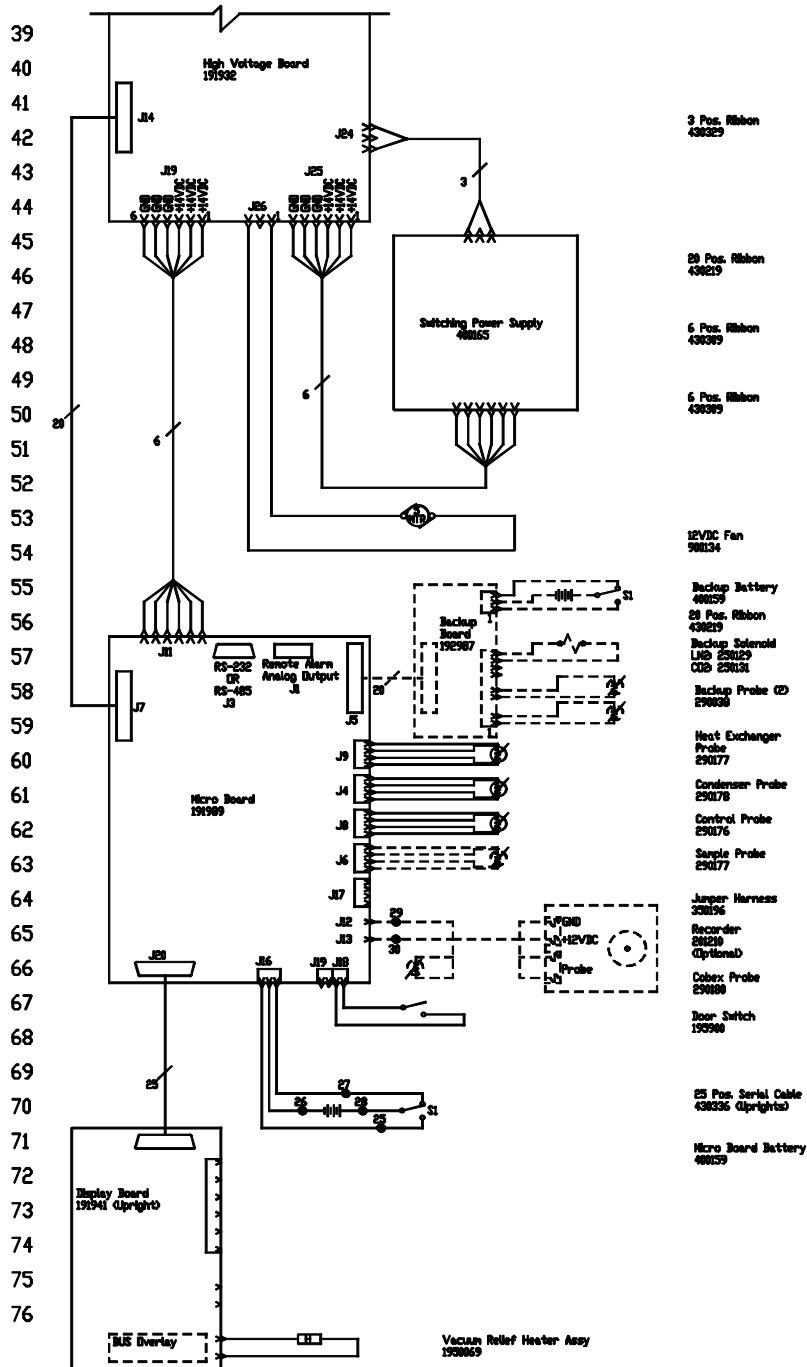


Electrical Schematic  
Model:  
120V, 60Hz, -86°C  
ULT Upright Freezers

5603-70-1-0 REV 3  
Page 1 of 3



**Section 10**  
Electrical Schematics



Electrical Schematic  
Model:  
120V, 60Hz, -86°C  
ULT Upright Freezers

5603-70-1-0 REV 3  
Page 2 of 3

77  
78  
79  
80  
81  
82  
83  
84  
85  
86  
87  
88  
89  
90  
91  
92  
93  
94  
95  
96  
97  
98  
99  
100  
101  
102  
103  
104  
105  
106  
107

| WIRE # | COLOR   | GAUGE | WIRE # | COLOR   | GAUGE |
|--------|---------|-------|--------|---------|-------|
| 1      | BLK     | 14    | 29     | BLK     | 22    |
| 2      | RED     | 14    | 30     | RED     | 14    |
| 3      | GRN/YEL | 14    | 31     | GRN/YEL | 14    |
| 4      | RED     | 14    | 32     | BLK     | 14    |
| 5      | BLK     | 14    | 33     | BLU     | 14    |
| 6      | BLK     | 18    | 34     | BLK     | 14    |
| 7      | BLK     | 18    | 35     | BLU     | 14    |
| 8      | BLK     | 18    | 36     | BLK     | 14    |
| 9      | BLK     | 18    | 37     | BLU     | 14    |
| 10     | -       | -     | 38     | BLK     | 14    |
| 11     | -       | -     | 39     | BLU     | 14    |
| 12     | BLK     | 18    | 40     | BLK     | 14    |
| 13     | BLK     | 18    | 41     | BLU     | 14    |
| 14     | GRN/YEL | 18    | 42     | BLK     | 14    |
| 15     | GRN/YEL | 18    | 43     | BLU     | 14    |
| 16     | -       | -     | 44     | -       | -     |
| 17     | GRN/YEL | 14    | 45     | -       | -     |
| 18     | GRN/YEL | 14    | 46     | -       | -     |
| 19     | -       | -     | 47     | -       | -     |
| 20     | -       | -     | 48     | -       | -     |
| 21     | -       | -     | 49     | -       | -     |
| 22     | -       | -     | 50     | -       | -     |
| 23     | -       | -     | 51     | -       | -     |
| 24     | -       | -     | 52     | YEL/WHI | 14    |
| 25     | WHI     | 18    | 53     | YEL     | 14    |
| 26     | BLK     | 18    | 54     | BLU/WHI | 14    |
| 27     | BLK     | 18    | 55     | BLU     | 14    |
| 28     | RED     | 18    |        |         |       |

