



# Ultra Low Temperature Freezers

**Revco RLE series, Forma 89000  
series, HERAFreeze HLE series and  
Thermo Scientific TLE series**

## **Installation and Operation**

328398H01

Rev. C

August 2017

Visit us online to register your warranty  
[www.thermofisher.com/labwarranty](http://www.thermofisher.com/labwarranty)



**thermo scientific**

**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.

© 2017 Thermo Fisher Scientific Inc. All rights reserved.

# Table of Contents

<b>Models</b> .....	<b>1</b>
<b>Safety Precautions</b> .....	<b>2</b>
<b>Unpacking</b> .....	<b>4</b>
<b>Packing List</b> .....	<b>5</b>
<b>General Recommendations</b> .....	<b>6</b>
Temperature Monitoring .....	6
General Usage .....	6
Initial Loading.....	6
Battery Door Opening / Closing .....	7
<b>Operating Standards</b> .....	<b>8</b>
Electrical Specifications.....	8
<b>Installation</b> .....	<b>9</b>
Location .....	9
Wiring .....	9
Leveling .....	10
Backup System (Optional) .....	10
Super Insulated Cabinet Construction.....	10
Door Operation .....	10
Pressure Equalization Port .....	12
Installing the Remote Alarm Connector .....	12
Intended Use .....	13
<b>Start Up</b> .....	<b>14</b>
Initial Start Up.....	14
Standby Mode .....	21
<b>Operation</b> .....	<b>22</b>
Operation Overview .....	22
Home Screen .....	22
Settings.....	23
Event Log.....	30
Alarms.....	31
Users.....	32
Reports .....	35
Chart.....	36
<b>Health Status and Alarm Management</b> .....	<b>40</b>
Health Status Overview .....	40
Notifications / Cautions .....	42
Alarms / Warning.....	43
<b>Backup System (Optional)</b> .....	<b>45</b>
CO <sub>2</sub> and LN <sub>2</sub> Precautions .....	45

## Table of Contents

Installation.....	46
Start Up .....	47
Operation.....	47
<b>Chart Recorders (Optional).....</b>	<b>48</b>
Set Up and Operation .....	48
Changing Chart Paper .....	49
Calibration Adjustment .....	49
<b>Maintenance and Troubleshooting.....</b>	<b>50</b>
Cleaning the Condenser .....	50
Cleaning the Condenser Filter.....	50
Gasket Maintenance .....	50
Defrosting the Freezer.....	51
Battery Maintenance .....	51
<b>Warranty .....</b>	<b>52</b>
<b>Warranty (International) .....</b>	<b>53</b>
<b>Appendix A: Alarm Summary.....</b>	<b>54</b>
<b>Appendix B: Event Log Detail.....</b>	<b>60</b>

# 1 Models

<b>Brand - Model</b>	<b>Size (xxx)</b>	<b>Voltage (*)</b>
Forma - 89xxx86*	300 / 400 / 500 / 600	A / D / V
Thermo Scientific –TLExxx86*	300 / 400 / 500 / 600	A / D / V
HERAfreeze –HLExxx86*	300 / 400 / 500 / 600	A / D / V
Revco -RLExxx86*	300 / 400 / 500 / 600	A / D / V

**Note** *Energy Star is not applicable to models of size 300 and 400.*

## 2 Safety Precautions

In this manual, the following symbols and conventions are used:



This symbol used alone indicates important operating instructions which reduce the risk of injury or poor performance of the unit.



**CAUTION:** This symbol, in the context of a CAUTION, indicates a potentially hazardous situation which if not avoided could result in minor to moderate injury or damage to the equipment.



**WARNING:** This symbol indicates potentially hazardous situations which, if not avoided, could result in serious injury or death.



**WARNING:** This symbol indicates situations where dangerous voltages exist and potential for electrical shock is present.



The snowflake symbol indicates extreme low temperatures and high risk of frostbite. Do not touch bare metal or samples with unprotected body parts.



This symbol indicates a need to use gloves during the indicated procedures. If performing decontamination procedures, use chemically resistant gloves. Use insulated gloves for handling samples and when using liquid nitrogen.



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 this 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 OEM exact replacement equipment or parts. Before use, confirm that the product has not been altered in any way.



**WARNING:** Your unit must be properly grounded in conformity with national and local electrical codes. Never connect the unit to overloaded power sources.



**WARNING:** Disconnect the unit from all power sources before cleaning, troubleshooting, or performing other maintenance on the product or its controls.



