

# **Thermo** Forma

**Model:  
916/917**

Ultra Low Temperature  
Upright Freezer  
17 cu. ft. capacity

Operating and Maintenance Manual

**Manual No. 7050916 Rev. 7**

**CAUTION!** All internal adjustments and maintenance must be performed by qualified service personnel.



**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.

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

**MANUAL NO. 7050916**

--	20775/SI-8506	6/3/02	Updated to 40°C Pepi thermostat from 43°C Klixon	ccs
7	20762/FR-1606	5/20/02	Added 'Normally Open' description to expansion tank solenoid valve	ccs
	20845/FR-1606	5/22/02	Updated refrigerant charge on schematics (oil return issue)	ccs
--	20684/FR-1596	3/28/02	Added sample pull-down curve after specs	ccs
6	20644/FR-1557	3/8/02	Added two expansion tanks to refig system	ccs
5	20251/FR-1526	2/19/02	Cordset connector	aks
4	20522/FR-1557	12/7/01	Updated 917 schematic to 12 FLA	ccs
3	20400/FR-1557	11/16/01	Added expansion tank with solenoid and relay timer (Release 11)	ccs
--	19363/FR-1429	10/18/00	New thermostat hot condenser temp reaches 43°C	ccs
2	19404/FR-1441	9/22/00	New temperature control board	ccs
1	19154/FR-1421	9/1/00	Updated parts and refrigeration drawings (dryer part number)	ccs
0	--	6/12/00	New Manual	aks
<b>REV</b>	<b>ECR/ECN</b>	<b>DATE</b>	<b>DESCRIPTION</b>	<b>By</b>

## Do You Need Information or Assistance on Thermo Forma Products?

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

1-740-373-4763	Direct
1-888-213-1790	Toll Free, U.S. and Canada
1-740-373-4189	FAX
<a href="http://www.thermoforma.com">http://www.thermoforma.com</a>	Internet Worldwide Web Home Page
<a href="mailto:service@thermoforma.com">service@thermoforma.com</a>	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 Forma products.

Whatever Thermo Forma 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 Forma  
Millcreek Road, PO Box 649  
Marietta, OH 45750

International customers, please contact your local Thermo Forma distributor.

# General Safety Notes used in this Manual

---



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

Ce symbole attire l'attention de l'utilisateur sur des instructions importantes de fonctionnement et/ou d'entretien. Il peut être utilisé seul ou avec d'autres symboles de sécurité. Lire attentivement le texte d'accompagnement.

Wichtige Betriebs- und/oder Wartungshinweise. Lesen Sie den nachfolgenden Text sorgfältig.

Importante instrucciones de operacion y/o mantenimiento. Lea el texto acompanante cuidadosamente.

---

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



Ce symbole attire l'attention de l'utilisateur sur des risques électriques potentiels. Seules des personnes qualifiées doivent appliquer les instructions et les procédures associées à ce symbole.

Gefahr von Stromschlägen. Nur qualifizierte Personen sollten die Tätigkeiten ausführen, die mit diesem Symbol bezeichnet sind.

Potencial de riesgos electricos. Solo personas das capacitadas deben ejecutar los procedimientos asociadas con este simbolo.

---

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



Risques potentiels liés à l'énergie. L'équipement en entretien ou en maintenance doit être éteint et mis sous clé pour éviter des blessures possibles.

Geräte, an denen Wartungs- oder Servicearbeiten durchgeführt werden, müssen abgeschaltet und abgeschlossen werden, um Verletzungen zu vermeiden.

El equipo recibiendo servicio o mantenimiento debe ser apagado y asegurado para prevenir danos.



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

Présence de surface(s) chaude(s) pouvant causer des brûlures sur la peau non protégée, ou sur des matières pouvant être endommagées par des températures élevées.

Heiße Oberfläche(n) können ungeschützter Haut Verbrennungen zufügen oder Schäden an Materialien verursachen, die nicht hitzebeständig sind.

Superficies calientes que pueden causar quemaduras a piel sin proteccion o a materiales que pueden estar danados por elevadas temperaturas.

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

## Table of Contents

### Section 1 - Installation and Start-Up

1.1 Unloading and Moving the Freezer	1-1
1.2 Getting to Know Your Freezer	1-1
1.3 Environmental Conditions	1-3
1.4 Installing the Wall Bumpers	1-3
1.5 Installing the Shelves	1-3
1.6 Location	1-3
1.7 Connecting the Alarm Battery	1-3
1.8 Remote Alarm Contacts	1-3
1.9 Optional Temperature Recorder	1-4
a. Connecting recorder battery	1-4
b. Installing the chart paper	1-4
1.10 Attaching the Power Cord	1-4
1.11 Power Switch (mains disconnect)	1-4
1.12 Water-cooled Condenser	1-4
1.13 Electrical Requirement and Connection	1-4
1.14 Factory Settings	1-4
1.15 Start Up and Loading	1-4

### Section 2 - Operation . . . . .2-1

2.1 Changing the Control Temperature Set Point	2-1
2.2 Changing the Over Temperature Alarm Set Point	2-1
2.3 Alarms	2-2
2.4 Silencing the Over Temperature Alarm	2-2

### Section 3 - Calibration . . . . .3-1

3.1 Temperature Display Calibration	3-1
3.2 Recorder Calibration	3-1
a. Changing the recorder range	3-1
b. Calibrating the chart recorder	3-1

### Section 4 - Routine Maintenance . . . . .4-1

4.1 Cleaning the Cabinet Exterior	4-1
4.2 Cleaning the Air Filter	4-1
4.3 Cleaning the Condenser	4-1
a. Cleaning the Water-cooled Condenser	4-1
4.4 Defrosting the Chamber	4-1
4.5 Cleaning the Door Gasket	4-1
4.6 Cleaning the Vacuum Relief Port	4-1
4.7 Replacing the Battery	5-1
4.8 Preparing the Unit for Storage	5-1

### Section 5 - Service . . . . .5-1

5.1 Servicing the Refrigeration System	5-1
5.2 Troubleshooting Chart	5-2

### Section 6 - Specifications . . . . .6-1

### Section 7 - Parts List . . . . .7-1

### Section 8 - Refrigeration Schematics . . . . .8-1

### Section 9 - Electrical Schematics . . . . .9-1

### Section 10 - Warranty . . . . .10-1

## Section 1 - Installation and Start-Up

### 1.1 Unloading and Moving the Freezer

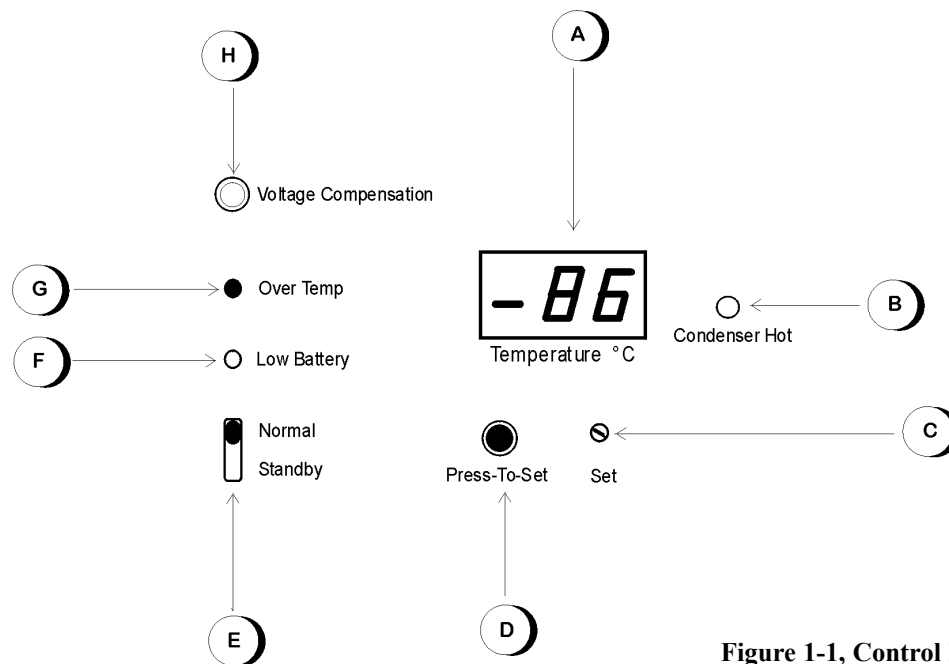
To remove the freezer from the pallet, use the 7/16" wrench to remove all the bolts securing the shipping bracket to the pallet.



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

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.

### 1.2 Getting to Know Your Freezer



**Figure 1-1, Control Panel**

**A** LED display - Shows actual chamber temperature and indicates the control temperature set point when the Press-To-Set button is pressed.

**B** Condenser Hot indicator - Lights when the thermostat on the condenser reaches approximately 40°C, indicating typically a dirty air filter, a clogged condenser, a fan failure or high ambient temperature.

**C** Set Point Adjustment screw - Used to set the control temperature set point.

**D** Press-To-Set button - Press to display control temperature set point.

The freezer can be easily pushed to the desired approved location, described in Section 1.6. If necessary, the doorstop may be removed to allow the door to swing fully open to move the unit through tight openings. When the freezer is in position, set the front caster brakes.



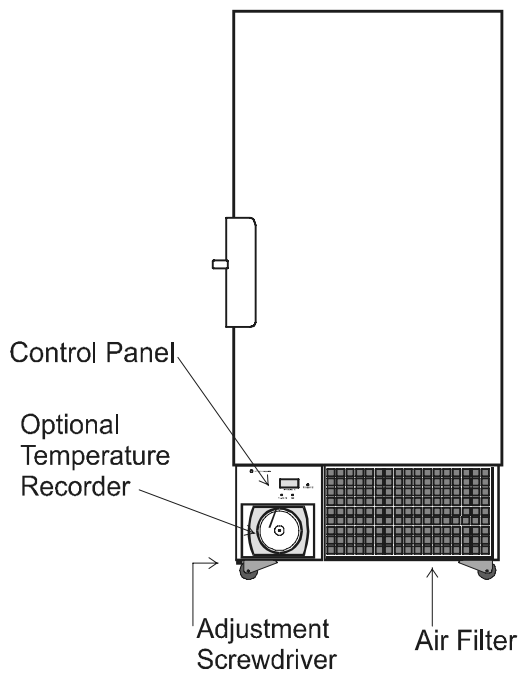
**The freezer must not be moved with the product load inside.**

**E** Normal/Standby switch - Silences the audible alarm.

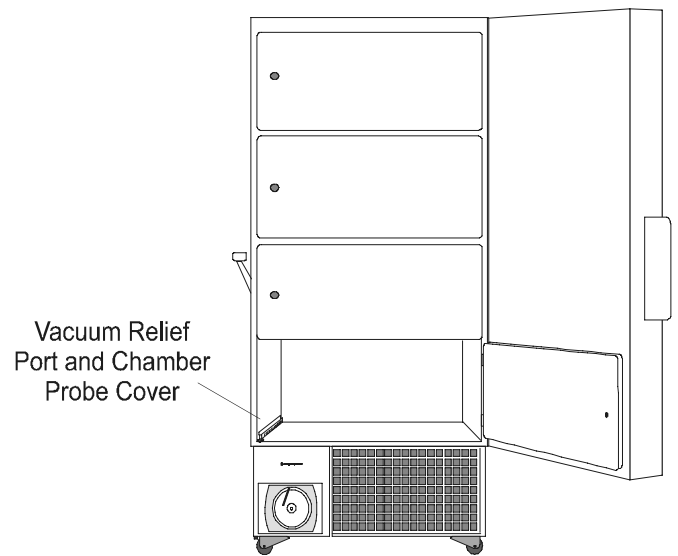
**F** Low Battery indicator - Flashes when the battery needs to be replaced.

**G** Over Temp indicator - Flashes when the chamber temperature rises above the Over Temperature Alarm setting.

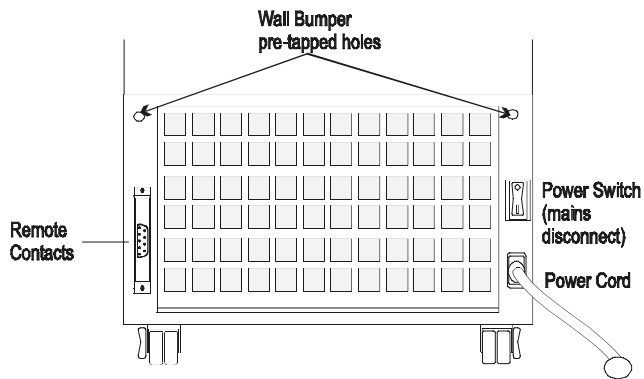
**H** Voltage Compensation indicator - Indicates when the incoming electrical power is being automatically adjusted to ensure that the compressor operates within specification.