UPRIGHT MODELS:  
13 CU.FT. (25V) - VUR24086A  
17 CU.FT. (25V) - VUR32086A  
23 CU.FT. (25V) - VUR40086A

RS-232 SPECIFICATION

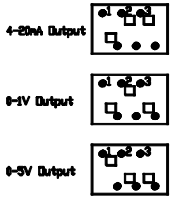
BAUD = 1200  
PARITY = N  
BITS = 8  
STOP BITS = 2

| RS-232 PIN |     |
|------------|-----|
| PIN 2      | TXD |
| PIN 3      | RXD |
| PIN 5      | GNB |

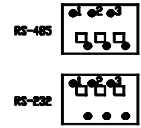
SENSE CONTACTS/ANALOG OUTPUT  
PIN 1 Analog Output +  
PIN 2 Analog Output -  
PIN 3 Not Connected  
PIN 4 Not Connected  
PIN 5 Normally Closed  
PIN 6 Common  
PIN 7 Normally Open

CONTACT RATING 1A @ 30V  
CONTACTS IN ALARM STATE

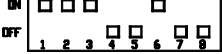
SV1 Settings for Analog Output



SV2 Settings for Communication Output



SV3 SETTINGS



- Expansion solenoid on 17 & 23 cuft uprights only.
- Door switches shown in open position.
- Battery switch shown in the OFF position.
- Circuit breaker switch shown in OFF position.
- Options and accessories shown in dashed lines.

| REV | NO.    | DATE     | BY  | APP'D | DESCRIPTION OF REVISION              |
|-----|--------|----------|-----|-------|--------------------------------------|
| 3   | 4275   | 10-10-17 | END | YFI   | UPDATED "UPRIGHT MODEL" DETAILS      |
| 2   | FR-666 | 10-10-17 | END | CVB   | 1200V BACKUP BEARD TO 120V/7         |
| 1   | FR-666 | 10-07-16 | END | SHS   | 250V BIP CIRCUIT WAS 250V/6          |
| 0   | FR-670 | 10-05-14 | END | END   | RELEASED FOR PRODUCTION / RELEASE 11 |

THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE RELEASED TO OTHERS FOR ANY PURPOSE, NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC



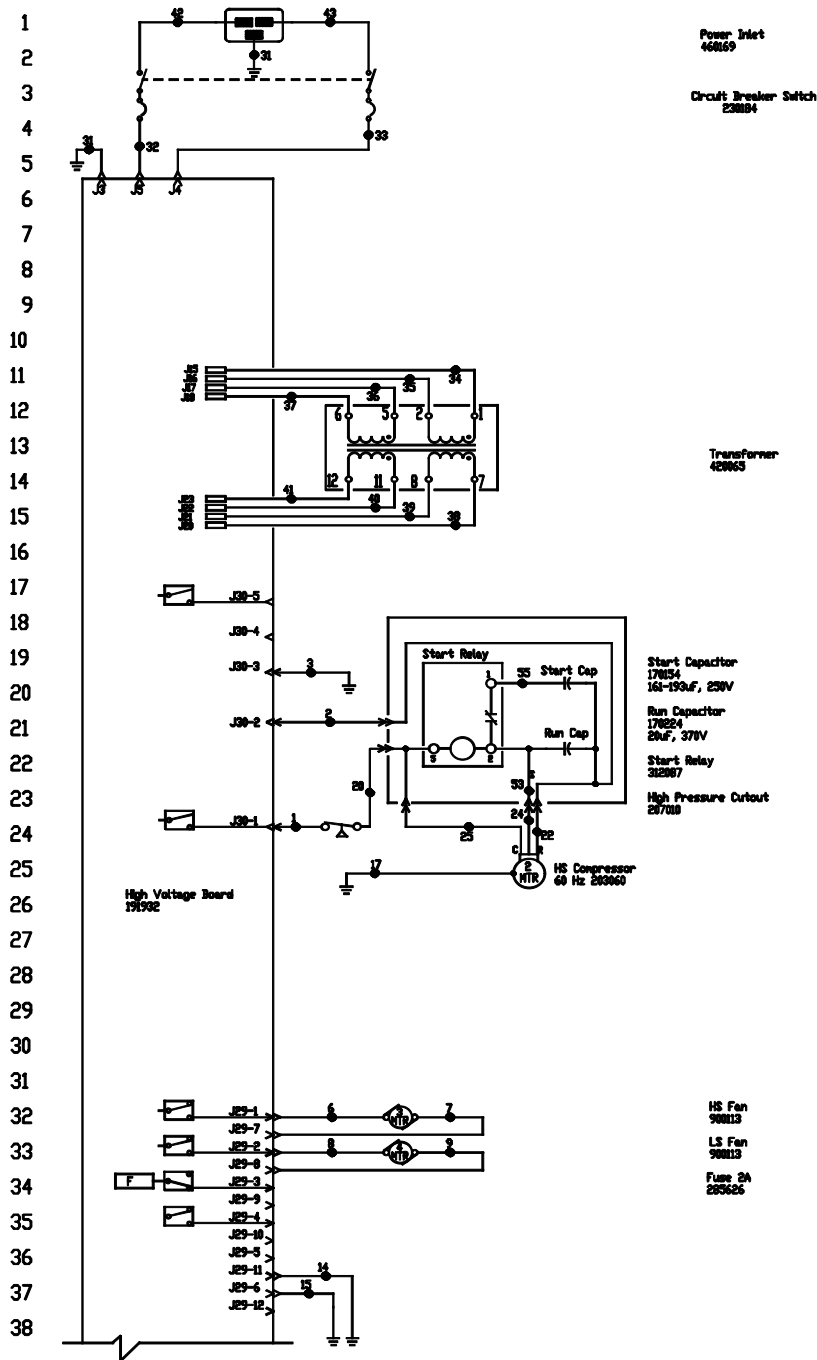
|                 |          |          |          |
|-----------------|----------|----------|----------|
| MODEL/PART NAME | UNIT     | DATE     | SCALE    |
| FR-666          | ULT      | 10-10-17 | 1:1      |
| PRINT COLOR     | DATE     | SCALE    |          |
| BLACK           | 10-10-17 | 1:1      |          |
| AWG/IN          | AWG/IN   | AWG/IN   | AWG/IN   |
| 14/0.075        | 18/0.090 | 22/0.054 | 24/0.047 |

