



-30°C Laboratory and Plasma Freezers V and W

Installation and Operation

325099H12 Rev. B

October 2014

Visit us online to register your warranty
www.thermoscientific.com/labwarranty

IMPORTANT Read this instruction manual. Failure to follow the instructions in this manual can result in damage to the unit, injury to operating personnel, and poor equipment performance.

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

Material in this manual is for informational 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 from or related to the use of this manual.

© 2014 Thermo Fisher Scientific Inc. All rights reserved.

Table of Contents

Introduction	1
Specifications	2
Intended Use.....	2
Safety Precautions	3
Unpacking	4
Temperature Monitoring	4
Installation	5
Location.....	5
Leveling	5
Wiring	5
Drawers.....	5
Door Seal	6
Solid Doors.....	6
Final Checks	6
Control Panel	7
Control Panel Features.....	7
Display Functions	9
Programming Functions.....	10
Service Parameters.....	11
Operation	12
Temperature Settings	12
Start Up	12
Defrost	13
Automatic Defrost	13
Alarm Systems	14
Operating the Alarm	14
Local and Remote Alarms	14
Installing a Remote Alarm (Optional)	15
Alarm Test	15
Chart Recorders	17
Set Up and Operation.....	17
Changing Chart Paper	18
Power Supply	18
Calibration Adjustment.....	19
Drawers	20
Removing the Drawers.....	20
Reinstalling the Drawers	21
Changing Drawer Position.....	21
Cleaning	22
Cleaning the Drawers.....	22
Cleaning the Condenser	22
Troubleshooting	23
Warranty Statement	24

1 Introduction

This manual provides installation and operation instructions for laboratory and plasma freezers with preset temperature setpoints of -30°C.

The control system, standard on all models, includes:

- Key-operated power and alarm switch
- Preset temperature setpoint
- Digital temperature display with 0.1°C resolution
- Graphic temperature display
- Audible and visual power failure indicators
- Alarm silence, ringback, and automatic reset functions
- SureTemp™ alarm system test

Other standard features include:

- Keyed door locks
- Remote alarm contacts
- CFC-free refrigerant
- CFC-free foamed in-place urethane insulation
- Quiet, hermetically sealed refrigeration compressors

1.1 Specifications

Environmental Operating Conditions

Pollution Degree 2
 Installation Category II
 Altitude 2000m MSL (mean sea level)
 Humidity max 80% - non-condensing
 Voltage Tolerance $\pm 10\%$
 Ambient Temperature 15-32°C (59-90°F)
 Product Usage Indoor Use Only

Power Requirements

Model Number	Voltage/Frequency/Phase	Amperage	Plug Type	Power Module Plug
430	V: 230/50/1 W: 230/50/1	12	CEE 7/7 BS 1363	IEE C19
1230	V: 230/50/1 W: 230/50/1	12	CEE 7/7 BS 1363	IEE C19
2330	V: 230/50/1 W: 230/50/1	12	CEE 7/7 BS 1363	IEE C19
3030	V: 230/50/1 W: 230/50/1	12	CEE 7/7 BS 1363	IEE C19
5030	V: 230/50/1 W: 230/50/1	12	CEE 7/7 BS 1363	IEE C19

Intended Use

Be sure to read this section to determine whether your freezer model is intended for use as a medical device.



Models with the following prefixes **are** intended for use as medical devices:

FFPF, UFP, PPL, UFP

Models with the following prefixes **are not** intended for use as medical devices and **are not** intended for in vitro diagnostic tests:

FFGL, ULT, JLE, IUF

2 Safety Precautions

In this manual and on labels attached to this product, the words **WARNING** and **CAUTION** mean the following:



WARNING : potential danger of electric shock which, if not avoided, could result in serious injury or death.



CAUTION : a situation which, if not avoided, may result in minor or moderate injury, impaired performance, or damage to the equipment.

Before installing, using or maintaining this product, please be sure to read this manual and product warning labels carefully. Failure to follow these instructions may cause this product to malfunction, which could result in injury or damage.

Below are important safety precautions that apply to this product:

- Use this product only in the way described in the product literature and in this manual. Before using it, verify that this product is suitable for its intended use. If the equipment is used in a manner not specified by the manufacturer, the protection provided by the equipment may be impaired.
- Do not modify system components, especially the controller. Use Thermo Scientific exact replacement equipment or parts. Before use, confirm that the product has not been altered in any way.
- Your unit must be properly grounded in conformity with national and local electrical codes. Never connect the unit to overloaded power sources.
- Disconnect the unit from all power sources before cleaning, troubleshooting, or performing other maintenance on the product or its controls.