**WARNING:** “Caution, risk of fire”. This unit is charged with hydrocarbon refrigerants.

### **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 to the manufacturer without written authorization. When submitting a claim for shipping damage, request that the carrier inspect the shipping container and equipment.

## 4 Packing List

Inside the freezer cabinet is a bag containing:

- This manual
- A handle lock key
- A CD with user's manuals, including translated versions
- Certificates of conformance and calibration
- A remote alarm contact connector
- Posts for rear spacing

If you have ordered a field-installed chart recorder, the bag will also contain:

- Recorder installation instructions
- Extra inkless paper

If you have ordered a backup system, the cabinet will also contain:

- A hose assembly
- English and metric connectors

If specified on the order, the bag may also include:

- A QC temperature graph and test log
- Calibration information

If you have ordered the Proximity Access Card Option, the cards will be in a bag attached to the front of the freezer.

## 5 General Recommendations

### 5.1 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.

### 5.2 General Usage

This refrigeration system is designed to maintain ultra-low temperatures with safety in an ambient environment within 15°C to 32°C (59°F to 90°F), only when the freezer is used for storage.



**WARNING:** This unit is not a “rapid-freeze” device. Freezing large quantities of liquid, or high-water content items, will temporarily increase the chamber temperature and will cause the compressors to operate for a prolonged time period.

Avoid opening the door for extended time periods since chamber temperature air will escape rapidly. Also, keep the inner doors closed as much as possible. When room air, which is higher in humidity, replaces chamber air, frost may develop in the chamber more rapidly.

### 5.3 Initial Loading

**Allow the freezer to operate at the desired temperature for a minimum of 12 hours before loading.**

Load the freezer one shelf at a time, beginning with the top shelf. After loading each shelf, allow the freezer to recover to the desired set point before loading the next shelf. Repeat this process until the freezer is fully loaded.



**CAUTION:** Failure to follow these procedures or overloading the unit may cause undue stress on the compressors or jeopardize user product safety.

## 5.4 Battery Door Opening / Closing

To open the grille door, pull the door from the top right corner as shown in the figure below.

To close the grille door, push the door against frame to hold latch in position.



**Figure 1.** Door Opening

## 6 Operating Standards

The freezers described in this manual are classified for use as stationary equipment in a Pollution Degree 2 and Overvoltage Category II environment.

These units are designed to operate under the following environmental conditions:

- Indoor use
- Altitude up to 2000m
- Maximum relative humidity 60% for temperatures within 15°C to 32°C (59°F to 90°F).
- Main supply voltage fluctuations not to exceed  $\pm 10\%$  of the nominal voltage.

### 6.1 Electrical Specifications

The last character in the model number listed on the dataplate identifies the electrical specifications for your unit. Specific unit current rating is listed on the dataplate.

The voltage types are A, D and V as specified in the following table:

**Table 1. Revco RLE series, Forma 89000 series, HERAfreeze HLE series and Thermo Scientific TLE series Electrical Specifications**

Model	Voltage	Frequency	Current
300A	115 V	60 Hz	12.6 A
300D	208-230 V	60 Hz	5.6 A
300V	230 V	50 Hz	5.9 A
400A	115 V	60 Hz	13.0 A
400D	208-230 V	60 Hz	5.9 A
400V	230 V	50 Hz	5.3 A
500A	115 V	60 Hz	12.5 A
500D	208-230 V	60 Hz	5.0 A
500V	230 V	50 Hz	5.9 A
600A	115 V	60 Hz	12.7 A
600D	208-230 V	60 Hz	6.2 A
600V	230 V	50 Hz	6.1 A

## 7 Installation



**WARNING:** Do not exceed the electrical rating printed on the data plate located on the lower left side of the unit.

### 7.1 Location

Install the unit in a level area free from vibration with a minimum of 8 inch (20 cm) of space on the top and sides, 6 inch (15 cm) in back. Refer to Section 7.3 for further instructions on leveling cabinets. Allow enough clearance so that door can swing open at least 85°.

The rear spacing posts provided with the freezer can be used to ensure proper clearance. To install the spacing posts, screw them into the back in the rear deck area.

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 15°C to 32°C (59°F to 90°F).

### 7.2 Wiring



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



**CAUTION:** 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 the unit to overloaded power lines.



**CAUTION:** Do not position the unit in a way that impedes access to the disconnecting device or circuit breaker in the back of the unit.



**CAUTION:** Always connect the freezer to a dedicated (separate) circuit. Each freezer is equipped with a service cord and plug designed to connect it to a power outlet which delivers the correct voltage. Supply voltage must be within  $\pm 10\%$  of the freezer rated voltage.