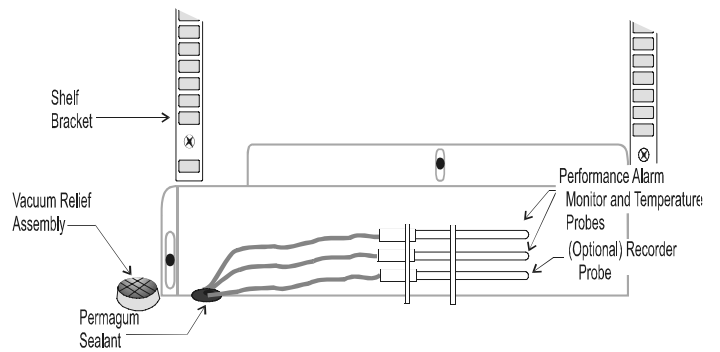
**Figure 1-2**  
**Freezer Exterior, front view**



**Figure 1-4, Vacuum Relief Port and Chamber Probe and Chamber Probe**



**Figure 1-3**  
**Freezer Base, rear view**



**Figure 1-5, Vacuum Relief Port and Chamber Probe Assembly (without cover)**

### 1.3 Environmental Conditions

The ULT Freezers are designed to be electrically safe in the following environmental conditions:

- Indoors
- Altitude: Up to 2,000 meters
- Temperature: 5°C to 43°C
- Humidity: 80% RH at or below 31°C, decreasing linearly to 50% RH at 40°C
- Mains Supply Fluctuations: ± 10% of nominal.
- Installation Category II <sup>1</sup>
- Pollution Degree 2 <sup>2</sup>
- Class of Equipment I
- Climatic Class T (Tropical)<sup>3</sup>

<sup>1</sup> 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.

<sup>2</sup> 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.

<sup>3</sup> Class T (Tropical) means that the freezers are electrically safe in a 43°C ambient.

### 1.4 Installing the Wall Bumpers (Refer to Figure 1-3)

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 1-3 for the locations of the pre-tapped holes.

### 1.5 Installing 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.

### 1.6 Location

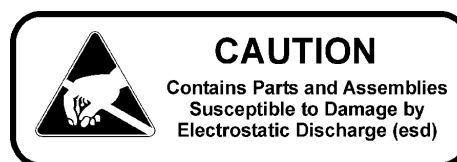
Locate the freezer on a firm, level surface in an area with an ambient temperature between 18°C and 32°C. Provide ample room to reach the mains disconnect switch (power switch) located on the rear of the freezer.



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

### 1.7 Connecting the Alarm Battery

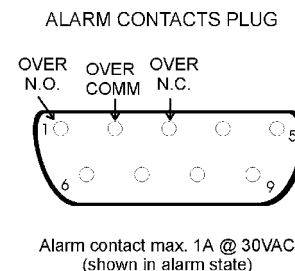
**NOTE:** The battery must be connected upon start-up so that the alarm will activate during an Overtemperature condition.



1. To gain access to the battery, remove the grill on the front of the freezer. The grill is attached to the freezer by friction plugs on each of the four corners and is readily pulled off. The battery is the rectangle fastened by Velcro to the compressor compartment frame and has a connector with red and black wires. Another red and black wire set with connector is secured to the wiring harness passing through the area of the freezer. Join the two wire connectors and replace the grill.
2. When the battery is connected, the OVERTEMP light will flash and the alarm will sound. Turn the NORMAL/STANDBY switch to the STANDBY position. This will silence the alarm. The alarm will remain silent until the unit is below alarm setpoint. The alarm setpoint is factory set for -65°C. If another setpoint is required, see Section 2.2.

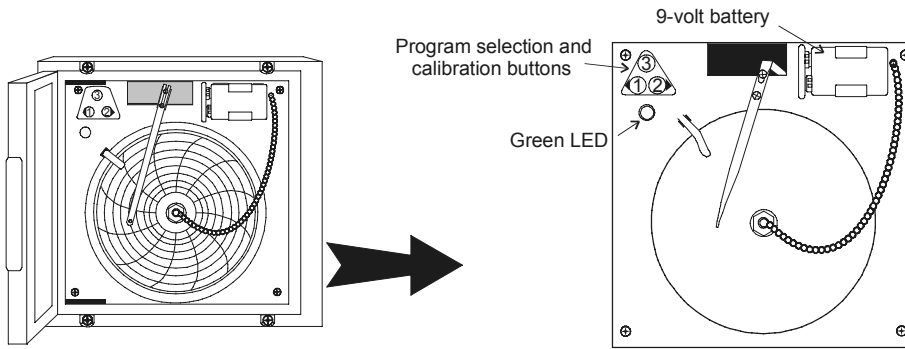
### 1.8 Remote Alarm Contacts

Remote alarm contacts are located on the rear of the unit. See Figure 1-6 for pin description, shown in the alarm state. The Thermo Forma part number for the plug required is 195482. Maximum rating for this plug is 1.0A @ 30VAC.



**Figure 1-6**  
**Remote Alarm Contact**





**Figure 1-8**  
Chart Recorder, detail view

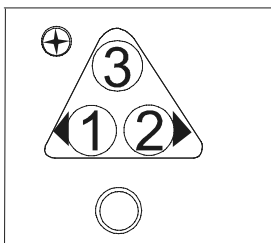
**1.9 Optional Temperature Recorder**

**a. Connecting recorder battery**

Open the glass door of the recorder and connect the 9-volt battery. The green light on the recorder will come on.

**b. Installing the chart paper**

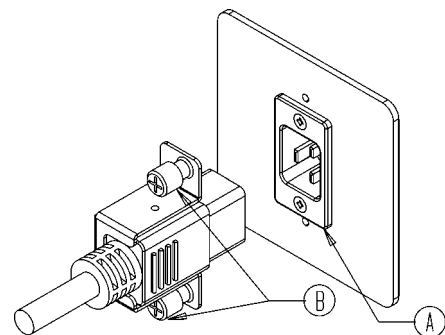
1. Open the glass door of the recorder and press button #3 until the pen begins to move outward.
2. Unscrew the knob at the center of the chart and remove the paper.
3. Install the new chart paper, position the paper to the correct time line and replace the knob.
4. Remove the cap from the felt pen and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



**Figure 1-9**  
Recorder Buttons

**1.10 Attaching the Power Cord**

1. Insert the power cord into the power outlet module (A).
2. Tighten screws (B) on the power cord retainer.



**Figure 1-10**  
Power Cord Assembly

**1.11 Power Switch (mains disconnect)**

The power switch is located on the rear of the unit, directly above the power cord. The power switch is also a circuit breaker that protects the entire unit.

**1.12 Water-cooled Condenser**

An accessory available for your freezer is the water-cooled condenser. Details for the use of this condenser follow:

**Water Supply**

- Connections:
  - Inlet - 1/2" compression
  - Outlet - 1/2" compression
- Maximum water pressure: 150 PSIG
- Water usage:
  - City water - 65°F (18°C), 1 GPM (3.4 liters)/HP
  - Tower water - 85°F (29°C), 3GPM (11.4 liters)/HP

**Water Strainer (P/N 780268)**

Included in the parts bag within the freezer is a water strainer recommended to be installed in the water inlet line. The connection size is 1/2" FPT.

**1.13 Electrical Requirement and Connection**

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.

**1.14 Factory Settings**

- Temperature: -80°C
- High Temperature Alarm: -65°C

If you wish to change any of these settings, see the appropriate section(s) in Section 2.

**1.15 Start Up and Loading**

- Turn the freezer on and allow it to run empty over night.
- When the empty freezer has stabilized over night at the control temperature set point, load the chamber with pre-frozen product.



**Make sure the power cord connection is completely seated.**



**The freezer was designed for the storage of pre-frozen product only. The addition of warm product may cause a temporary rise in the cabinet temperature.**

## Section 2 - Operation

### 2.1 Changing the Control Temperature Set Point

1. Remove the screwdriver on the front left corner of the control panel.
2. Press and hold in the PRESS-TO-SET key on the control panel.
3. Using the screwdriver, turn the SET screw until the desired temperature is displayed. Clockwise lowers the temperature and counter clockwise raises the temperature. The approved operating temperature range for the unit is  $-50^{\circ}\text{C}$  to  $-86^{\circ}\text{C}$ .

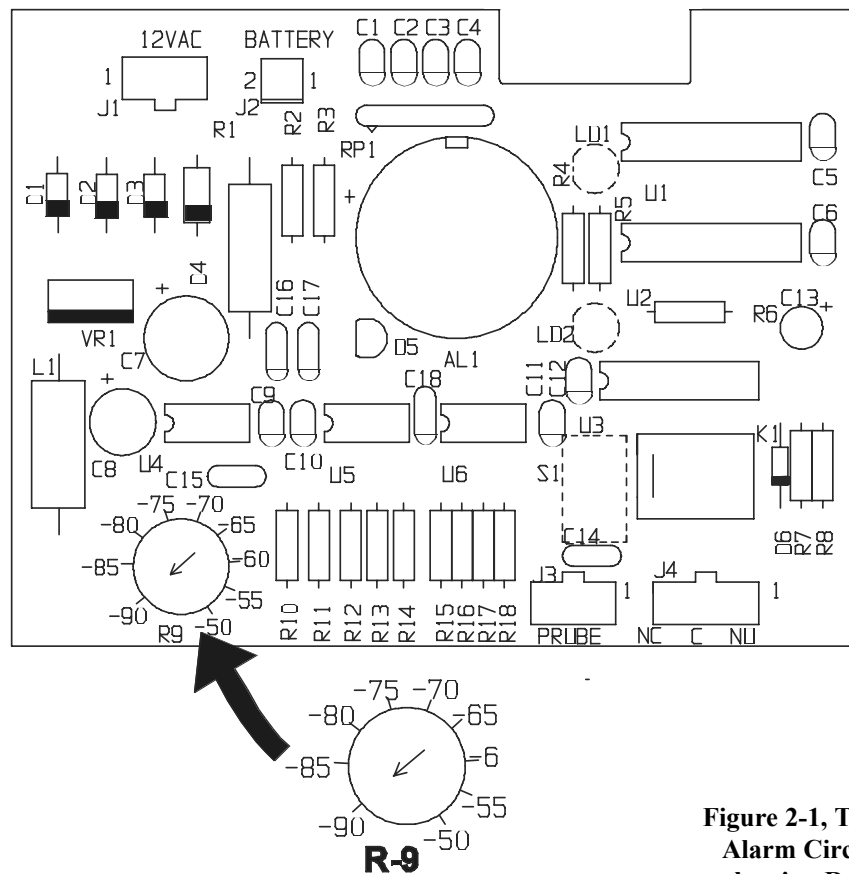


**The unit should NEVER be set to operate below  $-86^{\circ}\text{C}$ .**

### 2.2 Changing the Over Temperature Alarm Set Point

The following procedure sets the alarm for a precise temperature during initial set up. If a precise temperature is not required, set R9 to the desired temperature.

1. Set the unit to operate at the desired alarm point (see Section 2.1) and allow the unit to operate until it reaches the set point.
2. Remove the top two screws securing the control panel and loosen the bottom two screws. Carefully remove the control panel.
3. Remove the six screws securing the cover of the box attached to the inside of the control panel to expose the alarm board.
4. Turn the switch on the control panel to the NORMAL position.
5. If the unit alarms, slowly adjust R9 clockwise (warmer) until the alarm turns off. If the unit is not in alarm, slowly adjust R9 counterclockwise (colder) until the alarm turns on.
6. The alarm is now set for the desired temperature. Reinstall the box cover and secure with screws. Carefully install the control panel to its proper location and secure with screws.
7. Set the unit to the desired control temperature. See Section 2.1.



**Figure 2-1, Temperature Alarm Circuit Board showing R-9 Control**

## 2.3 Alarms

Alarm Description	Visual	Audible	Cause	Action Required
Over Temperature	Flashing (red)	On	<ul style="list-style-type: none"> <li>· Unit is above alarm setpoint</li> <li>· Addition of excessive product load</li> </ul>	<ul style="list-style-type: none"> <li>· Check freezer operation.</li> </ul>
Condenser Hot	On (red)*	---	<ul style="list-style-type: none"> <li>· Filter and/or condenser dirty.</li> <li>· Operating freezer in greater than 40°C ambient.</li> </ul>	<ul style="list-style-type: none"> <li>· See section 4.2 and 4.3 for cleaning filter and condenser.</li> </ul>
Low Battery	Flashing (red)	---	<ul style="list-style-type: none"> <li>· Rechargeable battery not connected.</li> <li>· Rechargeable battery needs replaced.</li> </ul>	<ul style="list-style-type: none"> <li>· See Section 1.6.</li> <li>· Replace battery. See section 4.7 for battery replacement instructions. See spare parts list for battery part number.</li> </ul>
Voltage Compensation	On (green)	---	<ul style="list-style-type: none"> <li>· Incoming voltage too high or too low</li> </ul>	<ul style="list-style-type: none"> <li>· Check if line voltage is within operating range. See 6 - Electrical. If within specification no action required.</li> </ul>

## 2.4 Silencing the Over Temperature Alarm:

- Move the switch from the NORMAL position to the STANDBY position. When the alarm condition has been corrected, the alarm will sound. The switch must then be returned to the NORMAL position.

**NOTE:** When the problem has been corrected and the alarm condition no longer exists an alarm will sound reminding the operator to press *Standby* to exit the STANDBY mode.