WARNING: Do not store flammable materials in this unit..

3 Unpacking

At delivery, examine the exterior for physical damage while the carrier's representative is present. If exterior damage is present, carefully unpack and inspect the unit and all accessories for damage.

If there is no exterior damage, unpack and inspect the equipment within five days of delivery. If you find any damage, keep the packing materials and immediately report the damage to the carrier. *Do not return goods without written authorization.* When submitting a claim for shipping damage, request that the carrier inspect the shipping container and equipment.

4 Temperature Monitoring



IMPORTANT NOTE: We recommend the use of a redundant and independent temperature monitoring system so that the freezer can be monitored continuously for performance commensurate with the value of product stored.

5 Installation



CAUTION : Improper operation of the equipment could result in dangerous conditions. Follow all instructions and operate within design limits noted on the dataplate.

5.1 Location

Install the unit in a level area free from vibration with a minimum of 6 inches of space on the sides and rear and 12 inches at the top. Undercounter units require 1 inch of clearance on top.

Do not position the equipment in direct sunlight or near heating diffusers, radiators, or other sources of heat. The ambient temperature range at the location must be 59 to 90°F (15 to 32°C).

5.2 Leveling

The unit must be level. If the unit is out of level, you may need to shim the corners with thin sheets of metal.

5.3 Wiring

Before connecting your freezer to a power source, be sure to check the dataplate for correct voltage. “V” models use plug type CEE 7/7; “W” models use plug type BS1363. Wiring diagrams are attached to the back of the cabinet.



CAUTION : Connect the equipment to the correct power source. Incorrect voltage can result in severe damage to the equipment.



WARNING : For personal safety and trouble-free operation, this unit must be properly grounded before it is used. Failure to ground the equipment may cause personal injury or damage to the equipment. Always conform to the National Electrical Code and local codes. Do not connect unit to already overloaded power lines.

Always connect the equipment to a dedicated (separate) circuit. Electrical codes require fuse or circuit breaker protection for branch circuit conductors. Use time delay fuses for #12 AWG circuits.

5.4 Drawers

Plasma freezers come standard with drawers. Additional drawers are offered as available options.

5.5 Door Seal

Door seal integrity is critical for freezers. A loose fitting gasket allows moist air to be drawn into the cabinet, resulting in quicker frost buildup on the evaporator coil, longer running time, poor temperature maintenance, and increased operation cost.

To check the door seal, complete the following steps:

1. Open the door.
2. Insert a strip of paper (a couple of inches wide) between the door gasket and the cabinet flange and close the door.
3. Slowly pull the paper strip from the outside. You should feel some resistance.
4. Repeat this test at 4-inch intervals around the door. If the door does not seal properly, replace the gasket.

5.6 Solid Doors

Solid doors stay open if opened 90 degrees. Solid door spring tension cannot be adjusted.

5.7 Final Checks

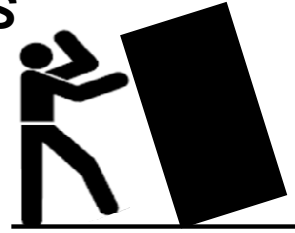
Before start up, complete the following steps:

1. Make sure that the unit is free of all wood or cardboard shipping materials, both inside and outside.
2. Verify that the unit is connected to a dedicated circuit.



WARNING: SAFETY INSTRUCTIONS

- REFRIGERATOR/FREEZER MUST BE SECURED BY THE ANTI-TIP BRACKET SUPPLIED
- UNLESS PROPERLY INSTALLED, REFRIGERATOR/FREEZER COULD TIP WHEN SHELVES/ DRAWERS ARE LOADED. INJURY AND DAMAGE TO EQUIPMENT AND CONTENTS MAY RESULT FROM REFRIGERATOR/ FREEZER TIPPING
- THIS REFRIGERATOR/FREEZER HAS BEEN DESIGNED TO MEET ALL RECOGNIZED INDUSTRY TIP STANDARDS FOR ALL NORMAL CONDITIONS.



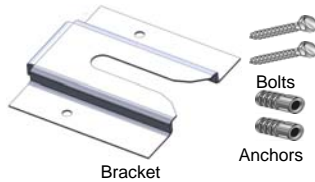
INSTALLATION INSTRUCTIONS:

Installation instructions are provided for wood and concrete floors. Any other type of construction may require special installation techniques as deemed necessary to provide adequate fastening of the Anti-Tip bracket to the floor. [For installation on floors other than wood or concrete, please contact technical support.](#)

The use of this bracket does not prevent the tipping of the Refrigerator/Freezer when not properly installed.