**CAUTION:** Never remove or disable the grounding prong from the service cord plug. If the prong is removed, the warranty is invalidated.

### 7.3 Leveling

Make sure that the floor is level. The unit must be level both front to back and side to side.

300 and 400 box capacity models are equipped with one or two leveling legs on the right hand side. These may be used to help prevent the unit from shifting during a door opening.

Be certain to lock the brakes for units equipped with casters.

### 7.4 Backup System (Optional)

If you are using a CO<sub>2</sub> or LN<sub>2</sub> backup system, refer to Section 11 for installation and operation instructions.

### 7.5 Super Insulated Cabinet Construction

In all models, the cabinet walls have a vacuum insulation core encapsulated by a sealed film laminate.



**CAUTION:** Never drill holes in or near the cabinet walls. Drilling could damage the insulation and make the unit inoperable.

### 7.6 Door Operation

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

Features include:

- One-hand operation
- A front-accessible lock
- Hasps for a standard padlock to provide additional security. Length of the shackle must be between 3/4 inch (1.9 cm) and 1 1/2 inch (3.8 cm).
- Durable construction for reliable operation and safe product storage.
- Door ramp alignment feature
- Optional controlled access to the freezer with Proximity Access cards.



**CAUTION:** When moving the freezer, always grasp cabinet surfaces; never pull the freezer by the latch handle.

#### 7.6.1 Opening the Door

For freezers with the Proximity Access Card option:

1. Remove the padlock if installed.

2. To unlock the door, pass the card in front of the freezer below the LCD display.
3. Grasp the latch handle and pull it toward yourself until the latch disengages from the cabinet strike.
4. Keep pulling by the latch handle to open the main door.

For freezers without the Access Card option:

1. Remove the padlock if installed.
2. Grasp the latch handle and pull it toward yourself until the latch disengages from the cabinet strike.
3. Keep pulling by the latch handle to open the main door.

### 7.6.2 Opening the Door During a Power Outage

In case of power outage and a unit that has the Proximity Access Card option, you may use a 9 volt battery to activate the system. To access the 9 volt terminal, remove the USB cover and locate the battery terminals.

Once the terminals are exposed, open the door by holding the 9 volt battery against the terminals and pass a valid proximity card below the display area. Once the door is open, remove the 9 volt battery.

**Note** *The terminals are polarized therefore orient the 9 volt battery properly.*

### 7.6.3 Closing the Door

**Note** *The latch does not self-engage automatically when you close the door. You must rotate the latch into the open position first.*

1. Grasp the latch handle (preferably with your left hand) and pull it toward yourself, rotating the latch into the open position.
2. Move the freezer door into the closed position and gently push the handle away from you, making sure that the latch engages fully with the cabinet strike.
3. Keep applying gentle pressure to the latch handle until the latch is securely in closed position.
4. Insert the key and rotate counterclockwise to lock.
5. Replace the padlock as required.

## 7.7 Pressure Equalization Port

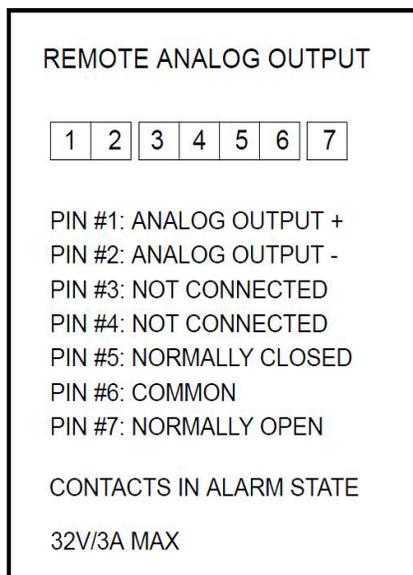
When an upright ultra-low temperature freezer door is opened, room temperature air rushes into the storage compartment. When the door is closed, the fixed volume of air is cooled rapidly. Pressure drops below atmospheric pressure, resulting in a substantial vacuum. Re-entry into the cabinet is impossible until internal pressures are returned to atmospheric pressure. Without a pressure equalization mechanism, it can take, in extreme cases, several hours before the door can easily be reopened.

All upright models feature a port that provides vacuum relief after door openings. The pressure equalization port is located in the door behind the eye-level panel on the front of the freezer. Although the port is designed to self-defrost, excessive frost accumulation on the inner door could eventually restrict air flow. Therefore you should periodically inspect the inner door and brush away any loose frost using a stiff nylon brush.

## 7.8 Installing the Remote Alarm Connector

The remote alarm contacts are located on the back of the freezer above and to the left of the power switch. After installing the wiring from the remote alarm to the connector, install the connector to the freezer micro-board.