***If the unit is left in the STANDBY mode, none of the alarms listed will sound, alarm lights only will turn on.***

### Section 3 - Calibration



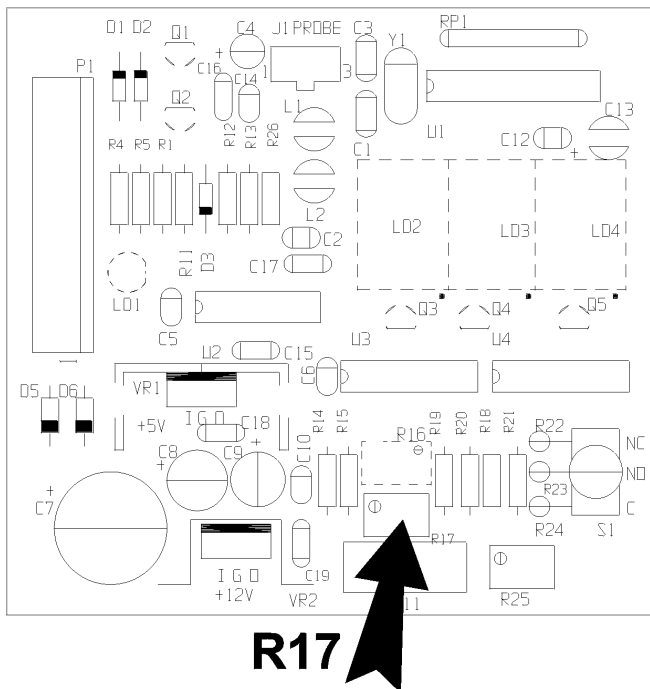
**Calibration must be performed when the unit is at operating temperature.**

**Required equipment:** Accurate low temperature remote bulb thermometer or thermocouple of known accuracy.

#### 3.1 Temperature Display Calibration

Place a measuring device near the probe cover. See Figures 1-4 and 1-5.

1. Allow the unit to stabilize at the operating temperature.
2. Remove the top two screws securing the control panel and loosen the bottom screws.
3. Carefully remove the control panel.
4. Remove the six screws securing the cover of box attached to the inside of the control panel. Remove the cover.
5. Adjust R17 until the control panel display matches the independent measuring device. Several turns of R17 may be required to achieve the desired temperature.
6. Reinstall cover to the box and secure.
7. Check the control temperature set point. Set point may have been altered during the calibration procedure.



**Figure 3-1, Temperature Control Board, R-17 Location**

#### 3.2 Optional Recorder Calibration

##### a. Changing the recorder 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.

NOTE: Changing ranges may require an offset calibration as outlined in Section 3.2.b.

Program No.	Range
Program 1	-40°C to 30°C
Program 2	0°C to 60°C
Program 3	-100°C to 38°C
Program 4	-5°C to 50°C
Program 5	0°C to 100°C
Program 6	-100°C to -200°C
Program 7	-115°C to 50°C
Program 8	-10°C to 70°C

**Table 3-1  
Recorder Range Chart**

##### b. Calibrating 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 from Thermo Forma. Refer to Parts List, Section 7.

# Thermo Forma

## PREVENTIVE MAINTENANCE

### SUVA Freezers

Your Thermo Forma 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.

Thermo Forma has 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 setpoint 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.

•Millcreek Road, Box 649 •Marietta, Ohio 45750 USA •740-373-4763  
•USA and Canada 888-213-1790 •Telefax: 740-373-4189 •email: [service@thermoforma.com](mailto:service@thermoforma.com)

### Preventive Maintenance for 900 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 probe location 4.5	Check and clean probe and cover, gaskets, hinges and doors of ice and snow.	<input checked="" type="checkbox"/>		
			<i>More frequent cleaning may be required, depending on use and environmental conditions.</i>	
4.2	Check air filter. Clean or replace as needed	<input checked="" type="checkbox"/>		
1.7	Check alarm back-up battery.	<input checked="" type="checkbox"/>		** Replace
--	Check condenser fan motor for unusual motor noise or vibration.		<input checked="" type="checkbox"/>	
--	* Verify and document calibration, at the minimum, annually.		<input checked="" type="checkbox"/>	
4.3	* 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.

## Section 4 - Routine Maintenance

### 4.1 Cleaning the 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.

### 4.2 Cleaning the Air Filter (minimum of twice a year\*)

1. Locate the grille on the right side of control panel. See Figure 1-2. Grasp the corner of the grille and gently pull to remove.
2. Remove the filter material and wash, using water and a mild detergent.
3. Dry by pressing between two towels.
4. Install the filter back into the grille and attach the grille.

\* Depending upon environmental conditions, the filter may need to be cleaned or replaced more frequently. If the filter becomes torn or excessively dirty, a replacement can be purchased from Thermo Forma. See the exploded parts list, Section 7, for filter part number. A filter kit (set of 5) part number 195517 is also available.

### 4.3 Cleaning the Condenser (minimum of twice a year\*)

1. Locate the grille on the right side of control panel. See Figure 1-2. Grasp the corner of the grille and gently pull to remove.
2. Using a vacuum cleaner, exercising care to not damage the condenser fins, clean the condenser.
3. Install the grille.

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

#### a. Cleaning 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.



**Do not use liquids that are corrosive to stainless steel or the brazing material (copper or nickel). Do not leave the unit on stand-by after cleaning.**

### 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.

### 4.4 Defrosting 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 the Alarm switch to Standby, or disconnect the battery.
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. Allow the freezer to operate empty overnight before reloading the product.

### 4.5 Cleaning the Door Gasket (minimum monthly\*)

Using a soft cloth, remove any frost build-up from the gasket and door(s).

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

### 4.6 Cleaning the Vacuum Relief Port (minimum monthly\*)

Using a soft cloth, remove any frost build-up from the vacuum relief, located in the front left corner of the chamber. See Figure 1-4 and 1-5.



**The vacuum relief port contains a small heating element. If the freezer is not disconnected from the electrical supply or turned off at the power switch, the heating element will continue to operate and may cause injury to personnel cleaning the freezer chamber.**

\*The vacuum relief port may need to be cleaned more frequently if dirt or excessive frost buildup prevents the door from closing properly.

#### 4.7 Replacing the Battery



***For a consistent and dependable charge, replace the battery every 2 years. Replacement batteries must be rechargeable and are available from Thermo Forma. 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.***

1. Turn off the power switch and unplug the unit from the AC power supply
2. To gain access to the battery, remove the grill on the front of the freezer. The grill is attached to the freezer by friction plugs on each of the four corners and is readily pulled off. The battery is the rectangle fastened by Velcro to the compressor compartment frame and has a connector with red and black wires. Another red and black wire set with connector is secured to the wiring harness passing through this area of the freezer.
3. Disconnect the two wire connectors, replace the battery pack and secure with the Velcro.
4. Reconnect the battery and replace the grill.
5. Plug unit in and turn power switch on.

#### 4.8 Preparing the Unit for Storage

Defrost the unit as described in Section 4.4. Disconnect the battery. This will prepare the unit for storage.

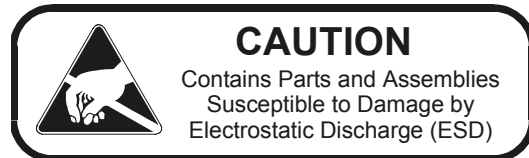


***The battery must be disconnected to prevent the battery from becoming completely discharged during storage.***

## Section 5 - Service



***Servicing must only be performed by service personnel who are qualified to repair cascade refrigeration systems. Always use standard safety practices when servicing the equipment.***



### 5.1 Servicing the Refrigeration System

Before opening the refrigeration system, use the troubleshooting chart to check out the electrical system. Electrical schematics and refrigeration drawings with parts are included with this manual.

Refer to the troubleshooting chart on the following page.



## 5.2 Troubleshooting Chart

Symptom	Possible Cause
No Temperature Display	<ul style="list-style-type: none"> <li>• Power line cord disconnected or not properly installed</li> <li>• External power circuit breaker tripped/open</li> <li>• Main power switch OFF</li> <li>• 24 volt transformer fuses open.</li> </ul>
Chamber Temperature Deviates from Set Point	<ul style="list-style-type: none"> <li>• Too much warm product added</li> <li>• Door open too long</li> <li>• Inadequate air circulation</li> <li>• Calibration</li> <li>• Dirty condenser</li> <li>• High ambient temperature</li> </ul>
Too Much Frost Build-Up	<ul style="list-style-type: none"> <li>• Door not properly sealed.</li> </ul>
Freezer not Being Refrigerated (unit is receiving power)	<ul style="list-style-type: none"> <li>• Compressor thermal overload open</li> <li>• Loss of refrigerant in either system</li> <li>• Defective compressor(s)</li> <li>• Defective temp control</li> <li>• Defective high pressure cut-off</li> </ul>
Display Problems in General	<ul style="list-style-type: none"> <li>• Defective temp control board.</li> </ul>
Condenser Hot Light On	<ul style="list-style-type: none"> <li>• Clogged air filter</li> <li>• Clogged condenser</li> <li>• Fan failure</li> </ul>

---

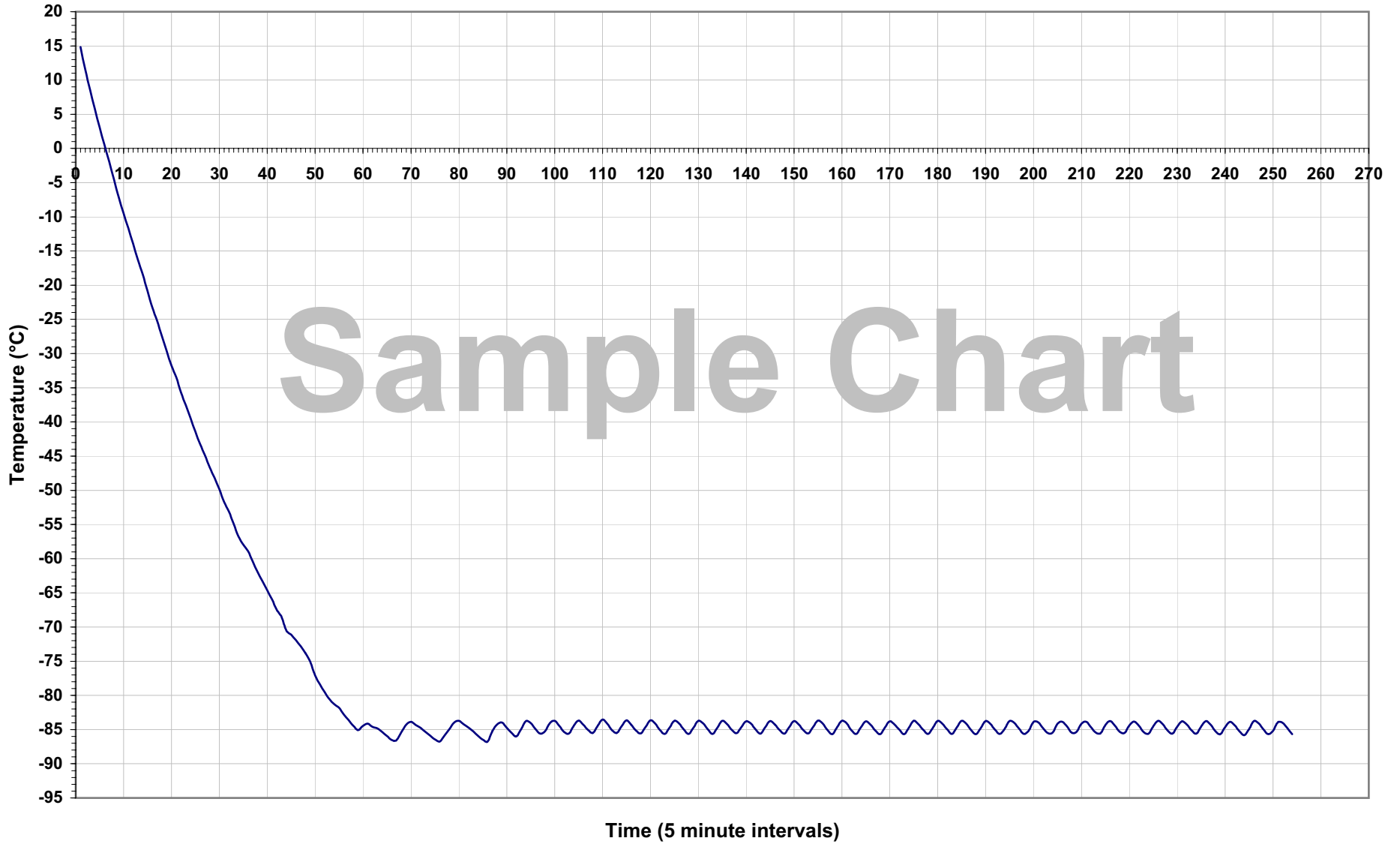
## Section 6 - Specifications

**MODEL 916/917**

<b>Temperature Range</b>	-50°C (-58°F) to -86°C (-123°F)		
<b>Exterior Dimensions</b>	33.3" W x 77.9" H x 37.0" F-B (84.6cm x 197.9cm x 94.0cm)		
	Add 3.0" (7.6cm) to W for latch/hinge; add 6.5" (16.5cm) to F-B for control panel/wall spacer.		
<b>Interior Dimensions</b>	23.0" W x 51.5" H x 25.3" F-B (58.4cm x 130.8cm x 64.3cm)		
<b>Capacity</b>	17 cu. ft. (489.9 liters)		
<b>Refrigeration</b>	Cascade system, (2) hermetically-sealed compressors		
<b>Insulation</b>	Non-CFC, foamed-in-place urethane; 5.0" (12.7cm) cabinet; 4.5" (11.4cm) door		
<b>Electrical</b>	<b>916</b> 120VAC, 1 PH, 60 Hz, 16 FLA Operating Range, 108VAC - 130VAC		
	<b>917</b> 230VAC, 1 PH, 50/60 Hz, 12.0 FLA Operating Range: 208VAC - 240VAC		
<b>Breaker Requirements</b>	<b>916</b> , 20 Amp, 120V <b>917</b> , 15 Amp, 230V		
<b>Automatic Voltage</b>	<b>Low:</b>		
	Cut In: 110V	Cut Out: 115V	Volts Boost: 10
	Cut In: 210V	Cut Out: 220V	Volts Boost: 18
	<b>High:</b>		
	Cut In: 125V	Cut Out: 120V	Volts Buck: 10
	Cut In: 235V	Cut Out: 225V	Volts Buck: 18
<b>Shipping Weight</b>	<b>Ocean:</b> 1025 lbs. (464.91kg) <b>Air/Container:</b> 785 lbs. (356.1kg) <b>Motor:</b> 785 lbs. (356.1kg)		