Materials Supplied

1. Anti-Tip Bracket (1)
2. 5/16" Lag Bolt (2)
3. Lag Screw Anchor (2), for concrete installation only
4. Bracket location template



Tools Required

Wood Floor

- Flashlight
- Tape Measure
- Drill
- 15/64" (6mm) Drill Bit
- 1/2" (13mm) Wrench
- 3/4" (19mm) Wrench

Concrete Floor

- Flashlight
- Tape Measure
- Hammer Drill
- 1/2" (13mm) Masonry Bit
- 1/2" (13mm) Wrench
- 3/4" (19mm) Wrench

Step 1

Locating the Bracket

- Determine where you want the centerline of the refrigerator/freezer to be
- Place the included template on the floor lined up with the centerline of the refrigerator/freezer and keep 6"-12" between the wall and the back of the unit
- On the floor, mark the location of Hole #1 & Hole #2 (also Hole #3 & Hole #4 for 50ft³ & 75ft³ models).

Step 2

Anti-Tip Bracket Installation

Wood Construction

- Drill 15/64" (6mm) pilot holes in locations marked in step 1
- Place bracket on floor aligned with holes
- Use supplied lag bolts to attach bracket to floor

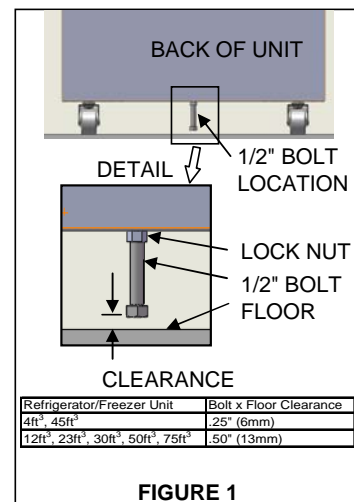
Concrete Construction

- Drill 1/2" (13mm) holes in locations marked in step 1 with masonry bit
- Slide Lag Screw Anchors into holes to be flush with floor surface
- Place bracket on floor aligned with holes
- Use supplied lag bolts to attach bracket to floor

Step 3

Adjusting Bolt in Refrigerator/Freezer

- Locate 1/2" bolt attached to bottom of cabinet
- Unscrew 1/2" bolt until there is the required clearance between floor and head of bolt as shown in Figure 1
- Tighten lock nut against bottom of unit



Step 4

Refrigerator/Freezer Positioning

- Line up 1/2" bolt installed in Step 3 with anti-tip bracket
- Roll or slide Refrigerator/Freezer into position until bolt stops against bracket
- Lock the casters

Step 5

Checking the Installation

- Complete the installation of the Refrigerator/Freezer per the installation instructions provided with the product.
- Check to see if the Anti-Tip bracket is installed properly by shining light under cabinet and confirming bolt in cabinet is secured by bracket on floor