Electrical Schematic  
Model:  
120V, 60Hz, -86°C  
ULT Upright Freezers

5603-70-1-0 REV 3  
Page 3 of 3

**Section 10**  
Electrical Schematics

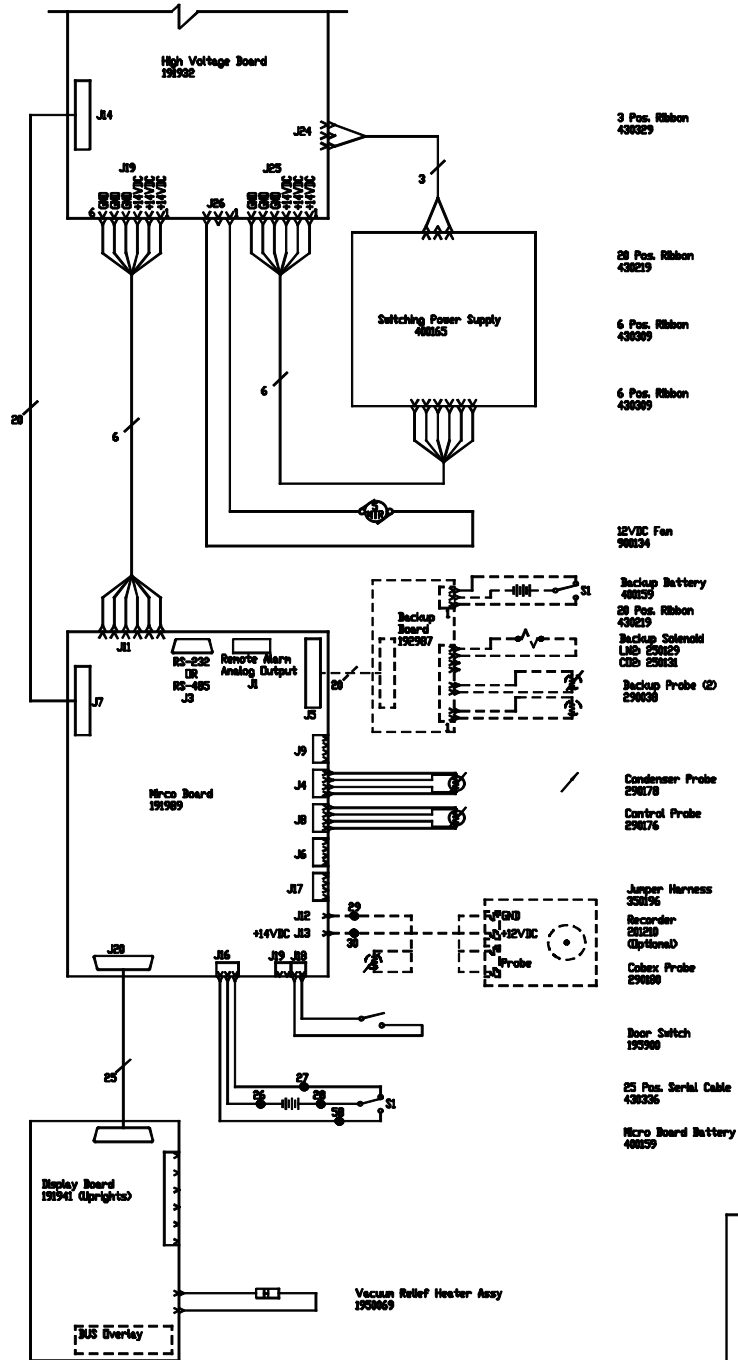
**POWER CONNECTION**  
120V, 16, 2V, 60HZ, 100 FLA  
OPERATING RANGE 100-132V



Electrical Schematic  
Model:  
120V, 60Hz, -40°C  
ULT Upright Freezers

5722-70-1-0 REV 4  
Page 1 of 3

39  
40  
41  
42  
43  
44  
45  
46  
47  
48  
49  
50  
51  
52  
53  
54  
55  
56  
57  
58  
59  
60  
61  
62  
63  
64  
65  
66  
67  
68  
69  
70  
71  
72  
73  
74  
75  
76



Electrical Schematic  
Model:  
120V, 60Hz, -40°C  
ULT Upright Freezers

5722-70-1-0 REV 4  
Page 2 of 3

| WIRE # | COLOR   | GAUGE | WIRE # | COLOR   | GAUGE |
|--------|---------|-------|--------|---------|-------|
| 1      | BLK     | 14    | 29     | BLK     | 22    |
| 2      | RED     | 14    | 30     | RED     | 14    |
| 3      | GRN/YEL | 14    | 31     | GRN/YEL | 14    |
| 4      |         |       | 32     | BLK     | 14    |
| 5      |         |       | 33     | BLU     | 14    |
| 6      | BLK     | 18    | 34     | BLU     | 14    |
| 7      | BLK     | 18    | 35     | BLU     | 14    |
| 8      | BLK     | 18    | 36     | BLK     | 14    |
| 9      | BLK     | 18    | 37     | BLU     | 14    |
| 10     |         |       | 38     | BLK     | 14    |
| 11     |         |       | 39     | BLU     | 14    |
| 12     |         |       | 40     | BLK     | 14    |
| 13     |         |       | 41     | BLU     | 14    |
| 14     | GRN/YEL | 18    | 42     | BLK     | 14    |
| 15     | GRN/YEL | 18    | 43     | BLU     | 14    |
| 16     |         |       | 44     | -       | -     |
| 17     | GRN/YEL | 14    | 45     | -       | -     |
| 18     |         |       | 46     | -       | -     |
| 19     |         |       | 47     | -       | -     |
| 20     | BLK     | 14    | 48     | -       | -     |
| 21     |         |       | 49     | -       | -     |
| 22     |         |       | 50     | -       | -     |
| 23     |         |       | 51     | -       | -     |
| 24     | YEL     | 14    | 52     | -       | -     |
| 25     | BLK     | 14    | 53     | YEL     | 14    |
| 26     | BLK     | 18    | 54     | -       | -     |
| 27     | BLK     | 18    | 55     | BLU     | 14    |
| 28     | RED     | 18    | 56     | -       | -     |
|        |         |       | 57     | -       | -     |
|        |         |       | 58     | WHT     | 18    |