**Thermo Forma - Representative Pulldown and Cycling Curve**

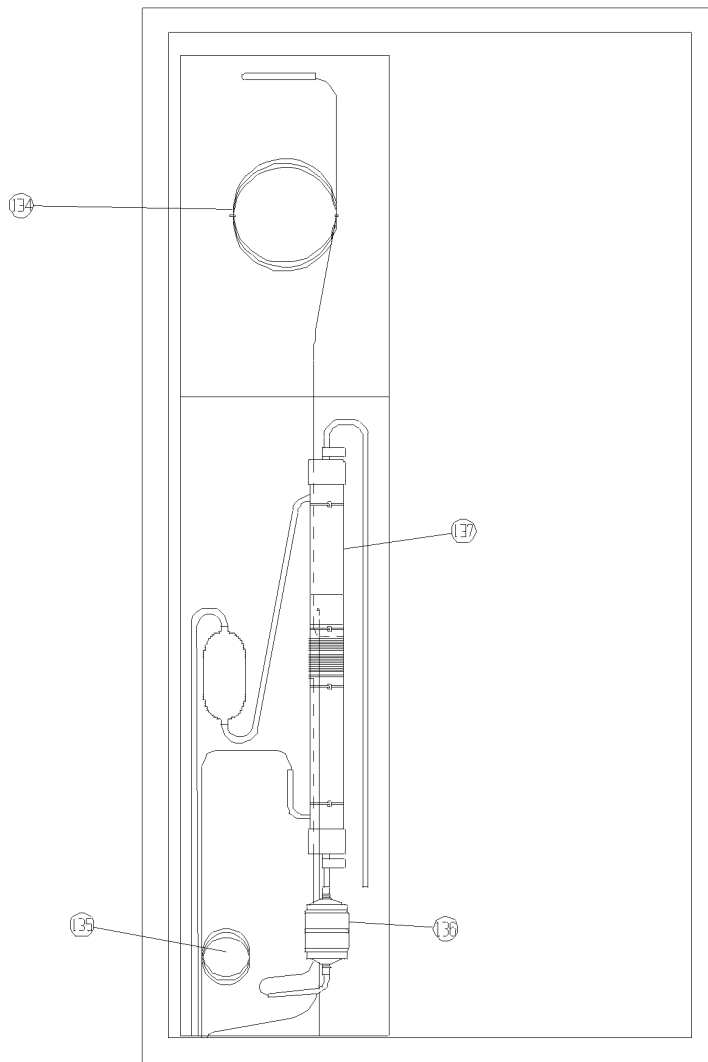
for 17 Ft.<sup>3</sup> -86°C Freezers



Sample Chart

6-2

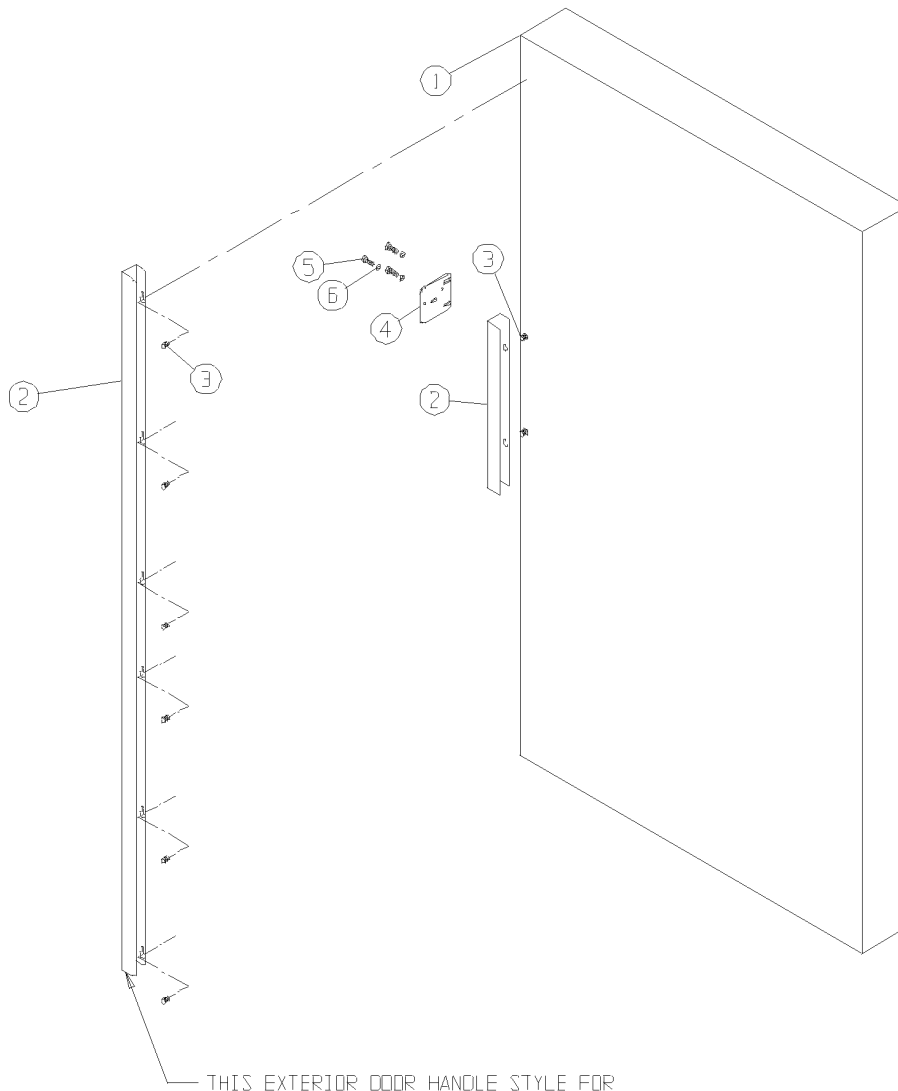
134	227928	Low stage cap. tube
135	227927	High stage cap. tube
136	209016	Dryer
137	211028	Heat exchanger



HEAT EXCHANGER ASSEMBLY

8516-206-0-0 REV. D  
Page 1 of 1

1	195570	Door (13 & 17 cuft)	18	195170	Interior door (4 door 13&17 cuft)
	195649	Door (23 cuft)		195171	Interior door (4 door 23 cuft)
	195650	Door (28 cuft)		195233	Interior door (5 door 13&17 cuft)
2	189934	Handle		195234	Interior door (5 door 23 cuft)
	195548	Handle (28 cuft)		195511	Interior door (4 door 28 cuft)
3	510022	#10-32 x 1/2 Hex head screw	19	285658	Knob (13, 17 & 23 cf)
4	121062	Can latch strike		120400	Knob (28 cf)
5	20003	1/4-20 x 3/4 Bolt	20	23043	Washer
6	23062	1/4 Star washer	21	195169	Latch
7	24042	#8-32 x 1/2 F screw	22	515083	Spacer
8	510303	Can latch	23	23044	Shoulder washer
9	189192	.063 Shim	24	23021	Flat washer
	189288	.125 Shim	25	23080	Spring washer
10	22115	#6-32 Screw	26	590008	#8-32 x 7/8 screw
11	23020	#6 Washer	27	23023	1/4 Flat washer
12	23015	#6-32 Cap nut	28	23033	1/4 Int. lockwasher
13	420308	Door gasket (13&17 cuft)	29	510305	Washer
	420309	Door gasket (23 cuft)	30	121032	Latch
	103104	Door gasket (28 cuft)	31	23057	Washer
14	116092	Door hinge	32	20058	1/4 x 3/4 Screw
15	189287	Door stop (23 & 28 cuft only)			
16	201120	Recorder			
17	245231	Recorder pen			



THIS EXTERIOR DOOR HANDLE STYLE FOR  
28 CF SINGLE DOOR FREEZER ONLY

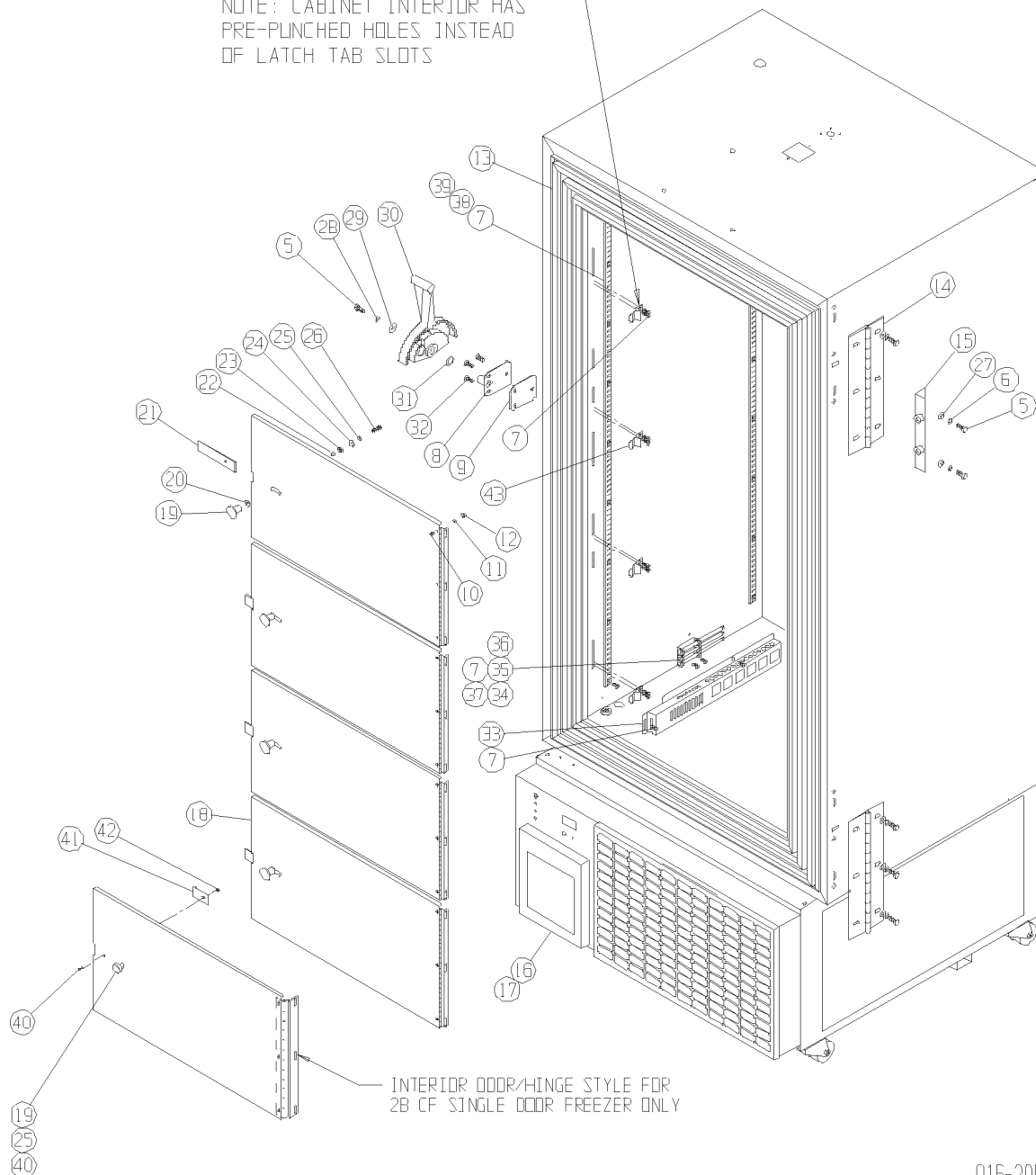
SINGLE DOOR CABINET ASSEMBLY

916-200-0-D REV. 4  
Page 1 of 2

- 33 189532 Probe Guard
- 34 195176 Probe mount
- 35 114020 Recorder probe grommet
- 36 30073 Control/alarm probe grommet
- 37 290165 Control/alarm probe
- 38 500177 Pilaster (900 only)
- 39 500175 Pilaster clip (flat) (900 only)
- 500184 Pilaster clip (tab) (900 only)
- 40 22051 8-32 x 1/4 SS PHP screw
- 41 195602 Interior door latch tab
- 42 23010 8-32 SS hex nut
- 43 195614 Interior door latch

- NOT SHOWN:
- B30011 Shelf (13 & 17 cuft)
  - B30025 Shelf (23 cuft)
  - B30028 Shelf (28 cuft)
  - 430263 Line cord (N. American 120V/20A)
  - 430264 Line cord (N. American 230V/15A)
  - 430270 Line cord (Continental Europe)
  - 430272 Line cord (Italy)
  - 430269 Line cord (Great Britain)
  - 430271 Line cord (India)
  - 430273 Line cord (Israel)
  - 430268 Line cord (Australia)