6 Control Panel

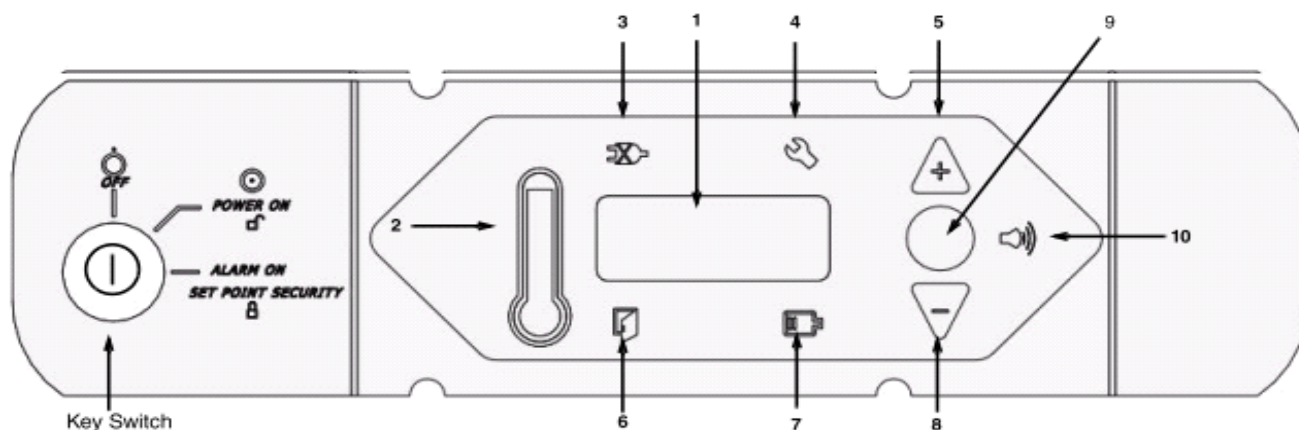


Figure 1. Laboratory Freezer Control Panel

6.1 Control Panel Features

The control panel is located on the top right side of your freezer. You can use the three pushbuttons (#5, #8, and #9 in Figure 1) to change the temperature display (#1) or to adjust temperature and alarm setpoints. The thermometer display (#2) provides a quick visual indicator of current cabinet temperature and alarm conditions.

1. Main temperature display — during normal operation, shows cabinet temperature in degrees Celsius, as measured by the sensor inside the cabinet. You can use the buttons to display other values such as setpoints and highest and lowest recorded temperatures. The number in the main display flashes when the value can be modified.
2. Thermometer — shows cabinet temperature and alarm conditions. There are 10 horizontal bars: 9 are displayed during normal operation, the tenth (top) bar indicates a warm alarm condition. The number of bars illuminated indicates approximate cabinet temperature. Depending on alarm settings, 4 or 5 bars illuminated indicate that the cabinet is at setpoint. For example, suppose that the cabinet temperature setpoint is -20°C and that the warm and cold alarm setpoints are -16°C and -24°C . Then the number of bars illuminated indicates cabinet temperature as shown in Table 1

Table 1. Thermometer display on control panel (setpoint -30°C)

Bars displayed	Temperature	Bars displayed	Temperature
bulb only	-34°C (cold alarm)	6 bars	-28°C
1 bar	-33	7 bars	-27
2 bars	-32	8 bars	-26 (warm alarm)
3 bars	-31	9 bars	-25
4 bars	-30 (setpoint)	10 bars	-24
5 bars	-19		

When cabinet temperature exceeds the warm alarm setpoint, the top bar of the thermometer flashes. When temperature is lower than the cold alarm setpoint, the bulb flashes. When you are in programming mode (described in Table 3) the thermometer shows the setpoint value you are changing.

3. Power failure — illuminated when the main power supply is interrupted. In this case the audible alarm also sounds.
4. Service required — illuminated when the controller is in service programming mode or when simulated warm or cold alarm conditions are failing to occur during an alarm test.
5. Increase — used to increase setpoint values in programming mode and for various display functions.
6. Door ajar — illuminated when the freezer door is open, the alarm is activated, and the key switch is turned to the alarm position).
7. Battery low — illuminated when the backup battery is low.
8. Decrease — pushbutton used to decrease setpoint values in programming mode and for various display functions.
9. Scan — pushbutton used to change the main display and for various other functions.
10. Audible alarm — illuminates during warm and cold alarm conditions.

For full descriptions of display, programming, and service functions, refer to the following tables.

6.2 Display Functions

Function	Meaning	Sequence	Display
Normal operation	Default display while freezer is running	—	Temperature display and control panel thermometer icon show cabinet temperature.
Cold excursion	Show coldest cabinet temperature since last startup or reset	Press ▼	Display shows coldest logged temperature while button is pressed.
Warm excursion	Show warmest cabinet temperature since last startup or reset	Press ▲	Display shows warmest logged temperature while button is pressed.
Mute	Silence audible alarm	Press ● (the Scan button between ▼ and ▽)	Display and thermometer show cabinet temperature, alarm icon continues to flash.
Reset	Return to default display after excursion or alarm condition	Press ▲ and ▼ simultaneously, hold for five seconds	Excursion values are reset; temperature display shows cabinet temperature.
SureTemp Alarm test	Tests alarm system by warming/cooling probe surface; key switch must be in alarm mode	Press ▲ and ● simultaneously	Display and thermometer icon show simulated cabinet temperature, alarms flash and sound as appropriate. Alarms clear when test is completed.

Table 2. Control Panel Display Functions

6.3 Programming Functions

You can enter programming mode by pressing and holding for 5 seconds. Pressing repeatedly scrolls through the available setpoint functions: cold alarm and warm alarm.