MODELS  
13 CUFT. (23V) - VWR2400A  
17 CUFT. (23V) - VWR2800A

RS-232 SPECIFICATION

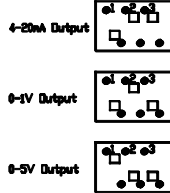
BAUD = 1200  
PARITY = N  
BITS = 8  
STOP BITS = 2

RS-232 PINS  
PIN 2 TXD  
PIN 3 RXD  
PIN 5 GND

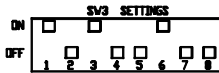
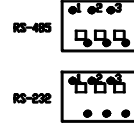
REMOTE CONTACTS/ANALOG OUTPUT  
PIN 1 Analog Output +  
PIN 2 Analog Output -  
PIN 3 Not Connected  
PIN 4 Not Connected  
PIN 5 Normally Closed  
PIN 6 Common  
PIN 7 Normally Open

CONTACT RATING 1A @ 30V  
CONTACTS IN ALARM STATE

SV1 Settings for Analog Output



SV2 Settings for Communication Output



1. Schematic represents single models.
2. Door switches shown in open position.
3. Battery switch shown in the OFF position.
4. Circuit breaker switch shown in OFF position.
5. Options and accessories shown in dashed lines.

| REV | DATE           | BY  | CHK | APP | DESCRIPTION OF REVISION              |
|-----|----------------|-----|-----|-----|--------------------------------------|
| 4   | 07/06 02-13-17 | KOB | TYI |     | UPDATED "MODELS" DETAILS             |
| 3   | 07-06/08-14-17 | KD  | CKG | KOB | SHW7 JACKUP MOVED TO SHW7            |
| 2   | 07-06/08-12-16 | CKG | SMS | DRP | SHW7 USE 200A                        |
| 1   | 06/06 02-07-15 | GLS | SMS | DRP | USE 60 PINS BLD TO REG. AS BULK CAFE |
| 0   | 07-07/08-12-14 | RV  | CKG | DRP | RELEASED FOR PROD / REL. II          |

Electrical Schematic  
Model:  
120V, 60Hz, -40°C  
ULT Upright Freezers

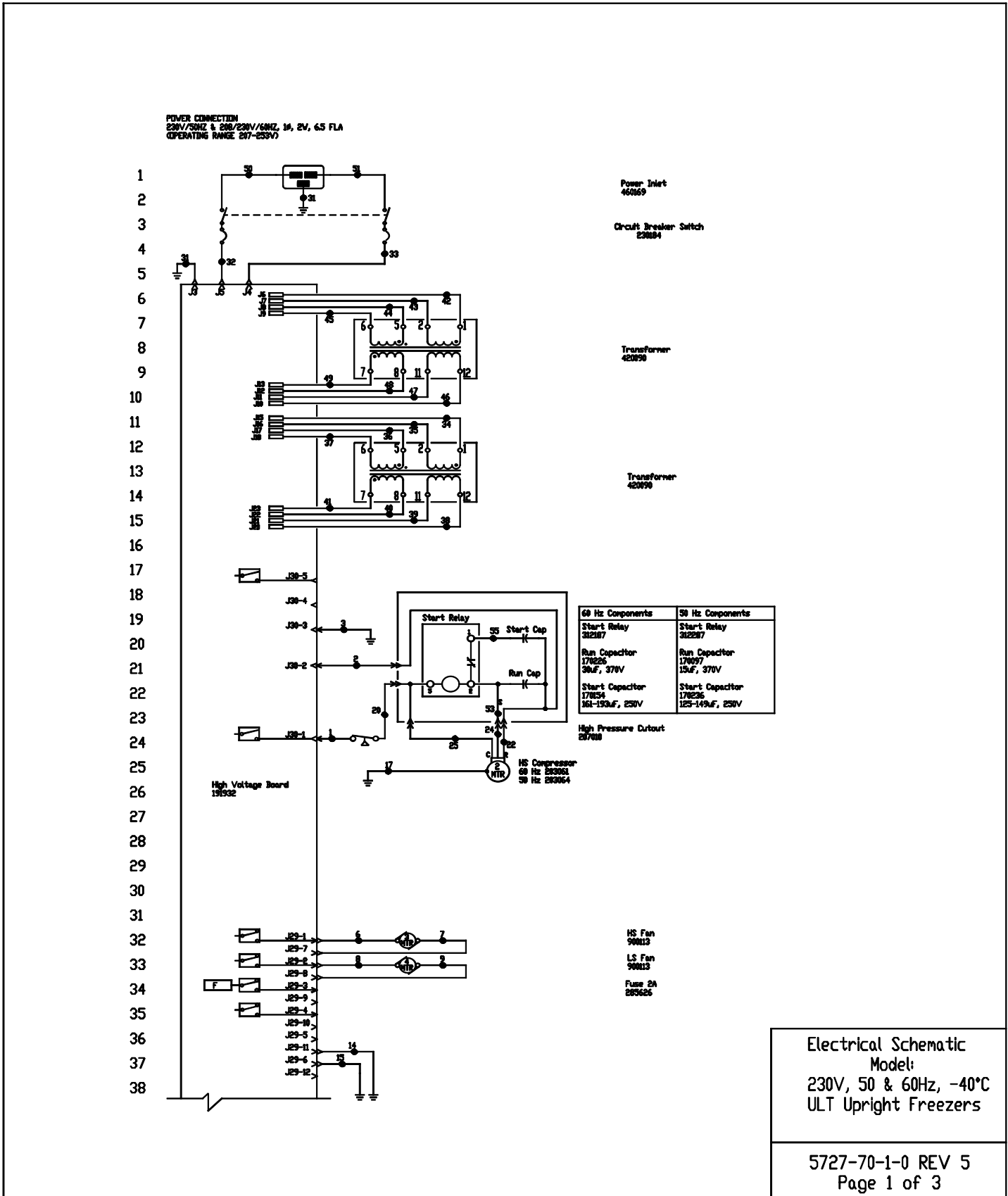


THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC.