The pin configuration is shown in Figure 2 below.



**Figure 2.** Remote Alarm Pin Configuration

The contacts will trip in the event of a power outage, high temperature alarm, low temperature alarm or door ajar alarm.

## 7.9 Intended Use

The -86°C freezer (refer to Section 1 for the specific model series) described in this manual are high performance units for professional use. These products are intended for use as cold storage in research use and as a general purpose laboratory freezer, storing samples or inventory at operating temperatures between -50°C and -80°C.

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

This unit is not intended for use in classified hazardous locations, nor to be used for the storage of flammable inventory.

## 8 Start Up

### 8.1 Initial Start Up

To start up the freezer, complete the following steps:

1. Plug the freezer into the power outlet.
2. Turn the power switch in back of the freezer, on the bottom right, to the ON position.
3. Once the freezer is powered up, the Thermo Scientific logo is displayed on the front screen. If this is the first time the unit is being turned on, an initial setup must be completed. Tap the Start Setup button to initiate the setup.



**Figure 3.** Main Screen

The first step of the setup is to select the language. This screen allows you to specify the preferred display language. Once the language is selected, tap the Next button.



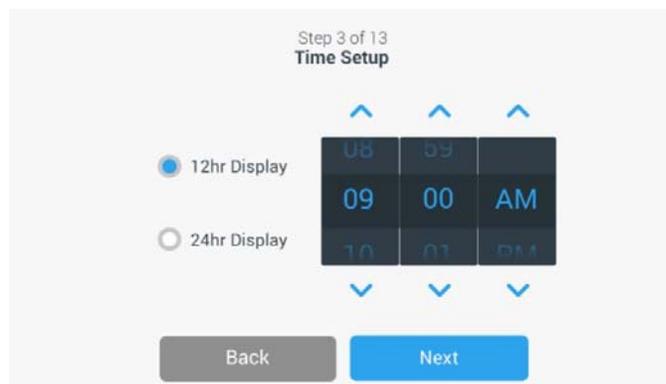
**Figure 4.** Language Selection Screen

The next screen allows you to select the date and date format. After setting the date, tap the Next button.



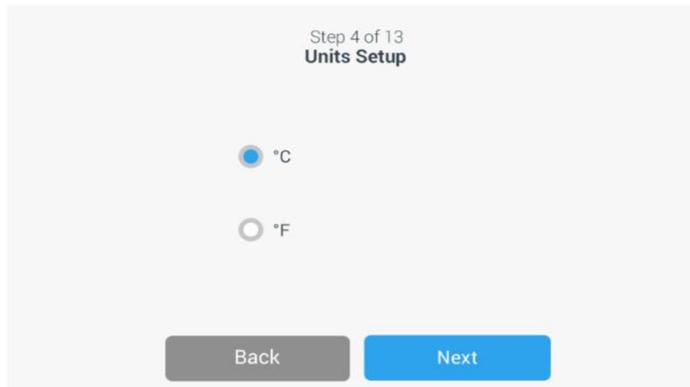
**Figure 5.** Date Setup Screen

The next screen allows you to specify the time and time format. After setting the time, tap the Next button.



**Figure 6.** Time Setup Screen

The next screen allows you to specify the temperature unit. After selecting the unit, tap the Next button.



**Figure 7.** Units Setup Screen

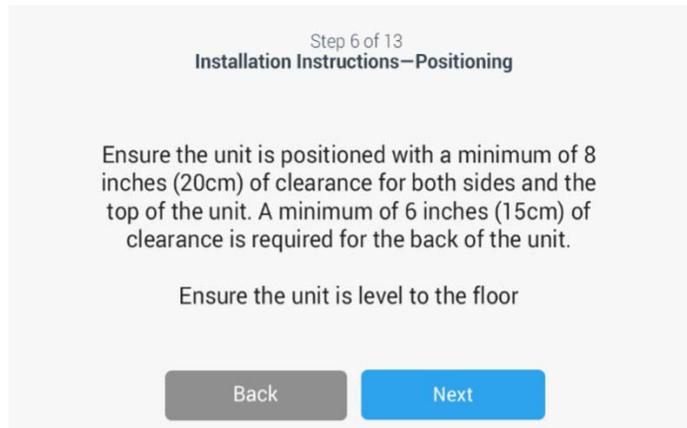
The next 3 screens provide installation instructions.

The first screen provides information regarding power source and ambient temperature conditions. Tap the Next button to continue through the installation instructions.