Function	Programming Sequence
Adjust temperature control setpoint	Enter programming mode by pressing ● and holding for 5 seconds. On release, the current temperature setpoint value flashes in the temperature display; use ▼ and ▲ to adjust it. Press ● to confirm the new value. The display automatically returns to normal operating mode 30 seconds after the last key entry or after scrolling through all available functions.
Adjust cold alarm setpoint	Enter programming mode by pressing ● and holding for 5 seconds. On release, the current cold alarm setpoint value flashes in the temperature display; use ▼ and ▲ to adjust it. The display automatically returns to normal operating mode 30 seconds after the last key entry or after scrolling through both available functions.
Adjust warm alarm setpoint	Enter programming mode and press ● repeatedly until the top of the thermometer is illuminated. The current warm alarm setpoint value then flashes in the temperature display; use ▼ and ▲ to adjust it. The display automatically returns to normal operating mode 30 seconds after the last key entry or after scrolling through both available functions.

Table 3. Setpoint Programming Functions

6.4 Service Parameters

You can access service parameters by entering programming mode with the controller key in the Power On position, then pressing \odot for an additional 5 seconds. On release of the button, the display will go blank, then display “SEr” with the service wrench icon illuminated. Then the *firmware checksum* (read-only) will be displayed for about 4 seconds. Pressing \odot repeatedly scrolls through the available service functions. While you are in service mode, the wrench icon is illuminated. For any flashing parameter you can use ∇ and Δ to adjust the value.

Parameter	Display	Notes
1. Offset	Value in main display, single bar illuminated in thermometer	Center air temperature calibration. Default value is 0 (maximum + or - 10.0).
2. Line voltage	Err	Not available for this model.
3. Network address	nEt (2 sec.); Adr (2 sec.); then value	Can only be modified by RS-485 communications software.
4. Defrost probe temperature	dEF (2 sec.); Prb (2 sec.); then value	Display only.

Table 4. Service Parameters

7 Operation

7.1 Temperature Settings

The factory default temperature setting is -30°C for automatic defrost freezers, including plasma freezers.

To change the factory temperature settings, refer to the instructions in Section 6.3.

7.2 Start Up

To start up the freezer, complete the following steps:

1. Plug in the power cord.
2. All freezers except undercounter models have a double pole circuit breaker switch located next to the power inlet. Make sure that is in the ON (“1”) position.
3. Insert the silver colored key in the switch on the control panel and turn the power on, turning the key switch to position 1. The compressor and the evaporator fans should start immediately.
4. Rotate the power switch to the ALARM ON position when the temperature drops below the warm alarm setpoint.
5. If desired, lock the cabinet door using the gold colored key. Place duplicate key copies in a safe place.

All controls should now be fully operational, the alarm activated, and all visual indicators active.



CAUTION : Do not operate strong radio emission sources such as walkie-talkies within 3 feet of the freezer. EMI and RFI can affect the performance of the control systems.

8 Defrost

8.1 Automatic Defrost

The defrosting process on all -30°C freezers initiates automatically in response to a built-in timer.

All models are set for one defrost cycle every six hours. The defrost cycle is 20 minutes. The cycle terminates automatically if during defrost the evaporator coil temperature exceeds 15°C.

9 Alarm Systems

9.1 Operating the Alarm

The alarm system is designed to provide visual and audible warning signals for both power failure and rise in temperature. The alarm is equipped with a battery backup.

Default cold and warm alarm values are -40°C and -20°C . These values may be adjusted, following the instructions in Section 6.3.

The alarm system is activated only when the key switch is turned to the Alarm On position. The audible warning signal sounds when there is a power failure or temperature alarm condition, or when the door is ajar for more than 3 minutes.

The Mute function (pressing the ● button) allows you to turn off the audio warning without turning off the visual indicators.

To turn off and reset flashing visual alarms, press ▲ and ▼ simultaneously.

There is also a ringback function after approximately 6 minutes if any alarm condition remains active.

9.2 Local and Remote Alarms

Freezer units can have either a factory-installed local alarm or an optional user-installed remote alarm. Operating and testing procedures are the same for both types of alarm.

The maximum distance between a freezer and a remote alarm depends on the wire gauge used. Refer to Table 5 below.

Wire Gauge	Total Wire Length (feet)	Distance to Alarm 1/2 Wire Length (feet)
20	530	265
18	840	420
16	1,330	665
14	2,120	1,060
12	3,370	1,685

Table 5. Wire Gauges and Distance to Remote Alarm

9.3 Installing a Remote Alarm (Optional)

Remote alarm terminals are located at the rear of the machine compartment. The terminals are: Common, Open on Fail (Normally Closed), and Close on Fail (Normally Open).

