

Model 8600 Series

-86C Ultra Low Temperature Freezer

Operating and Maintenance Manual 7048602 Rev. 0



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 1-3) terminates in an interconnect jack (Figure 1-5) behind the base front cover. (May be located differently in chests. See Section 1.) 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. s

Model	Capacity in Cubic Feet	Voltage
8602	13	230
8603	13	120
8604	17	120
8605	17	230
8606	23	230
8607	28	230
8656	23	120
Double Door Units		
8690	23	120
8691	13	230
8692	13	120
8693	17	120
8694	17	230
8695	23	230

MANUAL NUMBER 7048602

0	24020/FR-2145	1/17/10	Release 5 - Relay enclosure	ccs
REV	ECR/ECN	DATE	DESCRIPTION	By



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. s

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

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

©2010 Thermo Fisher Scientific. 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.



Hot surface(s) present which may cause burns to unprotected skin, or to materials which may be damaged by elevated temperatures.



Marking of electrical and electronic equipment, which applies to electrical and electronic equipment falling under the Directive 2002/96/EC (WEEE) and the equipment that has been put on the market after 13 August 2005.



This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96/EC. It is marked with the WEEE symbol. Thermo Fisher Scientific has contracted with one or more recycling/disposal companies in each EU Member State European Country, and this product should be disposed of or recycled through them. Further information on Thermo's compliance with this directive, the recyclers in your country and information on Thermo products will be available at www.thermo.com.

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

Do You Need Information or Assistance on Thermo Scientific Products?

If you do, please contact us 8:00 a.m. to 6:00 p.m. (Eastern Time) at:

1-740-373-4763

Direct

1-800-438-4851

Toll Free, U.S. and Canada

1-877-213-8051

FAX

<http://www.thermoscientific.com>

Internet Worldwide Web Home Page

service.led.marietta@thermofisher.com

Service E-Mail Address

Our **Sales Support** staff 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.

Our **Service Support** staff can supply technical information about proper setup, operation or troubleshooting of your equipment. We can fill your needs for spare or replacement parts or provide you with on-site service. We can also provide you with a quotation on our Extended Warranty for your Thermo Scientific products.

Whatever Thermo Scientific 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 direct factory trained technicians or a qualified service organization for on-the-spot repair. If your service need is covered by the warranty, we will arrange for the unit to be repaired at our expense and to your satisfaction.

Regardless of your needs, our professional telephone technicians are available to assist you Monday through Friday from 8:00 a.m. to 6:00 p.m. Eastern Time. Please contact us by telephone or fax. If you wish to write, our mailing address is:

Thermo Fisher Scientific
401 Millcreek Road, Box 649
Marietta, OH 45750

International customers, please contact your local Thermo Scientific distributor.

Table of Contents

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

Section 4	Alarms	4-1
	High Stage System Failure Alarm	4-2
	Error Messages	4-3
Section 5	Maintenance	5-1
	Clean Cabinet Exterior	5-1
	Clean Air Filter	5-1
	Clean the Condenser	5-1
	Clean the Water-cooled Condenser	5-2
	Clean the Door Gasket	5-3
	Defrost the Chamber	5-3
	Vacuum Relief Port	5-4
	Replace the Battery(s)	5-5
	Prepare the Unit for Storage	5-6
Section 6	Factory Installed Options	6-1
	BUS-Back-Up System	6-1
	Install Injection Assembly, Vent Stack and Solenoid	6-1
	Install the Temperature Probe	6-3
	Connect the Probe/Solenoid Harness	6-4
	BUS Control Panel	6-5
	Test the BUS	6-6
	Clean the Vent Stack	6-6
	Set the Optional BUS Set Point	6-6
	Disconnect the Fitting Assembly & Transfer Hose	6-7
	Chart Recorder	6-7
	Install the Chart Paper	6-7
	Change Recorder Temperature Range	6-8
	Calibrate the Chart Recorder	6-8
	Five Inner Door Option	6-9
	Water-cooled Condenser	6-9
	Datalogger	6-9
Section 7	Specifications	7-1
Appendix	Handling Liquid Nitrogen	A-1
	Handling Liquid CO₂	B-1
	First Aid	C-1

Section 1 Installation and Start-Up

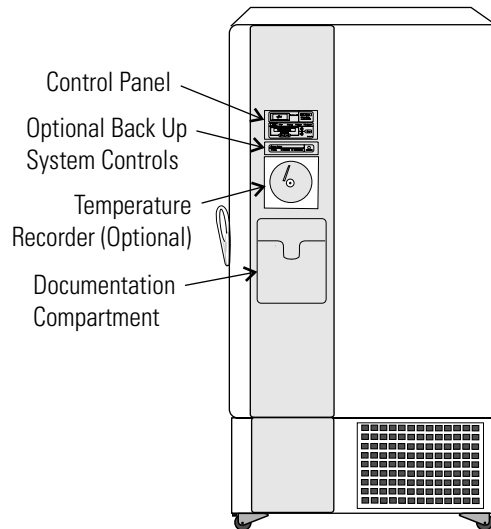


Figure 1-1. Front View

- Control Panel - keypad, displays and indicators.
- BUS (Optional Back Up System) panel.
- Optional temperature recorder - 7 day, one pen or datalogger.
- Documentation compartment - storage of user's manual and other documentation.

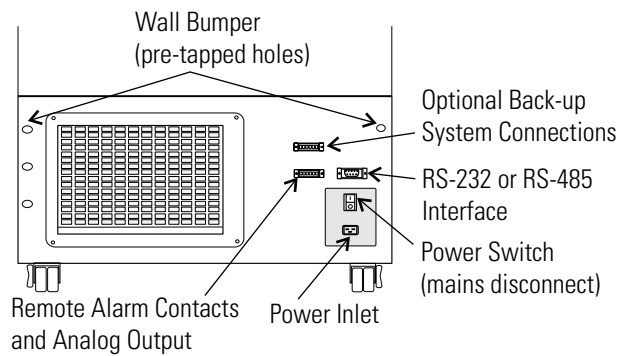


Figure 1-2. Rear View

- 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).

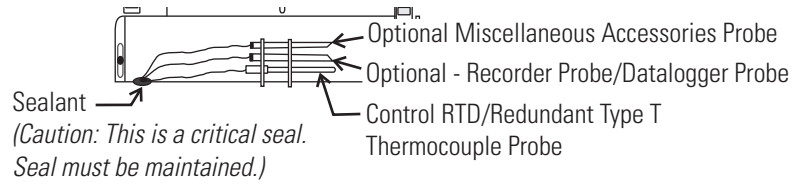


Figure 1-3. Chamber Probe

Figures 1-3 & 1-4

- Vacuum relief port - pressure equalization port
- Probe cover - houses control, optional recorder, datalogger or 1535 alarm probes

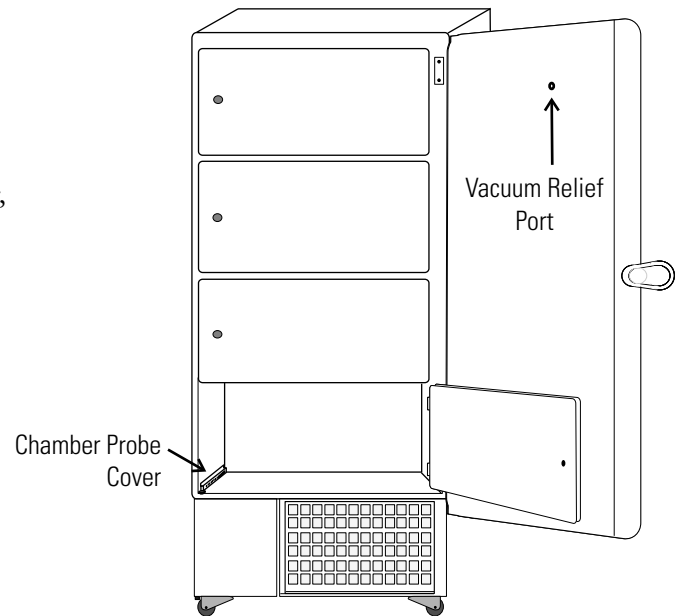


Figure 1-4. Vacuum Relief and Probe Cover Location

Figure 1-5

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

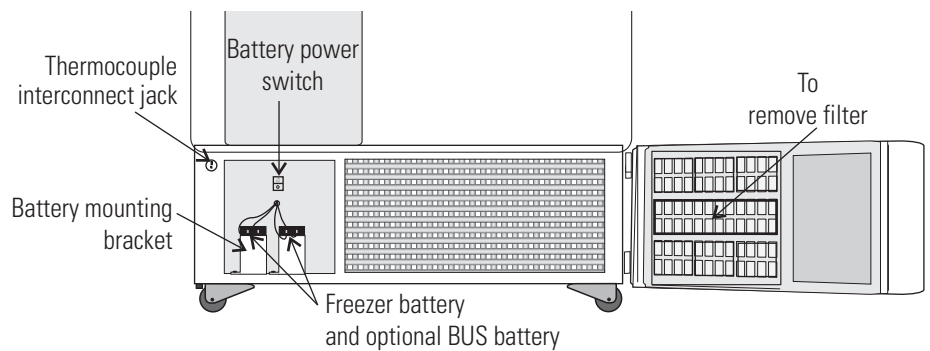


Figure 1-5. Battery(s) Location and Switch

Control Panel Keys, Displays & Indicators

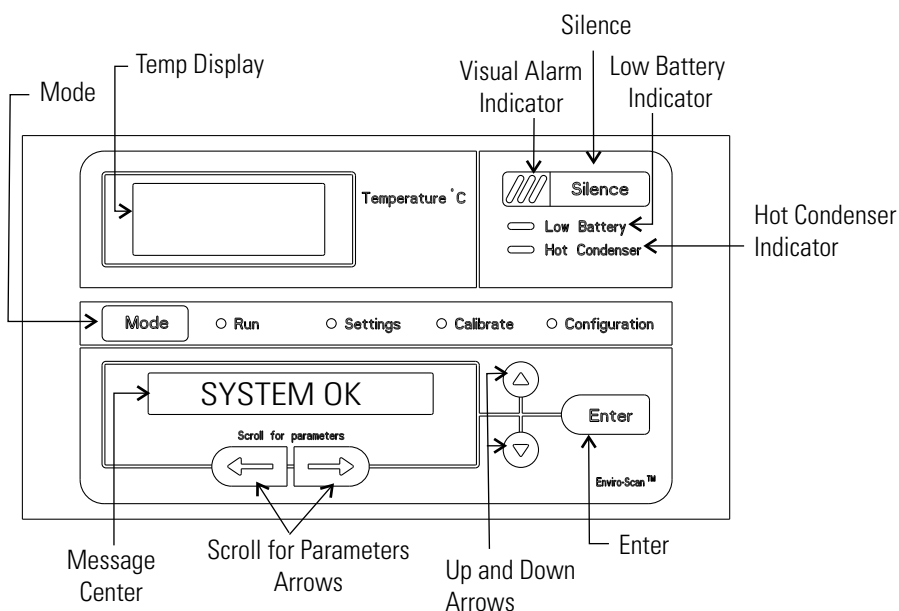


Figure 1-6. Control Panel

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 a cabinet alarm condition.

Silence - 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 - Stores the value into computer memory.

Keypad Operation

Model 8600 Series freezer has four basic modes which allow freezer setup: Run, Settings, Calibrate and Configuration.

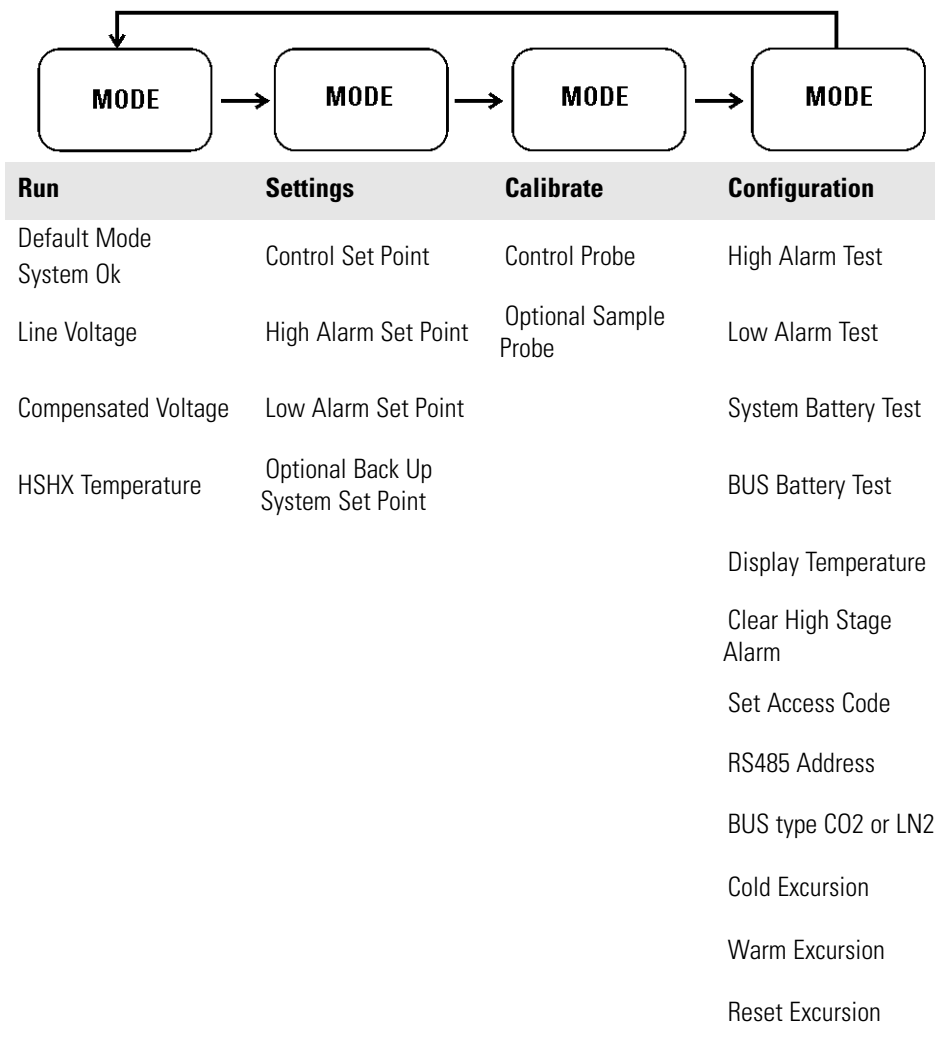
Run is the default mode for the freezer during normal operation.

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.



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 all changed values.

Keypad Operation (continued)

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

Silence 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 Section 4 - Alarms.

Install the Freezer

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

To remove the freezer from the pallet, use the 7/16" 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 in the following section. 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 The freezer must not be moved with the product load inside. s

Choose the Location

Locate the freezer on a firm, level surface in an area with an ambient temperature between 18°C and 32°C. Provide ample room to reach the mains disconnect switch (power switch) located on the rear of the freezer.

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 the front of the freezer for door opening. s

Install the Wall Bumpers

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

Qty	Stock #	Description	Purpose
2	510016	1/4-20 x 5-1/2" Bolt	Wall Bumper
2	380520	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 locations 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. s

RS-232 Communications

The Model 8600 Series freezer has a data communications interface. The factory default setting is RS-232.

The wiring identification for the interface is shown in Figure 1-7. 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.

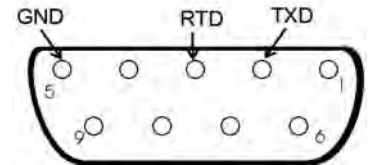


Figure 1-7. Wiring

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

Data format:

Baud1200
 Data bits8 (7 bit ASCII with leading zero)
 Start bits1
 Stop bits1
 Paritynone

RS-232 Communications (continued)

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)

Note The RS-232 is not compatible with Model 1535 Monitor/Alarm System. s

Remote Alarm Contacts and Analog Output

Model 8600 Series freezer has remote alarm contacts and analog output. See Figure 1-2 for the location of the remote alarm contacts. The remote alarm connector is shipped 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. 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 or microboard failures. Figure 1-8 shows the remote contacts in alarm state.

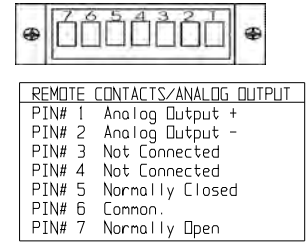


Figure 1-8. Remote Contacts in Alarm

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 1-9 for output specifications.

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

Figure 1-9. Output 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.

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

Connect the Unit to Electrical Power

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.

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.

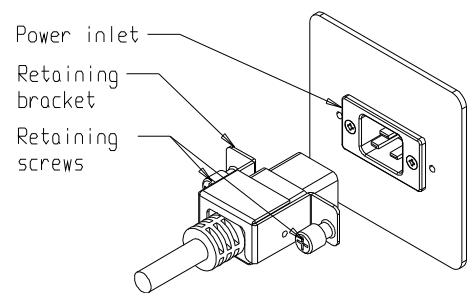


Figure 1-10. Secure Power Cord

Connect the Unit to Electrical Power (cont.)

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 indicator may illuminate.

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. s

Freezer Start-Up

With the freezer properly installed and connected to power, system set points can be entered. The following set points can be entered in Settings mode: Control temperature, high temperature alarm set point, low temperature alarm set point, and (optional) BUS set point. Default settings are shown in Table 1-1.

Table 1-1. Default Settings

Control Set Point	-80°C
High Temperature Alarm	-70°C
Low temperature alarm	-90°C
Optional BUS Set Point	-60°C

Caution If the set point is changed and the low temperature and high temperature alarms are set 10° from the set point, the alarm set points will adjust automatically to maintain a distance of at least 10° from set point. s

Set the Operating Temperature

All Model 8600 Series freezers have an operating temperature range of -50°C to -86°C, depending on ambient temperature. The freezer is shipped from the factory with a temperature set point of -80°C. To change the operating temperature set point:

1. Press the Mode key until the Settings indicator lights.
2. Press the right arrow until “SET PT = -XX” is displayed in the 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.

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 the right arrow until “HI ALM = -XX” is displayed in the 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.

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

Note At initial start-up, the high temperature alarm is disabled until the cabinet reaches set point or 12 hours elapse. s

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. s

Access Code

An access code of 000 is required to access the Settings, Calibrate or Configuration modes. If the access code is not at the default 000, a code must be entered to leave RUN mode. See Section 3 to modify the access code.

Run Mode

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

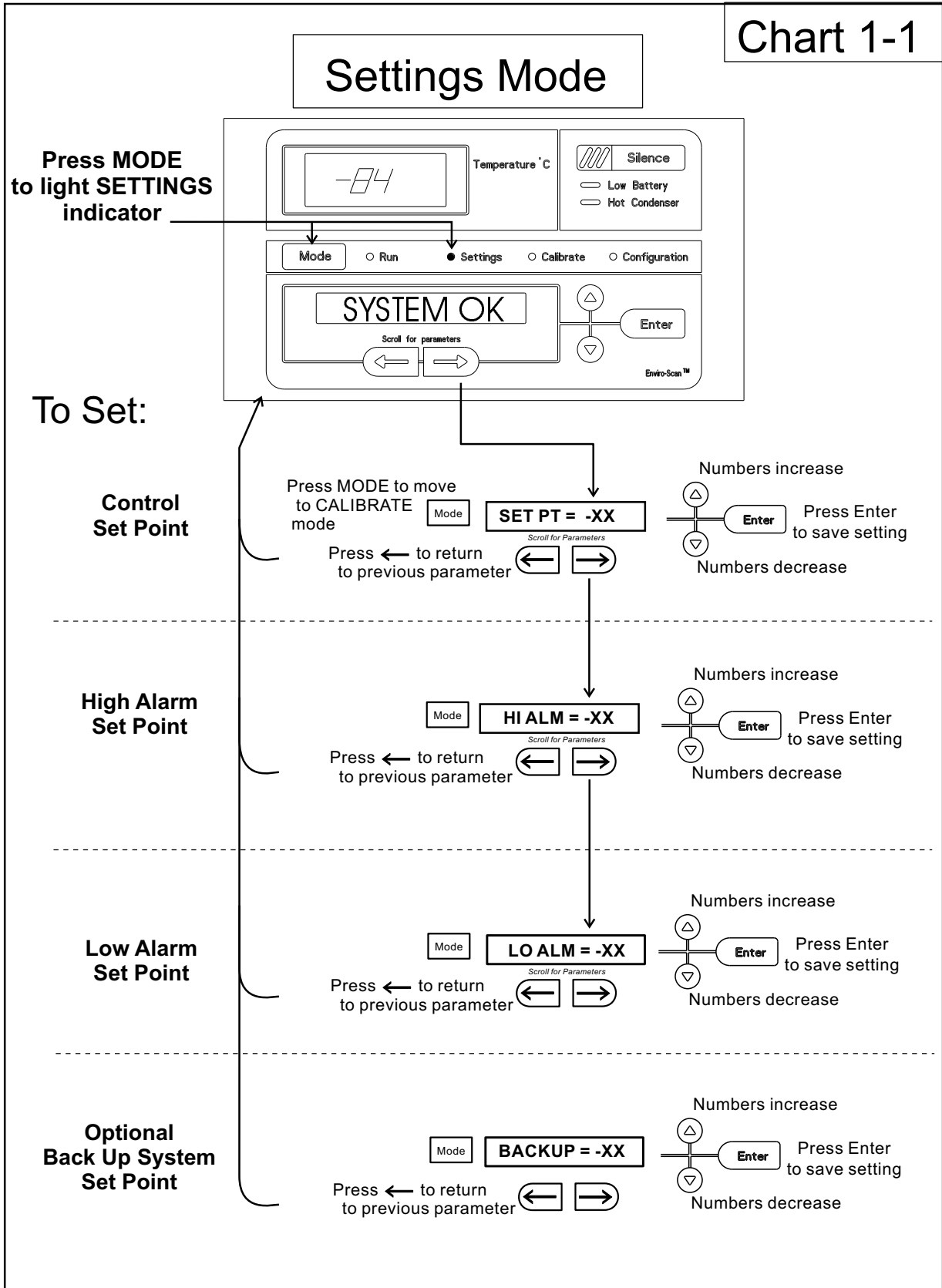
LINE VOLTAGE

COMPENSATED VOLTAGE

HSHX TEMPERATURE (heat exchanger temperature)

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 1-1



Section 2 Calibrate

After the freezer has stabilized, the control or sample probe may require calibration. 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.

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

Calibrate the Control Probe

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

1. Press the Mode key until the Calibrate indicator lights.
2. Press the right arrow until “CONT T = -XX.X” appears in the 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.

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 2-1 for calibration process functions.