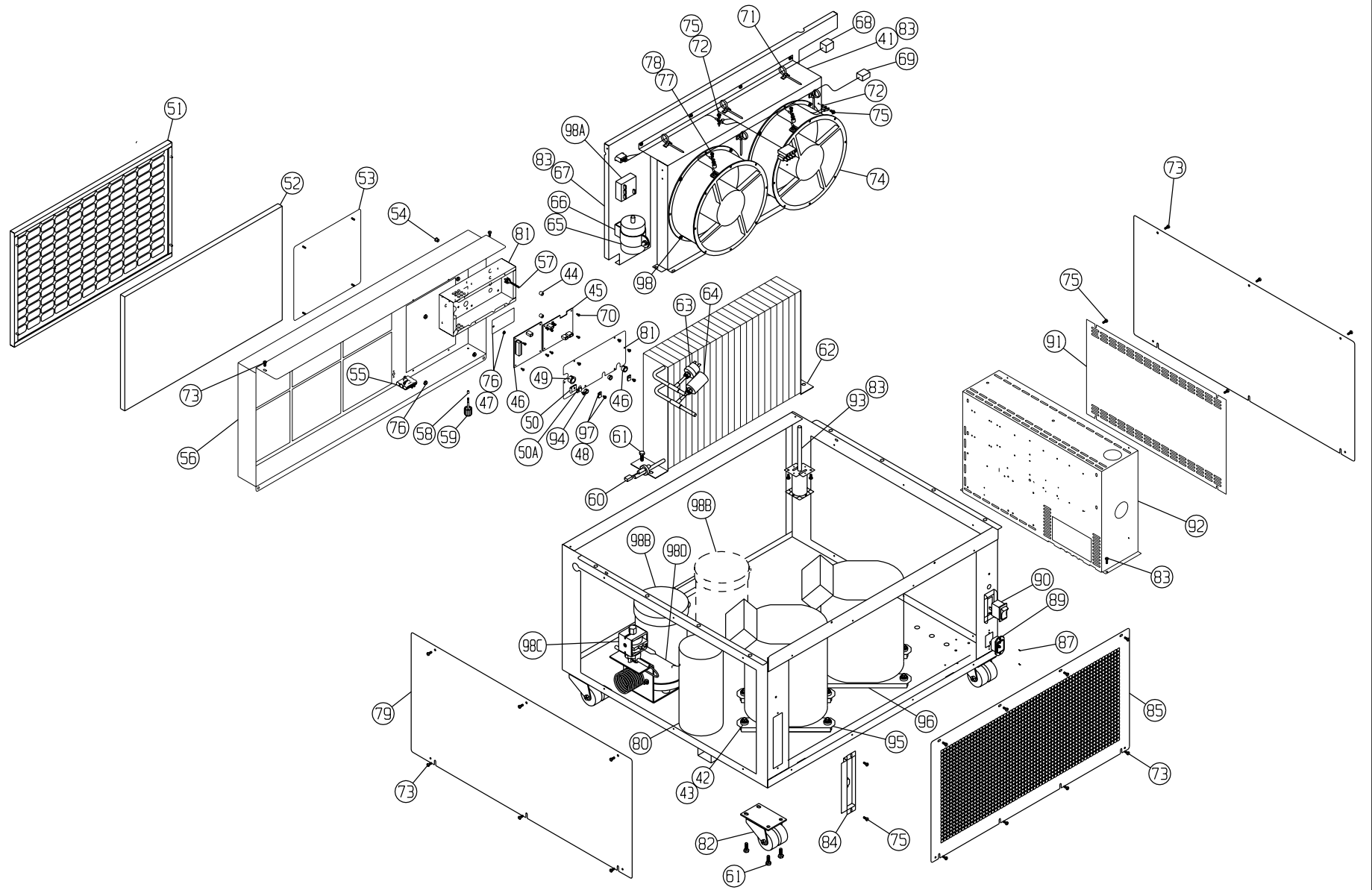
INTERIOR DOOR LATCH FOR 28 CF FREEZERS ONLY  
 NOTE: CABINET INTERIOR HAS PRE-PUNCHED HOLES INSTEAD OF LATCH TAB SLOTS



SINGLE DOOR CABINET ASSEMBLY

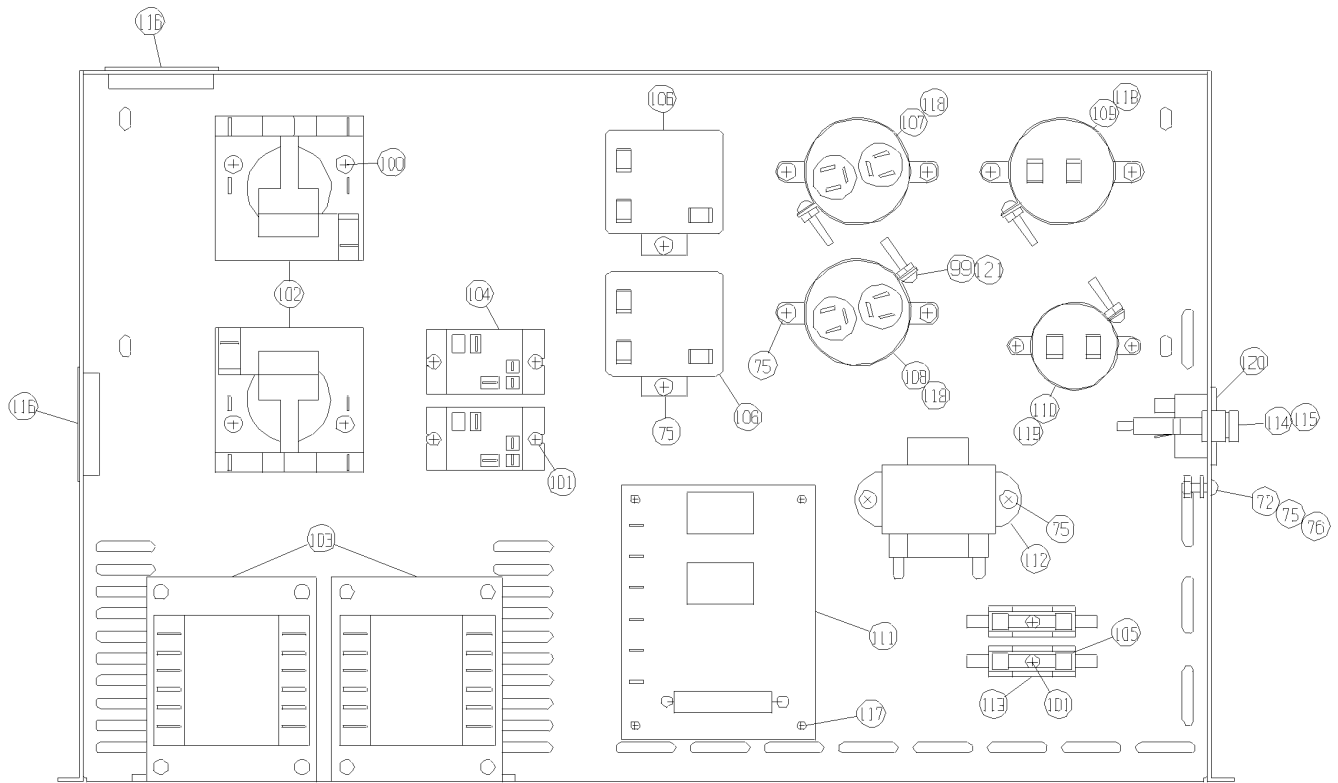


7-5



BASE ASSEMBLY



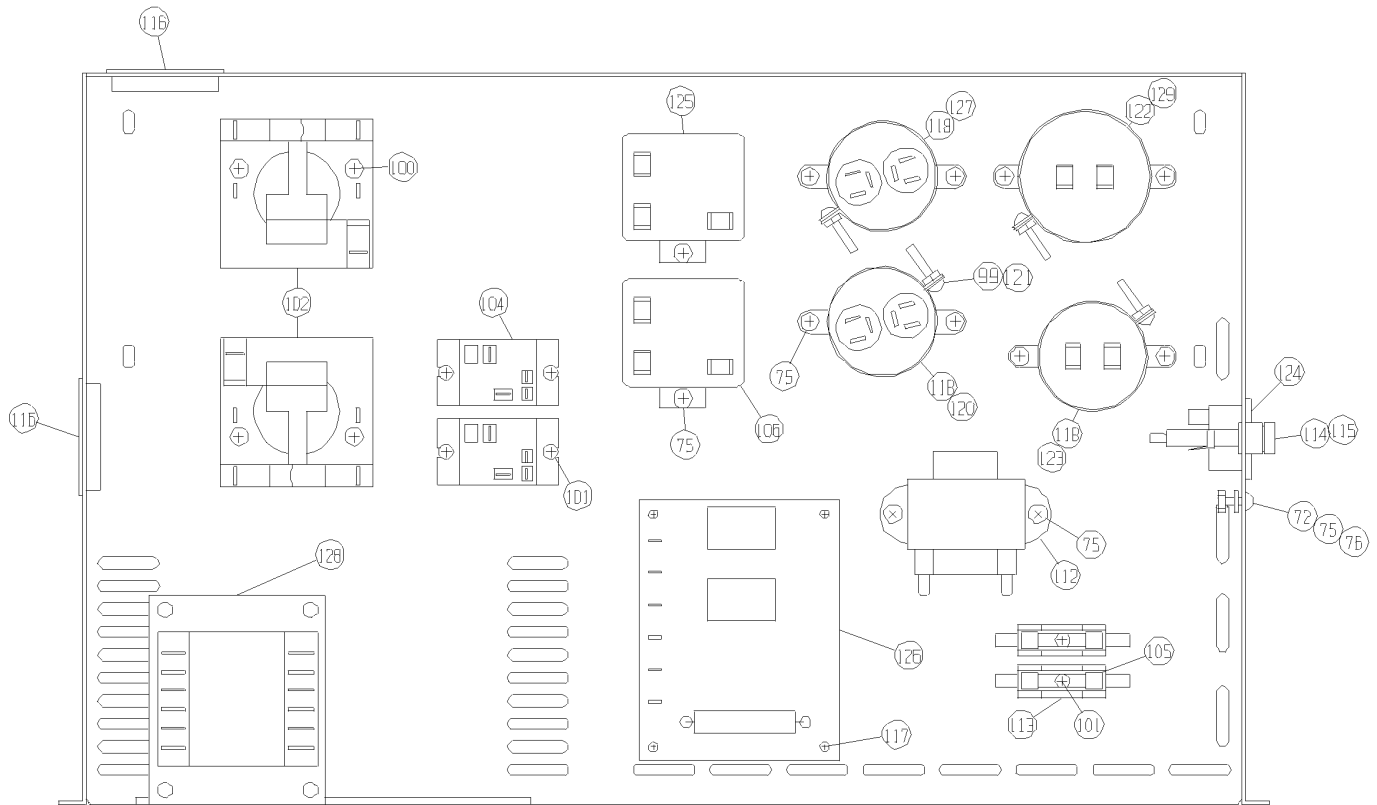


230 Volt relay enclosure

72	23059	#8 Ext. lockwasher	110	170159	Low stage start capacitor (N/A -40°C)
75	590020	#8-32 x 3/8 sens screw	111	191628	Boost/buck board
76	23002	#8-32 keps nut	112	420101	24V transformer
99	23001	#6-32 keps nut	113	230039	Fuse holder
100	22053	#8-32 x 1/2 screw	114	285632	Fuse holder
101	22049	#6-32 x 3/8 screw	115	230115	.15A slo blow fuse
102	300073	Boost & buck relay	116	330001	Snap bushing
103	420066	Transformer	117	59007	#4-40 x .375 screw
104	300261	Relay (Qty. 1, -40°C)	118	170091	1.75 dia. bracket
105	230110	1A slo blow fuse	119	170095	1.438 dia. bracket
106	300319	High & low stage start relay (H.S. only, -40°C)	120	460024	Outlet
107	170145	High stage run capacitor	121	22050	#6-32 x 1/2 screw
108	170149	Low stage run capacitor (N/A -40°C)			
109	170158	High stage start capacitor			