1. The remote alarm system has two keyhole slots on the back to hang the alarm system on the wall. Insert two screws, no longer than a #12 truss-head type and spaced two inches apart, into a wall and mount the alarm.
2. Make the following connections:
 - a. Connect the common terminal on the cabinet switch to the purple wire on the alarm.
 - b. Connect the normally closed terminal on the cabinet to the black (“open on fail”) wire on the alarm. This connection gives an alarm when the switch contacts open.
3. Plug the alarm system service cord into an electrical outlet. This alarm is designed for 115V/60 Hz, 115V/50 Hz, or 100V/50 Hz operation.



Note : If you want the alarm signal to sound when the switch contacts **close**, connect the normally open terminal on the cabinet to the red/white (“close on fail”) wire on the alarm. The purple and red/white wires must be tied together in this application.

The wiring diagram is attached to the inside of the alarm back cover.

9.4 Alarm Test

Your freezer is equipped with a SureTemp™ testing system which automatically tests the alarm probe and electronics. This may eliminate the need for other methods of warming or cooling the probe, such as ice baths.

Theory of Operation

During the alarm test, the temperature sensor is artificially heated and by a tiny, built-in thermoelectric heating unit which simulates warm conditions. The electronic control module notes the sensor temperature changes and the control panel displays these changes.

While this alarm testing procedure is very accurate and reliable, the temperature of the refrigerated space does not change during the alarm test.

Alarm Test Procedure



Note : This test automatically advances through all steps and stops.

1. Verify that the key position is in the Alarm On mode, and that the current warm and cold alarm setpoints are within normal ranges (the warm and cold simulations may not work if the setpoints are set to extreme values).
2. To start the alarm test, press Δ and \odot simultaneously. During the test the main display and thermometer bulb will indicate simulated (not actual) cabinet temperature.
3. When simulated temperature exceeds the warm alarm setpoint, the alarm sounds and the alarm icon on the control panel illuminates (#10 in Figure 1).
4. The temperature display begins to drop. After a few seconds, the temperature in the display is back in the operating range.
5. The alarm stops. The temperature on the display drops until the cold alarm sounds.
6. The test is now complete but the alarm continues to sound until the temperature on the display is back in the operating range.

If the simulated alarm conditions do not occur during the first five minutes of the alarm test, the service (wrench) icon illuminates and the test is terminated. You can also terminate the test immediately by turning the key switch to the second (Power On) position. When during the alarm test, the temperature display does not change or the service icon illuminates, check the sensor connections.

After an alarm test has terminated, there is a 10-minute delay before the test can be run again.

10 Chart Recorders

Panel-mounted six inch recorders are standard and factory-installed on all models except for undercounter models, for which free-standing recorders are provided.

Recorder operation begins when the system is powered on.

10.1 Set Up and Operation

To prepare the recorder to function properly, complete the following steps:

1. Open the recorder door to access the recorder.
2. Connect the nine volt DC battery located at the recorder's upper right corner. This battery provides back-up power.
3. Install clean chart paper (refer to Section 10.2 below).
4. Close the recorder door.



Note : The recorder may not respond until the system reaches temperatures within the recorder's range.

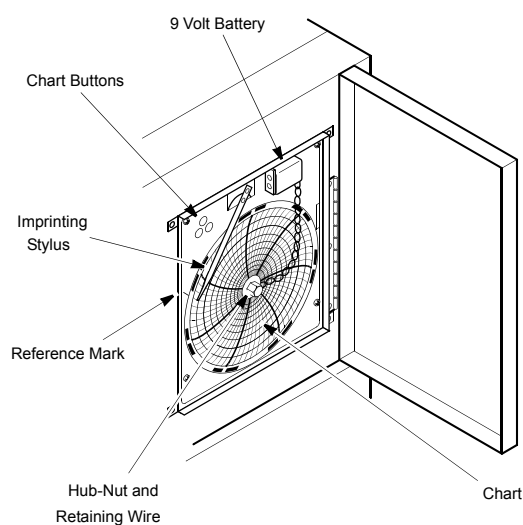


Figure 2. Six Inch Chart Recorder

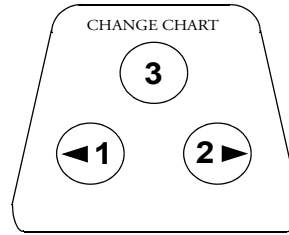


Figure 3. Pressure Sensitive Chart Buttons

10.2 Changing Chart Paper

To change the chart paper, complete the following steps:

1. Locate the pressure sensitive buttons at the front, upper left of the recorder panel.
2. Press and hold the change chart button (#3 on the upper left of the panel) for one second. The pen will move off the scale.
3. Unscrew the center nut, remove the old chart paper, and install new chart paper. Carefully align the day and time with the reference mark on the recorder panel (a small groove on the left side of the panel, shown in Figure 2).
4. Replace the center nut and hand tighten. Press the change chart button again (#3) to resume temperature recording.