**ThermoFisher SCIENTIFIC**  
BEE 649, MARICITA, CHD 4570

|                                      |  |
|--------------------------------------|--|
| MODEL/PART NAME                      | RESH-CHD, -SP UPRIGHT FREEZERS, 120V, 60Hz |
| 3/46 TITLE UNIT SCHEMATIC            |  |
| 3/46 RV                              | CKG KDG APPS                               |
| DATE                                 | 02-13-14                                   |
| SCALE                                | NONE                                       |
| PAPER COLOR                          |  |
| TOLERANCE UNLESS OTHERWISE SPECIFIED | DRAWING NUMBER                             |
| ANGLES: DECIMAL, 20±0.2, 30±0.2      | 5722-70-1                                  |
|                                      | SEC: D                                     |

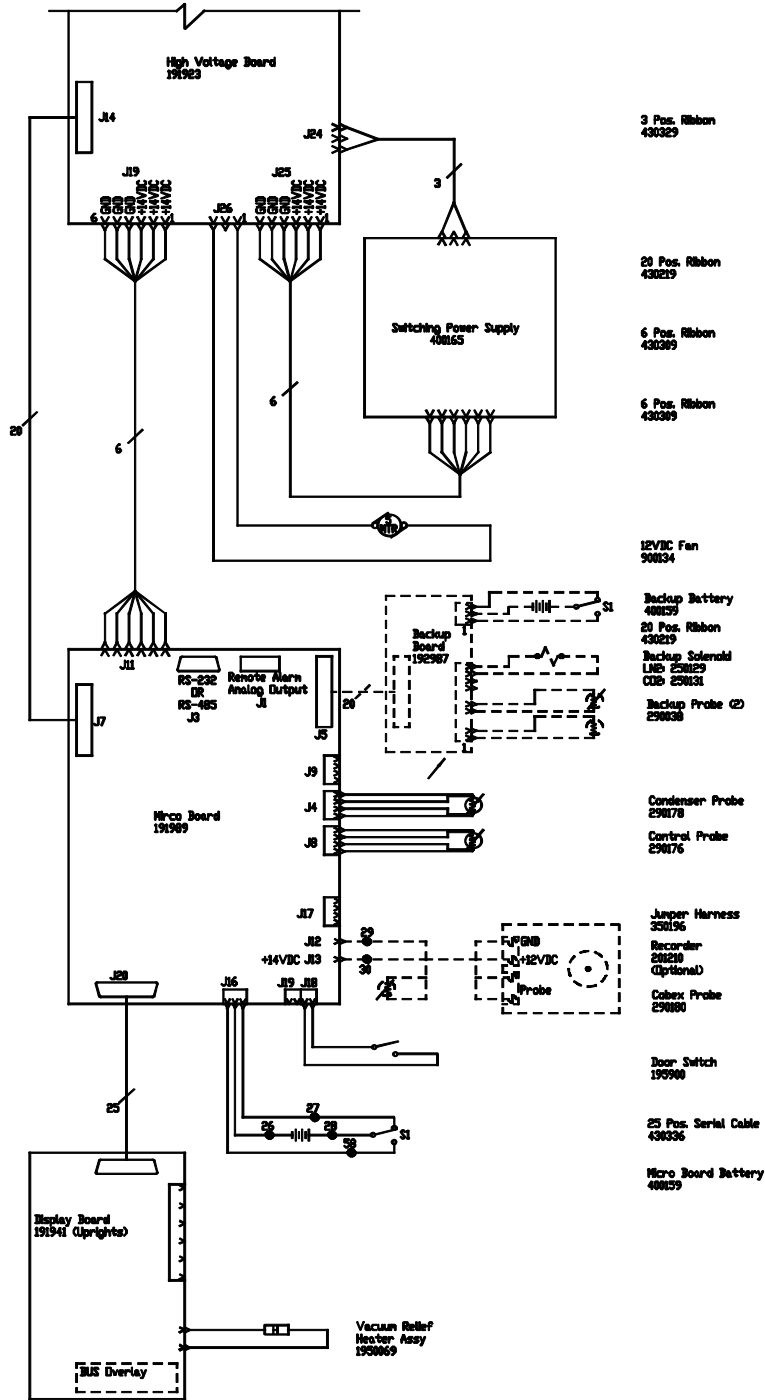
5722-70-1-0 REV 4  
Page 3 of 3



Electrical Schematic  
Model:  
230V, 50 & 60Hz, -40°C  
ULT Upright Freezers

**Section 10**  
Electrical Schematics

39  
40  
41  
42  
43  
44  
45  
46  
47  
48  
49  
50  
51  
52  
53  
54  
55  
56  
57  
58  
59  
60  
61  
62  
63  
64  
65  
66  
67  
68  
69  
70  
71  
72  
73  
74  
75  
76