Relay Assembly

916-204-0-D REV. 3  
Page 1 of 2

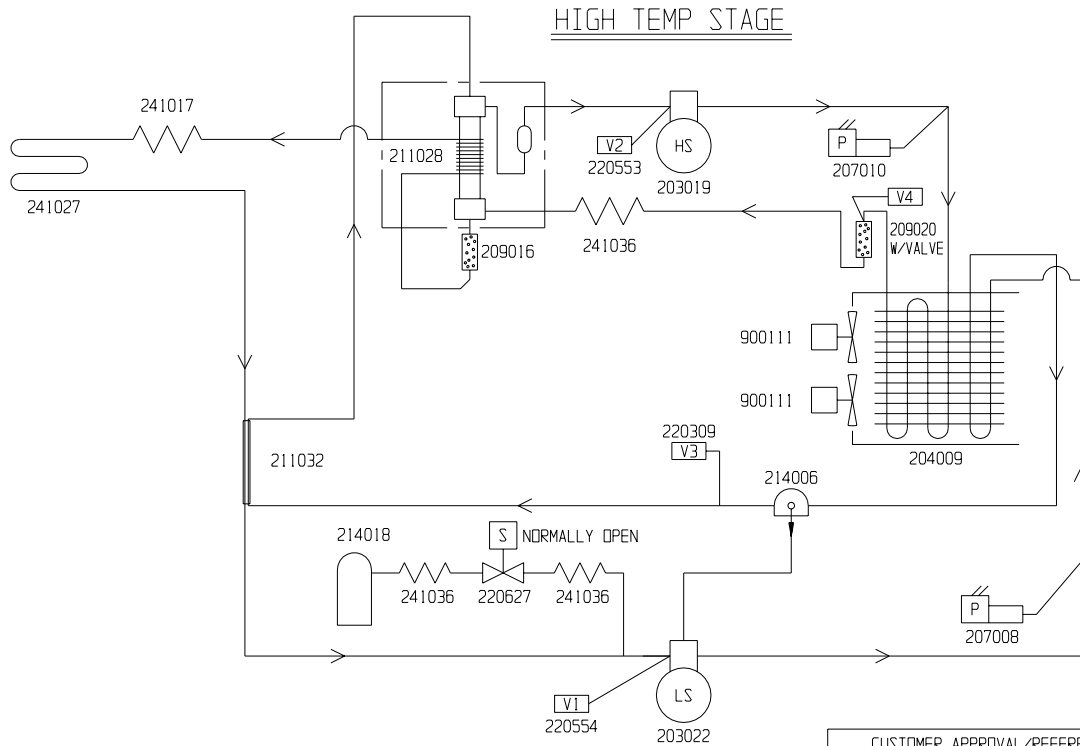


120 Volt relay enclosure

72	23059	#8 Ext. lockwasher	117	59007	#4-40 x 3/8 screw
75	590020	#8-32 x 3/8 sens screw	118	170091	1.75 dia. bracket
76	23002	#8-32 keps nut	119	170095	1.438 dia. bracket
99	23001	#6 keps nut	120	170160	Low stage run capacitor (N/A -40°C)
100	22053	#8-32 x 1/2 screw	121	24032	#8-32 x 3/8 F screw
101	22049	#6-32 x 3/8 screw	122	170157	2.5 dia. bracket
102	300073	Boost & buck relay	123	170010	Low stage start capacitor (N/A -40°C)
104	300261	Relay (Qty. 1, -40°C)	124	460024	Outlet
105	230110	1A slo blow fuse	125	300323	High Stage Start relay
106	300319	Low stage start relay (N/A -40°C)	126	190576	Boost/buck board
112	420101	24V transformer	127	170108	High stage run capacitor
113	230039	Fuse holder	128	420065	Transformer
114	285632	Fuse holder	129	170155	High stage start capacitor
115	230115	.15A slo blo fuse			
116	330001	Snap bushing			

Relay Assembly

916-204-0-D REV. 3  
Page 2 of 2



CASCADE REFRIGERATION

HIGH TEMPERATURE STAGE REFRIGERANT:

17 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:

17 CUFT UNITS: R-290 0.9 OZ. (26gr) MAX. ; VAC TO 3 PSIG  
 R-508B 15 OZ (425gr) ± 1/4 OZ. (7gr) OR  
 3 PSIG TO 134 PSIG ± 3 PSIG

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

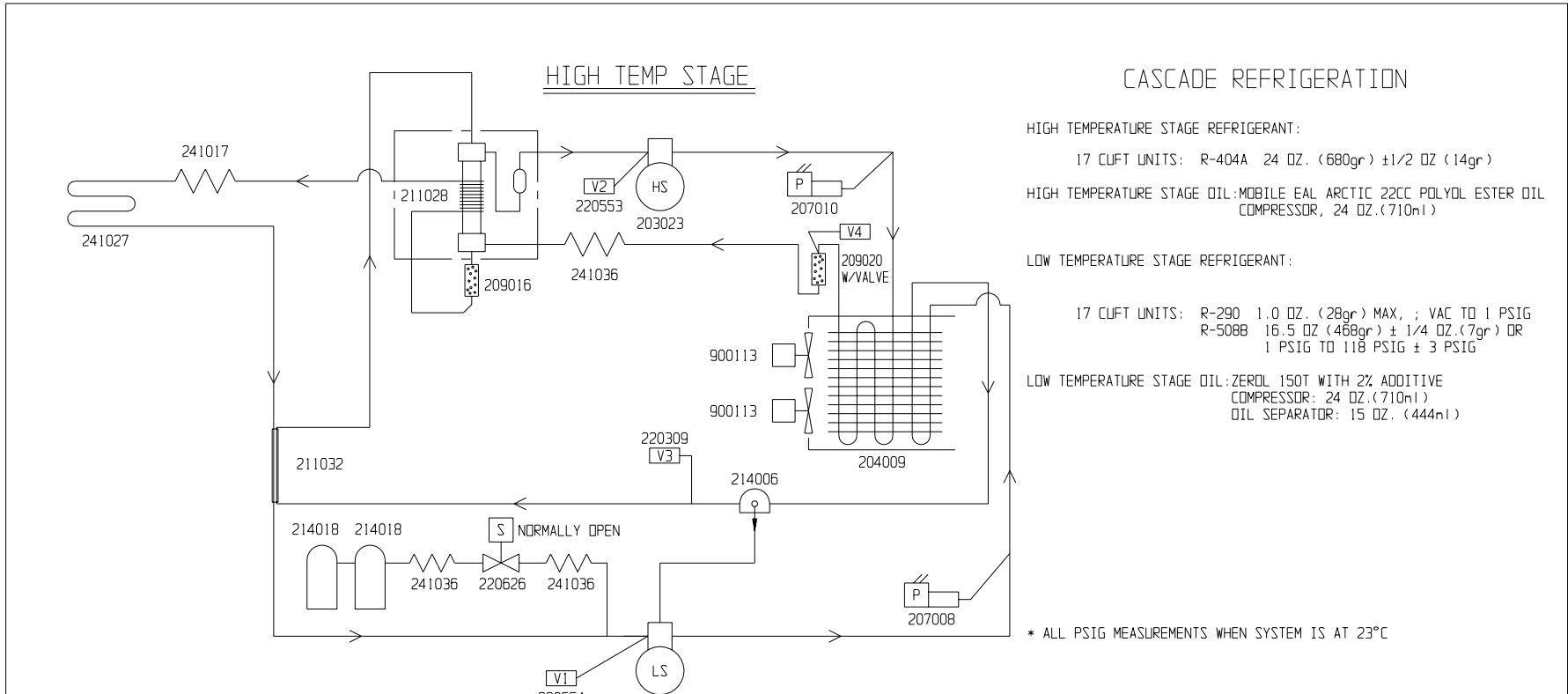
\* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C

LOW TEMP STAGE

1. COMMON TO: 717, 917, 982, 8517, 8582, 5467 & 5417 17 CU. FT. FREEZERS

CUSTOMER APPROVAL/REFERENCE						
APPROVED BY	3	FR-1616	05-21-02	DAS	KDG	REV. LOW STAGE REFRIG. CHARGES
APPROVING FIRM	2	FR-1606	05-08-02	RDS	KDG	NOTED SOLENOID AS NORMALLY OPEN
DATE OF APPROVAL	1	FR-1557	02-25-02	AFC	KDG	ADDED 241036 CAP. TUBE TO 220627
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 FORMA						
REV	ECR NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION
	0	FR-1557	11-12-01	LDC	KDG	RELEASED FOR PRODUCTION, REL. 11
DATE	04-21-98	DWN	LDC	CAD	LDC	APPD LON SCALE NONE
CUSTOMER						
JOB TITLE -86°C 17 CU. FT. UPRIGHT FREEZERS (230 V)						
DWG TITLE REFRIGERATION SCHEMATIC						
LOCATION		JOB NUMBER		DRAWING NUMBER		
				8517-91-0-B		

**Thermo Forma**  
 BOX 649, MARIETTA, OHIO 45750



CASCADE REFRIGERATION

HIGH TEMPERATURE STAGE REFRIGERANT:  
 17 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:  
 17 CUFT UNITS: R-290 1.0 OZ. (28gr) MAX, ; VAC TO 1 PSIG  
 R-508B 16.5 OZ (468gr) ± 1/4 OZ.(7gr) OR  
 1 PSIG TO 118 PSIG ± 3 PSIG

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

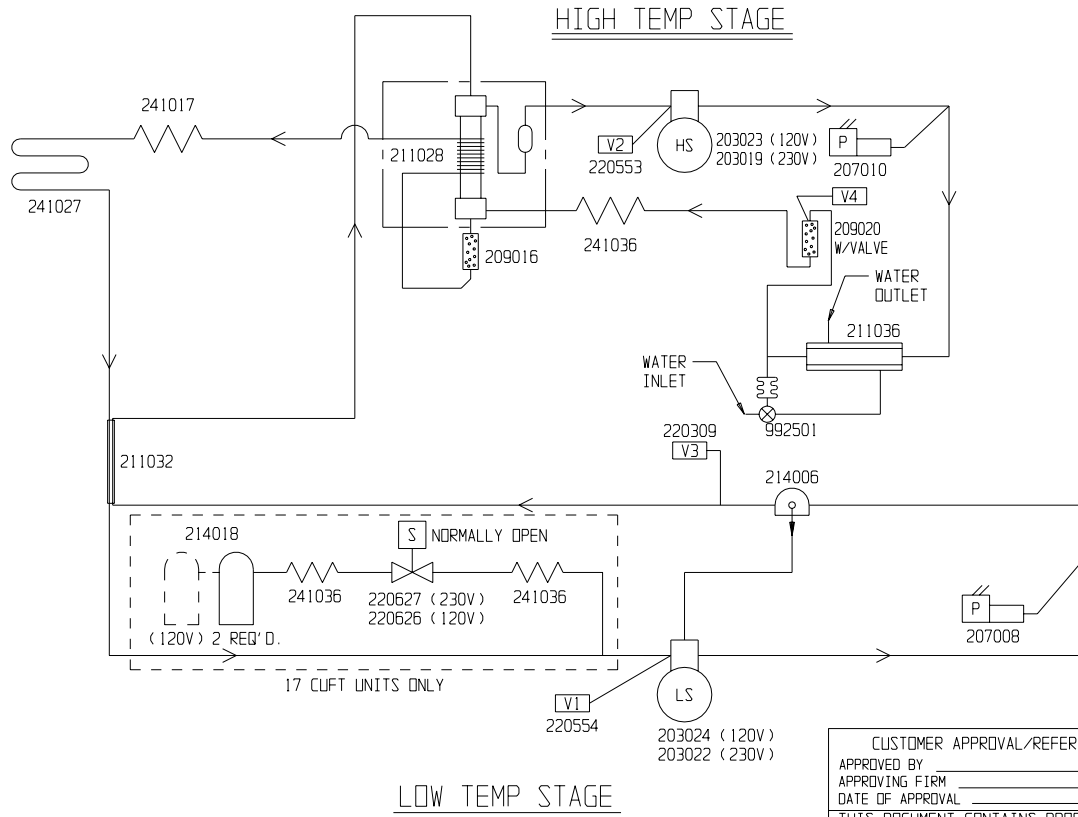
\* ALL PSIG MEASUREMENTS WHEN SYSTEM IS AT 23°C

LOW TEMP STAGE

1. COMMON TO: 716, 916, 981, 8516, 8581, 5466 & 5416 17 CU. FT. FREEZERS

CUSTOMER APPROVAL/REFERENCE							
APPROVED BY	_____						
APPROVING FIRM	_____						
DATE OF APPROVAL	2	FR-1616	05-21-02	DAS	KDG	LON	REV. LOW STAGE REFRIG. CHARGES
	1	FR-1606	05-08-02	RDS	KDG	LON	NOTED SOLENOID AS NORMALLY OPEN
	0	FR-1557	02-25-02	LDC	KDG	LON	RELEASED FOR PRODUCTION, REL. 11
REV	ECR NO.	DATE	BY	CAO	APPO	DESCRIPTION OF REVISION	
		04-21-98	DWN	LDC	CAO	LDC	APPO LON SCALE NDNE
CUSTOMER							
JOB TITLE -86°C 17 CU. FT. UPRIGHT FREEZERS (120 V)							
DWG TITLE REFRIGERATION SCHEMATIC							
LOCATION		JOB NUMBER		DRAWING NUMBER			
				8516-91-0-B			

**Thermo Forma**  
 BOX 649, MARIETTA, OHIO 45750



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)

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 OZ (369gr) ± 1/4 OZ. (.7gr) DR  
 20" TO 147 PSIG ± 3 PSIG

17 CUFT UNITS: R-290 1.0 OZ. (28gr) MAX, ; VAC TO 1 PSIG  
 (120 VOLTS) R-508B 16.5 OZ (468gr) ± 1/4 OZ. (.7gr) DR  
 1 PSIG TO 118 PSIG ± 3 PSIG

17 CUFT UNITS: R-290 0.9 OZ. (26gr) MAX, ; VAC TO 3 PSIG  
 (230 VOLTS) R-508B 15 OZ (425gr) ± 1/4 OZ. (.7gr) DR  
 3 PSIG TO 134 PSIG ± 3 PSIG