10.3 Power Supply

The recorder normally uses AC power when the system is operating. If AC power fails, the LED indicator on the recorder flashes to alert you to a power failure. The recorder continues sensing cabinet temperature and the chart continues turning for approximately 24 hours with back-up power provided by the nine-volt battery.

The LED indicator glows continuously when main power is functioning and the battery is charged.

When the battery is low, the LED flashes to indicate that the battery needs to be changed.

10.4 Calibration Adjustment

This recorder has been accurately calibrated at the factory and retains calibration even during power interruptions. If required, however, adjustments can be made as follows:

1. Run the unit continuously at the control setpoint temperature. Continue steady operation for at least two hours to provide adequate time for recorder response.
2. Measure cabinet center solution temperature with a calibrated temperature monitor.
3. Compare the recorder temperature to the solution temperature. If necessary, adjust the recorder by pressing the left (#1) or right (#2) chart buttons for five seconds.



Note : The stylus does not begin to move until the button is held for five seconds.

11 Drawers

11.1 Removing the Drawers

Larger models

To remove the drawers in 23 ft³ and larger models, complete the following steps (refer to Figure 4):

1. Pull the drawer toward you until the slides are fully extended.
2. Lift the back of the drawer to disengage the mounting tabs from the slots on the slides.



Note : The drawers fit snugly between the slides. Push the back of the drawer from underneath to remove the drawer.

3. Raise the back of the drawer almost to a vertical position and disengage the front mounting clips from the slides.

5 and 12 Cu. Ft. models

To remove the drawers on 5 and 12 cu.ft. models, depress the hooks located on both sides of the drawers (about 1/3 of the way back) and slide the drawers up and out.

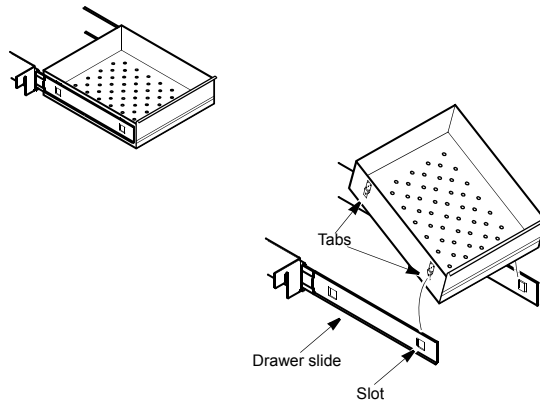


Figure 4. Drawer Removal (Upright Models)

11.2 Reinstalling the Drawers

To reinstall the drawers in 23 cu.ft. models and larger, complete the following steps (refer to Figure 4):

1. Pull both drawer slides toward you until the slides are fully extended.
2. Position the drawer between the slides and with the back facing end raised at a 45 degree angle, insert the front mounting clips into the slots on front of the slides.
3. Push the back of the drawer down between the slides and insert the drawer tabs into the back slots.



Note : The drawers fit snugly between the slides. Push on the back of the drawer from the inside to insert the drawer tabs completely into the slots. Make sure both drawer tabs are aligned with the slots on the slides before pushing the drawer down between the slides.

11.3 Changing Drawer Position

The drawer slides are adjustable to higher and lower positions in the cabinet. You can position these slides in the vertical slots which are spaced at one-inch intervals.

Drawer slides have a small wire safety clip at the front pilaster which prevents the slides from falling when the drawers are removed. To change the position of the drawer slides, complete the following steps:

1. Locate the safety clip.
2. Slip a small screwdriver under the bottom of the wire clip and pry the clip out toward the inside of the freezer.
3. Lift up the slide at the front. The slide is free to move from the front pilaster.
4. The drawer slide must be removed from the rear pilaster at approximately a 45 degree angle toward the center of the cabinet.
5. Pull the slide toward the front of the cabinet.
6. Reposition the slide to the desired vertical slot.
7. Replace the safety clip.



CAUTION : Drawer slides do *not* require lubrication. Additional lubricant could impede movement of the drawers when the lubricant is cold.

12 Cleaning

12.1 Cleaning the Drawers

To clean the drawers, use a solution of water and a mild detergent. Rinse the drawers and wipe them dry with a soft cloth.