Temperature Stabilization Periods

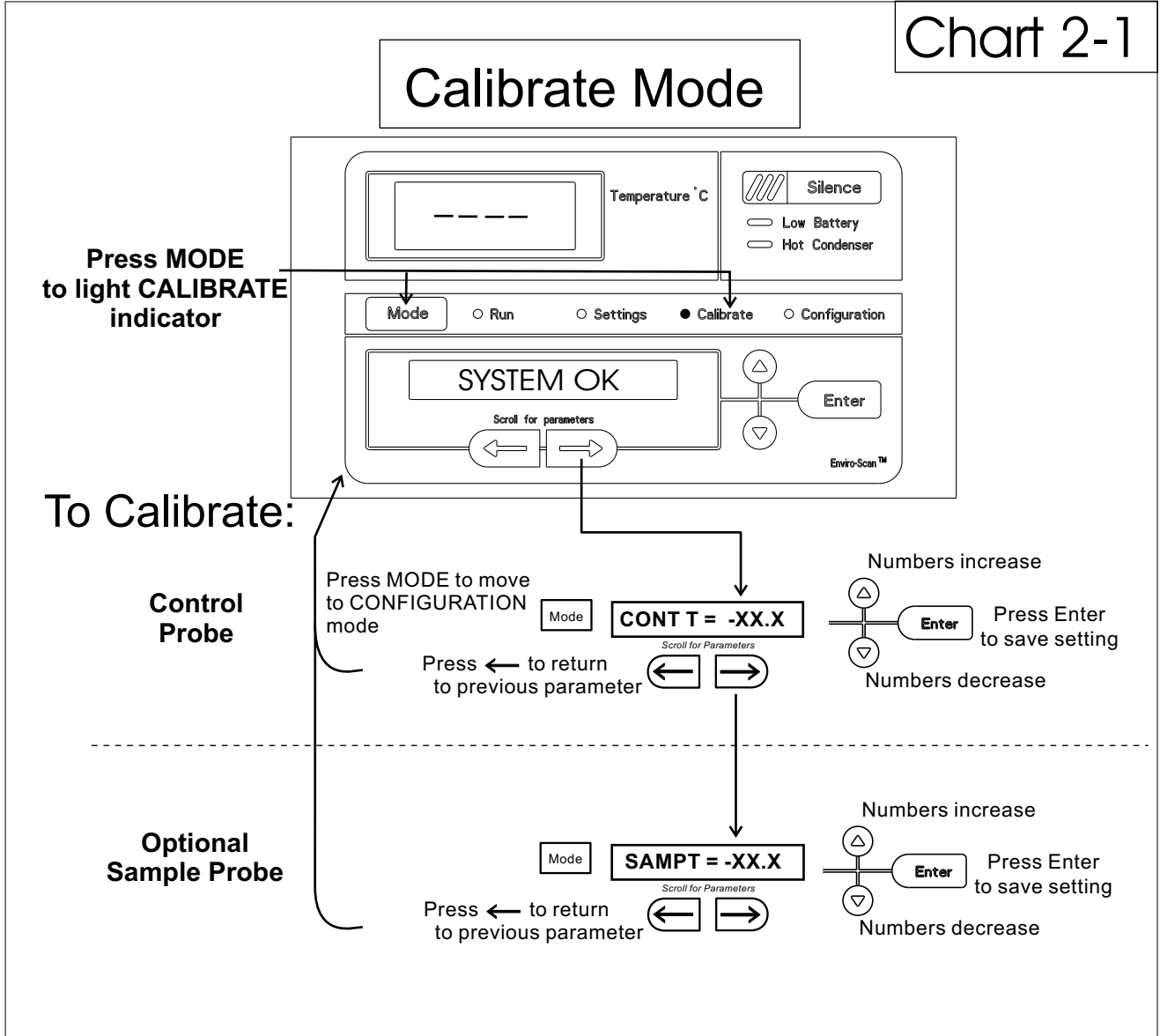
Startup - Allow 12 hours for the temperature in the cabinet to stabilize before proceeding.

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

Note During calibration, the temperature display is not available. s

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

Chart 2-1



Section 3 Configuration

Configuration mode is used for testing and custom setup of the freezer. The configuration functions 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 that the high alarm activates if the freezer temperature equals or exceeds 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 begins to increase until the high alarm set point has been reached. The audible alarm sounds and the alarm indicator flashes. Press the Silence key to silence the alarm.

Low Alarm Test

The low alarm test is used to verify that the low alarm activates if the freezer temperature equals or decreased to 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 begins to decrease until the low alarm set point has been reached. The audible alarm sounds and the alarm indicator flashes. Press the Silence key to silence the alarm.

System Battery Test

To test the freezer battery charge:

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. If a test fails, the audible alarm sounds, the alarm indicator and Low Battery indicator light. Press the Silence key and the alarm indicator will go off. The Low Battery light stays on until a future battery test is performed and passed.

BUS Battery Test

To test the charge of the back-up system (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. If a test is failed, the audible alarm sounds, the alarm indicator and the Low Battery indicator light. Press the Silence key. The audible alarm and alarm indicator will go off. The Low Battery light stays on. If this test fails, battery replacement is recommended.

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 occur, it may be 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 an 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 displays ACC CODE = 000. Press the up/down arrow until the desired access code is displayed (000 - 999). Press the left/right arrow 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. s

5. Press Enter to save the setting
6. Press the Mode key until the Run indicator lights. A 3-digit Access Code can be entered to avoid unauthorized personnel from changing the set points, calibration, or configuration. A setting of 000 bypasses the access code. The factory setting is 000.

RS485 Address

If the freezer is configured for RS485 communications, it must have a unique identification address. This address is set through 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/down arrow to select the appropriate address for the freezer (1 - 24).
5. Press Enter to save.

Back-Up System Type

This function, 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.

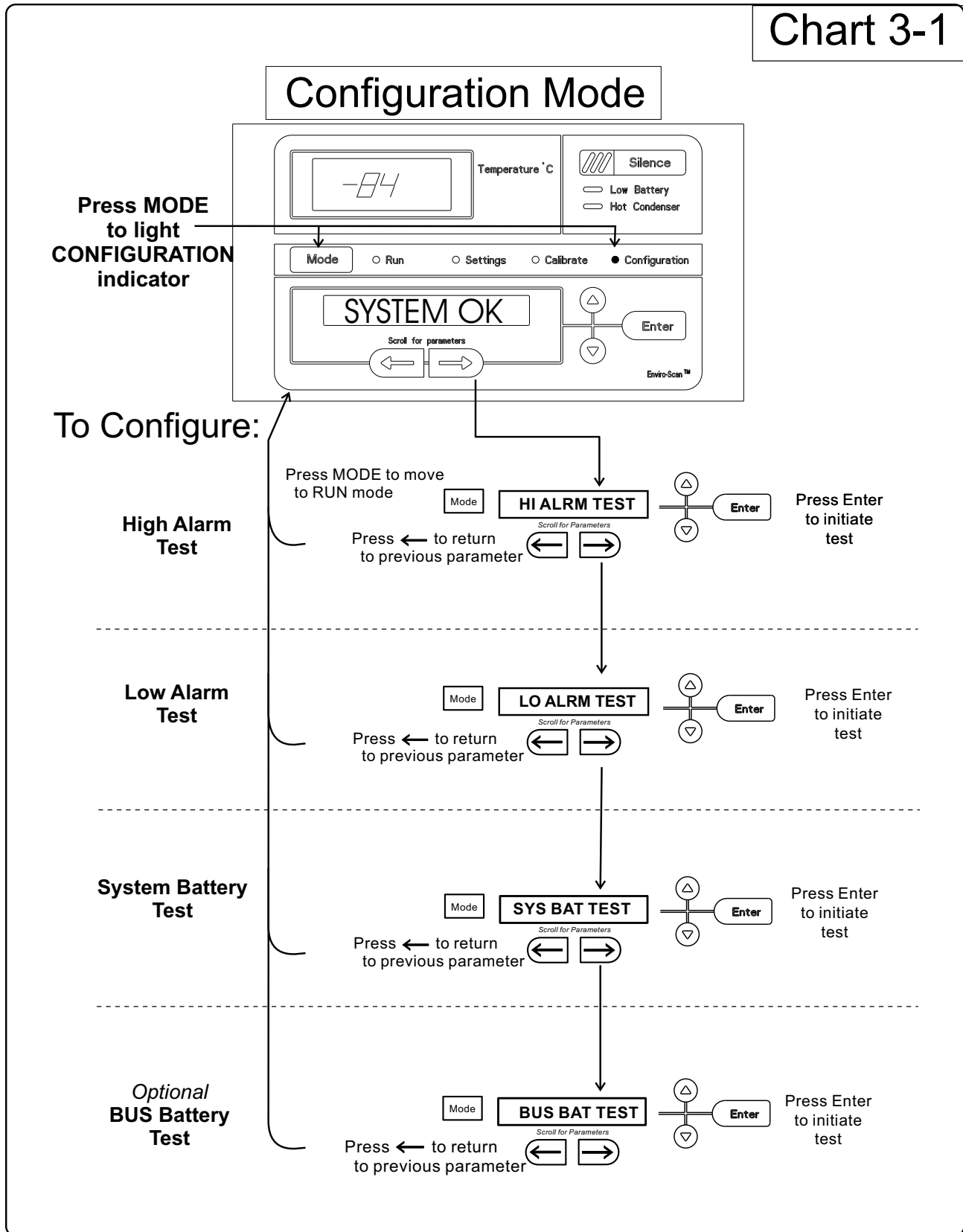
Warm Excursion

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

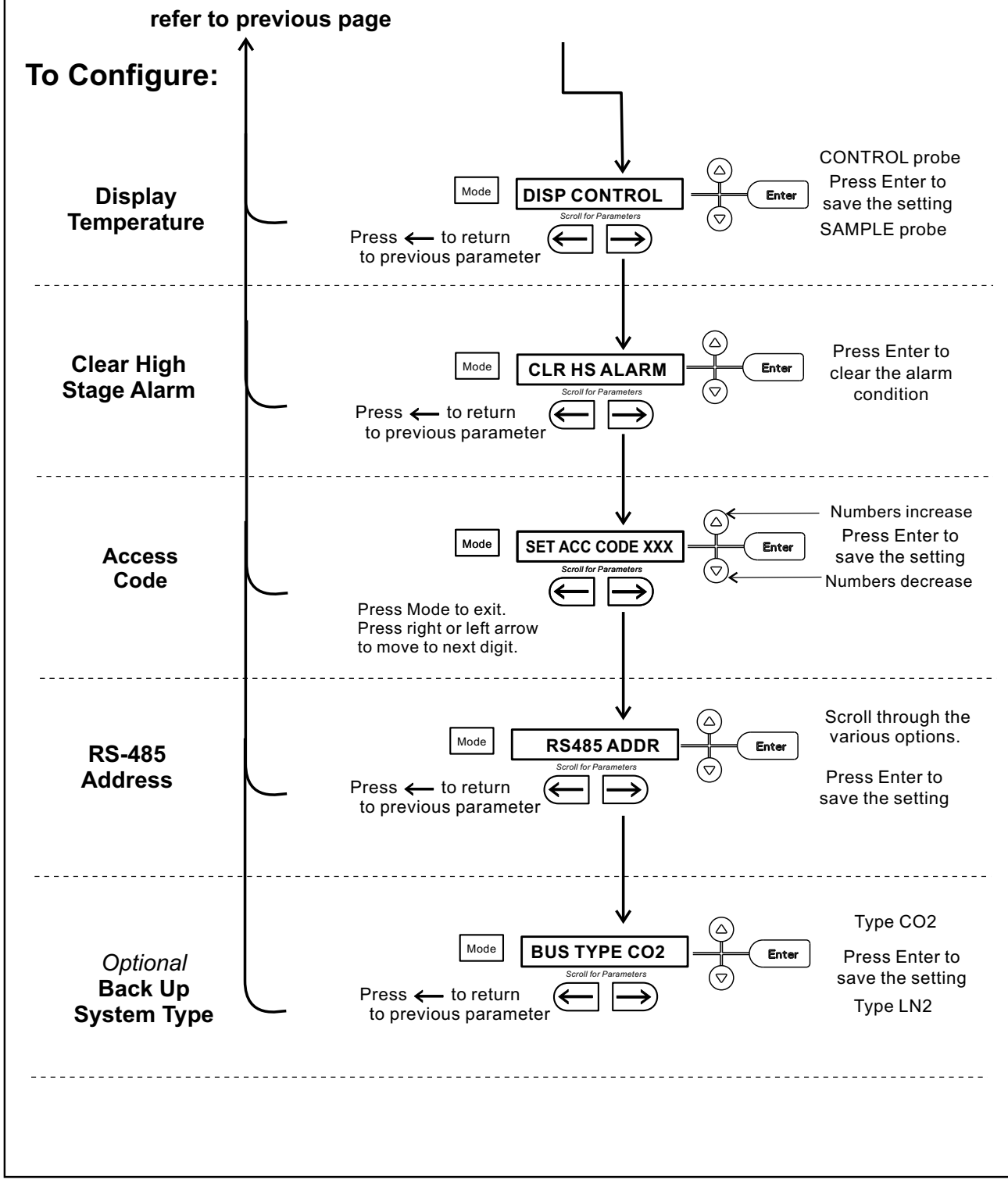
Reset Excursion

This function resets the cold and warm excursions.

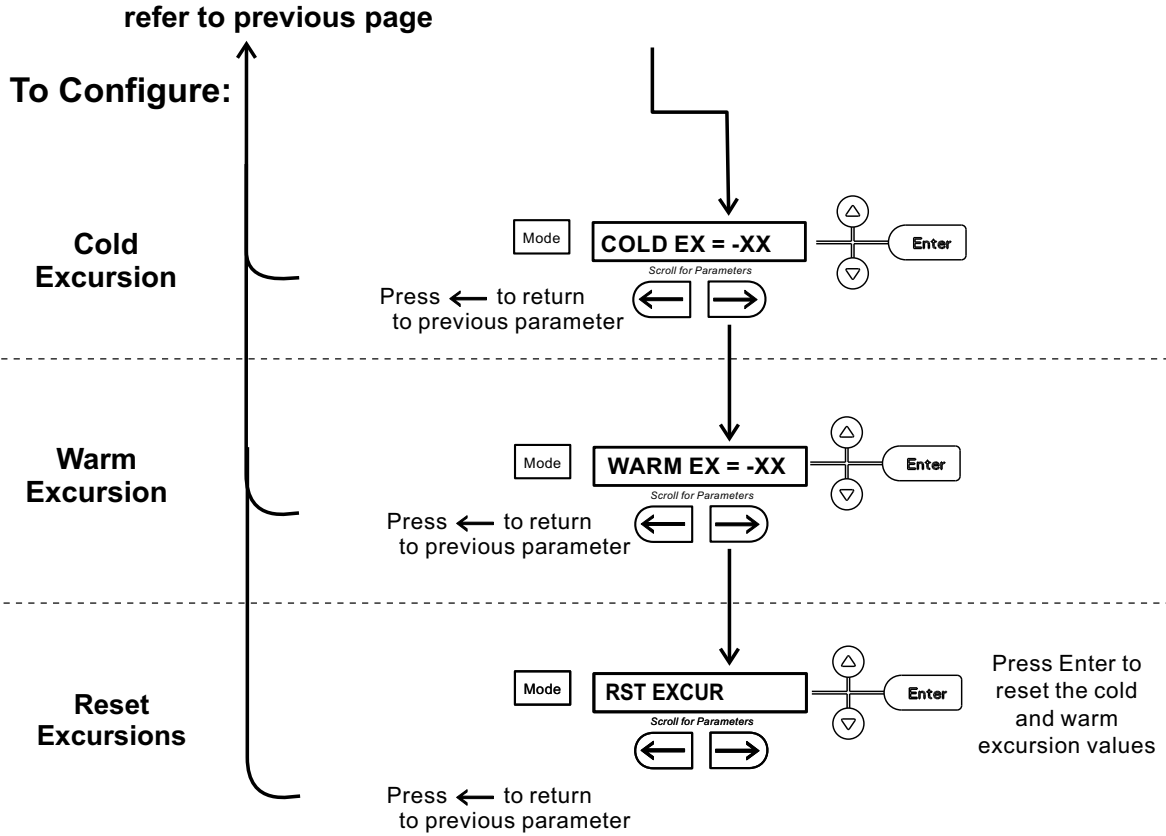
Chart 3-1



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



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



Section 4 Alarms

Model 8600 Series freezer alarm system is shown in the table below. When an alarm is active, the message appears in the LED message center. Press the Silence key to disable 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.

Table 4-1. Alarms

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	CNT PRB FLT	1 min.	15 min.	Yes
Heat Exchanger Probe Failure	HSHX PRB FLT	1 min.	15 min.	No
Condenser Probe	COND PRB FLT	1 min.	15 min.	No
Sample Probe Failure (optional)	SMPL PRB FLT	1 min.	15 min.	No
High Stage System Failure	HS SYST FAIL	1 min.	15 min.	YES
Condenser Hot Condition	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 +30 seconds.

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

High Stage System Failure Alarm

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 Silence during multiple alarms causes all active alarms to be silenced and to ringback 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

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>Name: Improper model selected.</p> <p>Description: Indicates that DIP SW3 has not selected a proper model or can't be accessed properly.</p> <p>Response: 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>Name: Voltage/Frequency failure</p> <p>Description: 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 (<180VAC < 270 for 230VAC unit / <85VAC < 160 for 115VAC unit); or the frequency measured over 10 cycles was not within a tolerable range (55 Hz < Freq < 70 Hz for 60 Hz units / 40 Hz < Freq < 55 Hz for 50 Hz units)</p> <p>Response: 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 4

Alarms

Error (cont.)	High End Message	Notes
Er02	"CNT PRB FLT"	<p>Name: Control (Cabinet) Sensor Failure</p> <p>Description: This condition indicates that the control sensor has failed to produce a valid reading for ≥ 12 consecutive reads (~60 seconds).</p> <p>Response: 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>Name: Heat Exchange Sensor Failure</p> <p>Description: This condition indicates that the heat exchange sensor has failed to produce a valid reading for ≥ 12 consecutive reads (~60 seconds).</p> <p>Response: The display will show "Er03" only when the button sequence to read the heat exchange sensor is depressed.</p>
Er05	N/A	<p>Name: Display Firmware Integrity Failure</p> <p>Description: The display firmware has failed to pass its CRC CCITT checksum integrity test.</p> <p>Response: 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>Name: Micro Firmware Integrity Failure</p> <p>Description: The micro firmware has failed to pass its CRC CCITT checksum integrity test.</p> <p>Response: 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>Name: Micro Fail - CS5521 SPI Failure / UISR Failure</p> <p>Description: 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>Response: 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"> 1. Remote alarm contacts will become active. 2. Buzzer will annunciate audibly and will have a ringback of 15 minutes. 3. "Seven segment" display will show "Er07". 4. The system will have 10 minute staging between the high stage compressor and the low stage compressor activation. 5. The system will go to bottom out temperatures.

Error (cont.)	High End Message	Notes
Er09	N/A	<p>Name: Stuck Button</p> <p>Description: This condition indicates that the display board has a stuck button.</p> <p>Response: The Er09 will show on the display periodically.</p>
Er11	"COND PRB FLT"	<p>Name: Condenser Probe Sensor Failure</p> <p>Description: This condition indicates that the condenser probe sensor has failed to produce a valid reading for ≥ 12 consecutive reads (~60 seconds).</p> <p>Response: The display shows "Er11".</p>
N/A	"SMPL PRB FLT"	<p>Name: Sample Probe Sensor Failure</p> <p>Description: This condition indicates that the sample probe sensor has failed to produce a valid reading for ≥ 12 consecutive reads (~60 seconds).</p> <p>Response: The message center shows "SMPL PRB FLT".</p>
dErr	N/A	<p>This is a general display error in which the value being displayed can not be represented within the characters provided.</p>
(four dashes) ---- in display	N/A	<p>Name: Lost Communication</p> <p>Description: 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 5 Maintenance

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

Clean Cabinet Exterior

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.

Clean Air Filter (minimum 4 x/year*)

1. Open the front lower door by grasping the bottom left corner.
2. Locate the grille on the door. See Figure 1-5. 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. Order part number 760203.

Clean the Condenser (minimum yearly*)

1. Open the front lower door by grasping the bottom left corner. See Figure 1-5.
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.

Clean the Water-cooled Condenser

The water-cooled condenser can be cleaned-in-place by using the CIP procedure. Cleaning solutions can be used, depending on type of deposits or build-up to be removed.

Caution Do not use liquids that are corrosive to stainless steel or the brazing material (copper or nickel).

CIP (Clean-In-Place) Procedure

1. Disconnect the unit from the water supply.
2. Drain the unit.
3. Rinse with fresh water and drain the unit again.
4. Fill with fresh water.
5. Add cleaning agent (solution and concentration dependent on deposits or build-up).
6. Circulate cleaning solution (if feasible).
7. Drain the cleaning solution.
8. Add and circulate a passivating liquid for corrosion inhibition of plate surfaces.
9. Drain this liquid.
10. Rinse with fresh water and drain.
11. Reconnect the water supply and fill the unit.
12. Return to service.

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 5-8.
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 (⏻).
10. Allow the freezer to operate empty overnight before reloading product.

Clean the Door Gasket (minimum monthly*)

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 door from closing properly.

Vacuum Relief Port

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 buildup on the inner doors.

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. The unit is designed with a “vacuum relief port” that allows the pressure to be equalized.

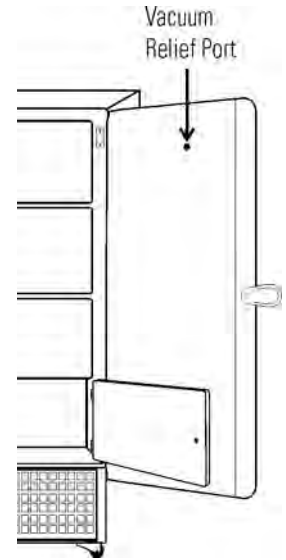


Figure 5-4. Location

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

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

Cabinets with the vacuum relief port operating normally, (i.e. vacuum relief port not iced over) requires 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 with door unlatched. The vacuum could release, resulting in a door opening and product loss. s

Vacuum Relief Port (continued)

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. Clear the vacuum relief tube completely free of ice to prevent rapid ice formation.

Factors that can affect the performance of the vacuum relief port include high ambient temperature, high humidity conditions and frequent door openings. Perform maintenance weekly, or as needed.

Note 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. s

Replace the Battery(s)

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 with three bolts. See Figure 5-2.
2. Directly above the battery(s) is the battery power switch. Turn the battery power switch to the Off position.
3. Disconnect the battery connections.
4. Remove the tape securing the battery.
5. 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.

Caution 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 Thermo. 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. s

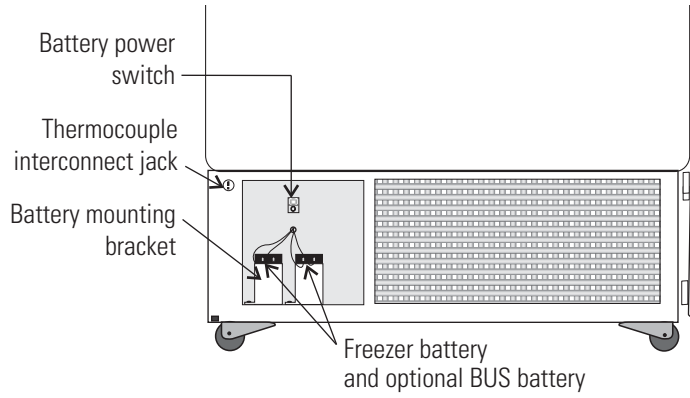


Figure 5-2. Battery Location

Prepare the Unit for Storage

Defrost the unit as described previously. This will prepare the unit for storage. Turn off the battery power switch. Turn off the freezer power switch. Disconnect power to the battery(s) and to the freezer.

PREVENTIVE MAINTENANCE

Freezers

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

The following is a condensed list of preventive maintenance requirements. See the specified section of the instruction 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 us at the number below.

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

Tips:

- Fill an upright by starting at the bottom near the probe and add racks to one shelf at a time. Allow freezer to recover to set point between shelves.
- Fill a chest by starting at the left side near the probe. Filling with room temperature racks will result in a long pull-down time.
- Fill unit with frozen product to help overall performance; frozen water jugs, for example.
- Always make certain the vacuum relief port is free of frost and ice, to allow for timely re-entry into the freezer after a door opening.