**Figure 8.** Installation Instructions (Power and Temperature) Screen

The second screen provides information regarding unit positioning, spacing and leveling. Tap the Next button after reviewing.



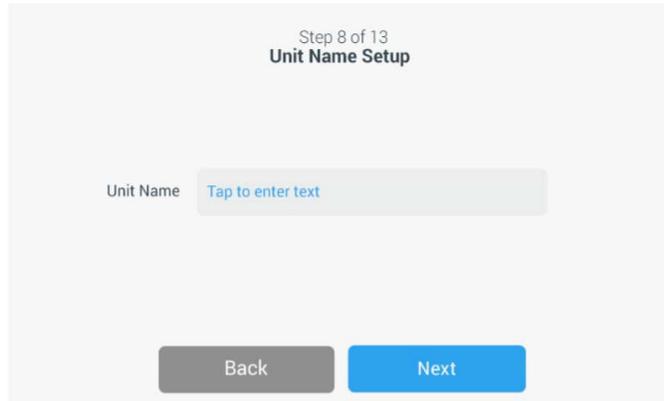
**Figure 9.** Installation Instructions (Positioning) Screen

The third screen provides information regarding initial freezer loading. Tap the Next button after reviewing.



**Figure 10.** Installation Instructions (Before Usage) Screen

The next screen allows you to identify an individual freezer by specifying a Unit Name. After entering a name, tap the Next button.



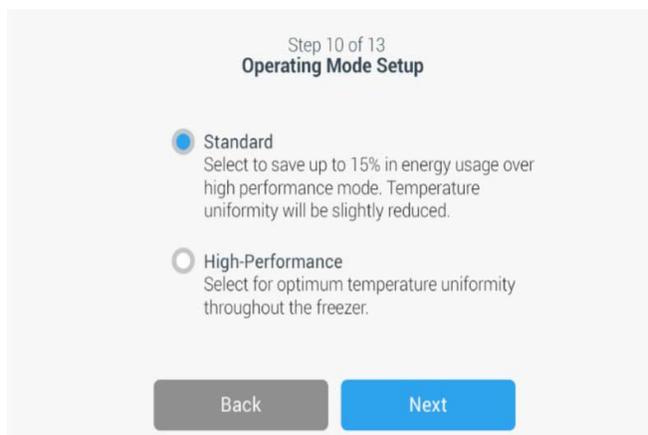
**Figure 11.** Unit Name Setup Screen

The next screen allows you to specify the temperature, warm and cold alarm setpoints. After setting the temperatures, tap the Next button.



**Figure 12.** Setpoints and Alarms Screen

This screen allows you to select the Operating Mode. Once the operating mode has been selected, tap the Next button.



**Figure 13.** Operating Mode Setup Screen

This screen allows you to specify the Access Mode.



**Figure 14.** Access Mode Setup Screen

If Secured Access is selected, at least one administrative account will need to be created. This can be done in step 12.

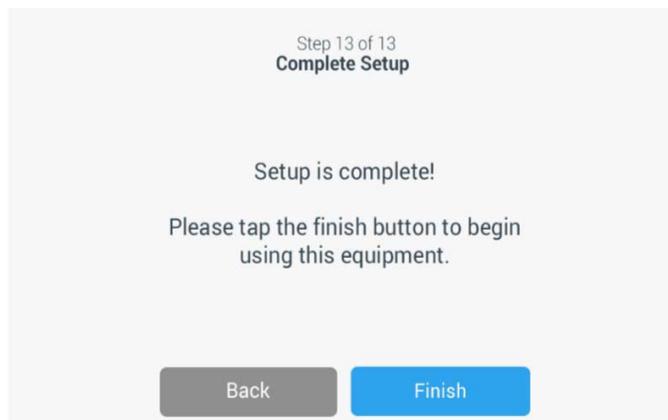
This screen allows you to enter the details of the administrator. All the fields marked with a red asterisk have to be filled. Once the details are entered, tap the Next button to proceed to step 13.

The figure consists of three sequential screenshots of the 'Step 12 of 13 Access Mode Setup' screen. Each screenshot shows a different set of form fields that must be filled out. The fields are marked with a red asterisk (\*). The 'Next' button is highlighted in blue, and the 'Back' button is greyed out.

- Screenshot 1:** Shows the first four fields: First Name\*, Last Name\*, Email\*, and User Name\*. The 'Next' button is visible at the bottom right.
- Screenshot 2:** Shows the next four fields: Email\*, User Name\*, Password\*, and Confirm Password\*. The 'Next' button is visible at the bottom right.
- Screenshot 3:** Shows the final four fields: Confirm Password\*, Phone: Office\* (with a checkmark icon), User Role\* (with a dropdown arrow), and Access ID\* (with an 'X' icon). The 'Next' button is visible at the bottom right.