For instructions on removing and reinstalling drawers, refer to Section 10.

12.2 Cleaning the Condenser



CAUTION : Condensers should be cleaned at least every six months. In heavy traffic areas, condensers load with dirt more quickly. Failure to keep the condenser clean can result in equipment warm-up or erratic temperatures.



WARNING : Be sure to disconnect the unit from main power before cleaning the condenser.



CAUTION : Never clean near condensers with your fingers. Some surfaces are sharp.

In all models, the condenser is located in the top machine compartment. To clean the condenser:

1. Disconnect the power.
2. Remove the top front grill.
3. Use a vacuum cleaner with hose and brush attachments to clean the front face of the finned surface.
4. Clean up any loose dust and replace the front grill.
5. Reconnect the power.

13 Troubleshooting



WARNING : Troubleshooting procedures involve working with high voltages which can cause injury or death. Troubleshooting should only be performed by trained personnel.

This section is a guide to troubleshooting equipment problems.

Problem	Cause	Solution
Unit does not operate or Power Failure Indicator is on.	Power supply	<p>Check that the cord is securely plugged in.</p> <p>All models except undercounter models have a double pole circuit breaker located next to the power inlet. Make sure that it is in the ON (“1”) position. Try cycling the switch to OFF (“0”) then ON (“1”).</p> <p>Plug another appliance into the outlet to see if it is live.</p> <p>If the outlet is dead, check the circuit breaker or fuses.</p>
Temperature fluctuates.	Cold control	Make sure that the cold control is set correctly. Refer to Section 6.
	Condenser	Make sure the condenser is clean. Refer to Section 12.2 on page 22.
Unit warms up.	Door is open.	Make sure the door is completely closed.
	Warm product recently loaded in unit.	Allow ample time to recover from loading warm product.
	Power supply	Check for proper voltage to the unit. If there is no voltage to the unit, call an electrician.
	Compressor	<p>If the compressor is not running, check if the unit has a power failure alarm. If the power failure alarm light is on, have an electrician check for proper voltage to the unit.</p> <p>If the compressor is running, open the door and look through the slotted air intake in the bottom of the evaporator cover to see if icing is present on the evaporator. If icing is present and there is no air flow behind evaporator, call technical service for assistance. The evaporator fans may be inoperative.</p> <p>If the compressor is running and there is air flow behind the evaporator, contact an authorized service provider or call the technical support hot line for assistance.</p>

Table -6. Troubleshooting Procedures

14 Warranty Statement

Domestic • 24 Months Full Warranty Parts and Labor
International Warranty • 24 Months Full Warranty for Parts Only

During the first twenty four (24) months from shipment, Thermo Fisher Scientific Inc, through its authorized Dealer or service organizations, will at its option and expense repair or replace any part found to be non-conforming in material or workmanship. Thermo Fisher Scientific Inc reserves the right to use replacement parts, which are used or reconditioned. Replacement or repaired parts will be warranted for only the unexpired portion of the original warranty.

This warranty does not apply to damage caused by (i) accident, misuse, fire, flood or acts of God; (ii) failure to properly install, operate or maintain the products in accordance with the printed instructions provided, (iii) causes external to the products such as, but not limited to, power failure or electrical power surges, (iv) improper storage and handling of the products, (v) use of the products in combination with equipment or software not supplied by Thermo Fisher; or (vi) installation, maintenance, repair, service, relocation or alteration of the products by any person other than Thermo Fisher or its authorized representative. To obtain proper warranty service, you must contact the nearest authorized service center or Dealer. Thermo Fisher Scientific, Inc's own shipping records showing date of shipment shall be conclusive in establishing the warranty period. At Thermo Fisher's option, all non-conforming parts must be returned to Thermo Fisher postage paid and replacement parts are shipped FOB Thermo Fisher's location.

Limitation of Liability

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 FISHER DOES NOT WARRANT THAT THE PRODUCTS ARE ERROR-FREE OR WILL ACCOMPLISH ANY PARTICULAR RESULT.

THERMO FISHER SHALL NOT BE LIABLE FOR ANY INDIRECT OR CONSEQUENTIAL DAMAGES INCLUDING, WITHOUT LIMITATION, DAMAGES TO LOST PROFITS OR LOSS OF PRODUCTS.

Thermo Fisher Scientific Inc.

275 Aiken Road
Asheville, NC 28804
United States

www.thermofisher.com

Thermo
SCIENTIFIC

325099H12 Rev. B