- Millcreek Road, Box 649 • Marietta, Ohio 45750 USA • 740-373-4763
- USA and Canada 888-213-1790 • Telefax: 740-373-4189 • email: services.controlenv@thermo.com

Preventive Maintenance for Model 8600 Series Freezers

Refer to Manual Section	Action	Monthly	Yearly	Every 2 Years
--	Verify ambient temperature, <90°F	<input checked="" type="checkbox"/>		
--	* Adjust door handle for firm latching, as needed	<input checked="" type="checkbox"/>		
Figure 1-4 for probe location	Check and clean probe cover, vacuum relief port, gaskets, hinges, and inner doors of ice and snow	<input checked="" type="checkbox"/>		
5		<i>More frequent cleaning may be required, depending on use and environmental conditions</i>		
5	Check air filter. Clean or replace as needed	<input checked="" type="checkbox"/>		
1, 3	Check alarm back-up battery	<input checked="" type="checkbox"/>		** Replace
--	Check condenser fan motor for unusual motor noise or vibration		<input checked="" type="checkbox"/>	
2	* Verify and document calibration, at the minimum, annually		<input checked="" type="checkbox"/>	
5	* Clean condenser compartment and wipe off condenser		<input checked="" type="checkbox"/>	

* Qualified service technicians only

** Dispose of properly, according to all state and federal regulations

To minimize ice build-up inside of freezer:

- Locate the freezer away from drafts or heating/cooling vents
- Keep the number of door openings to a minimum
- Minimize the length of time door is open
- Make sure door latches securely after opening

Section 6 Factory Installed Options

Details for the factory installed options available, or already installed, are listed below.

Back-Up System (BUS) - P/N 1950445, 1950447

Caution 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. s

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.

Install Injection Assembly, Vent Stack and Solenoid

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.

Injection Assembly, Vent Stack and Solenoid (cont.)

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 Permagem.
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 CO₂ supply cylinder, it must be equipped with a siphon tube. s

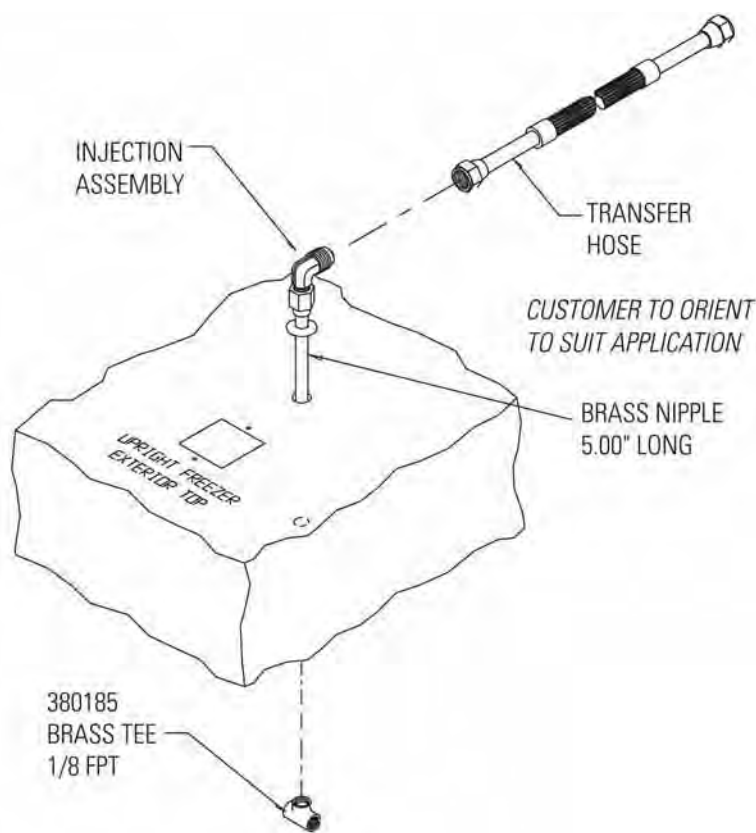


Figure 6-1. Injection Assembly

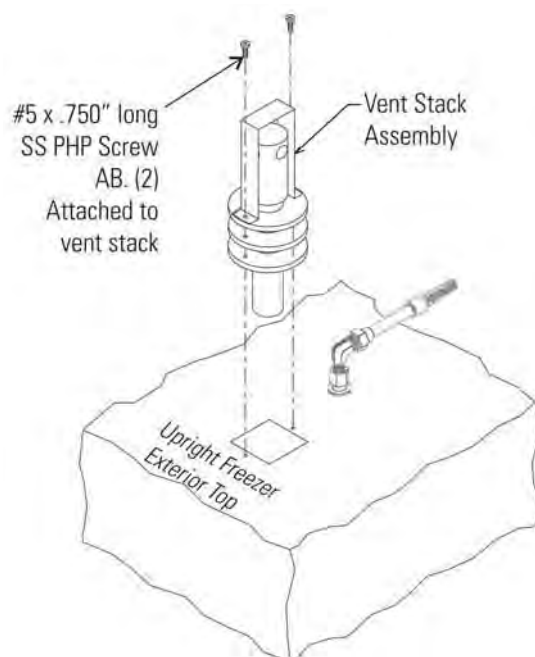


Figure 6-2. Vent Stack

Install the Temperature Probe

1. Locate the 0.500" pre-punched hole in the upper left hand 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 6-4).
2. As shown in Figure 6-3, 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 6-4 shows the back-up probe mounted on the interior left side wall of the freezer.

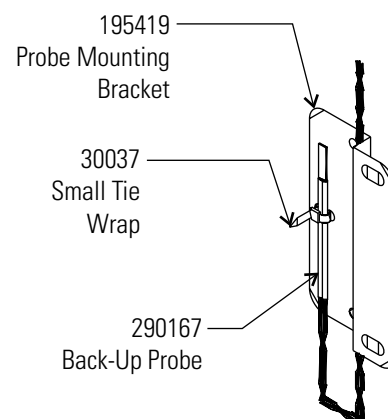


Figure 6-3. Secure Probe

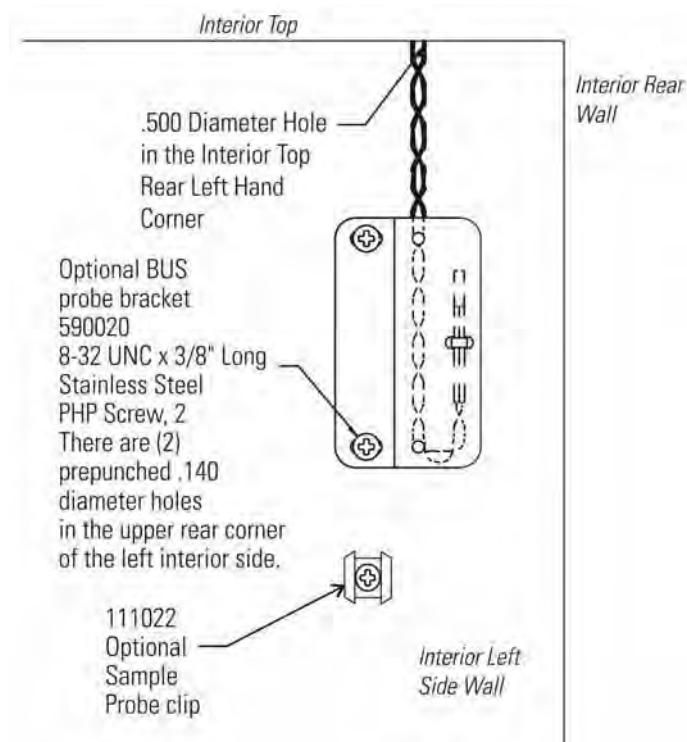


Figure 6-4. Mounted

Connect the Probe/Solenoid Harness

1. Remove the four screws on the freezer back panel and use them to mount the tie wrap anchors as shown in Figure 6-5. 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.

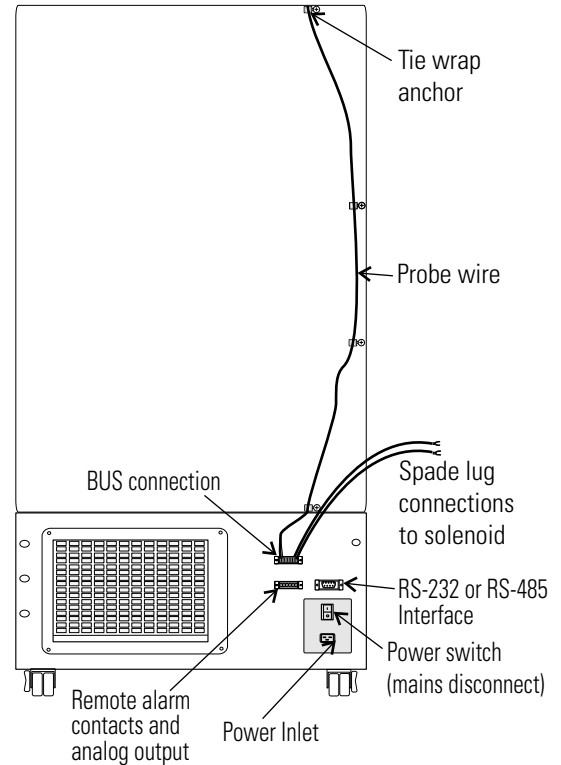


Figure 6-5. Routing

4. Connect power to the freezer. Turn the freezer On, with battery switch Off.
 - a. The Solenoid Engaged light on the BUS control panel will illuminate (no injection occurs). This light stays on until the unit is below BUS setpoint.
 - b. The Low Battery indicator may also illuminate.
5. Turn the battery switch to Standby mode (⏻) to charge both batteries.

BUS Control Panel

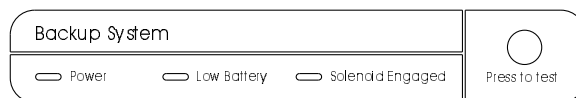


Figure 6-6. BUS Control Panel

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 suppresses oxygen levels and may cause suffocation if area is not well ventilated. Refer to Appendix A for the proper handling of liquid LN₂. s

Caution Make sure the pressure relief valve on any LN₂ tank is adjusted to 30 PSI maximum blow-off. s

Warning Carbon dioxide gas suppresses oxygen levels and may cause suffocation if area is not well ventilated. Refer to Appendix B for the proper handling of liquid CO₂. s

Power - indicates the unit has AC power.

Low Battery - battery charge is low. The battery needs replaced or recharged.

Solenoid Engaged - BUS has opened the solenoid so it can inject gas (CO₂ or LN₂).

Press-To-Test - Activates the solenoid and injects LN₂ or CO₂ into the freezer chamber as long as the button is depressed. The solenoid engaged indicator should light. If the Low Battery indicator lights during the test, replace the BUS battery.

Note Solenoid will not engage if door is open. s

Set the Optional BUS Set Point

The optional back up system is designed to inject CO₂ or LN₂ 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 the next or previous parameter.

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

Caution Changing the operating temperature set point can affect the BUS set point. The BUS set point will self-adjust to maintain a temperature of at least 10°C above the operating temperature set point. s

Caution The BUS set point cannot be set colder than the high temperature alarm set point. (See Section 1.) If the back-up system is installed with CO₂, then -65°C is the coldest BUS set point that can be used (if cabinet set point is -75°C or colder). s

Test the BUS

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 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.

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 & 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

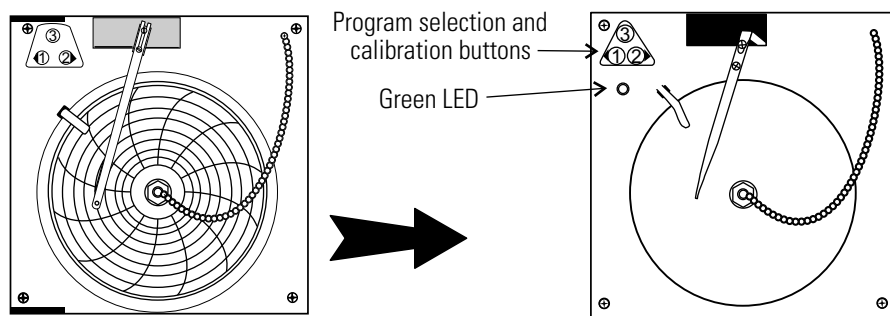


Figure 6-7. Recorder Buttons Location

Install the Chart Paper

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 6-8.
Buttons

Change Recorder Temperature Range

The chart recorder contains eight temperature ranges and is factory-programmed for the freezer.

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.

Program	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

Calibrate the Chart Recorder

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 1-4).
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 may be purchased. Refer to exploded parts drawings. s

Datalogger

Dataloggers and ELPRO evaluation software provide monitoring and documentation of temperature and alarm conditions. The dataloggers have a memory capacity of 64,000 measured values or data points. Temperature is measured, stored and displayed. Alarm conditions are recorded. Optional evaluation software permits data to be downloaded to a PC. A variety of statistical information is provided through calculations, analysis, graphs and printed reports. Refer to the ELPRO documentation for operating instructions for the datalogger.

Water-cooled Condenser

The water-cooled condenser (P/N 195964 [13cf], 195965 [17, 23, 28 cf], 195967 [12, 17, 20 cf chests]) is a factory installed option and requires a qualified technician at freezer installation. The installation should include proper adjustment of the regulating valve, which controls the discharge pressure. Specifications for this option are displayed in Table 6-1.

Table 6-1. Specifications

Water Source	Tower	City
Water Pressure	Not to exceed 150 psig	
Water Temperature Range	Not to exceed 29.4C (85F)	
Inlet Connection	0.5" compression	
Outlet Connection	0.5" compression	
Flow Rate Required	3.0 gallons (11.4 liters) per minute	1.0 gallon (3.8 liters) per minute
Drain Required	No (return line is required)	Yes

Five Inner Door Option

The five inner door option (P/N 189405 [13cf], 189406 [17cf], 189407 [23cf] & 195652 [28cf]) is factory installed. The freezer is converted to accommodate four adjustable specimen shelves with the fifth "shelf" as the bottom of the freezer chamber.

Section 7 Specifications

Model	8602	8603	8604	8605	8606
Temperature Range	-50°C (-58°F) to -86°C (-123°F) in an 18C to 32C* (64.4F to 89.6F) ambient				
Exterior Dimensions	33.3"W x 77.8"H x 31.0" 84.6 x 197.6 x 78.7 cm	33.3"W x 77.8"H x 31.0" 84.6 x 197.6 x 78.7 cm	33.3"W x 77.8"H x 37.0" 84.6 x 197.6 x 94.0 cm	33.3"W x 77.8"H x 37.0" 84.6 x 197.6 x 94.0 cm	40.8"W x 77.8"H x 37.0" 103.6 x 197.6 x 94.0 cm
Interior Dimensions	23.0"W x 51.5"H x 19.3" 58.4 x 130.8 x 49.0 cm	23.0"W x 51.5"H x 19.3" 58.4 x 130.8 x 49.0 cm	23.0"W x 51.5"H x 25.3" 58.4 x 130.8 x 64.3 cm	23.0"W x 51.5"H x 25.3" 58.4 x 130.8 x 64.3 cm	30.6"W x 51.5"H x 25.3" 77.7 x 130.8 x 64.3 cm
Capacity	13.0 cu. ft. (368.1 liters)	13.0 cu. ft. (368.1 liters)	17.3 cu. ft. (489.9 liters)	17.3 cu. ft. (489.9 liters)	23.0 cu. ft. (651.3 liters)
Refrigeration	Two 1 HP (2545 BTUH each)				
Insulation	Non-CFC, foamed-in-place urethane: 5.0" (12.7 cm) cabinet; 4.5" (11.4 cm) door				
Electrical	230V, 50/60Hz, 12.0 FLA Operating Range: 208VAC-240VAC	120V, 60 Hz, 16.0 FLA Operating Range: 108-130V	120V, 60 Hz, 16.0 FLA Operating Range: 108-130V	230V, 50/60Hz, 12.0 FLA Operating Range: 208VAC-240VAC	230V, 50/60Hz, 12.0 FLA Operating Range: 208VAC-240VAC
Breaker Requirements	15 amp, 230V, Dedicated Circuit, 15 Amp Time Delay Breaker	20 amp, 120V, Dedicated Circuit, 20 amp Time Delay Breaker	20 amp, 120V, Dedicated Circuit, 20 amp Time Delay Breaker	15 amp, 230V, Dedicated Circuit, 15 Amp Time Delay Breaker	15 amp, 230V, Dedicated Circuit, 15 Amp Time Delay Breaker
Shipping Weight Motor	712 lbs. (323.0 kg)	712 lbs. (323.0 kg)	795 lbs. (360.6 kg)	795 lbs. (360.6 kg)	900 lbs. (408.2 kg)

* Compressors may not cycle off with cabinet running at -86C in a 32C ambient.

Section 7
Specifications

Model	8607	8656	8690	8691	8692
Temperature Range	-50°C (-58°F) to -86°C (-123°F) in an 18C to 32C* (64.4F to 89.6F) ambient				
Exterior Dimensions	46.8"W x 77.8"H x 37.0" 118.9x197.6x94.0cm	40.8"W x 77.8"H x 37.0" 103.6x197.6x94.0cm	40.8"W x 77.8"H x 37.0" 103.6 x 197.6 x 94.0 cm	33.3"W x 77.8"H x 31.0" 84.6 x 197.6 x 78.7 cm	33.3"W x 77.8"H x 31.0" 84.6 x 197.6 x 78.7 cm
Interior Dimensions	36.6"W x 51.5"H x 27.0" 93.0x130.8x68.6cm	30.6"W x 51.5"H x 25.3" 77.7x130.8x64.3cm	30.6"W x 51.5"H x 25.3" 77.7 x 130.8 x 64.3 cm	23.0"W x 51.5"H x 19.3" 58.4 x 130.8 x 49.0 cm	23.0"W x 51.5"H x 19.3" 58.4 x 130.8 x 49.0 cm
Capacity	28.0 cu. ft. (792.8 liters)	23.0 cu. ft. (651.3 liters)	23.0 cu. ft. (651.3 liters)	13.0 cu. ft. (368.1 liters)	13.0 cu. ft. (368.1 liters)
Refrigeration	Two 1 HP (2545 BTUH each)				
Insulation	Non-CFC, foamed-in-place urethane: 5.0" (12.7 cm) cabinet; 4.5" (11.4 cm) door				
Electrical	230V,50/60Hz, 12.0FLA Operating Range: 208VAC-240VAC	120V, 60 Hz, 16.0 FLA Operating Range: 108-130V	120V, 60 Hz, 16.0 FLA Operating Range: 108-130V	230V,50/60Hz, 12.0FLA Operating Range: 208VAC-240VAC	120V, 60 Hz, 16.0 FLA Operating Range: 108-130V
Breaker Requirements	15 amp, 230V, Dedicated Circuit, 15 Amp Time Delay Breaker	20 amp, 120V, Dedicated Circuit, 20 amp Time Delay Breaker	20 amp, 120V, Dedicated Circuit, 20 amp Time Delay Breaker	15 amp, 230V, Dedicated Circuit, 15 Amp Time Delay Breaker	20 amp, 120V, Dedicated Circuit, 20 amp Time Delay Breaker
Shipping Weight Motor	980 lbs. (444.5 kg)	900 lbs. (408.2 kg)	900 lbs. (408.2 kg)	712 lbs. (323.0 kg)	712 lbs. (323.0 kg)

Model	8693	8694	8695
Temperature Range	-50°C (-58°F) to -86°C (-123°F) in an 18C to 32C* (64.4F to 89.6F) ambient		
Exterior Dimensions	33.3"W x 77.8"H x 37.0" 84.6 x 197.6 x 94.0 cm	33.3"W x 77.8"H x 37.0" 84.6 x 197.6 x 94.0 cm	40.8"W x 77.8"H x 37.0" 103.6 x 197.6 x 94.0 cm
Interior Dimensions	23.0"W x 51.5"H x 25.3" 58.4 x 130.8 x 64.3 cm	23.0"W x 51.5"H x 25.3" 58.4 x 130.8 x 64.3 cm	30.6"W x 51.5"H x 25.3" 77.7 x 130.8 x 64.3 cm
Capacity	17.3 cu. ft. (489.9 liters)	17.3 cu. ft. (489.9 liters)	23.0 cu. ft. (651.3 liters)
Refrigeration	Two 1 HP (2545 BTUH each)		
Insulation	Non-CFC, foamed-in-place urethane: 5.0" (12.7 cm) cabinet; 4.5" (11.4 cm) door		
Electrical	120V, 60 Hz, 16.0 FLA Operating Range: 108-130VAC	230V, 50/60Hz, 12.0FLA Operating Range: 208VAC-240VAC	230V, 50/60Hz, 12.0FLA Operating Range: 208VAC-240VAC
Breaker Requirements	20 amp, 120V, Dedicated Circuit, 20 amp Time Delay Breaker	15 amp, 230V, Dedicated Circuit, 15 Amp Time Delay Breaker	15 amp, 230V, Dedicated Circuit, 15 Amp Time Delay Breaker
Shipping Weight Motor	795 lbs. (360.6 kg)	795 lbs. (360.6 kg)	900 lbs. (408.2 kg)

* Compressors may not cycle off with cabinet running at -86C in a 32C ambient.

Certifications

Declaration of Conformity is available from the factory

Safety Specifications

Indoor Use Only

Altitude - up to 2,000 meters

Temperature - 5°C to 43°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 $\pm 10\%$ of the nominal voltage

Installation Category II ¹

Pollution Degree 2 ²

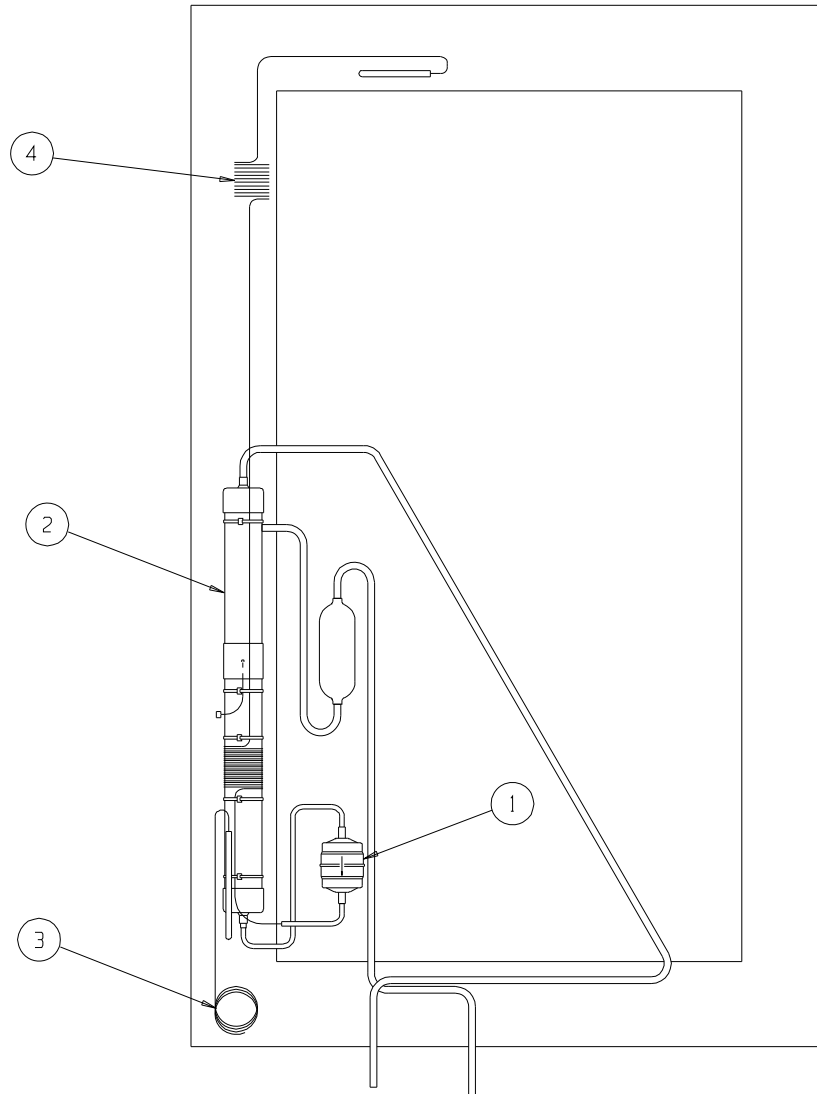
Class of Equipment I

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.