Electrical Schematic  
Model:  
230V, 50 & 60Hz, -40°C  
ULT Upright Freezers

5727-70-1-0 REV 5  
Page 2 of 3

| WIRE # | COLOR | GAUGE | WIRE # | COLOR   | GAUGE |
|--------|-------|-------|--------|---------|-------|
| 78     |       |       |        |         |       |
| 79     | 1     | 14    | 29     | BLK     | 22    |
|        | 2     | 14    | 30     | RED     | 22    |
|        | 3     | 14    | 31     | GRN/YEL | 14    |
| 80     | 4     |       | 32     | BLK     | 14    |
|        | 5     |       | 33     | BLU     | 14    |
|        | 6     | 18    | 34     | BLK     | 14    |
| 81     | 7     | 18    | 35     | BLU     | 14    |
|        | 8     | 18    | 36     | BLK     | 14    |
| 82     | 9     | 18    | 37     | BLU     | 14    |
|        | 10    | 18    | 38     | BLK     | 14    |
|        | 11    |       | 39     | BLU     | 14    |
| 83     | 12    |       | 40     | BLK     | 14    |
|        | 13    |       | 41     | BLU     | 14    |
| 84     | 14    | 18    | 42     | BLK     | 14    |
|        | 15    | 18    | 43     | BLU     | 14    |
| 85     | 16    |       | 44     | BLK     | 14    |
|        | 17    | 14    | 45     | BLU     | 14    |
|        | 18    |       | 46     | BLK     | 14    |
| 86     | 19    |       | 47     | BLU     | 14    |
|        | 20    | 14    | 48     | BLK     | 14    |
| 87     | 21    |       | 49     | BLU     | 14    |
|        | 22    | 18    | 50     | BLK     | 14    |
|        | 23    |       | 51     | BLU     | 14    |
| 88     | 24    | 14    | 52     | -       | -     |
|        | 25    | 14    | 53     | YEL     | 14    |
|        | 26    | 18    | 54     | -       | -     |
| 89     | 27    | 18    | 55     | BLU     | 14    |
|        | 28    | 18    | 56     | -       | -     |
| 90     |       |       | 57     | -       | -     |
| 91     |       |       | 58     | WHT     | 18    |

MODELS:  
13 CUFT. (230V) - VVR2404B, VVR2404V  
17 CUFT. (230V) - VVR3204B, VVR3204V

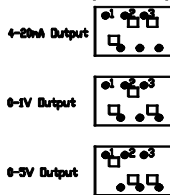
RS-232 SPECIFICATION

BAUD = 1200  
PARITY = N  
BITS = 8  
STOP BITS = 2

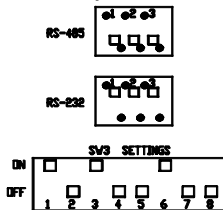
RS-232 PINS:  
PIN 2 TXD  
PIN 3 RXD  
PIN 5 GND

REMOTE CONTACTS/ANALOG OUTPUT  
PINS 1 Analog Output +  
PINS 2 Analog Output -  
PINS 3 Not Connected  
PINS 4 Not Connected  
PINS 5 Normally Closed  
PINS 6 Common  
PINS 7 Normally Open  
CONTACT RATING 1A @ 30V  
CONTACTS IN ALARM STATE

SV1 Settings for Analog Output



SV2 Settings for Communication Output



- Schematic represents upright models.
- Door switch shown in open position.
- Battery switch shown in the OFF position.
- Circuit breaker switch shown in OFF position.
- Options and accessories shown in dashed lines.

| REV | REV NO. | DATE     | BY  | CHK | APP'D | DESCRIPTION OF REVISION          |
|-----|---------|----------|-----|-----|-------|----------------------------------|
| 5   | 4/2/00  | 02-10-07 | WHD | TFI |       | UPDATED "MODEL#" DETAILS         |
| 4   | FR-2006 | 09-14-07 | ICJ | ICJ | END   | 210007 BACKUP BOARD TO 210007    |
| 3   | FR-2006 | 06-25-06 | END | SMS | IMP   | SHOWN WAS 2006, 2007 AND 2008    |
| 2   | FR-2004 | 04-25-04 | ICJ | SMS | IMP   | 2004 WAS 2004                    |
| 1   | FR-2004 | 04-25-04 | ICJ | SMS | IMP   | WIRE @ PIN 14 TO REMAIN BLT CURR |
| 0   | FR-2700 | 10-25-04 | WV  | ICJ | IMP   | RELEASED FOR PINS / REL II       |

Electrical Schematic  
Model:  
230V, 50 & 60Hz, -40°C  
ULT Upright Freezers



THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE, WITH OR WITHOUT WRITING, WITHOUT THE WRITTEN PERMISSION FROM THOMAS FISHER SCIENTIFIC.

**Thomas Fisher SCIENTIFIC**  
BOX 649, MARICCA, OHIO 45870

MATERIAL:  
PART COLOR:  
TOLERANCE UNLESS OTHERWISE SPECIFIED:  
ASSEMBLED BY: JSM  
REWORK: JSM

SWG TITLE: UNIT SCHEMATIC  
SWG NO: CAD ICG APP'D: DATED: 10-10-04 SCALE: NONE

SHAWING NUMBER: 5727-70-1  
SIZE: D

5727-70-1-0 REV 5  
Page 3 of 3



## VWR FREEZER WARRANTY - USA

The Warranty Period starts two weeks from the date your equipment is shipped from our facility. This allows for shipping time so the warranty will go into effect at approximately the same time your equipment is delivered. The warranty protection extends to any subsequent owner during the warranty period.

During the first two years of the warranty period, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at VWR's expense, labor included. The VWR ULT Freezers include an additional three year warranty on the compressors, parts only, F.O.B. factory. Installation and calibration is not covered by this warranty agreement. The Technical Services Department must be contacted for warranty determination and direction prior to any work being performed. Expendable items, i.e., glass, filters, pilot lights, light bulbs and door gaskets are excluded from this warranty.

Replacement or repair of component parts or equipment under this warranty shall not extend the warranty to either the equipment or to the component part beyond the original two year warranty period. VWR must give prior approval for the return of any components or equipment.

**THIS WARRANTY IS EXCLUSIVE AND IN LIEU OF ALL OTHER WARRANTIES, WHETHER WRITTEN, ORAL, OR IMPLIED. NO WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE SHALL APPLY. VWR shall not be liable for any indirect or consequential damages including, without limitation, damages relating to lost profits or loss of products.**

Your local VWR Sales Office is ready to help with comprehensive site preparation information before your equipment arrives. Printed instruction manuals carefully detail equipment installation, operation, and preventive maintenance.