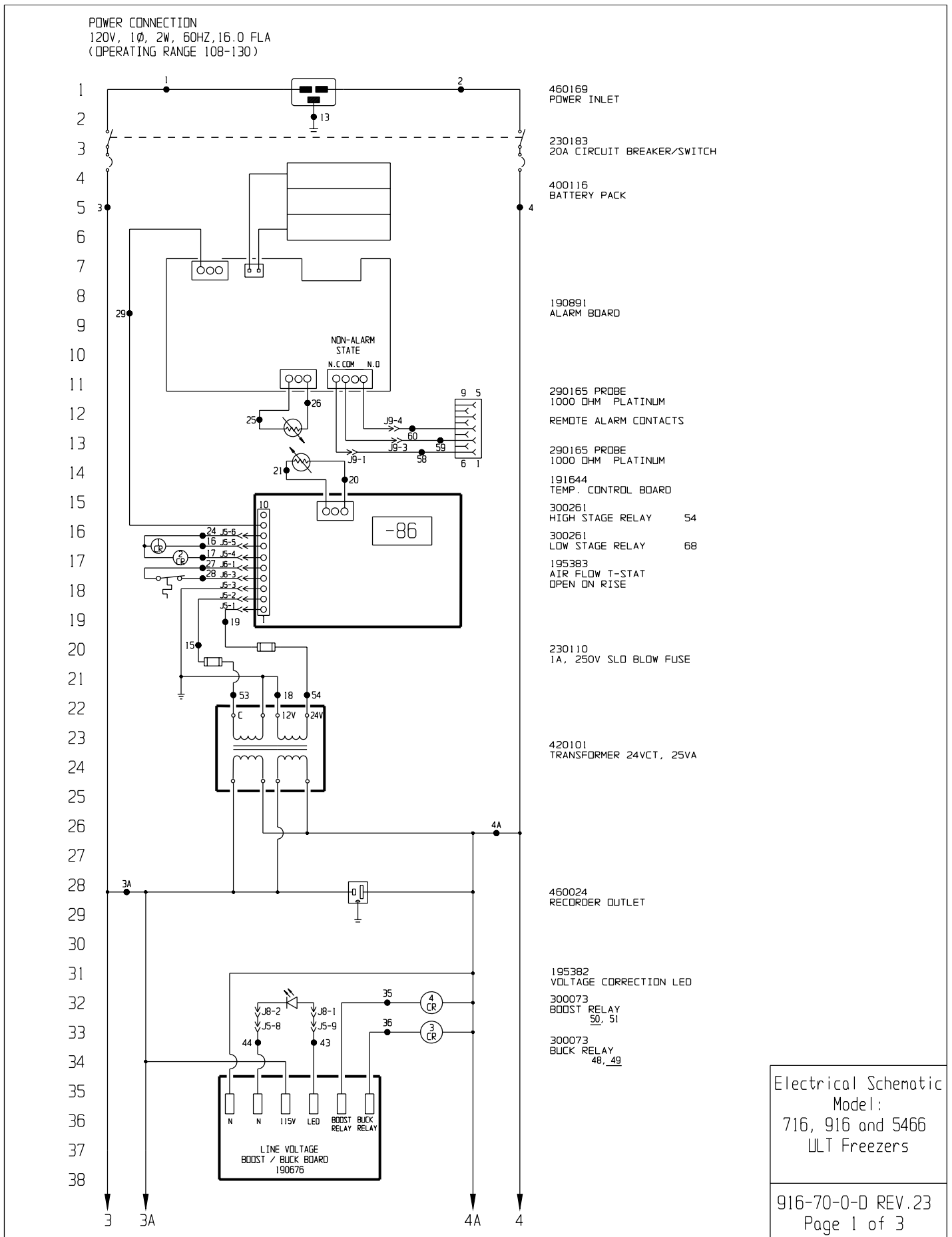
LOW TEMPERATURE STAGE OIL: ZEROL 150T WITH 2% ADDITIVE  
 COMPRESSOR: 24 OZ. (710ml)  
 OIL SEPARATOR: 15 OZ. (444ml)

\*\*\*SET WATER REGULATING VALVE FOR 170 PSIG

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

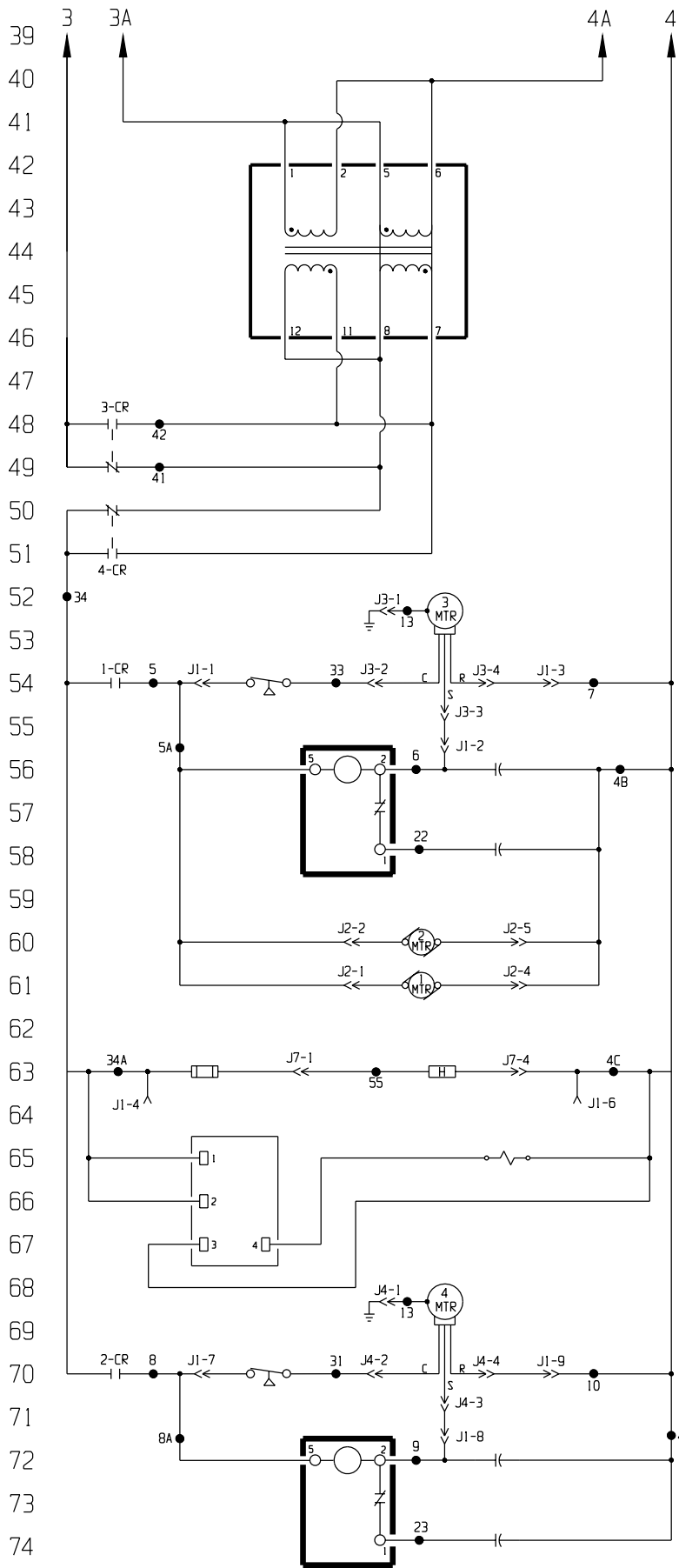
CUSTOMER APPROVAL/REFERENCE		6	FR-1616	05-21-02	DAS	KDG	LON	REV. LOW STAGE REFRIG. CHARGES
APPROVED BY	_____	5	FR-1606	05-08-02	RDS	KDG	LON	NOTED SOLENOID AS NORMALLY OPEN
APPROVING FIRM	_____	4	FR-1557	02-25-02	LDC	KDG	LON	ADDED EXPANSION TANK 17 CF 120V
DATE OF APPROVAL	_____	3	FR-1557	11-12-01	LDC	KDG	LON	ADDED EXPANSION TANK 17 CF 230V
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 FORMA								
REV	ECR NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION		
		12-10-99	DWN	CAW	CAW	APPD	CAW	SCALE NONE
CUSTOMER								
JOB TITLE -86°C 13 & 17 CU. FT. UPRIGHT H2O COOLED FREEZERS								
DWG TITLE REFRIGERATION SCHEMATIC								
LOCATION			JOB NUMBER			DRAWING NUMBER		
						195145-90-0-B		

**Thermo Forma**  
 BOX 649, MARIETTA, OHIO 45750



Electrical Schematic  
Model:  
716, 916 and 5466  
ULT Freezers

916-70-0-D REV.23  
Page 1 of 3



420065  
TRANSFORMER

203023  
HIGH STAGE COMPRESSOR  
120V

207010  
HIGH PRESSURE CUTOUT

300323  
START RELAY

170108  
RUN CAPACITOR  
35UF, 370V

170155  
START CAPACITOR  
189-227UF, 250V

900113  
TUBEAXIAL FAN

900113  
TUBEAXIAL FAN

230115  
0.15A SLO-BLOW FUSE

195153  
VACUUM RELIEF PORT HEATER  
12.5W @ 120VAC

220626  
SOLENOID VALVE

300359  
8 HR. TIMER/RELAY

203024  
LOW STAGE COMPRESSOR  
120V

207008  
HIGH PRESSURE CUTOUT

300319  
START RELAY

170160  
RUN CAPACITOR  
40UF, 370V

170010  
START CAPACITOR  
88-108UF, 250V

Electrical Schematic  
Model:  
716, 916 and 5466  
ULT Freezers

916-70-0-D REV.23  
Page 2 of 3

77  
78  
79  
80  
81  
82  
83  
84  
85  
86  
87  
88  
89  
90  
91  
92  
93  
94  
95  
96  
97  
98  
99  
100

WIRE REFERENCE CHART

WIRE #	GAUGE	COLOR
1	14	BRN
2	14	BLU
3	14	BLK
3A	14	BLK
4	14	BLU
4A	18	BLU
4B	18	BLU
4C	18	BLU
4D	18	BLU
5	14	BRN
5A	18	BRN
6	18	RED
7	14	YEL
8	14	BLK
8A	18	BLK
9	18	PUR
10	14	ORG
13	14	GRN/YEL
15	22	RED
16	22	YEL
17	22	ORG
18	22	BLK
19	22	WHT
20	22	RED
21	22	WHT
22	18	GRY
23	18	GRY
24	22	BLU
25	22	RED

WIRE REFERENCE CHART


WIRE #	GAUGE	COLOR
26	22	WHT
27	22	BLK
28	22	BLK
29	22	BLK
31	14	BLK
33	14	BRN
34	14	BRN
34A	18	BRN
35	18	ORG
36	18	YEL
41	14	RED
42	14	GRY
43	18	RED
44	18	BLU
48	18	BRN
53	22	RED
54	22	WHT
55	18	BRN
58	24	RED
59	24	BLK
60	24	WHT

REMOTE CONTACTS	
PIN# 1	NORMALLY CLOSED
PIN# 2	COMMON
PIN# 3	NORMALLY OPEN

CONTACT RATING: 1A @ 30V

23	FR-1557	02-18-02	AT	KDG	CCS	ADDED TIMER & SOLENOID (REL. 11)
22	FR-1441	09-11-00	RSB	KDG	LON	CHANGE 190860 TEMP. BOARD TO 191644
21	FR-1402	05-24-00	AT	GLS	LON	REVISED ALARM BD PWR CONNECTION
20	FR-1284	04-28-99	AT	KDG	LON	REVISED MODELS IN TITLE BLOCK
19	N/A	11-03-98	AT	AT	LON	SUVA & EMC UPDATES
REV	ECN NO.	DATE	BY	CAD	APPO	DESCRIPTION OF REVISION

Electrical Schematic  
Model:  
716, 916 and 5466  
ULT Freezers



**ATTENTION**  
OBSERVE PRECAUTIONS  
ELECTROSTATIC  
SENSITIVE DEVICES

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 FORMA

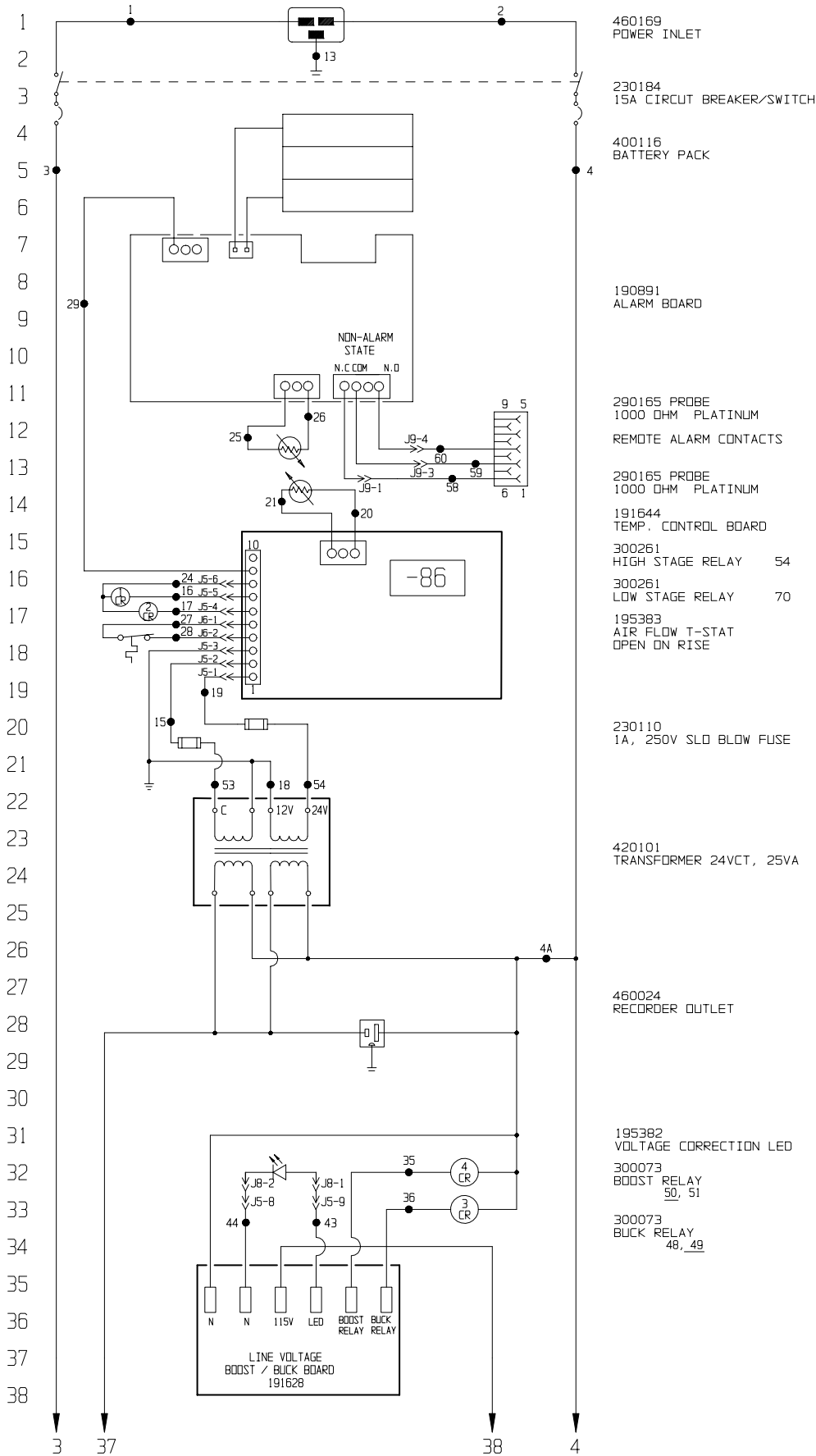
**Thermo Forma**  
BOX 649, HARRISBURG, OHIO 45750