REV	ECN NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION
0	N/A	12-04-02	PDK	PDK	LDN	RELEASED FOR PRODUCTION
1	FR-1943	06-13-07	CRM	KDG	LDN	CHG. 211039/41 HEAT EXCHG. TO 211050
2	FR-2055	07-29-09	NWM	KDG	CCS	CHANGED 209016 DRYER TO 209017

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



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, MARIETTA, OHIO 45750

MODEL/PART NAME: 8600 UP-RIGHT FREEZER

DWG TITLE: HEAT EXCHANGER ASSEMBLY

DWN: PDK CAD: PDK APPD: LDN DATE: 12-04-02 SCALE: NTS

MATERIAL: N/A

PAINT COLOR: N/A

TOLERANCE UNLESS OTHERWISE SPECIFIED

ANGLES: DECIMAL: .XX=±
.XXX=±

DRAWING NUMBER

8602-205-1

SIZE

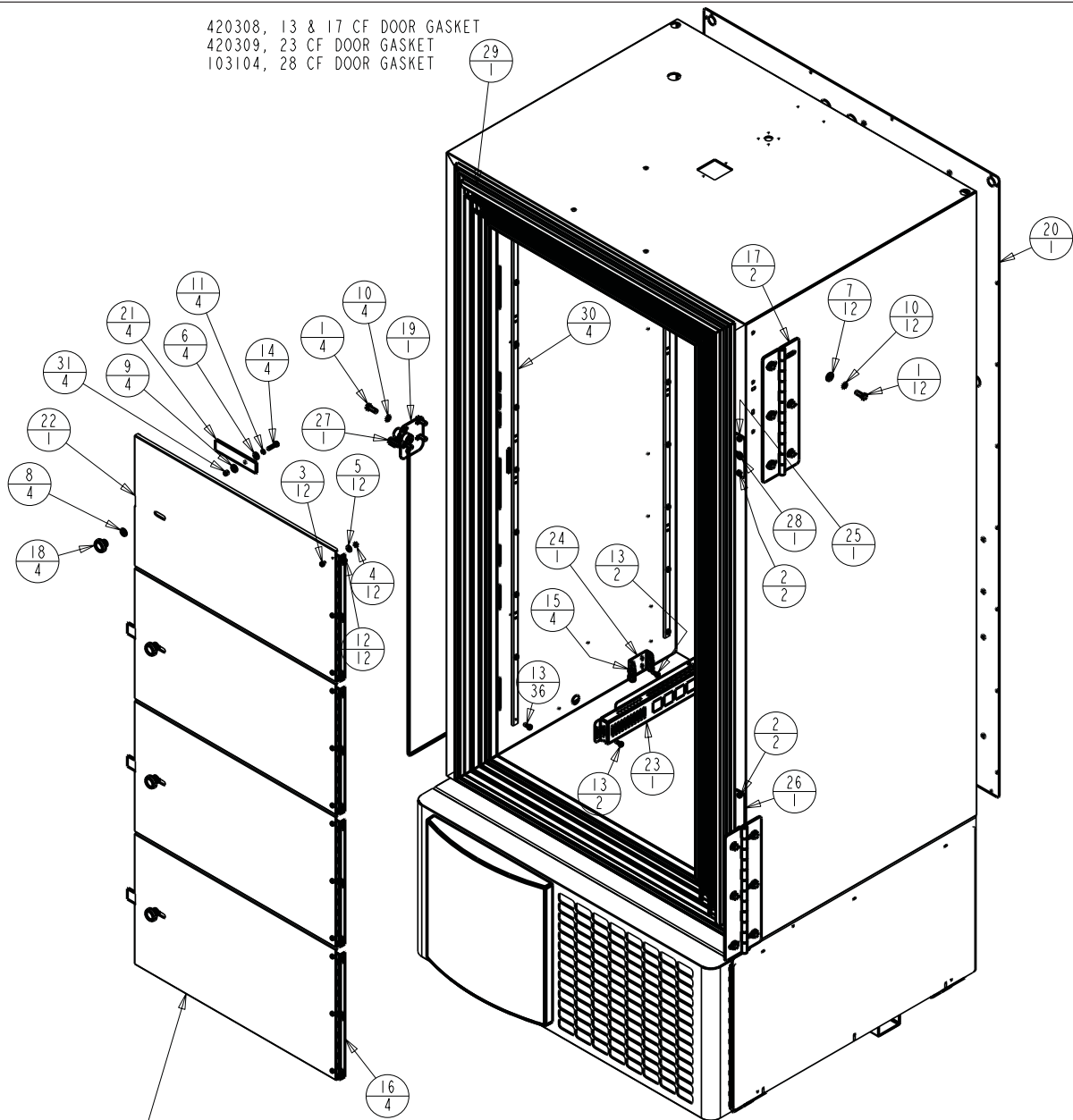
B

Exploded Drawing
for
Upright Freezer
Heat Exchanger Assy

8602-205-1-D REV.2

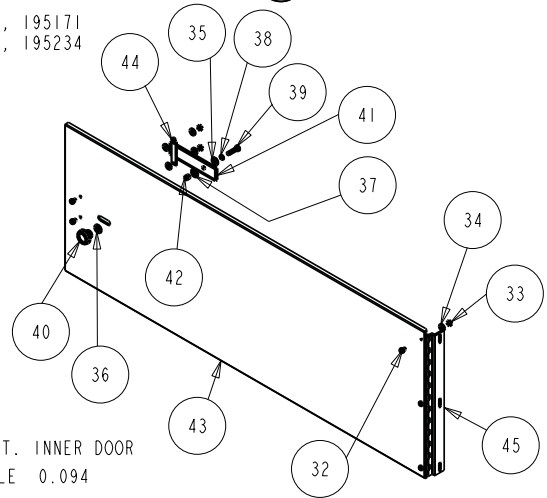
Page 1 of 1

420308, 13 & 17 CF DOOR GASKET
 420309, 23 CF DOOR GASKET
 103104, 28 CF DOOR GASKET



116077, USE ON INNER DOOR 195170, 195171
 116069, USE ON INNER DOOR 195233, 195234

195170, 13 & 17 CF INNER DOOR "4 DOOR"
 195171, 23 CF INNER DOOR "4 DOOR"
 195233, 13 & 17 CF INNER DOOR "5 DOOR"
 195234, 23 CF INNER DOOR "5 DOOR"



28 CU. FT. INNER DOOR
 SCALE 0.094

Exploded Parts
 Drawing
 Model 8600 Series
 Upright Freezer
 8602-200-1-B Rev. 5
 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	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	59008	#8-32 X 7/8 SS PHP SCREW
15	114020	5/8" X 1/2" ID GROMMET
16	116077	FRONT PANEL HINGE
17	116092	EXTERIOR FREEZER DOOR HINGE
18	120400	BLACK PLASTIC KNOB
19	121069	FREEZER CAM LATCH STRIKE
20	189921	EXTERIOR BACK 13 & 17
21	195169	LATCH TAB
22	195170	13/17 CU. FT. INNER DOOR
23	195866	PROBE GUARD
24	195867	PROBE MOUNT
25	195874	CABINET CABLE COVER PLATE
26	195879	CABINET CABLE BLANK COVER PLATE
27	195900	SINGLE DOOR SWITCH ASSY.
28	330010	1/2" SPLIT SNAP BUSHING
29	420308	13 & 17 CU. FT. SINGLE DOOR FRAME GASKET
30	500177	PILSATER 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

REV	ECN NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION
3	FR-1789	06-02-05	DHG	DHG	LDN	REMOVED VRP IN TOP OF CABINET, DMHVRP
4	FR-2055	05-15-09	GJG	SAG	LDN	CHANGED 28 CU. FT. INNER DOOR
5	SI-10308	08-11-10	CAC	LDC	CCS	285659 KNOB TO 120400

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

ThermoFisher
SCIENTIFIC

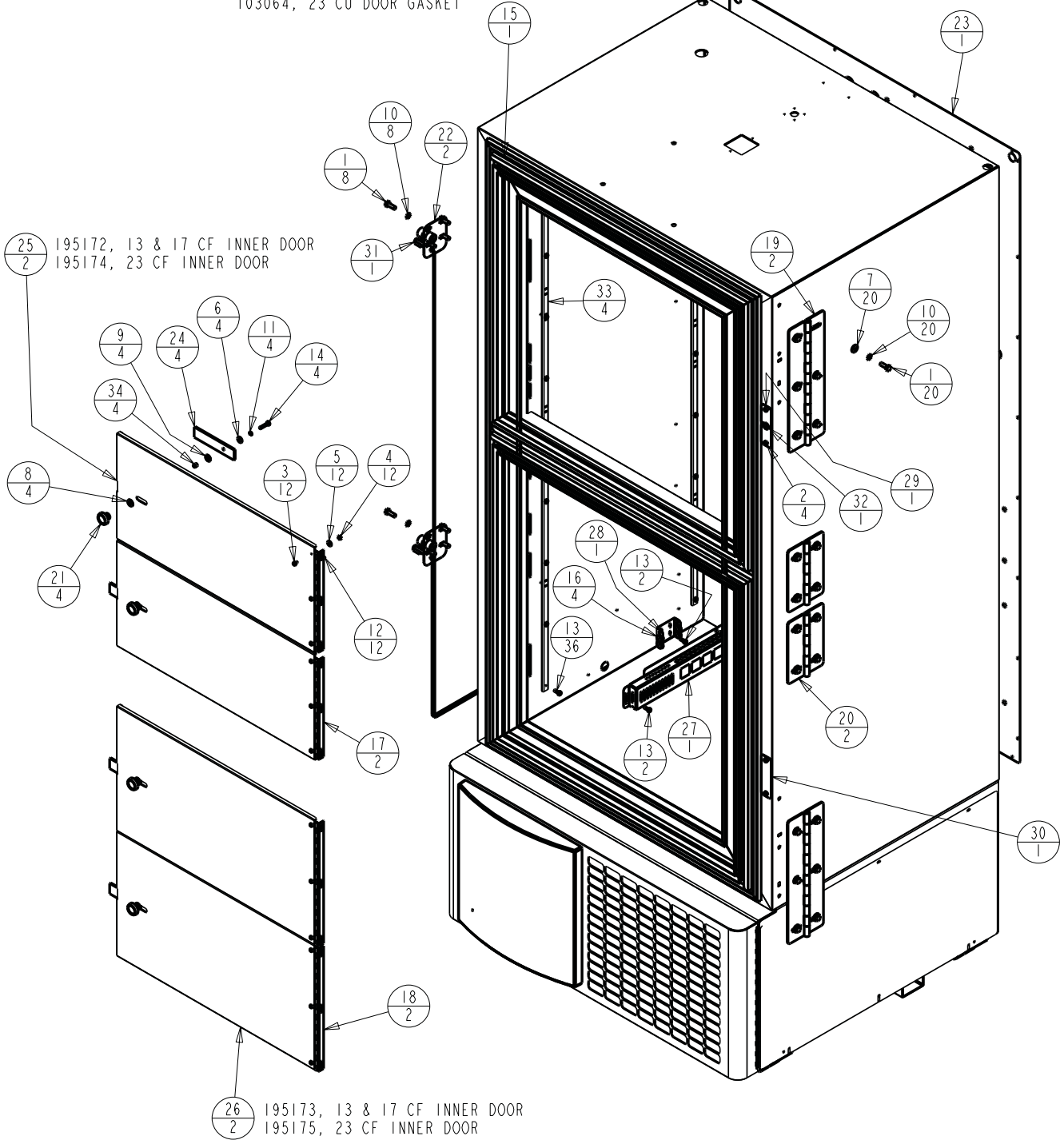
BOX 649, MARIETTA, OHIO 45750

MODEL/PART NAME: 8600 UP-RIGHT FREEZER					
DWG TITLE: 8602 UP-RIGHT FREEZER ASSEMBLY					
DWN: PDK	CAD: PDK	APPD: MAH	DATE: 10-30-02	SCALE: 0.094	
MATERIAL: N/A					
PAINT: N/A					
TOLERANCE UNLESS OTHERWISE SPECIFIED			DRAWING NUMBER	SIZE	
ANGLES: DECIMAL: .XX± .xxx±			8602-200-1	B	

**Exploded Parts
Drawing
Model 8600 Series
Upright Freezer**

**8602-200-1-B Rev. 5
Page 2 of 2**

103063, 13 & 17 CF DOOR GASKET
 103064, 23 CU DOOR GASKET



Exploded Parts
 Drawing
 Model 8600 Series
 Upright Freezer
 8602-200-2-D Rev. 6
 Page 1 of 2

REV	ECN NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION
4	SI-9595	05-15-06	DHG	KDG	LDN	CHG. TO COMMON SWITCH ASSY 195900
5	SI-10308	08-12-10	CAC	LDC	LDN	285659 KNOB TO 120400
6	FR-1999	09-01-10	CAS	KDG	CCS	REMOVED 180312 PLASTIC, OBS. 4/08

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	59008	#8-32 X 7/8 SS PHP SCREW
15	103063	DOUBLE DOOR FRAME GASKET
16	114020	5/8" X 1/2" ID GROMMET
17	116069	FRONT PANEL HINGE
18	116077	FRONT PANEL HINGE
19	116092	EXTERIOR FREEZER DOOR HINGE
20	116093	EXTERIOR FREEZER DOOR HINGE
21	120400	BLACK PLASTIC KNOB
22	121069	FREEZER CAM LATCH STRIKE
23	189921	EXTERIOR BACK 13 & 17
24	195169	LATCH TAB
25	195172	13/17 CU. FT. INNER DOOR, TOP
26	195173	23 CU. FT. INNER DOOR, TOP
27	195866	PROBE GUARD
28	195867	PROBE MOUNT
29	195874	CABINET CABLE COVER PLATE
30	195879	CABINET CABLE BLANK COVER PLATE
31	195900	TOP DOOR SWITCH ASSY.
32	330010	1/2" SPLIT SNAP BUSHING
33	500177	PILSATER STRIPS
34	515083	1/4 DIA. X 1/4L SS SPACER

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

MODEL/PART NAME: 8600 UP-RIGHT FREEZER				
DWG TITLE: 8691 UP-RIGHT FREEZER ASSEMBLY				
DWN: PDK	CAD: PDK	APPD: MAH	DATE: 10-30-02	SCALE: 0.094
MATERIAL: N/A				
PAINT: N/A				
TOLERANCE UNLESS OTHERWISE SPECIFIED			DRAWING NUMBER	SIZE
ANGLES: DECIMAL: .XX± .xxx±			8602-200-2	B

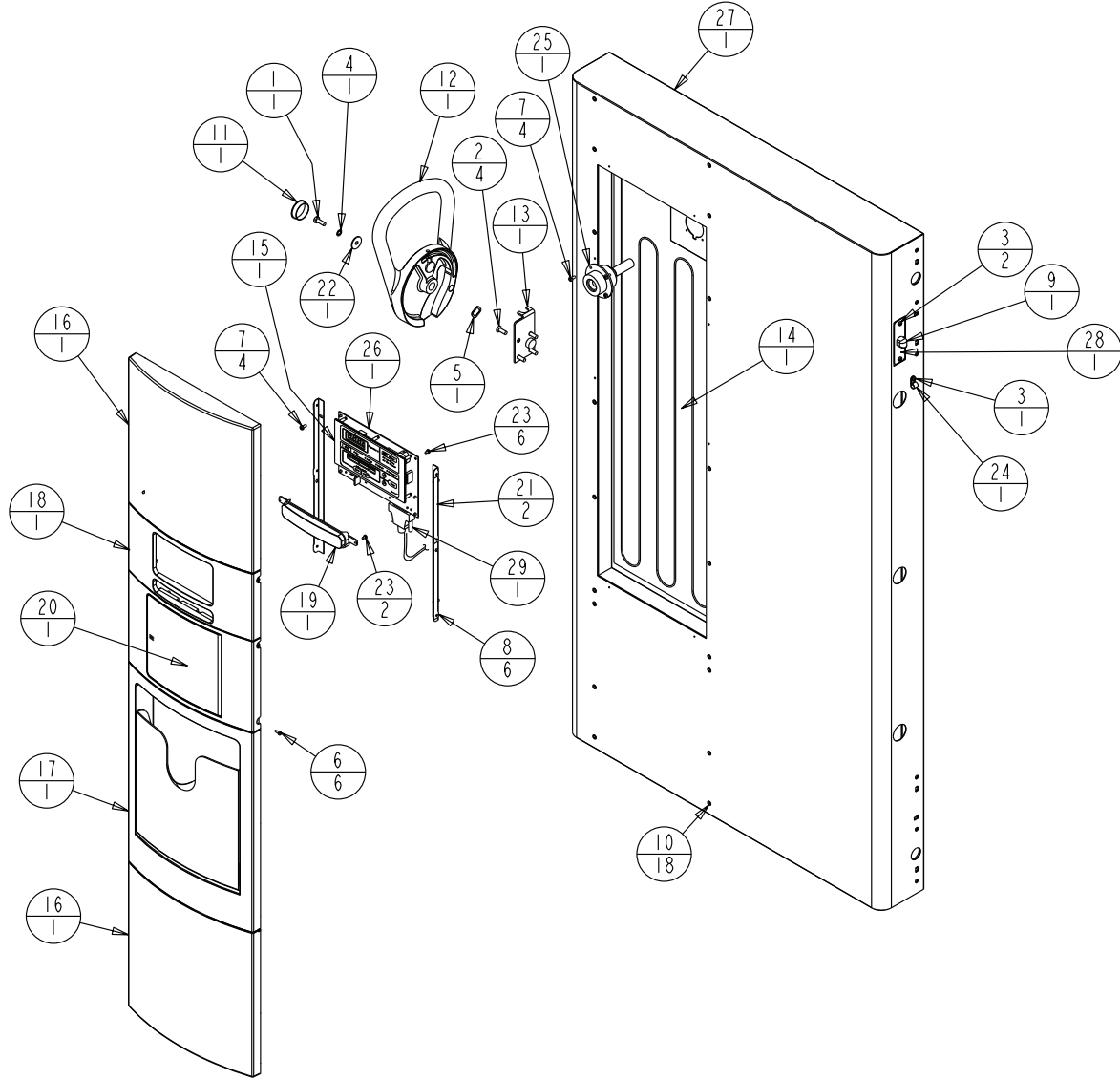
ThermoFisher
SCIENTIFIC

BOX 649, MARIETTA, OHIO 45750

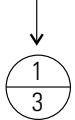
Exploded Parts
Drawing
Model 8600 Series
Upright Freezer

8602-200-2-D Rev. 6
Page 2 of 2

195771, 23 CU. FT. UPRIGHT FREEZER DOOR "8600"
 195772, 28 CU. FT. UPRIGHT FREEZER DOOR "8600"



Part Number



Quantity


Upright Freezer
 Door
 Assembly
 Single Door

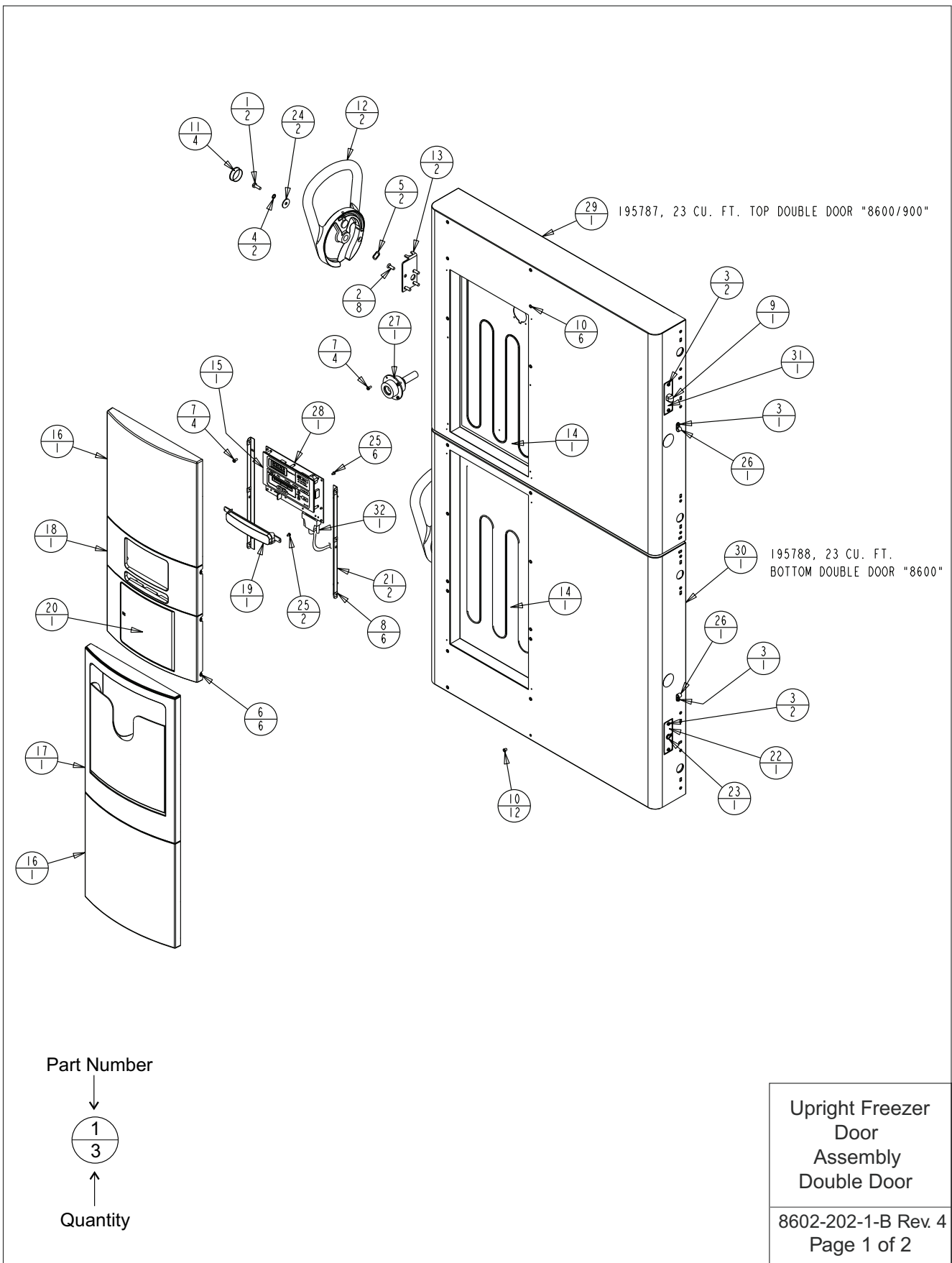
8602-201-1-B Rev. 4
 Page 1 of 2

REV	ECN NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION
0	N/A	10-30-02	PDK	PDK	MAH	RELEASED FOR PRODUCTION
1	FR-1673	06-06-03	DHG	PDK	LDN	REVISED CONTROL PANEL FASTENERS
2	FR-1776	03-08-04	AT	pdk	LDN	CORRECTED CONTROL PANEL STK NUMBER
3	FR-1789	11-18-04	RSB	KDG	LDN	ADDED VACUUM RELIEF PORT
4	FR-2145	12-28-10	HCE	KDG	CCS	DISPLAY 191802 TO 191941, ROHS ENCL.

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	111028	TINNERMAN TUBULAR SPEED CLIP
11	117038	1-3/8" DIA. THERMO WHITE HOLE PLUG
12	121068	121068 FINISHED HANDLE/LATCH ASSEMBLY
13	121075	CAM LATCH MOUNT
14	132115	HEATER, 5W, 14VDC
15	140367	CONTROL PANEL ASSEMBLY
16	180301	THERMO CONTROL CENTER BLANK PANEL
17	180304	THERMO DATA STORAGE POCKET ASSEMBLY
18	180305	CONTROL CENTER DISPLAY BEZEL
19	180306	THERMO BACK-UP SYSTEM BLANK PANEL
20	180308	THERMO CONTROL CENTER RECORDER BLANK
21	195837	MOUNTING ANGLE FOR 180305
22	510305	1" OD FLAT WASHER
23	590027	#6-32 X 1/4 SS PHP EXT SEMS SCREW
24	600085	5/16 NYLON CABLE CLAMP
25	1950069	HEATED VACUUM RELIEF PORT
26	191941_	FREEZER DISPLAY BOARD
27	195770	13 & 17 CU. FT. UPRIGHT FREEZER DOOR "8600"
28	195830	UPRIGHT DOOR WIREWAY COVER PLATE
29	430336	15 FT, RS-232 CABLE 25 POS.