If equipment service is required, please call the Technical Services Office at 1-800-438-4851 (USA and Canada) or 1-740-373-4763. We're ready to answer your questions on equipment warranty, operation, maintenance, service, and special applications. Outside the USA, contact your local distributor for warranty information.



Rev. 9 12/15

## **VWR FREEZER WARRANTY - INTERNATIONAL (EXCLUDING EUROPE)**

The Warranty Period starts two months from the date your equipment is shipped from our facility. This allows for shipping time so the warranty will go into effect at approximately the same time your equipment is delivered. The warranty protection extends to any subsequent owner during the warranty period.

During the first two years of the warranty period, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at VWR's expense, labor excluded. The VWR ULT Freezers include an additional three year warranty on the compressors, parts only, F.O.B. factory. Installation and calibration is not covered by this warranty agreement. The Technical Services Department must be contacted for warranty determination and direction prior to any work being performed. Expendable items, i.e., glass, filters, pilot lights, light bulbs and door gaskets are excluded from this warranty.

Replacement or repair of component parts or equipment under this warranty shall not extend the warranty to either the equipment or to the component part beyond the original two year warranty period. VWR must give prior approval for the return of any components or equipment.

**THIS WARRANTY IS EXCLUSIVE AND IN LIEU OF ALL OTHER WARRANTIES, WHETHER WRITTEN, ORAL, OR IMPLIED. NO WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE SHALL APPLY. VWR shall not be liable for any indirect or consequential damages including, without limitation, damages relating to lost profits or loss of products.**

Your local VWR Sales Office is ready to help with comprehensive site preparation information before your equipment arrives. Printed instruction manuals carefully detail equipment installation, operation, and preventive maintenance.

Outside the USA, for information or technical assistance, please contact your local VWR representative, or visit [vwr.com](http://vwr.com) for warranty information.



Rev. 9 12/15

## **VWR FREEZER WARRANTY - INTERNATIONAL (EUROPE ONLY)**

VWR warrants that this product will be free from defects in material and workmanship for a period of two (2) years from date of delivery. If a defect is present, VWR will, at its option and cost, repair, replace, or refund the purchase price of this product to the customer, provided it is returned during the warranty period. The VWR ULT Freezers include an additional three year warranty on the compressors, parts only, F.O.B. factory.

Installation and calibration is not covered by this warranty agreement. This warranty does not apply if the product has been damaged by accident, abuse, misuse, or misapplication, or from ordinary wear and tear. Expendable items, i.e., glass, filters, pilot lights, light bulbs and door gaskets are excluded from this warranty. If the required maintenance and inspection services are not performed according to the manuals and any local regulations, such warranty turns invalid, except to the extent, the defect of the product is not due to such non performance.

Items being returned must be insured by the customer against possible damage or loss. This warranty shall be limited to the aforementioned remedies. IT IS EXPRESSLY AGREED THAT THIS WARRANTY WILL BE IN LIEU OF ALL WARRANTIES OF FITNESS AND IN LIEU OF THE WARRANTY OF MERCHANTABILITY.



Rev. 9 12/15

# Appendix A Handling Liquid Nitrogen



**Warning** Contact of liquid nitrogen or cold gas with the skin or eyes may cause serious freezing (frostbite) injury. ▲

## Handle liquid nitrogen carefully.

The extremely low temperature can freeze human flesh very rapidly. When spilled on a surface the liquid tends to cover it completely and intimately, cooling a large area. The gas issuing from the liquid is also extremely cold. Delicate tissue, such as that of the eyes, can be damaged by an exposure to the cold gas which would be too brief to affect the skin of the hands or face.

## Never allow any unprotected part of your body to touch objects cooled by liquid nitrogen.

Such objects may stick fast to the skin and tear the flesh when you attempt to free yourself. Use tongs to withdraw objects immersed in the liquid, and handle the object carefully.

## Wear protective clothing.

Protect your eyes with a face shield or safety goggles (safety glasses without side shields do not give adequate protection). Always wear gloves when handling anything that is, or may have been, in immediate contact with liquid nitrogen. Insulated gloves are recommended, but heavy leather gloves may also be used. The gloves should fit loosely, so that they can be thrown off quickly if liquid should splash into them. When handling liquid in open containers, it is advisable to wear high-top shoes. Trousers (which should be cuffless if possible) should be worn outside the shoes.

## Introduction

The safe handling and use of liquid nitrogen in cryogenic refrigerators and dewar flasks is largely a matter of knowing the potential hazards and using common-sense procedures based on that knowledge. There are two important properties of liquid nitrogen that present potential hazards:

1. It is extremely cold. At atmospheric pressure, liquid nitrogen boils at -320°F (-196°C).
2. Very small amounts of liquid vaporize into large amounts of gas. One liter of liquid nitrogen becomes 24.6 cu. ft. (700l) of gas.

The safety precautions in this booklet must be followed to avoid potential injury or damage which could result from these two characteristics. Do not attempt to handle liquid nitrogen until you read and fully understand the potential hazards, their consequences, and the related safety precautions. Keep this booklet handy for ready reference and review.

**Note** Argon is an inert gas whose physical properties are very similar to those of nitrogen. The precautions and safe practices for the handling and use of liquid argon are the same as those for liquid nitrogen. ▲

### Use only containers designed for low temperature liquids.

Cryogenic containers are specifically designed and made of materials that can withstand the rapid changes and extreme temperature differences encountered in working with liquid nitrogen. Even these special containers should be filled SLOWLY to minimize the internal stresses that occur when any material is cooled. Excessive internal stresses can damage the container.

**Do not cover or plug the entrance opening of any liquid nitrogen refrigerator or dewar. Do not use any stopper or other device that would interfere with venting of gas.**

These cryogenic liquid containers are generally designed to operate with little or no internal pressure. Inadequate venting can result in excessive gas pressure which could damage or burst the container. Use only the loose-fitting necktube core supplied or one of the approved accessories for closing the necktube. Check the unit periodically to be sure that venting is not restricted by accumulated ice or frost.