MODEL/PART NAME: 716, 916 & 5466 UPRIGHT ULT FREEZERS			
DWG TITLE: ELECTRICAL SCHEMATIC			
OWN: AT	CAD: AT	APPO: CBL	DATE: 7-2-96
SCALE: NONE			
MATERIAL:			
PAINT COLOR:			
TOLERANCE UNLESS OTHERWISE SPECIFIED		DRAWING NUMBER	SIZE
ANGLES:	DECIMAL: .XX±	916-70-0	D

916-70-0-D REV.23  
Page 3 of 3



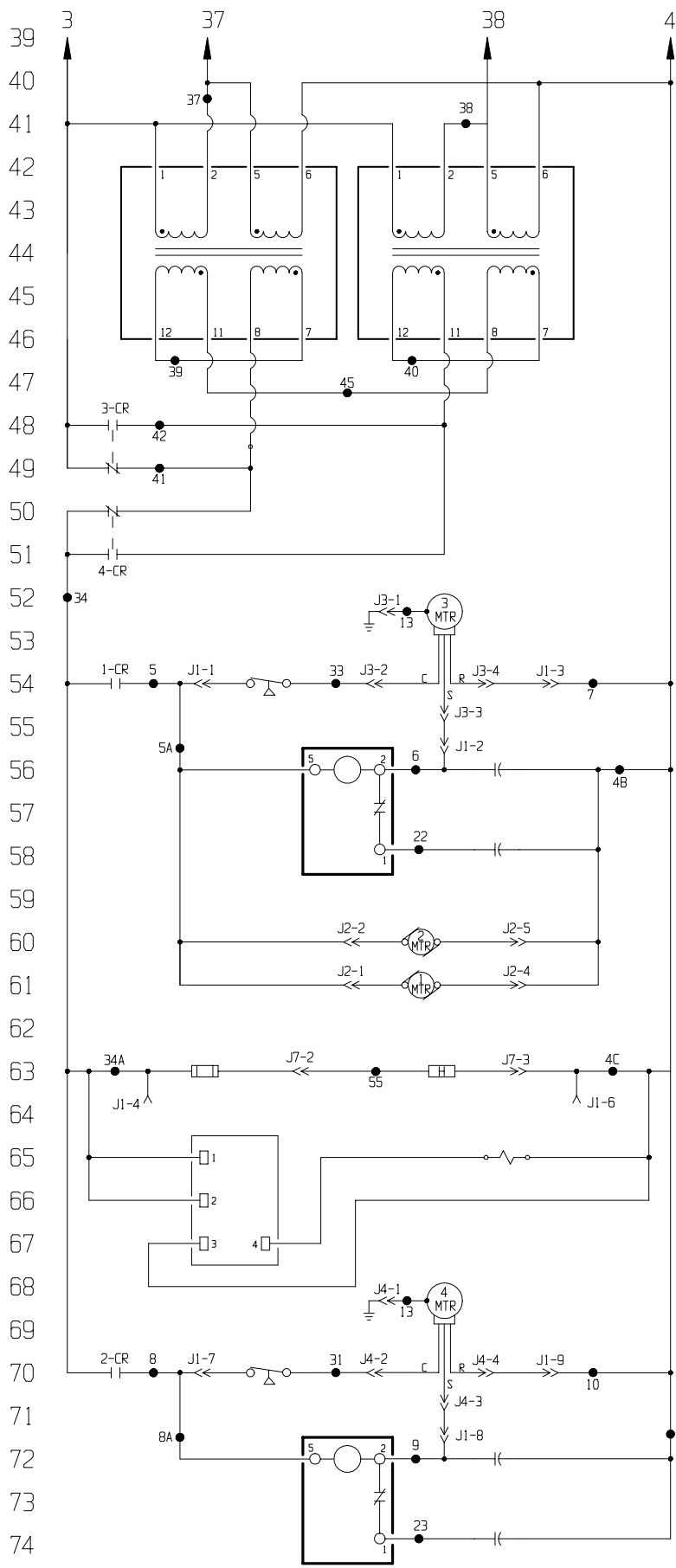
POWER CONNECTION  
 230V, 1 $\phi$ , 2W, 50/60HZ, 12.0 FLA  
 (OPERATING RANGE 208-240)



- 460169  
POWER INLET
- 230184  
15A CIRCUIT BREAKER/SWITCH
- 400116  
BATTERY PACK
- 190891  
ALARM BOARD
- 290165 PROBE  
1000 OHM PLATINUM
- REMOTE ALARM CONTACTS
- 290165 PROBE  
1000 OHM PLATINUM
- 191644  
TEMP. CONTROL BOARD
- 300261  
HIGH STAGE RELAY 54
- 300261  
LOW STAGE RELAY 70
- 195383  
AIR FLOW T-STAT  
OPEN ON RISE
- 230110  
1A, 250V SLD BLOW FUSE
- 420101  
TRANSFORMER 24VCT, 25VA
- 460024  
RECORDER OUTLET
- 195382  
VOLTAGE CORRECTION LED
- 300073  
BOOST RELAY  
50, 51
- 300073  
BUCK RELAY  
48, 49

Electrical Schematic  
 Model:  
 717, 723, 917, 923,  
 929, 5463 and 5467  
 Upright Freezer

917-70-0-D REV.30  
 Page 1 of 3



420066  
130VA TRANSFORMER, 2

203019  
HIGH STAGE COMPRESSOR  
230V

207010  
HIGH PRESSURE CUTOOUT

300319  
START RELAY

170145  
RUN CAPACITOR  
20UF, 440V

170158  
START CAPACITOR  
108-130UF, 250V

900111  
TUBEAXIAL FAN

900111  
TUBEAXIAL FAN

230115  
0.15A SLO-BLOW

195152  
VACUUM RELIEF PORT HEATER  
12.5W @ 230VAC

220627  
SOLENOID VALVE

300363  
8 HR. TIMER/RELAY

203022  
LOW STAGE COMPRESSOR  
230V

207008  
HIGH PRESSURE CUTOOUT

300319  
START RELAY

170149  
RUN CAPACITOR  
10UF, 370V

170159  
START CAPACITOR  
43-52UF, 250V

Electrical Schematic  
Model:  
717, 723, 917, 923,  
929, 5463 and 5467  
Upright Freezer

917-70-0-D REV.30  
Page 2 of 3

WIRE REFERENCE CHART

WIRE REFERENCE CHART

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

WIRE #	GAUGE	COLOR
1	14	BRN
2	14	BLU
3	14	BRN
4	14	BLU
4A	18	BLU
4B	18	BLU
4C	18	BLU
4D	18	BLU
5	14	BRN
5A	18	BRN
6	18	RED
7	14	YEL
8	14	BLK
8A	18	BLK
9	18	PUR
10	14	DRG
13	14	GRN/YEL
15	22	RED
16	22	YEL
17	22	DRG
18	22	BLK
19	22	WHT
20	22	RED
21	22	WHT
22	18	GRY
23	18	GRY
24	22	BLU
25	22	RED

WIRE #	GAUGE	COLOR
26	22	WHT
27	22	BLK
28	22	BLK
29	22	BLK
31	14	BLK
33	14	BRN
34	14	BRN
34A	18	BRN
35	18	DRG
36	18	YEL
37	18	BLK
38	18	BLK
39	14	DRG
40	14	DRG
41	14	RED
42	14	GRY
43	18	RED
44	18	BLU
45	14	YEL
48	18	BRN
53	22	RED
54	22	WHT
55	18	BRN
58	24	RED
59	24	BLK
60	24	WHT


REMOTE CONTACTS  
 PIN# 1 NORMALLY CLOSED  
 PIN# 2 COMMON  
 PIN# 3 NORMALLY OPEN

CONTACT RATING: 1A @ 30V

30	FR-1557	12-07-01	AT	KDG	AKS	CORRECT 10.0 FLA RATING TO 12.0 FLA
29	FR-1557	10-26-01	AT	KDG	LDN	ADDED TIMER & SOLENOID (REL. 11)
28	FR-1441	09-11-00	RSB	KDG	LDN	CHANGE 190860 TEMP. BOARD TO 191644
27	FR-1402	05-24-00	AT	GLS	LDN	REVISED ALARM BD PWR CONNECTION
26	FR-1357	01-31-00	AT	KDG	LDN	CHG. H.S. RUN CAPACITOR FROM 170097
REV	ECN NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION

Electrical Schematic  
 Model:  
 717, 723, 917, 923,  
 929, 5463 and 5467  
 Upright Freezer

**ATTENTION**  
 OBSERVE PRECAUTIONS  
 ELECTROSTATIC  
 SENSITIVE DEVICES



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 FORMA

**Thermo Forma**  
 BOX 649, WARRETTA, OHIO 45750

MODEL/PART NAME: 717, 723, 917, 923, 929, 5463 & 5467 UR FREEZERS

DWG TITLE: ELECTRICAL SCHEMATIC

DWN: AT	CAD: AT	APPD: CBL	DATE: 7-2-96	SCALE: NONE
---------	---------	-----------	--------------	-------------

MATERIAL:

PAINT COLOR:

TOLERANCE UNLESS OTHERWISE SPECIFIED	DRAWING NUMBER	SIZE
ANGLES: DECIMAL: .XX±	917-70-0	D

917-70-0-D REV.30  
 Page 3 of 3

## **THERMO FORMA 900 & 8500 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 year of the warranty period, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo Forma's expense, labor included. The 900 Series ULT Freezers include a second year warranty on the compressors, parts only, F.O.B. factory. The 8500 Series ULT Freezers include an additional four year warranty on the compressors, parts only, F.O.B. factory. Installation and calibration is not covered by this warranty agreement. The Thermo Forma Service 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, effective March 1, 2000, the foamed-in-place cabinet design carries a unit production lifetime warranty. Please contact your sales representative or Thermo Forma 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 one year warranty period. The Thermo Forma Service 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 Forma 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 Forma 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 Thermo Forma Service Office at 1-888-213-1790 (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.



## **THERMO FORMA 900 & 8500 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 Thermo Forma Service Dept.

During the first year of the warranty period, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo Forma's expense, labor excluded. The 900 Series ULT Freezers include a second year warranty on the compressors, parts only, F.O.B. factory. The 8500 Series ULT Freezers include an additional four year warranty on the compressors, parts only, F.O.B. factory. Installation and calibration is not covered by this warranty agreement. The Thermo Forma Service 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, effective March 1, 2000, the foamed-in-place cabinet design carries a unit production lifetime warranty. Please contact your sales representative or Thermo Forma 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 one year warranty period. The Thermo Forma Service 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 Forma 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 Forma 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 Forma (1-888-213-1790 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.



# Declaration of Conformity

**Manufacturer's Name:** Thermo Forma, Inc.

**Manufacturer's Address:** 401 Millcreek Road  
Marietta, Ohio 45750  
U.S.A.

**Product Description:** Laboratory Freezer

**Product Designations:** 917

**Year of Initial Marking (CE):** 1996

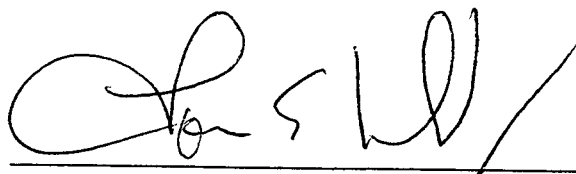
**Affected Units:** Release Level 11  
Release Level (REL#) shown on Serial Tag

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

**EMC:** 89/336/EEC  
**LVD:** 73/23/EEC

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

<b>EMC:</b>	<b>LVD:</b>
EN 61326-1:1997	EN 61010-1:1993
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



Louis E. Urschel, Jr.  
V. P. of Quality & Service

**Thermo** Forma

15 November 2001

**Thermo** Forma

**Millcreek Road, P.O. Box 649  
Marietta, Ohio 45750  
U.S.A.**

---

**Telephone: (740) 373-4763**

**Telefax: (740) 373-4189**