<p>THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSE WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC</p>  <p>BOX 649, MARIETTA, OHIO 45750</p>	MODEL/PART NAME: ULT UP-RIGHT SERIES FREEZER					<p>Upright Freezer Door Assembly Single Door</p>
	DWG TITLE: 8600 SINGLE DOOR BOM ASSEMBLY					
	DWN: PDK	CAD: PDK	APPD: MAH	DATE: 10-30-02	SCALE: 0.094	
	MATERIAL: N/A					
PAINT: N/A					<p>8602-201-1-B Rev. 4 Page 2 of 2</p>	
TOLERANCE UNLESS OTHERWISE SPECIFIED		DRAWING NUMBER		SIZE		
ANGLES: DECIMAL: .XX±		8602-201-1		B		
		.xxx±				



REV	ECN NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION
0	N/A	11-05-02	PDK	PDK	MAH	RELEASED FOR PRODUCTION
1	FR-1673	06-06-03	DHG	PDK	LDN	REVISED CONTROL PANEL FASTENERS
2	FR-1776	03-09-04	AT	pdk	LDN	CORRECTED CONTROL PANEL STK NUMBER
3	FR-1789	11-19-04	RSB	KDG	LDN	ADDED VACUUM RELIEF PORT
4	FR-2145	12-28-10	HCE	KDG	CCS	DISPLAY 191802 TO 191941, ROHS ENCL.

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	111028	TINNERMAN TUBULAR SPEED CLIP
11	117038	1-3/8" DIA. THERMO WHITE HOLE PLUG
12	121068	121068 FINISHED HANDLE/LATCH ASSEMBLY
13	121075	CAM LATCH MOUNT
14	132114	HEATER, 3W, 14VDC
15	140367	CONTROL PANEL ASSEMBLY
16	180301	THERMO CONTROL CENTER BLANK PANEL
17	180304	THERMO DATA STORAGE POCKET ASSEMBLY
18	180305	CONTROL CENTER DISPLAY BEZEL
19	180306	THERMO BACK-UP SYSTEM BLANK PANEL
20	180308	THERMO CONTROL CENTER RECORDER BLANK
21	195837	MOUNTING ANGLE FOR 180305
22	195874	CABINET CABLE COVER PLATE
23	330010	1/2" SPLIT SNAP BUSHING
24	510305	1" OD FLAT WASHER
25	590027	#6-32 X 1/4 SS PHP EXT SEMS SCREW
26	600085	5/16 NYLON CABLE CLAMP
27	1950069	HEATED VACUUM RELIEF PORT
28	191941_	FREEZER DISPLAY BOARD
29	195785	13/17 CF TOP DOUBLE DOOR "8600/900"
30	195786	13/17 CF BOTTOM DOUBLE DOOR "8600"
31	195830	UPRIGHT DOOR WIREWAY COVER PLATE
32	430336	15 FT, RS-232 CABLE 25 POS.

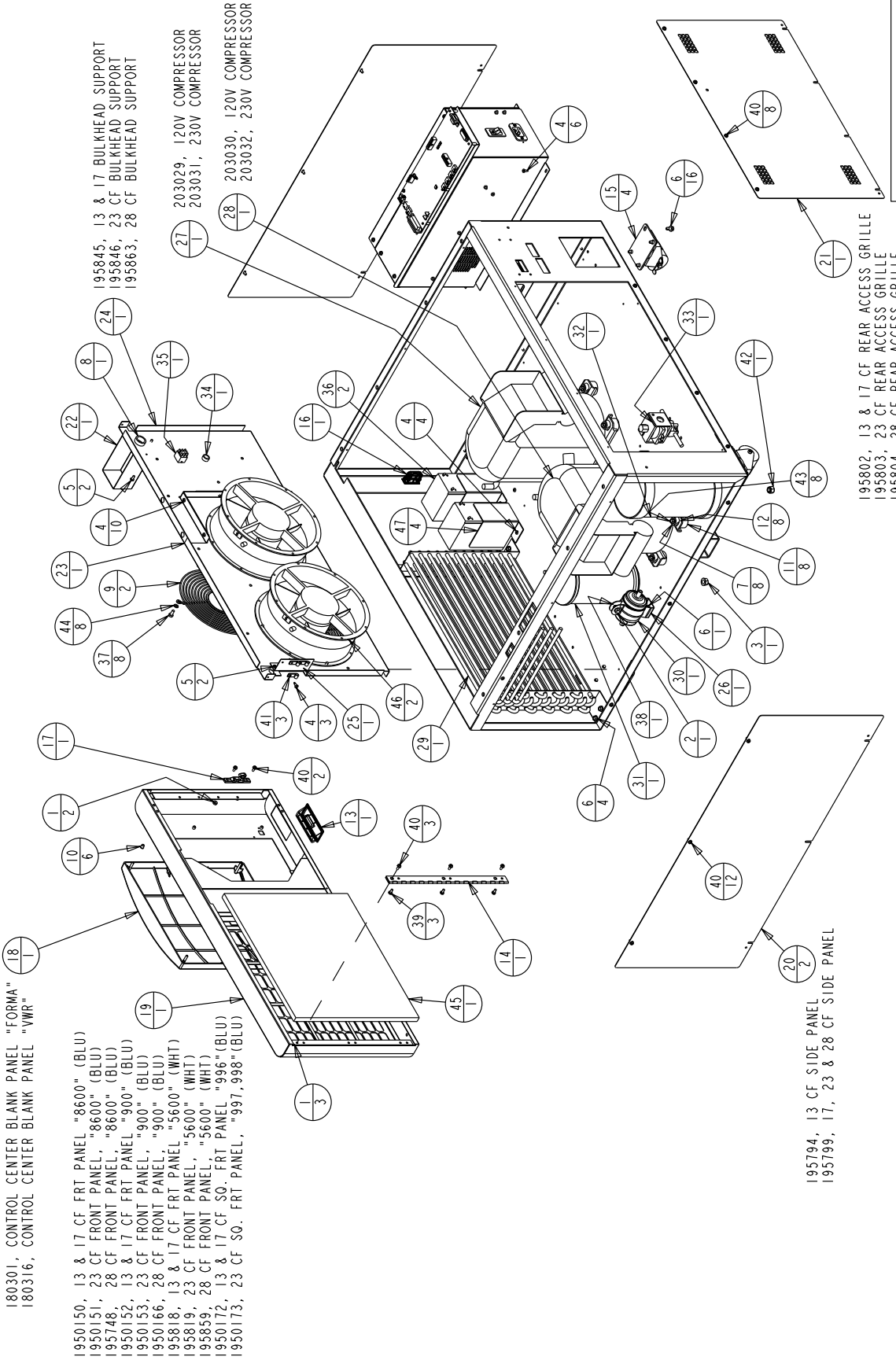
<p>THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSE WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC</p> <p>ThermoFisher SCIENTIFIC</p> <p>BOX 649, MARIETTA, OHIO 45750</p>	MODEL/PART NAME: ULT UP-RIGHT SERIES FREEZER					<p>Upright Freezer Door Assembly Double Door</p> <p>8602-202-1-B Rev. 4 Page 2 of 2</p>
	DWG TITLE: 8600 DOUBLE DOOR BOM ASSEMBLY					
	DWN: PDK	CAD: PDK	APPD: MAH	DATE: 11-05-02	SCALE: 0.094	
	MATERIAL: N/A					
	PAINT: N/A					
TOLERANCE UNLESS OTHERWISE SPECIFIED			DRAWING NUMBER		SIZE	
ANGLES: DECIMAL: .XX=± .xxx=±			8602-202-1		B	

180301, CONTROL CENTER BLANK PANEL "FORMA"
180316, CONTROL CENTER BLANK PANEL "VWR"

1950150, 13 & 17 CF FRT PANEL "8600" (BLU)
1950151, 23 CF FRONT PANEL, "8600" (BLU)
195748, 28 CF FRONT PANEL, "8600" (BLU)
1950152, 13 & 17 CF FRT PANEL "900" (BLU)
1950153, 23 CF FRONT PANEL, "900" (BLU)
1950166, 28 CF FRONT PANEL, "900" (BLU)
195818, 13 & 17 CF FRT PANEL "5600" (WHT)
195819, 23 CF FRONT PANEL, "5600" (WHT)
195859, 28 CF FRONT PANEL, "5600" (WHT)
1950172, 13 & 17 CF SO. FRT PANEL "996" (BLU)
1950173, 23 CF SO. FRT PANEL, "997,998" (BLU)

195845, 13 & 17 BULKHEAD SUPPORT
195846, 23 CF BULKHEAD SUPPORT
195863, 28 CF BULKHEAD SUPPORT

203029, 120V COMPRESSOR
203031, 230V COMPRESSOR
203030, 120V COMPRESSOR
203032, 230V COMPRESSOR



195794, 13 CF SIDE PANEL
195799, 17, 23 & 28 CF SIDE PANEL

195802, 13 & 17 CF REAR ACCESS GRILLE
195803, 23 CF REAR ACCESS GRILLE
195804, 28 CF REAR ACCESS GRILLE

Assembly Drawing
Upright Freezer
Base
8602-203-1-D Rev. 5
Page 1 of 2

REV	ECM NO.	DATE	BY	CAD APPD	DESCRIPTION OF REVISION
3	SI-9962	03-26-08	KDG	LDN	CHG. FORMA FRONT PANEL TO BLUE/VBL
4	FR-2004	07-01-08	KDG	LDN	ADDED FRONT PANEL FOR 5607
5	FR-2004	07-28-08	KDG	CCS	ADDED FRONT PANEL FOR 996.997.998

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	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	24049 1/4 ZP FLAT WASHER
8	30016 1" SNAP BUSHING
9	108020 10" WIRE FAN GUARD
10	111028 TINNERMAN TUBULAR SPEED CLIP
11	114033 COMPRESSOR MOUNTING FOOT
12	114034 COMPRESSOR MOUNTING SLEEVE
13	115032 BLACK ABS PLASTIC PULL
14	116115 FRONT PANEL HINGE
15	120011 DUAL WHEEL CASTER
16	121071 LATCH CATCH, PART OF 121071 ASSEMBLY
17	121071 LATCH KEEPER, PART OF 121071 ASSEMBLY
18	180301 THERMO CONTROL CENTER BLANK PANEL
19	195746 13&17 CU. FT. UR FRZ BASE FRONT PNL
20	195799 SIDE PANEL 17, 23 AND 28 CU. FT. UPRIGHT
21	195802 13/17 REAR ACCESS GRILLE
22	195829 MULLION/DOOR SWITCH WIRE COVER
23	195844 UR FRZ FAN BULKHEAD
24	195845 13&17 CU. FT. FAN BULKHEAD SUPPORT
25	195882 REFRIGERATION LINE SUPPORT BRACKET
26	200126 2" RIGID HANGER
27	203031 230V HIGH STAGE COMPRESSOR
28	203032 230V LOW STAGE COMPRESSOR
29	204009 REFRIGERATION CONDENSER
30	209020 LIQUID LINE FILTER DRYER WITH ACCESS PORT
31	214006 OIL SEPARATOR
32	214018 10.000" H X 5.000" DIA. EXPANSION TANK

BILL OF MATERIALS	
ITEM NO.	PART NO. PART DESCRIPTION
33	220626 120V - 50/60 HZ SOLENOID VALVE
34	330002 5/8" SNAP BUSHING
35	360248 MINI SNAP-IN POWER SWITCH
36	400159 SEALED LEAD ACID BATTERY - 12 VOLT - 7.2 Ah
37	510035 #12-24 X 1/2 SS HH CAP SCREW
38	550043 1/4-20 X 1" L ZP CARRIAGE BOLT
39	590020 #8-32 X 3/8 SS PHP EXT SEMS SCREW
40	590029 #8-32 X 3/8 SS PHP EXT SEMS SCREW W/PATCH
41	600080 1/4 ALUM CLAMP W/LINER
42	610053 3/8-16 NYLON INSERT LOCK NUT
43	680008 1/4-20 X 1-3/4 SELF TAPPING SCREW
44	730087 #12 SS EXT TOOTH LOCKWASHER
45	760203 AIR FILTER
46	900113 10" TUBEAXIAL FAN, 115V
47	1950074 BATTERY MOUNTING BRACKET

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

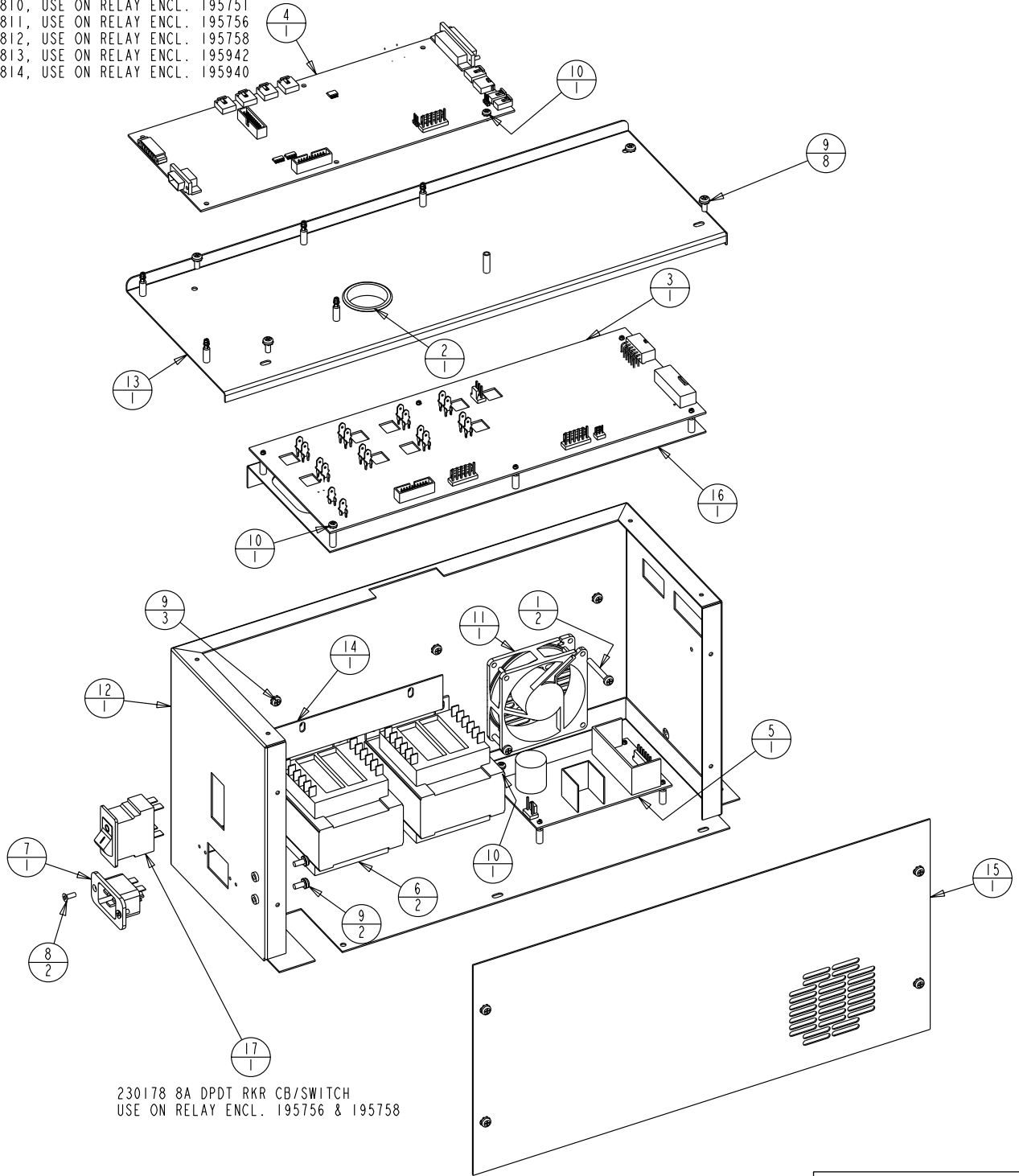
ThermoFisher
SCIENTIFIC

BOX 649, MARIETTA, OHIO 45750

MODEL/PART NAME: 8600 UP-RIGHT FREEZER	
DWG TITLE: 8605 UP-RIGHT FREEZER ASSEMBLY	
DWN: PDK	CAD: PDK
APPD: MAH	DATE: 10-30-02
SCALE: 0.094	
MATERIAL: N/A	
PAINT: N/A	
TOLERANCE UNLESS OTHERWISE SPECIFIED DECIMAL: .XX±	DRAWING NUMBER 8602-203-1
ANGLES: .XXX±	SIZE B

Assembly Drawing
Upright Freezer
Base
8602-203-1-D Rev. 5
Page 2 of 2

191804, USE ON RELAY ENCL. 195754, 195755
 191810, USE ON RELAY ENCL. 195751
 191811, USE ON RELAY ENCL. 195756
 191812, USE ON RELAY ENCL. 195758
 191813, USE ON RELAY ENCL. 195942
 191814, USE ON RELAY ENCL. 195940



230178 8A DPDT RKR CB/SWITCH
 USE ON RELAY ENCL. 195756 & 195758

1. COMMON TO: 195751, 195754, 195755, 195756, 195758, 195940 & 195942

Part Number

15
1

Quantity

Exploded View
 Drawing
 230V Relay Enclosure
 Assembly

8602-204-1-D Rev. 5
 Page 1 of 2

REV	ECN NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION
1	FR-1673	03-06-03	DHG	KDG	LDN	MADE COMMON TO 195940 & 195942
2	FR-1789	07-28-04	ADT	KDG	LDN	CHG. MICRO BOARD FOR VACUUM RELIEF
3	PIP-111	08-02-04	TJ	KDG	LDN	REMOVED 114031 GROMMET EDGING
4	FR-1806	08-23-04	JDL	KDG	LDN	SPECIFIED AMPERAGE OF CB SWITCHES
5	FR-2008	02-02-09	RTB	SAG	CCS	REFLECTS CHANGES MADE TO METALWORK

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	191658	HIGH VOLTAGE BOARD 230V
4	191804	MICRO BOARD (-86 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

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

MODEL/PART NAME: RELAY ENCLOSURE ASSEMBLY					
DWG TITLE: 230 VOLT RELAY ENCLOSURE ASSY (HIGH END)					
DWN: DHG	CAD: DHG	APPD: MAH	DATE: 07-26-01	SCALE: 0.250	
MATERIAL: -					
PAINT: N/A					
TOLERANCE UNLESS OTHERWISE SPECIFIED			DRAWING NUMBER		SIZE
ANGLES: DECIMAL: .XX=± .xxx=±			8602-204-1		B

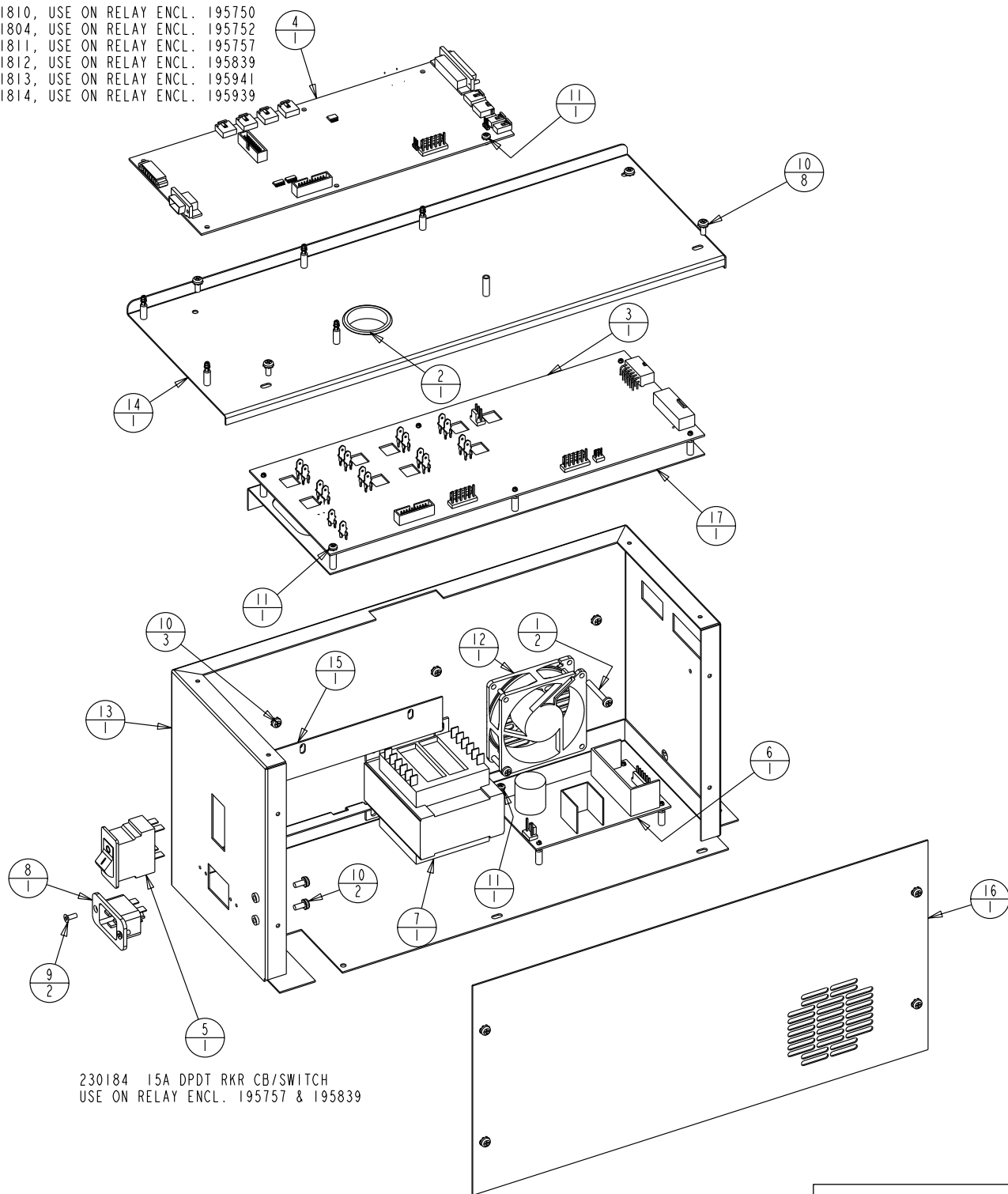
**Exploded View
Drawing
230V Relay Enclosure
Assembly**

8602-204-1-D Rev. 5
Page 2 of 2

**ThermoFisher
SCIENTIFIC**

BOX 649, MARIETTA, OHIO 45750

191810, USE ON RELAY ENCL. 195750
 191804, USE ON RELAY ENCL. 195752
 191811, USE ON RELAY ENCL. 195757
 191812, USE ON RELAY ENCL. 195839
 191813, USE ON RELAY ENCL. 195941
 191814, USE ON RELAY ENCL. 195939



230184 15A DPDT RKR CB/SWITCH
 USE ON RELAY ENCL. 195757 & 195839

1. COMMON TO: 195750, 195752, 195757, 195839, 195939 & 195941

Exploded View
 Drawing
 120V Relay Enclosure
 Assembly
 8602-204-2-D Rev. 5
 Page 1 of 2