**Figure 15.** Secured Access Mode Setup Screen

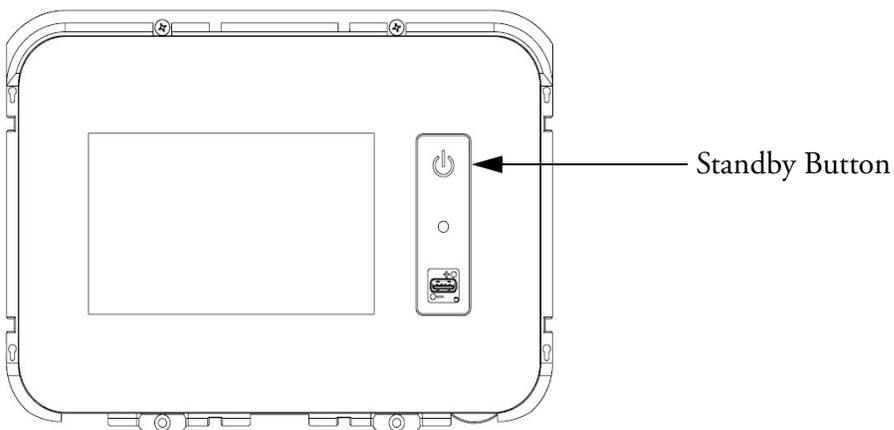
This screen indicates your initial setup is complete. Press the Finish button to complete initial setup or the Back button to make changes.



**Figure 16.** Complete Setup Screen

## 8.2 Standby Mode

There is a standby button on the front of the unit. When pressed and held for ~3 seconds the user will be prompted to confirm that the unit should be placed in standby mode. Upon confirmation, the unit will enter standby mode. **The refrigeration system will be shut down and the unit will not cool while in standby mode.** To exit standby mode and restore normal operation, the standby power button must be pressed and held for ~1 second.



**Figure 17.** Standby Button

## 9 Operation

### 9.1 Operation Overview

Once you have successfully completed the initial start up procedures, the freezer starts operating normally and the only actions required are:

- Setting the operating and alarm set points, refer to Section 9.3.1.
- Activating the CO<sub>2</sub> or LN<sub>2</sub> backup system if installed. For instructions on backup settings and activating the system, refer to Section 11.

### 9.2 Home Screen

The Home Screen below is the default screen.



**Figure 18.** Home Screen

The various options available on the home screen are:

- The vertical panel on the left is the navigation bar that provides access to all functions of the unit.
- The colored icon in the middle indicates the health of the unit. There are four icons to denote this:
  - Green heart with a check mark indicates operation is normal.
  - Yellow triangle with an exclamation mark indicates a notification.
  - Red bell indicates an alarm condition.
  - Red bell with a diagonal indicates the alarm has been snoozed.
- A login button to login into the system. Refer to Section 9.2.1 for more information.

- The setpoint temperature is displayed. This can be changed by tapping the Setpoint button at the bottom of the screen.
- The temperature chart can be viewed by tapping the Chart button at the bottom of the screen. Refer to Section 9.8 for more information.
- The “i” icon at the bottom right corner of the screen is the onboard help button. Tap this icon to display an onboard help box with text explaining all of the features available on that particular screen.
- The back button to navigate to the previous screen.

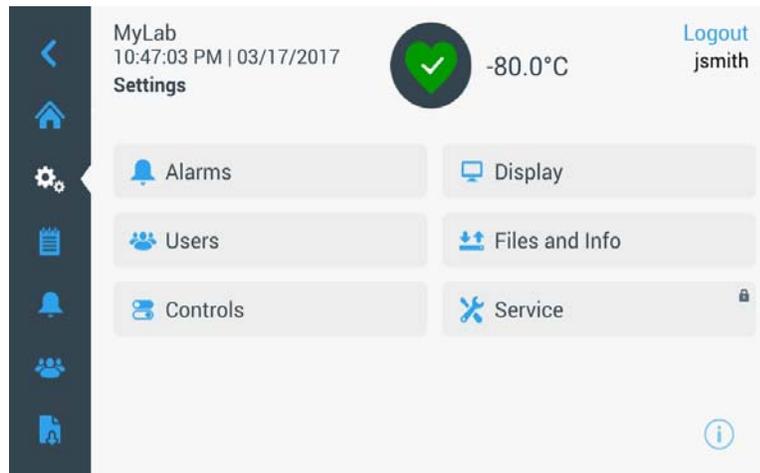
**9.2.1 User Login**

When the system is in Secured Access Mode, the user has to login by entering their username and password.