### Use proper transfer equipment.

Use a phase separator or special filling funnel to prevent splashing and spilling when transferring liquid nitrogen into or from a dewar or refrigerator. The top of the funnel should be partly covered to reduce splashing. Use only small, easily-handled dewars for pouring liquid. For the larger, heavier containers, use a cryogenic liquid withdrawal device to transfer liquid from one container to another. Be sure to follow instructions supplied with the withdrawal device. When liquid cylinders or other large storage containers are used for filling, follow the instructions supplied with those units and their accessories.

## Introduction (continued)

### Do not overfill containers.

Filling above the bottom of the necktube (or specified maximum level) can result in overflow and spillage of liquid when the necktube core or cover is placed in the opening.

### Never use hollow rods or tubes as dipsticks.

When a warm tube is inserted into liquid nitrogen, liquid will spout from the top of the tube due to gasification and rapid expansion of liquid inside the tube.



### **Warning** Nitrogen gas can cause suffocation without warning! ▲

### Store and use liquid nitrogen only in a well-ventilated place.

As the liquid evaporates, the resulting gas tends to displace the normal air from the area. In closed areas, excessive amounts of nitrogen gas reduce the concentration of oxygen and can result in asphyxiation. Because nitrogen gas is colorless, odorless and tasteless, it cannot be detected by the human senses and will be breathed as if it were air. Breathing an atmosphere that contains less than 18% oxygen can cause dizziness and quickly result in unconsciousness and death.

**Note** The cloudy vapor that appears when liquid nitrogen is exposed to the air is condensed moisture; not the gas itself. The issuing gas is invisible.

### Never dispose of liquid nitrogen in confined areas or places where others may enter.

Disposal of liquid nitrogen should be done outdoors in a safe place. Pour the liquid slowly on gravel or bare earth where it can evaporate without causing damage. Do not pour the liquid on pavement.

## Handling Liquid Carbon Dioxide



**Warning** High concentrations of CO<sub>2</sub> gas can cause asphyxiation! OSHA Standards specify that employee exposure to carbon dioxide in any eight-hour shift of a 40-hour work week shall not exceed the eight-hour time weighted average of 5000 PPM (0.5% CO<sub>2</sub>). The short term exposure limit for 15 minutes or less is 30,000 PPM (3% CO<sub>2</sub>). Carbon dioxide monitors are recommended for confined areas where concentrations of carbon dioxide gas can accumulate. ▲

### Store and use liquid CO<sub>2</sub> only in a well-ventilated place.

As the liquid evaporates, the resulting gas tends to displace the normal air from the area. In closed areas, excessive amounts of CO<sub>2</sub> gas reduce the concentration of oxygen and can result in asphyxiation. Because CO<sub>2</sub> gas is colorless, odorless and tasteless, it cannot be detected by the human senses and will be breathed as if it were air. Breathing an atmosphere that contains less than 18% oxygen can cause dizziness and quickly result in unconsciousness and death.

**Note** The cloudy vapor that appears when liquid CO<sub>2</sub> is exposed to the air is condensed moisture, not the gas itself. The issuing gas is invisible.

### Never dispose of liquid CO<sub>2</sub> in confined areas or places where others may enter.

Disposal of liquid CO<sub>2</sub> should be done outdoors in a safe place. Pour the liquid slowly on gravel or bare earth where it can evaporate without causing damage. Do not pour the liquid on pavement.

## First Aid

If a person seems to become dizzy or loses consciousness while working with liquid nitrogen or carbon dioxide, move to a well-ventilated area immediately. If breathing has stopped, apply artificial respiration. If breathing is difficult, give oxygen. Call a physician. Keep warm and at rest.

If exposed to liquid or cold gas, restore tissue to normal body temperature (98.6° F) as rapidly as possible, followed by protection of the injured tissue from further damage and infection. Remove or loosen clothing that may constrict blood circulation to the frozen area. Call a physician. Rapid warming of the affected part is best achieved by using water at 108° F. Under no circumstance should the water be over 112° F, nor should the frozen part be rubbed either before or after rewarming. The patient should neither smoke nor drink alcohol.

**FIELD INSTALLED ACCESSORIES**

| <b>NA VWR #</b> | <b>EU VWR #</b> | <b>DESCRIPTION</b>  | <b>-86°C Upright Freezers</b> | <b>-40°C Upright Freezers</b> | <b>3 cu ft (85L) Chest Freezers</b> | <b>Other Chest Freezers</b> |
|-----------------|-----------------|---|-------------------------------|-------------------------------|-------------------------------------|-----------------------------|
| 10185-060       | 471-1144        | CO2 BUS Kit VWR Collection Upright Freezers Field Installed Option                  | x                             | x                             | x                                   | x                           |
| 10185-062       | 471-1145        | LN2 BUS Kit VWR Collection Upright Freezers Field Installed Option                  | x                             | x                             | x                                   | x                           |
| 10185-064       | 471-1146        | Chart Recorder, Pen VWR Collection Upright Freezers *CI, Field Installed Option     | x                             | x                             |                                     |                             |
| 10185-072       | 471-1147        | Chart Recorder, Pen VWR Collection Chest Freezers *CI, Field Installed Option       |                               |                               | x                                   | x                           |
| 10185-068       | 471-1148        | Chart Recorder, Inkless VWR Collection Upright Freezers *CI, Field Installed Option | x                             | x                             |                                     |                             |
| 10185-076       | 471-1149        | Chart Recorder, Inkless VWR Collection Chart Freezers *CI, Field Installed Option   |                               |                               | x                                   | x                           |
| 10185-082       | 471-1150        | S.S. Shelf Kit 13 Cu Ft ULT   | x                             | x                             |                                     |                             |
| 10185-084       | 471-1151        | S.S. Shelf Kit 17.3 Cu Ft ULT   | x                             | x                             |                                     |                             |
| 10185-086       | 471-1152        | S.S. Shelf Kit 23 Cu Ft ULT   | x                             |                               |                                     |                             |
| 10185-088       | 471-1153        | S.S. Shelf Kit 28 Cu Ft ULT   | x                             |                               |                                     |                             |
| 89202-604       | 471-1154        | Replacement Back-up Battery   | x                             | x                             | x                                   | x                           |