REV	ECN NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION
1	FR-1673	03-06-03	DHG	KDG	KDG	MADE COMMON TO 195939 & 195941
2	FR-1789	07-28-04	ADT	KDG	LDN	CHG. MICRO BOARD FOR VACUUM RELIEF
3	PIP-111	08-02-04	TJ	KDG	LDN	REMOVED 114031 GROMMET EDGING
4	FR-1806	08-23-04	JDL	KDG	LDN	SPECIFIED AMPERAGE OF CB SWITCHES
5	FR-2008	02-02-09	RTB	SAG	CCS	REFLECTS CHANGES MADE TO METAL WORK

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	191680	HIGH VOLTAGE BOARD 120V
4	191810	MICRO BOARD (-86 LOW END)
5	230183	20A 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

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

ThermoFisher
SCIENTIFIC

BOX 649, MARIETTA, OHIO 45750

MODEL/PART NAME: RELAY ENCLOSURE ASSEMBLY

DWG TITLE: 120 VOLT RELAY ENCLOSURE ASSY (LOW END)

DWN: DHG CAD: DHG APPD: MAH DATE: 07-26-01 SCALE: 0.250

MATERIAL: -

PAINT: N/A

TOLERANCE UNLESS OTHERWISE SPECIFIED
ANGLES: DECIMAL: .XX±
.xxx±

DRAWING NUMBER

8602-204-2

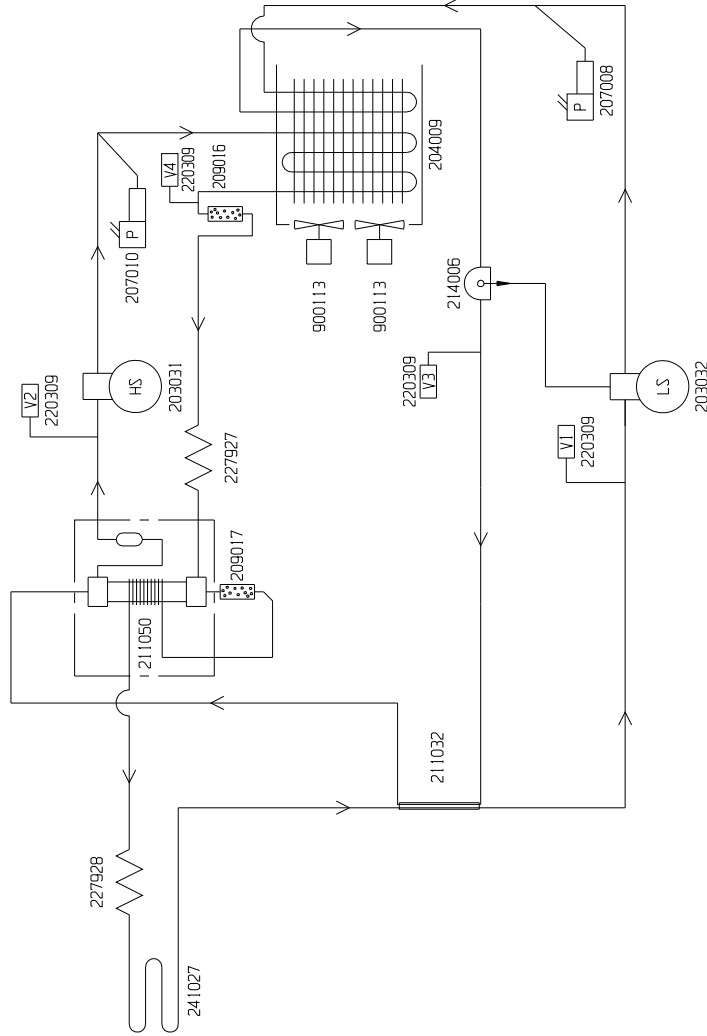
SIZE

B

Exploded View
Drawing
120V Relay Enclosure
Assembly

8602-204-2-D Rev. 5
Page 2 of 2

HIGH TEMP STAGE



LOW TEMP STAGE

CASCADE REFRIGERATION

HIGH TEMPERATURE STAGE REFRIGERANT:

13 CUFT UNITS: R-404A 24 OZ. (680g) ±1/2 OZ (14g)

HIGH TEMPERATURE STAGE OIL: MOBILE EAL ARCTIC 22CC POLYOL ESTER OIL COMPRESSOR; 24 OZ. (710ml)

LOW TEMPERATURE STAGE REFRIGERANT:

13 CUFT UNITS: R-290 0.5 OZ. (14g) MAX.; VAC TO 20" R-508B 13.5 OZ (382g) ± 1/4 OZ. (7g) OR 20" TO 166 PSIG ± 3 PSIG

LOW TEMPERATURE STAGE OIL: ZEROL 150T WITH 2% ADDITIVE COMPRESSOR; 24 OZ (710ml) 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

ThermoFisher
SCIENTIFIC
BOX 649, MARIETTA, OHIO 45750

1. COMMON TO: 702, 902, 991, 995, 8602, 8627, 8691, 5702 & 5602 13 CU. FT. FREEZERS (230 V)

10	FR-2080	10-10-10	NM	KDG	CCS	147	TO 166 PSIG, OLE TO DANFOSS DRIVER
9	FR-2080	07-27-09	NM	KDG	LDN	209017	WAS 209016, 13 TO 13.5 OZ
8	FR-2049	04-06-09	NM	SAG	LDN	209016	WAS 209020, ADD (1) 220309
7	FR-2004	07-25-08	KDG	KDG	LDN		ADDED COMMON TO 995
6	FR-1943	06-13-07	CRM	SAG	LDN	211050	WAS 211039
5	FR-1962	06-06-07	SUN	SAG	LDN	227928	WAS 241017, 227927 WAS 241036
4	FR-1867	07-26-05	RDS	DHG	LDN	REMOVED (2)	220553 ADDED (2), 220309
REV	ECR NO.	DATE	BY	CAO	APPD	DESCRIPTION OF REVISION	
DATE	05-14-02	LDWN	MAH	CAO	KOG	APPD	MAH SCALE NONE
CUSTOMER							
JOB TITLE	-86°C 13 CU. FT. UPRIGHT FREEZERS (230 V)						
DWG TITLE	REFRIGERATION SCHEMATIC						
LOCATION	JOB NUMBER	DRAWING NUMBER					
		8602-90-0-B					

CASCADE REFRIGERATION

HIGH TEMPERATURE STAGE REFRIGERANT:

13 CUFT UNITS: R-404A 24 OZ. (680g) ±1/2 OZ (14g)

HIGH TEMPERATURE STAGE OIL: MOBILE EAL ARCTIC 22CC POLYOL ESTER OIL COMPRESSOR, 24 OZ. (710ml)

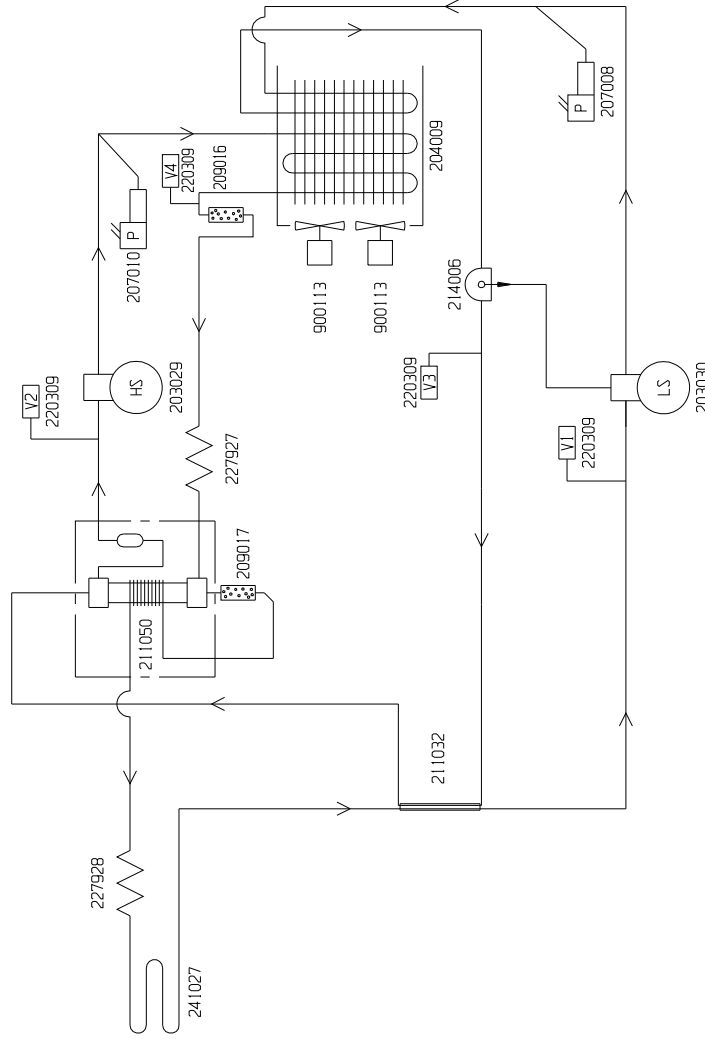
LOW TEMPERATURE STAGE REFRIGERANT:

13 CUFT UNITS: R-290 0.5 OZ. (14g) MAX. ; VAC TO 20" R-508B 13.5 OZ (382g) ± 1/4 OZ (.7g) DR 20" TO 166 PSIG ± 3 PSIG

LOW TEMPERATURE STAGE OIL: ZEROL 150T WITH 2% ADDITIVE COMPRESSOR: 24 OZ. (710ml) OIL SEPARATOR: 15 OZ. (444ml)

* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C

HIGH TEMP STAGE



LOW TEMP STAGE

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

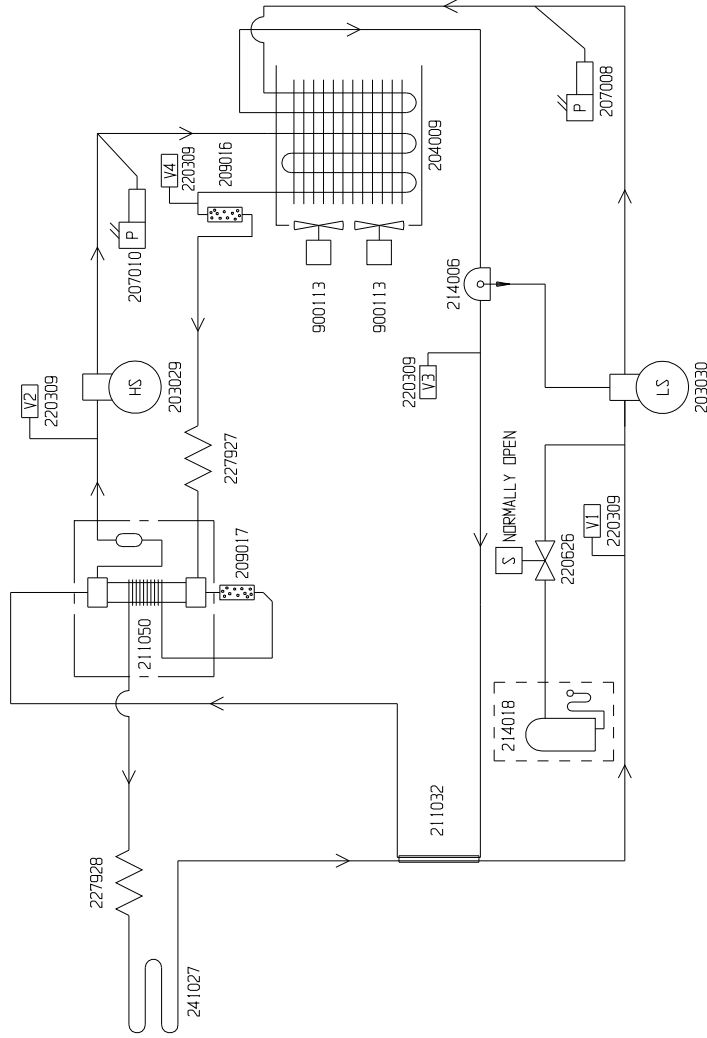
BOX 649, MARIETTA, OHIO 45750

1. COMMON TO: 703, 903, 992, 8603, 8692, 5703 & 5603 13 CU. FT. FREEZERS (120 V)

8	FR-2080	10-10-10	NM	KDG	CCS	147 TO 166 PSIG, DUE TO DANFOSSE DRIVER
7	FR-2080	07-27-09	NM	KDG	LDN	209017 WAS 209016, 13 TO 13.5 OZ
6	FR-2049	04-06-09	NM	SAG	LDN	209016 WAS 209020, ADD (1), 220309
5	FR-1943	06-13-07	CRM	SAG	LDN	211050 WAS 211039
4	FR-1962	06-06-07	SJN	SAG	LDN	227928 WAS 241017, 227927 WAS 241036
3	FR-1867	07-26-05	RDS	DHG	MAH	REMOVED (2) 220553 ADDED (2) 220309
2	FR-1729	08-27-03	RDS	KDG	MAH	220554 RETOLLOCK VALVE TO 220553
REV	ECR NO.	DATE	BY	CAI	APPD	DESCRIPTION OF REVISION
		05-14-02	DWN	MAH	CAJ	SCALE NONE
CUSTOMER						
JOB TITLE -66°C 13 CU. FT. UPRIGHT FREEZERS (120 V)						
DWG TITLE REFRIGERATION SCHEMATIC						
LOCATION					JOB NUMBER	DRAWING NUMBER
						8603-90-0-B

CASCADE REFRIGERATION

HIGH TEMP STAGE



HIGH TEMPERATURE STAGE REFRIGERANT:

17 CUFT UNITS: R-404A 24 OZ. (680g) ±1/2 OZ (14g)

HIGH TEMPERATURE STAGE OIL: MOBILE EAL ARCTIC 22CC POLYOL ESTER OIL COMPRESSOR, 24 OZ. (710ml)

LOW TEMPERATURE STAGE REFRIGERANT:

17 CUFT UNITS: R-290 0.8 OZ. (23g) MAX. ; VAC TO 2 PSIG
R-508B 14.4 OZ (408g) ± 1/4 OZ (7g) DR
2 PSIG TO 137 PSIG ± 3 PSIG

LOW TEMPERATURE STAGE OIL: ZEROL 150T WITH 2% ADDITIVE
COMPRESSOR: 24 OZ. (710ml)
OIL SEPARATOR: 15 OZ. (444ml)

* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C

LOW TEMP STAGE

THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR REPRODUCING OR TRANSMITTING IN ANY MANNER WITHOUT WRITTEN PERMISSION FROM THERMO FISHER SCIENTIFIC

ThermoFisher
SCIENTIFIC

BOX 668, MARLETTA, OHIO 45750

1. COMMENT TO: 704, 904, 993, 8604, 8693, 5704 & 5604 17 CU. FT. FREEZERS (120 V)

10	FR-2080	10-10-10	NWM	KDG	CCS	123 TO 137 PSIG, OLE TO DMFOSS DRYER				
9	FR-2080	07-27-09	NWM	KDG	LDN	209017 WAS 209016, 15 TO 14.4 OZ				
8	FR-2049	04-06-09	NWM	SAG	LDN	209016 WAS 209020, ADD (1) 220309				
7	FR-1943	06-13-07	CRM	SAG	LDN	211050 WAS 211039				
6	FR-1962	06-06-07	SJN	SAG	LDN	227928 WAS 241017, 227927 WAS 241036				
5	FR-1867	07-26-05	RDS	DHG	MAH	REMOVED (2) 226553 ADDED (2) 220309				
4	FR-1729	08-27-03	RDS	KDG	MAH	2220554 ROTILOCK VALVE TO 2220553				
REV	ECR NO.	DATE	BY	CAJ/APPO	DESCRIPTION OF REVISION					
		05-14-02	DWN	MAH	CAJ	KDG	APPO	MAH	SCALE	NONE

CUSTOMER

JOB TITLE -86°C 17 CU. FT. UPRIGHT FREEZERS (120 V)

DWG TITLE REFRIGERATION SCHEMATIC

LOCATION

JOB NUMBER

DRAWING NUMBER

8604-90-0-B

CASCADE REFRIGERATION

HIGH TEMPERATURE STAGE REFRIGERANT:

17 CUFT UNITS: R-404A 24 OZ. (680g) ±1/2 OZ (14g)

HIGH TEMPERATURE STAGE OIL: MOBILE EAL ARCTIC 22CC POLYOL ESTER OIL COMPRESSOR, 24 OZ. (710ml)

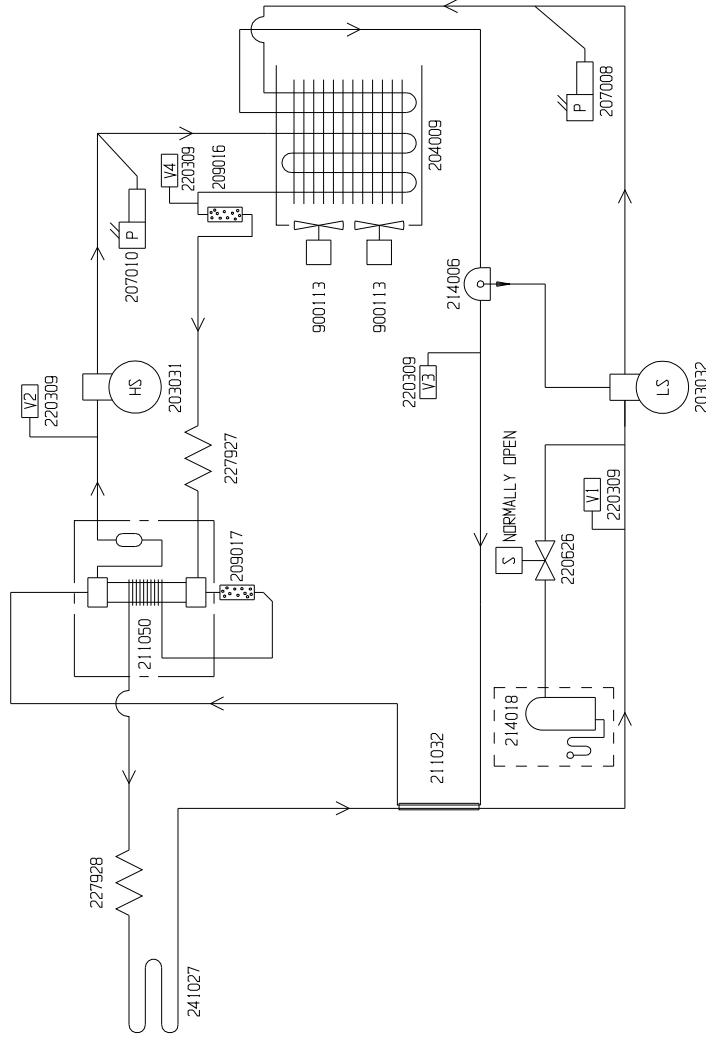
LOW TEMPERATURE STAGE REFRIGERANT:

17 CUFT UNITS: R-290 0.9 OZ. (26g) MAX. ; VAC TO 3 PSIG
R-508B 16 OZ (453g) ± 1/4 OZ (7g) OR
3 PSIG TO 152 PSIG ± 3 PSIG

LOW TEMPERATURE STAGE OIL: ZEROL 150T WITH 2% ADDITIVE
COMPRESSOR: 24 OZ. (710ml)
OIL SEPARATOR: 15 OZ. (444ml)

* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C

HIGH TEMP STAGE



LOW TEMP STAGE

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

ThermoFisher
SCIENTIFIC
BOX 649, MARLETTA, OHIO 45750

1. COMMON TD: 705, 905, 994, 997, 8605, 8694, 5705 & 5605 17 CU. FT. FREEZERS (230 V)

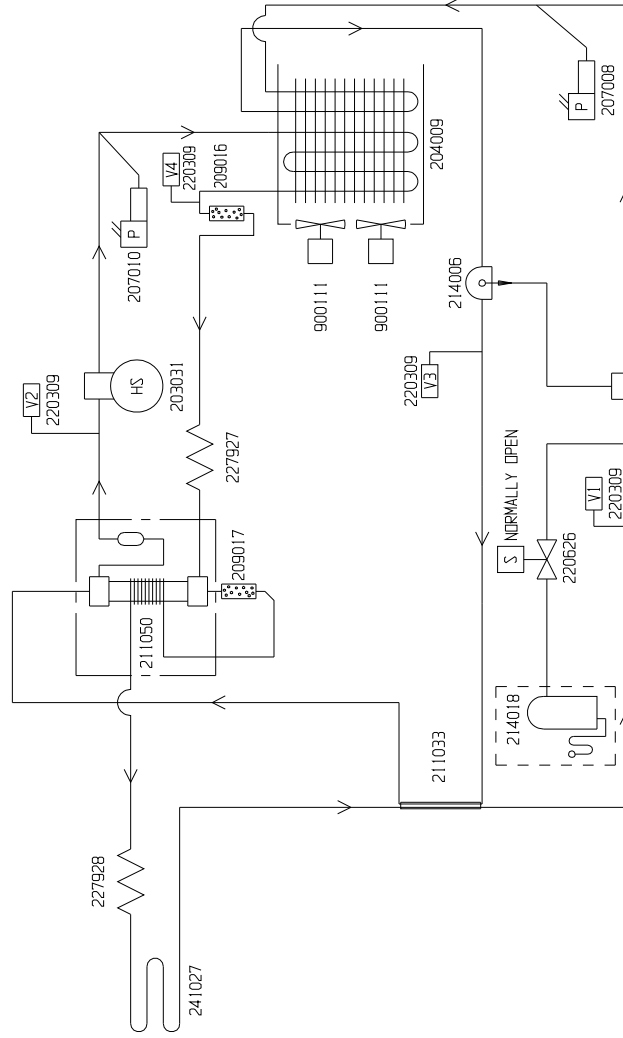
11	FR-2080	10-10-10	NM	KDG	CCS	134 TO 152 PSIG, OLE TO DAMPOSS DRYER
10	FR-2080	07-27-09	NM	KDG	LDN	209017 WAS 209016, 15 TO 16 OZ
9	FR-2049	04-06-09	NM	SAG	LDN	209016 WAS 209020, ADD (1) 220309
8	FR-2004	07-25-08	KDG	KDG	LDN	ADDED COMMON TO 997
7	FR-1943	06-13-07	ERM	SAG	LDN	211050 WAS 211039
6	FR-1962	06-06-07	SJN	SAG	LDN	227928 WAS 241017, 227927 WAS 241036
5	FR-1867	07-25-05	ROD	DHG	LDN	REMOVED (2) 220553 ADDED (2) 220309
REV	ECR NO.	DATE	BY	CA/APPD	DESCRIPTION OF REVISION	
	DATE	05-14-02	DWN	MAH	CAO KDG APPD MAH	SCALE NONE
CUSTOMER						
JOB TITLE -86°C 17 CU. FT. UPRIGHT FREEZERS (230 V)						
DWG TITLE REFRIGERATION SCHEMATIC						
LOCATION					JOB NUMBER	DRAWING NUMBER
					8605-90-0-B	

CASCADE REFRIGERATION

HIGH TEMP. STAGE REFRIGERANT: R-404A, 24 OZ. (680g) ± 1/2 OZ. (14g)
 HIGH TEMPERATURE STAGE OIL: MIBILE EAL ARCTIC 22CC POLYOL ESTER OIL COMPRESSOR, 24 OZ. (710ml)
 LOW TEMP. STAGE REFRIGERANT: R-290 1.0 OZ. (28g) MAX.; VAC TO 5 PSIG R-508B 18.3 OZ. (518g) ± 1/4 OZ. (7g) OR 5 PSIG TO 163 PSIG ± 3 PSIG
 LOW TEMPERATURE STAGE OIL: ZEROL 150T WITH 2% ADDITIVE COMPRESSOR, 24 OZ. (710ml) OIL SEPARATOR, 15 OZ. (444ml)

* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C

HIGH TEMP STAGE



LOW TEMP STAGE

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

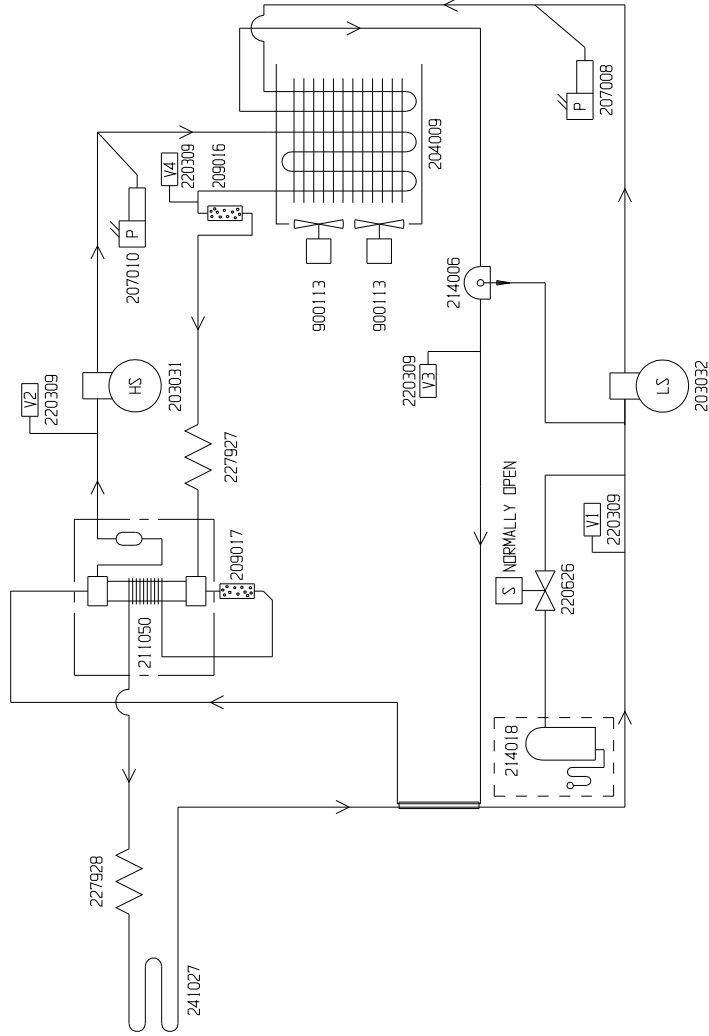
BOX 649, MARLETTA, OHIO 45750

1. COMMON TO: 706, 906, 995, 998, 8606, 8695, 5706 & 5606 23 CU.FT. FREEZERS (230 V)

11	FR-2080	10-10-10	NM	KDG	CCS	152 TO 163 PSIG, DUE TO DAMPSS DRIVER				
10	FR-2080	07-27-09	NM	KDG	LDN	209017 WAS 209016, 17.3 TO 18.3 OZ				
9	FR-2049	04-06-09	NM	SAG	LDN	209016 WAS 209020, ADD (1) 220309				
8	FR-2004	07-24-08	KDG	KDG	LDN	ADDED COMMON TO 998				
7	FR-1943	06-13-07	CRM	SAG	LDN	211050 WAS 211039				
6	FR-1962	06-06-07	SJN	SAG	LDN	227928 WAS 241017, 227927 WAS 241036				
5	FR-1867	07-26-05	RDS	DHG	LDN	REMOVED (2), 220553, ADDED (2) 220309				
REV	ECR NO.	DATE	BY	CAO	APPD	DESCRIPTION OF REVISION				
		05-14-02	DWN	MAH	CAO	KDG	APPD	MAH	SCALE	NONE
CUSTOMER										
JOB TITLE -86°C 23 CU. FT. UPRIGHT FREEZERS (230 V)										
DWG TITLE REFRIGERATION SCHEMATIC										
LOCATION										
JOB NUMBER										
DRAWING NUMBER										
8606-90-0-B										

CASCADE REFRIGERATION

HIGH TEMP STAGE



HIGH TEMPERATURE STAGE REFRIGERANT:

28 CUFT UNITS: R-404A 24 OZ. (680g) ± 1/2 OZ (14g)

HIGH TEMPERATURE STAGE OIL: MOBILE EAL ARCTIC 22CC POLYOL ESTER OIL COMPRESSOR, 24 OZ. (710ml)

LOW TEMPERATURE STAGE REFRIGERANT:

28 CUFT UNITS: R-290 1.0 OZ. (28g) MAX. ; VAC TO 4 PSIG
R-508B 18.7 OZ (529g) ± 1/4 OZ (7g) OR
4 PSIG TO 157 PSIG ± 3 PSIG

LOW TEMPERATURE STAGE OIL: ZEROL 150T WITH 2% ADDITIVE
COMPRESSOR: 24 OZ. (710ml)
OIL SEPARATOR: 15 OZ. (444ml)

* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C

11	FR-2080	10-10-10	NM	KDG	CCS	144	TO 157 PSIG, DUE TO DANFOSS DRYER
10	FR-2080	07-27-09	NM	KDG	LON	209017	WAS 209016, 17.7 TO 18.7 OZ
9	FR-2049	04-06-09	NM	SAG	LON	209016	WAS 209020, ADD (1) 220309
8	FR-2004	07-01-08	KDG	KDG	LON		ADDED COMMON TO 5607
7	FR-1943	06-13-07	CRM	SAG	LON		211050 WAS 211041
6	FR-1962	06-11-07	S	JN	SAG	LON	227928 WAS 241017, 227927 WAS 241036
5	FR-1867	07-26-05	ROS	DHG	LON		REMOVED (2) 220653 ADDED (2) 220309
REV	ECR NO.	DATE	BY	CAO	APPO	DESCRIPTION	OF REVISION
	DATE	05-14-02	DWN	MAH	CAO	KDG	APPO MAH SCALE NONE
CUSTOMER							
JOB TITLE -86°C 28 CU. FT. UPRIGHT FREEZERS (230 V)							
DWG TITLE REFRIGERATION SCHEMATIC							
LOCATION						JOB NUMBER	DRAWING NUMBER
						8607-90-0-B	

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, MARLETTA, OHIO 45750

1. COMMON TO: 8607, 907 & 5607 28 CU. FT. FREEZERS (230 V)

DRAWING NUMBER: 8656-90-0-B

REV	NO.	DATE	BY	APP'D	DESCRIPTION OF REVISION	
2	FR-1883	12-20-05	JAS	KDG	LON	CORRECTLY SHOWN EXP. TANK & VALVE
3	FR-1962	06-06-07	S.JN	SAG	LON	227928 WAS 241017. 227927 WAS 241036
4	FR-1943	06-13-07	CRM	SAG	LON	211050 WAS 211039
5	FR-2049	04-06-09	NMM	SAG	LON	209016 WAS 209020.ADD (1) 220309
6	FR-2080	07-27-09	NMM	KDG	LON	209017 WAS 209016. 15 TO 16 OZ
7	FR-2080	10-10-10	NMM	KDG	CCS	133 TO 145 PSIG. DUE TO DANFOSS DRIER

CASCADE REFRIGERATION

HIGH TEMPERATURE STAGE REFRIGERANT:

23 CUFT UNITS: R-404A 24 OZ. (680g) ± 1/2 OZ (14g)

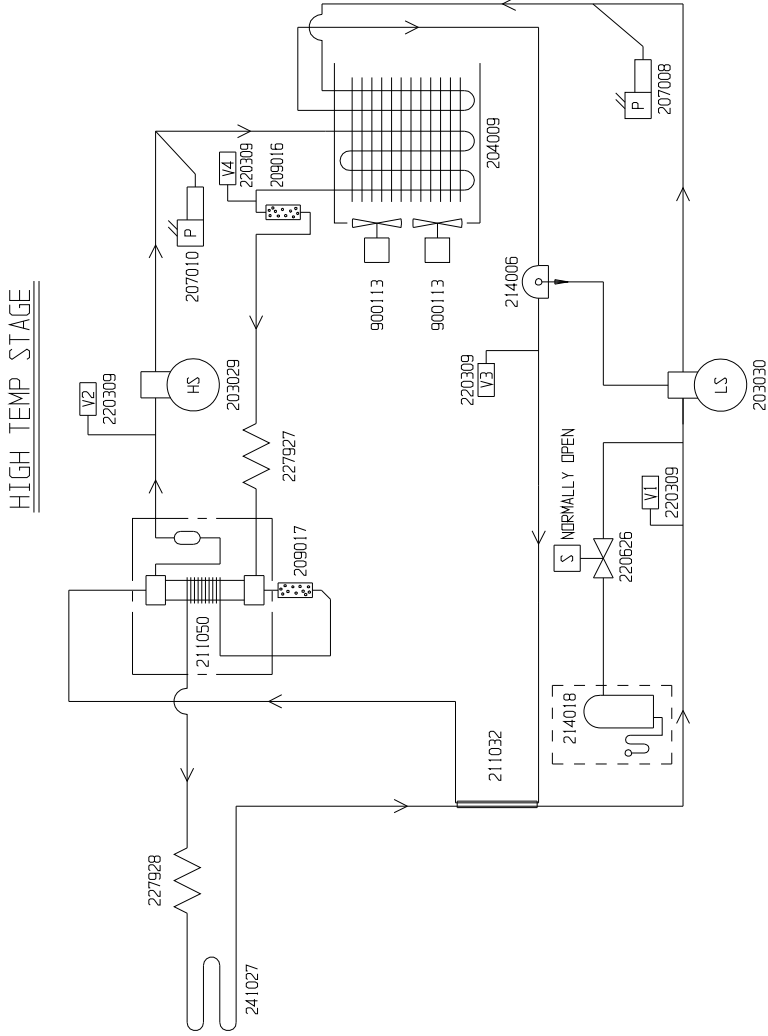
HIGH TEMPERATURE STAGE OIL: MOBILE EAL ARCTIC 22CC POLYOL ESTER OIL COMPRESSOR, 24 OZ. (710ml)

LOW TEMPERATURE STAGE REFRIGERANT:

23 CUFT UNITS: R-290 0.9 OZ. (26g) MAX. ; VAC TO 4 PSIG
R-5088 16 OZ (453g) ± 1/4 OZ. (7g) OR
4 PSIG TO 145 PSIG ± 3 PSIG

LOW TEMPERATURE STAGE OIL: ZEROL 150T WITH 2% ADDITIVE
COMPRESSOR: 24 OZ. (710ml)
OIL SEPARATOR: 15 OZ. (444ml)

* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C



LOW TEMP STAGE

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

BDX 649, MARIETTA, OHIO 45750

MODEL/PART NAME: -86° 23 CU. FT. UP-RIGHT FREEZERS (120 VOLT)	PAINT COLOR: N/A	DRAWING NUMBER: 8656-90-0	SIZE: B
DWG TITLE: REFRIGERATION SCHEMATIC	TOLERANCE UNLESS OTHERWISE SPECIFIED: .XX±		
DWN: ADT	CD: pdk	APPD: MAH	DATE: 03-09-04
SCALE: NONE	ANGLES: DECIMAL: .XXX±		
MATERIAL: N/A			

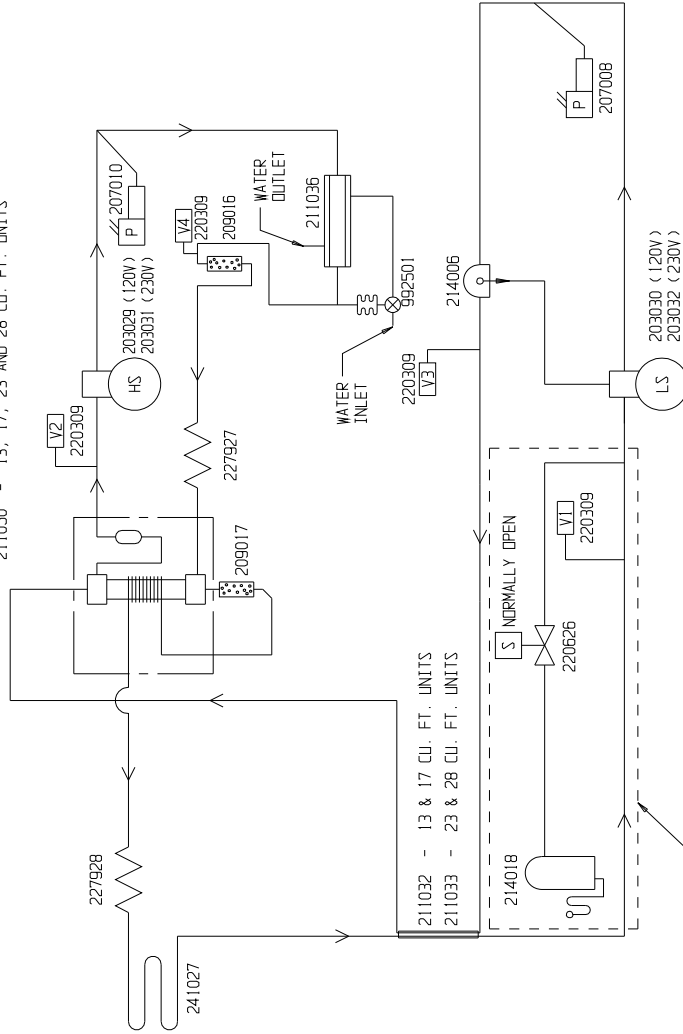
1. COMMON TO: 756, 956, 990, 8656, 8690, 5756 & 5656 23 CU. FT. FREEZERS (120 V)

DRAWING NUMBER: 195964-90-1-B

REV	NO.	DATE	BY	CAO	APPD	DESCRIPTION OF REVISION
3	FR-1963	06-13-07	SJM	SAG	LON	227928 WAS 241017, 227927 WAS 241028
4	FR-1943	06-13-07	CRM	SAG	LON	MADE 211050 STANDARD
5	FR-2049	04-06-09	NMM	KOG	LON	209016 WAS 209020, ADD (1) 220309
6	FR-2080	07-27-09	NMM	KOG	LON	209017 WAS 209016 IN CAB., CHG. CHARGES
7	FR-2080	10-10-10	NMM	KOG	CCS	UPDATED PRESSURES DUE TO DANFLOSS ORDER

HIGH TEMP STAGE

211050 - 13, 17, 23 AND 28 CU. FT. UNITS



17, 23 AND 28 CU. FT. UNITS ONLY

LOW TEMP STAGE

CASCADE REFRIGERATION

HIGH TEMPERATURE STAGE REFRIGERANT:

- 13 CUFT UNITS: R-404A 24 OZ. (680gr) ±1/2 OZ (14gr)
- 17 CUFT UNITS: R-404A 24 OZ. (680gr) ±1/2 OZ (14gr)
- 23 CUFT UNITS: R-404A 24 OZ. (680gr) ±1/2 OZ (14gr)
- 28 CUFT UNITS: R-404A 24 OZ. (680gr) ±1/2 OZ (14gr)

HIGH TEMPERATURE STAGE OIL: MOBILE EAL ARCTIC 22CC POLYOL ESTER OIL COMPRESSOR, 24 OZ. (710ml)

LOW TEMPERATURE STAGE REFRIGERANT:

- 13 CUFT UNITS: R-290 0.5 OZ. (14gr) MAX. ; VAC TO 20" R-508B 13.5 OZ. (382gr) ± 1/4 OZ. (.7gr) DR 20" TO 166 PSIG ± 3 PSIG
- 17 CUFT UNITS: R-290 0.8 OZ. (23gr) MAX. ; VAC TO 2 PSIG (120 VOLTS) R-508B 14.4 OZ. (406gr) ± 1/4 OZ. (.7gr) DR 2 PSIG TO 137 PSIG ± 3 PSIG
- 17 CUFT UNITS: R-290 0.9 OZ. (26gr) MAX. ; VAC TO 3 PSIG (230 VOLTS) R-508B 16 OZ. (453gr) ± 1/4 OZ. (.7gr) DR 3 PSIG TO 152 PSIG ± 3 PSIG
- 23 CUFT UNITS: R-290 1.0 OZ. (28gr) MAX. ; VAC TO 5 PSIG R-508B 18.3 OZ. (516gr) ± 1/4 OZ. (.7gr) DR 5 PSIG TO 163 PSIG ± 3 PSIG
- 28 CUFT UNITS: R-290 1.0 OZ. (28gr) MAX. ; VAC TO 4 PSIG R-508B 18.7 OZ. (529gr) ± 1/4 OZ. (.7gr) DR 4 PSIG TO 157 PSIG ± 3 PSIG

LOW TEMPERATURE STAGE OIL: ZEROL 150T WITH 2% ADDITIVE COMPRESSOR: 24 OZ. (710ml)

OIL SEPARATOR: 15 OZ. (444ml)

***SET WATER REGULATING VALVE FOR 170 PSIG

* 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

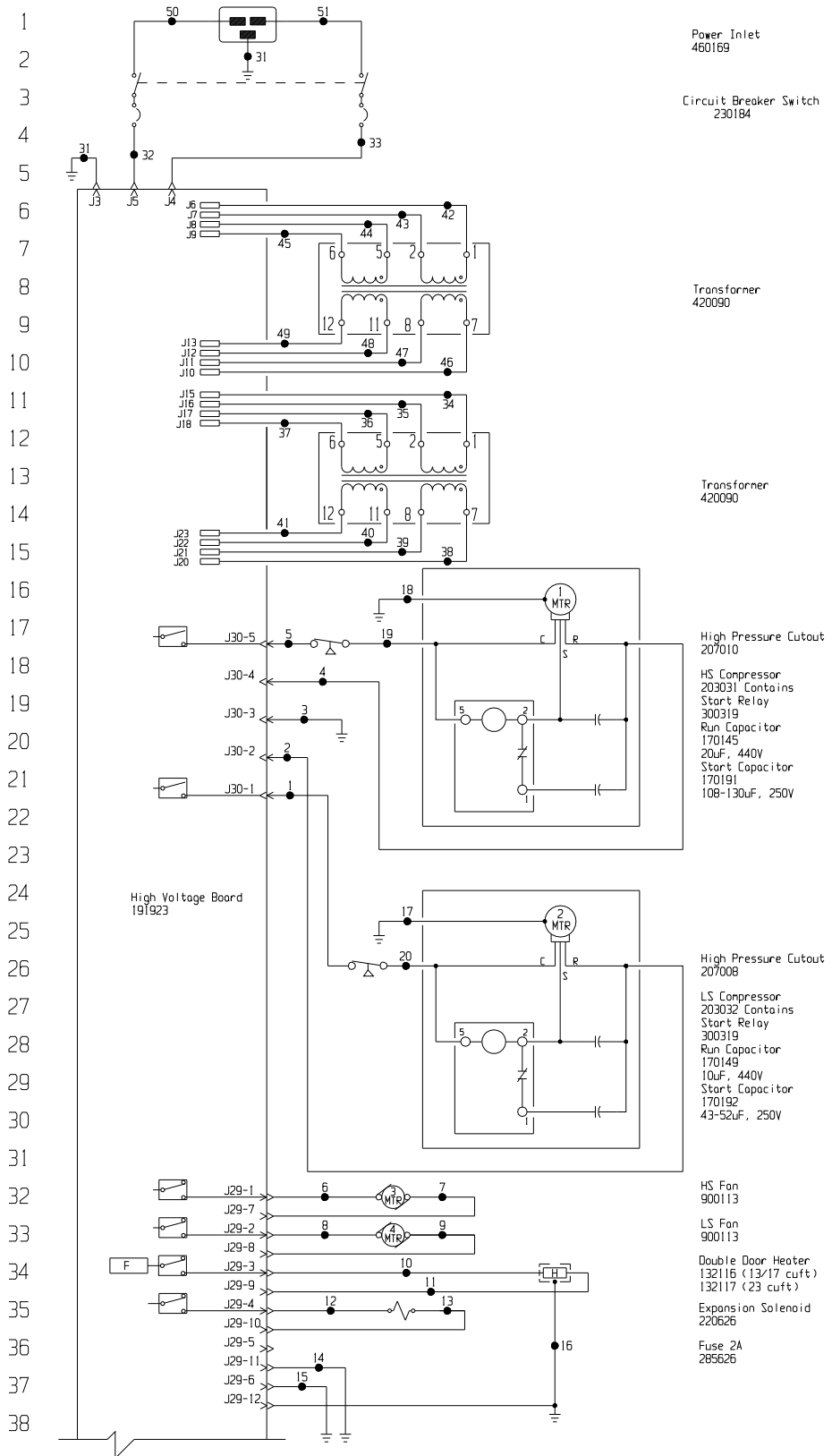
ThermoFisher
SCIENTIFIC

BOX 649, MARIETTA, OHIO 45750

COMMON TO: 195965

MODEL/PART NAME: -86°C 13, 17, 23 & 28 CU. FT. UPRIGHT H2O COOLED	DWG TITLE: FREEZER REFRIGERATION SCHEMATIC
DWN: DHG	CAO: DHG
APPD: MAH	DATE: 01-23-03
MATERIAL: N/A	SCALE: NONE
PAIN'T COLOR: N/A	
TOLERANCE UNLESS OTHERWISE SPECIFIED: .XX±	DRAWING NUMBER: 195964-90-1
ANGLES: DECIMAL: .XX±	SIZE: B

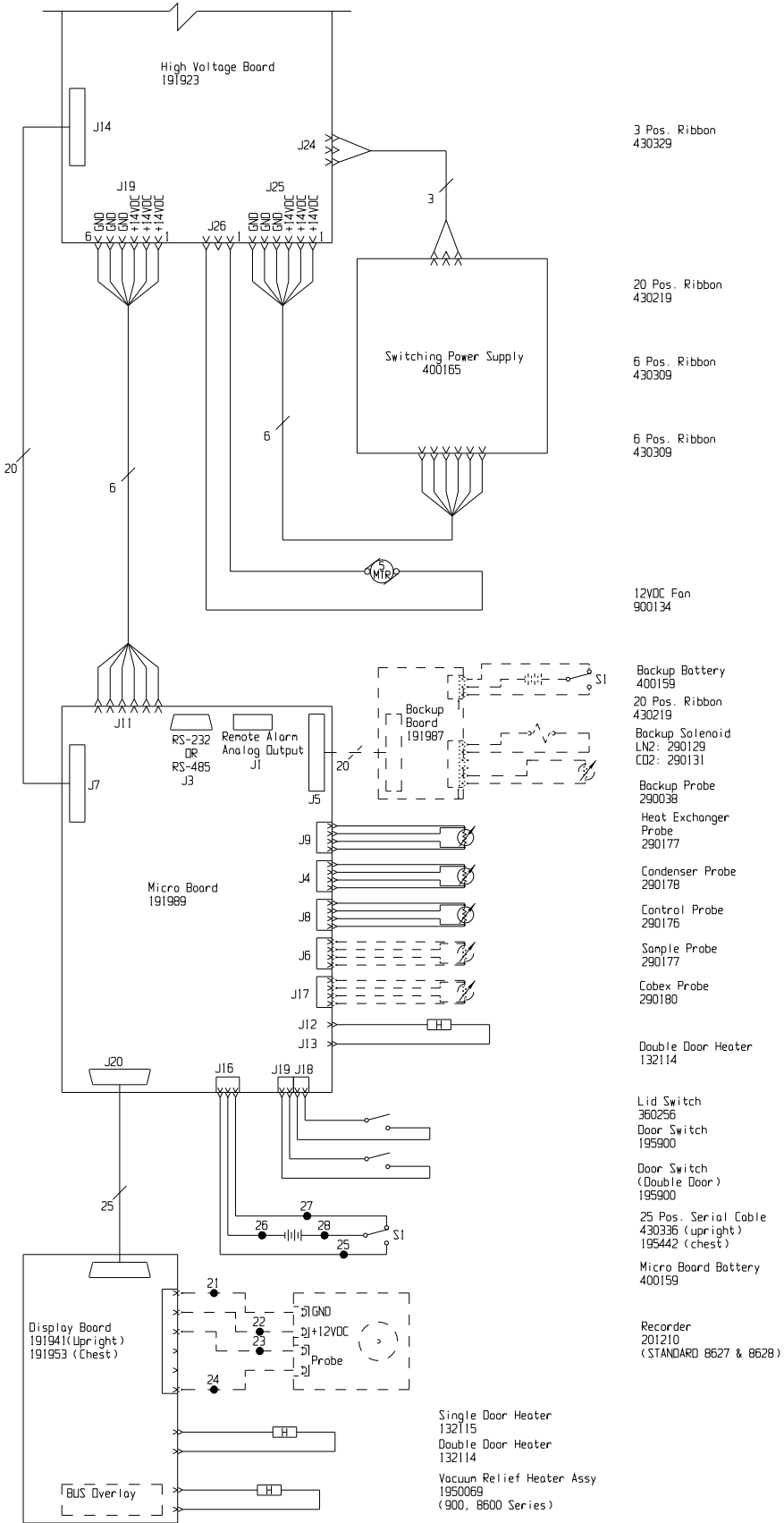
POWER CONNECTION
 230V, 1Ø, 2W, 50/60HZ, 12.0 FLA
 (OPERATING RANGE 208-240)