When the system is running in Full Access mode, the login feature is restricted to service technicians to access the Service screens.

**9.3 Settings**

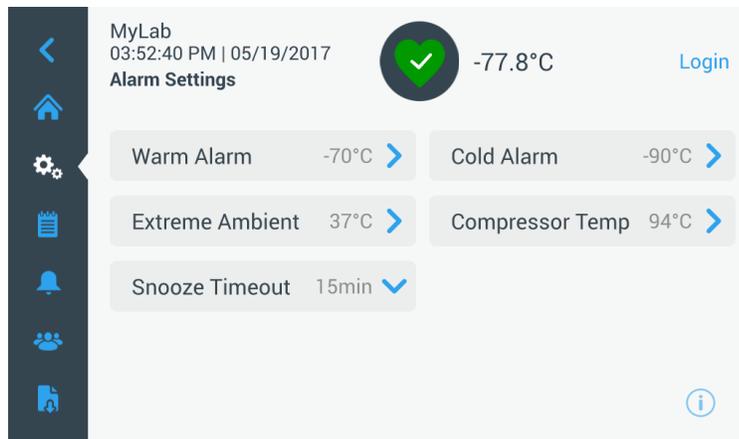
The second tab on the navigation panel is the Settings icon. The following screen will be displayed once the Settings icon is tapped:



**Figure 19.** Settings Screen

### 9.3.1 Alarms

The Alarm Settings screen provides the option to set the warm and cold alarm setpoints. Tap the Warm Alarm or Cold Alarm button and the Setpoints screen will be displayed permitting parameter adjustments. For more information, refer to Section 9.3.5.



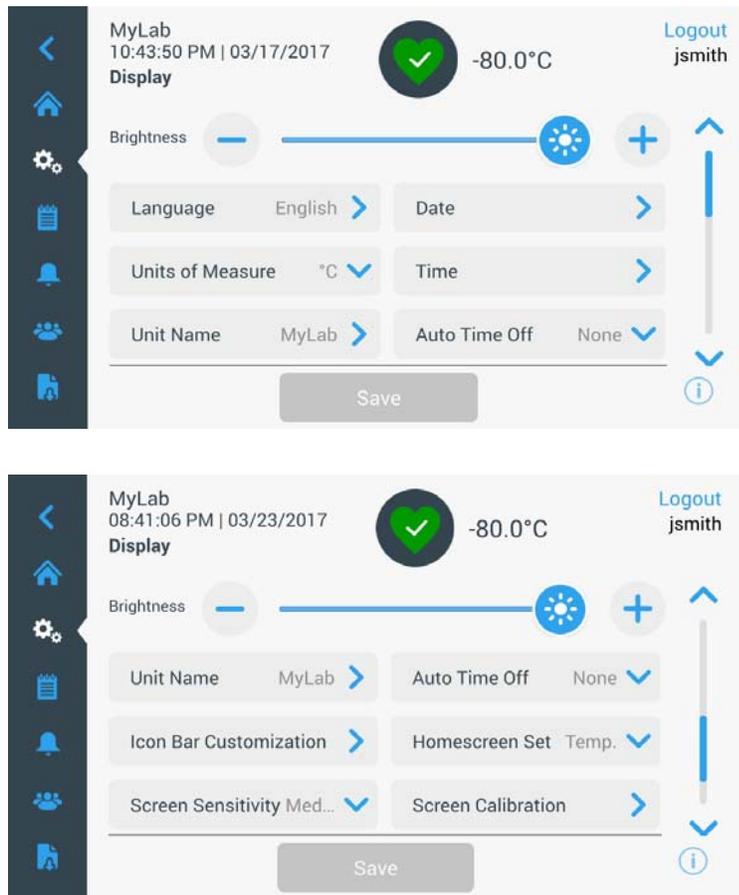
**Figure 20.** Alarm Settings Screen

- **Warm Alarm:** The range of the warm alarm temperature is  $-40^{\circ}\text{C}$  to within  $5^{\circ}\text{C}$  of setpoint.

**Note** *The warm alarm will be disabled for 12 hours from a warm start condition.*

- **Cold Alarm:** The range is  $-99^{\circ}\text{C}$  to within  $5^{\circ}\text{C}$  of setpoint.
- **Extreme Ambient:** This allows the user to set the extreme ambient alarm setpoint. The range is  $32^{\circ}\text{C}$  to  $40^{\circ}\text{C}$ . Default is  $37^{\circ}\text{C}$ .
- **Compressor Temp:** It is an alarm setpoint for the second stage compressor sump temperature. The range is  $70^{\circ}\text{C}$  to  $98^{\circ}\text{C}$ . Default is  $94^{\circ}\text{C}$ .
- **Snooze Timeout:** This sets the time to snooze the audible alarm for an active alarm.