Electrical Schematic
 for
 230V -86C Chest
 and Upright Freezers

8602-70-2-D REV. 0
 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



- 3 Pos. Ribbon 430329
- 20 Pos. Ribbon 430219
- 6 Pos. Ribbon 430309
- 6 Pos. Ribbon 430309
- 12VDC Fan 900134
- Backup Battery 400159
- 20 Pos. Ribbon 430219
- Backup Solenoid LN2: 290129 CO2: 290131
- Backup Probe 290038
- Heat Exchanger Probe 290177
- Condenser Probe 290178
- Control Probe 290176
- Sample Probe 290177
- Cobex Probe 290180
- Double Door Heater 132114
- Lid Switch 360256
- Door Switch 195900
- Door Switch (Double Door) 195900
- 25 Pos. Serial Cable 430336 (upright) 195442 (chest)
- Micro Board Battery 400159
- Recorder 201210 (STANDARD 8627 & 8628)
- Single Door Heater 132115
- Double Door Heater 132114
- Vacuum Relief Heater Assy 1950069 (900, 8600 Series)

Electrical Schematic for 230V -86C Chest and Upright Freezers

8602-70-2-D REV. 0
Page 2 of 3

WIRE #	COLOR	GAUGE	WIRE #	COLOR	GAUGE	
77	1	BLK	14	25	WHT	18
	2	RED	14	26	BLK	18
78	3	GRN/YEL	14	27	BLK	18
	4	RED	14	28	RED	18
79	5	BLK	14	29		
	6	BLK	18	30		
80	7	BLK	18	31	GRN/YEL	14
	8	BLK	18	32	BLK	14
81	9	BLK	18	33	BLU	14
	10	BLK	18	34	BLK	14
82	11	BLK	18	35	BLU	14
	12	BLK	18	36	BLK	14
83	13	BLK	18	37	BLU	14
	14	GRN/YEL	18	38	BLK	14
84	15	GRN/YEL	18	39	BLU	14
	16	GRN/YEL	18	40	BLK	14
85	17	GRN/YEL	14	41	BLU	14
	18	GRN/YEL	14	42	BLK	14
86	19	BLK	14	43	BLU	14
	20	BLK	14	44	BLK	14
87	21	BLK	18	45	BLU	14
	22	RED	18	46	BLK	14
88	23	BLK	18	47	BLU	14
	24	RED	18	48	BLK	14
				49	BLU	14
				50	BLK	14
				51	BLU	14

Uprights

SIZE	13CF	17CF	23CF	28CF
MODEL	8602	8605	8606	8607
	8691	8694	8695	-
	5602	5605	5606	5607

Chests

SIZE	12CF	17CF	20CF
MODEL	8612	5819	8620
	5612	5219	5620
	5812	5155	5820
	5211	5175	5220
	5152	-	5156
	5172	-	5176

RS-232 SPECIFICATION

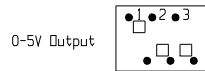
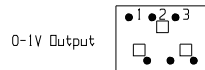
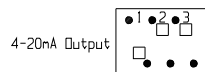
BAUD = 1200
 PARITY = N
 BITS = 8
 STOP BITS = 2

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

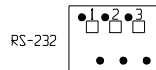
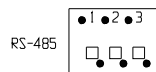
RS-232 PORT	
PIN# 2	TXD
PIN# 3	RXD
PIN# 5	GND

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

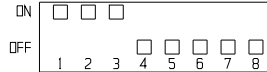
SW1 Settings for Analog Output



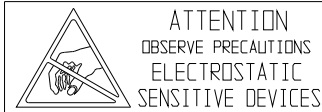
SW2 Settings for Communication Output



SW3 SETTINGS



1. Schematic represents single & double door upright models and chest models. All heaters are on uprights only. Expansion solenoid on 17, 23 & 28 uprights only.
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.



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, MARIETTA, OHIO 45750

REV	ENCL NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION
0	N/A	10-10-10	HCE	KDG	CCS	RELEASED FOR PRODUCTION / RELEASE 5

MODEL/PART NAME: 8600 SERIES, 230V, ULT FREEZERS
 DWG TITLE: UNIT SCHEMATIC
 DWG: HCE CAD: KDG APPD: LON DATE: 10-10-10 SCALE: NONE

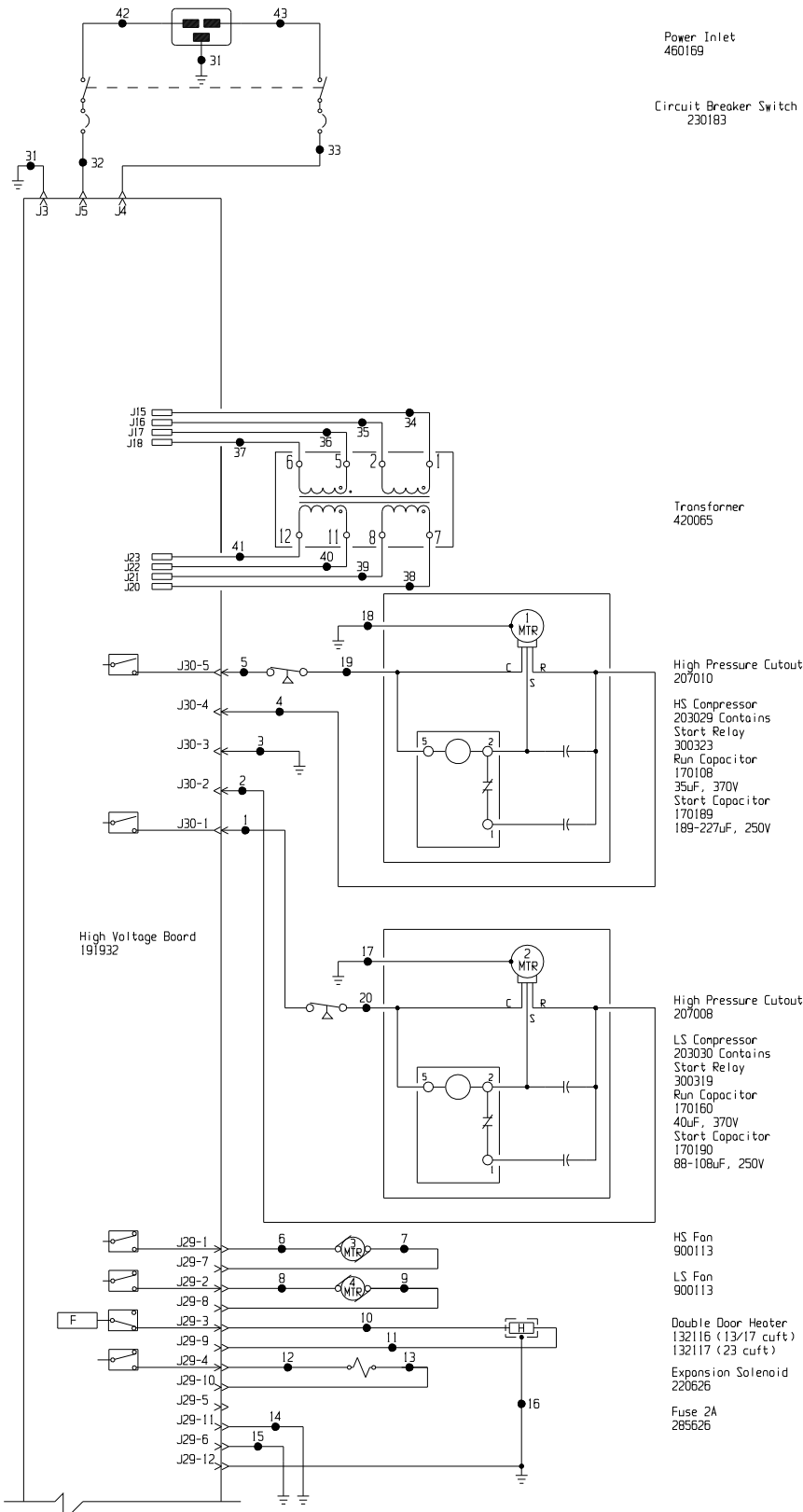
TOLERANCE UNLESS OTHERWISE SPECIFIED	DRAWING NUMBER	SIZE
ANGLES: DECIMAL: .XX± XXX±	8602-70-2	D

Electrical Schematic for
 230V -86C Chest
 and Upright Freezers

8602-70-2-D REV. 0
 Page 3 of 3

POWER CONNECTION
 120V, 1 ϕ , 2W, 60HZ, 16.0 FLA
 (OPERATING RANGE 108-130V)

1
2
3
4
5
6
7
8
9
10
11
12
13
14
15
16
17
18
19
20
21
22
23
24
25
26
27
28
29
30
31
32
33
34
35
36
37
38

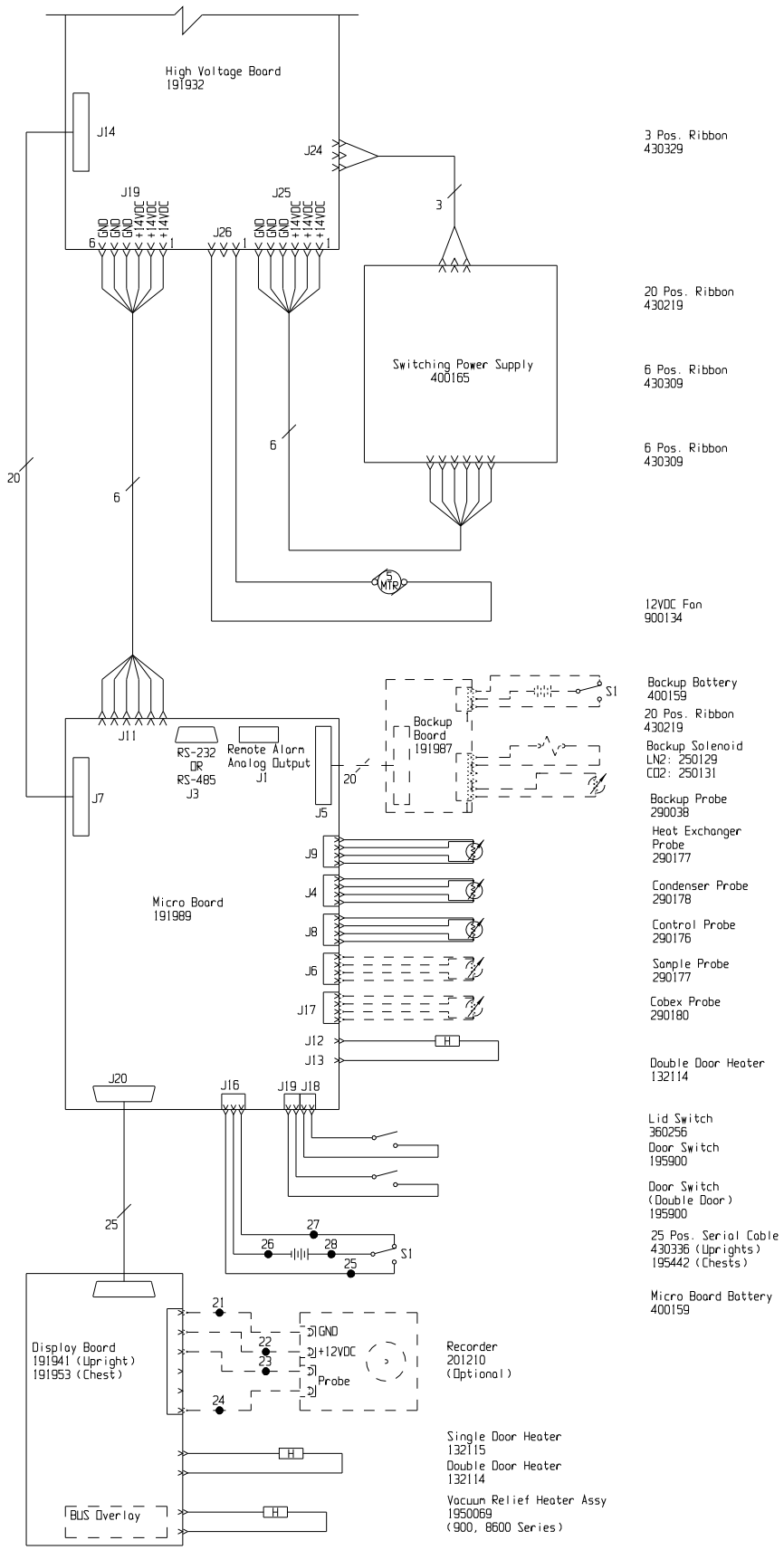


- Power Inlet
460169
- Circuit Breaker Switch
230183
- Transformer
420065
- High Pressure Cutout
207010
- HS Compressor
203029 Contains
Start Relay
300323
Run Capacitor
170108
35uF, 370V
Start Capacitor
170189
189-227uF, 250V
- High Pressure Cutout
207008
- LS Compressor
203030 Contains
Start Relay
300319
Run Capacitor
170160
40uF, 370V
Start Capacitor
170190
88-108uF, 250V
- HS Fan
900113
- LS Fan
900113
- Double Door Heater
13216 (13/17 cuft)
13217 (23 cuft)
- Expansion Solenoid
220526
- Fuse 2A
285626

Electrical Schematic
 for
 -86C 120V Chest
 and Upright Freezers

8603-70-2-D REV. 0
 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
for
-86C 120V Chest
and Upright Freezers

8603-70-2-D REV. 0
Page 2 of 3

WIRE #	COLOR	GAUGE	WIRE #	COLOR	GAUGE
1	BLK	14	25	WHT	18
2	RED	14	26	BLK	18
3	GRN/YEL	14	27	BLK	18
4	RED	14	28	RED	18
5	BLK	14	29		
6	BLK	18	30		
7	BLK	18	31	GRN/YEL	14
8	BLK	18	32	BLK	14
9	BLK	18	33	BLU	14
10	BLK	18	34	BLK	14
11	BLK	18	35	BLU	14
12	BLK	18	36	BLK	14
13	BLK	18	37	BLU	14
14	GRN/YEL	18	38	BLK	14
15	GRN/YEL	18	39	BLU	14
16	GRN/YEL	18	40	BLK	14
17	GRN/YEL	14	41	BLU	14
18	GRN/YEL	14	42	BLK	14
19	BLK	14	43	BLU	14
20	BLK	14			
21	BLK	18			
22	RED	18			
23	BLK	18			
24	RED	18			

Uprights

SIZE	13CF	17CF	23CF
MODEL	8603	8604	8656
	8692	8693	8690
	5603	5604	5656

Chests

SIZE	12CF	17CF	20CF
MODEL	8615	5818	8621
	5615	5218	5621
	5815	5154	5821
	5215	-	5221
	5153	-	5157

RS-232 SPECIFICATION

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

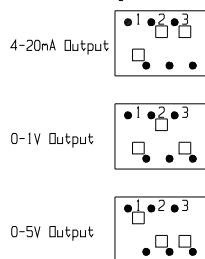
RS-232 PORT
PIN# 2 TXD
PIN# 3 RXD
PIN# 5 GND

REMOTE CONTACTS/ANALOG OUTPUT

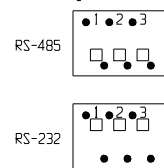
PIN# 1 Analog Output +
 PIN# 2 Analog Output -
 PIN# 3 Not Connected
 PIN# 4 Not Connected
 PIN# 5 Normally Closed
 PIN# 6 Common.
 PIN# 7 Normally Open

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

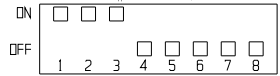
SW1 Settings for Analog Output



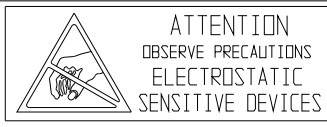
SW2 Settings for Communication Output



SW3 SETTINGS



1. Schematic represents single & double door upright models and chest models. All heater are for uprights only. Expansion solenoid on 17 & 23 cuft uprights only.
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.



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



REV	NO.	DATE	BY	APPD	DESCRIPTION OF REVISION	
0	N/A	10-10-10	HCE	KDG	CCS	RELEASED FOR PRODUCTION / RELEASE 5

MODEL/PART NAME: 8600 SERIES, 120V, ULT FREEZERS
 DWG TITLE: UNIT SCHEMATIC
 DWN: HCE CAD: KDG APPD: LON DATE: 10-10-10 SCALE: NONE
 MATERIAL:
 PAINT COLOR:
 TOLERANCE UNLESS OTHERWISE SPECIFIED: .XX±
 ANGLES: DECIMAL: .XXX±

Electrical Schematic for -86C 120V Chest and Upright Freezers
 8603-70-2-D REV. 0
 Page 3 of 3

THERMO FISHER SCIENTIFIC 900 & 8600 SERIES ULT FREEZER WARRANTY

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 Thermo's expense, labor included. The 900 Series ULT Freezers include an additional two year warranty on the compressors, parts only, F.O.B. factory. The 8600 Series 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.

In addition to the standard warranty, the foamed-in-place cabinet design carries a unit production lifetime warranty (foamed-in-place cabinet, evaporator and foamed-in-place door; parts only). Please contact your sales representative or Thermo for additional information.

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. The Technical Services Department 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.

Thermo shall not be liable for any indirect or consequential damages including, without limitation, damages relating to lost profits or loss of products.

Your local Thermo 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 your Technical Services Department 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. 4 4/09

THERMO FISHER SCIENTIFIC 900 & 8600 SERIES ULT FREEZER INTERNATIONAL DEALER WARRANTY

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. Dealers who stock our equipment are allowed an additional four months for delivery and installation, providing the warranty card is completed and returned to the Technical Services Department.

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 Thermo's expense, labor excluded. The 900 Series ULT Freezers include an additional two year warranty on the compressors, parts only, F.O.B. factory. The 8600 Series 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.

In addition to the standard warranty, the foamed-in-place cabinet design carries a unit production lifetime warranty (foamed-in-place cabinet, evaporator and foamed-in-place door; parts only). Please contact your sales representative or Thermo for additional information.

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. The Technical Services Department 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.

Thermo shall not be liable for any indirect or consequential damages including, without limitation, damages relating to lost profits or loss of products.

Your local Thermo 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 contact your local distributor or Thermo (1-800-438-4851 in 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. 4 4/09

Appendix A Handling Liquid Nitrogen

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

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 Because 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. s

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.

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! s

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. s

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.

Appendix B Handling Liquid CO₂

Warning High concentrations of CO₂ 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% CO₂). The short term exposure limit for 15 minutes or less is 30,000 PPM (3% CO₂). Carbon dioxide monitors are recommended for confined areas where concentrations of carbon dioxide gas can accumulate. s

Store and use liquid CO₂ 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 CO₂ gas reduce the concentration of oxygen and can result in asphyxiation. Because CO₂ 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 CO₂ is exposed to the air is condensed moisture; not the gas itself. The issuing gas is invisible. s

Never dispose of liquid CO₂ in confined areas or places where others may enter.

Disposal of liquid CO₂ 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.

Appendix C 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.

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Forma[®] Laboratory Freezer
Product Designations: 8602
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release Level [REL#] shown on Serial Tag)

This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


Dallas Kemper
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Forma[®] Laboratory Freezer
Product Designations: 8603
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release Level [REL#] shown on Serial Tag)

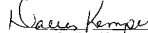
This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


Dallas Kemper
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Forma[®] Laboratory Freezer
Product Designations: 8604
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release Level [REL#] shown on Serial Tag)

This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


Dallas Kemper
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Laboratory Freezer
Product Designations: 8605
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release level [REL#] shown on Serial Tag)


This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


Dallas Kemper
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Forma® Laboratory Freezer
Product Designations: 8606
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release Level [REL#] shown on Serial Tag)

This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


Dallas Kemper
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Forma® Laboratory Freezer
Product Designations: 8607
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release Level [REL#] shown on Serial Tag)

This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


Dallas Kemper
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

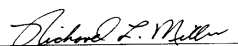
Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Blood Bank Freezer
Product Designations: 8627
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release level [REL#] shown on Serial Tag)

This product conforms to the following European Union Directive(s):

EMC: 89/336/EEC
MDD: 93/42/EEC

This product conforms to the following Harmonized, International and National Standards:

EMC:	MDD:
EN 61326-1:1997	EN 61010-1
EN 50081-1:92	Amendments 1 and 2
EN 50082-1:97	EN 60335-2-24 (applicable sections)
	CSA C22.2 No. 1010.1
	UL 471 (applicable sections)
	UL 61010A-1


Richard L. Miller, COE
Regulatory Compliance Manager

ThermoFisher
SCIENTIFIC

25 June 2007

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Forma® Laboratory Freezer
Product Designations: 8656
Year of Initial CE Marking: 2004
Affected Serial Numbers: Release 4
(Release Level [REL#] shown on Serial Tag)

This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


Dallas Kemper
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Laboratory Freezer
Product Designations: 8690
Year of Initial CE Marking: 2004
Affected Serial Numbers: Release 4
(Release level [REL#] shown on Serial Tag)

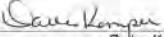
This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


04 November 2010
Dallas Kemper
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Forma® Laboratory Freezer
Product Designations: 8691
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release Level [REL#] shown on Serial Tag)

This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, international and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


03 August 2010
Dallas Kemper
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Forma® Laboratory Freezer
Product Designations: 8692
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release Level [REL#] shown on Serial Tag)

This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


03 August 2010
Dallas Kemper
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Forma® Laboratory Freezer
Product Designations: 8693
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release Level [REL#] shown on Serial Tag)

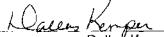
This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


03 August 2010
Dallas Kemper
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Forma® Laboratory Freezer
Product Designations: 8694
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release Level [REL#] shown on Serial Tag)

This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


Dallas Kemper 03 August 2010
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Declaration of Conformity

Manufacturer's Name: Thermo Fisher Scientific (Asheville) LLC
Manufacturer's Address: 401 Millcreek Road
Marietta, Ohio 45750
U.S.A.
Product Description: Forma® Laboratory Freezer
Product Designations: 8695
Year of Initial CE Marking: 2002
Affected Serial Numbers: Release 4
(Release Level [REL#] shown on Serial Tag)

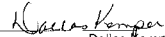
This product conforms to the following European Union Directive(s):

EMC: 2004/108/EC
LVD: 2006/95/EC

This product conforms to the following Harmonized, International and National Standards:

EMC:	LVD:
EN 61326-1:2006	EN 61010-1:2001
EN 61000-3-2	Amendments 1 and 2
EN 61000-3-3	EN 60335-2-89 (applicable sections)
	CSA C22.2 No. 61010-1
	CSA C22.2 No. 120 (applicable sections)
	UL 471 (applicable sections)
	UL 61010A-1 2 nd Edition

Declaration Date: 01 February 2009


Dallas Kemper 03 August 2010
Quality Engineer
FDA Official Correspondent for Regulatory Compliance

ThermoFisher
SCIENTIFIC

Thermo Fisher Scientific
Controlled Environment Equipment
401 Millcreek Road
Marietta, Ohio 45750
United States

www.thermofisher.com