**9.3.2 Display** From the Settings screen, pressing the Display button will show the Display screen. Various display settings can be adjusted.



**Figure 21.** Display Screen

- **Brightness:** Use the slide control or the +/- buttons to adjust the brightness level of the display.
- **Language:** To change the display language, tap the Language button and select the desired language.
- **Date:** To set the date and the date format, tap the Date button.
- **Units of Measure:** To change the displayed unit of measure, tap the Units of Measure button.
- **Time:** To set the time and time format, tap the Time button.
- **Unit Name:** To enter or change the unit name, tap the Unit Name button.

- **Auto Time Off:** Select the time range for when the LCD will automatically darken. The default option None, will leave the LCD constantly illuminated.
- **Icon Bar Customization:** This is used to customize the bottom three icons in the navigation bar.
- **Homescreen Set:** To select the default homescreen, tap the Homescreen Set button.
- **Screen Sensitivity:** Select to modify the touch sensitivity of the screen.
- **Screen Calibration:** Tap to run a screen auto calibration routine.

Select the Save button after making the necessary changes.

### 9.3.3 Users

Access Mode is used to change the access mode for the system (full or secured), add a user to the system, and to import and/or export a user database.

The Users screen can also be accessed by using the User Icon on the left navigation bar. For more information, refer to Section 9.6.



**Figure 22.** Users Screen

**9.3.4 Files and Info** The following screen is displayed when the Files and Info button is selected:

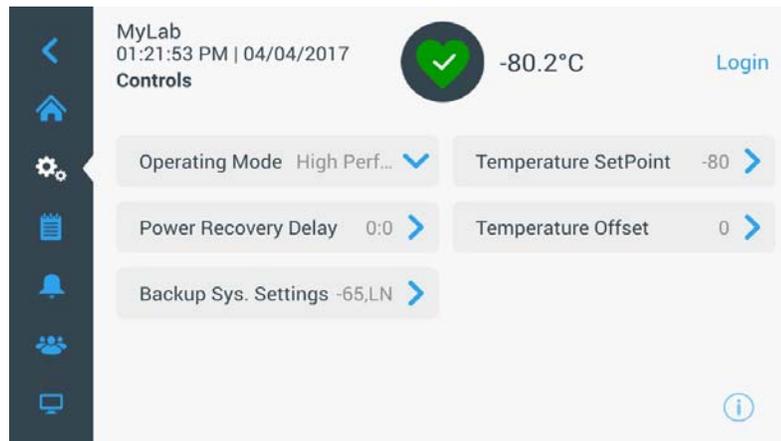


**Figure 23.** Files and Info Screen

- **Factory Reset:** Select to reset the settings, including temperature setpoint to factory defaults.
- **Configuration Files:** This is used to export or import configuration files. Configuration files may be uploaded to other freezers.
- **Reports:** This is used to export temperature and event data. A date range can be provided too. Refer to Section 9.7 for more information.
- **Contact Us:** Tap this button to view or modify the service contact information.

**9.3.5 Controls** Temperature setpoints can be set in the Controls screen. To select a temperature setpoint,

Select the Controls button to navigate to the Controls screen.



**Figure 24.** Controls Screen

- **Operating Mode:** The default setting is the Standard mode, which provides a balance between power consumption and peak variation performance. High Performance mode provides minimum temperature peak variation.

**Note** *Customers performing on-site temperature calibration may observe as much as a 2°C variation when an external probe is placed next to the freezer control probe. This variation is due to optimization of the control system to ensure temperature uniformity throughout the chamber.*

- **Power Recovery Delay:** Tap this button to set the time delay upon startup after power failure. Default is 0.
- **Temperature Offset:** This is used for calibration. Range is -10°C to +7°C. Default is 0.
- **Backup System Settings:** This screen allows the user to select the backup type and backup setpoint for units with a backup system installed. For more information, refer to Section 11.

- Tap the Temperature Setpoint button to display the Setpoints screen:



**Figure 25.** Setpoints Screen

- The setpoint and temperature alarm parameters may be adjusted by swiping the spin control up / down or pressing the up / down arrows.

**Note** *A setpoint change may automatically change the warm and / or cold alarm setpoints as well to prevent unnecessary alarms.*

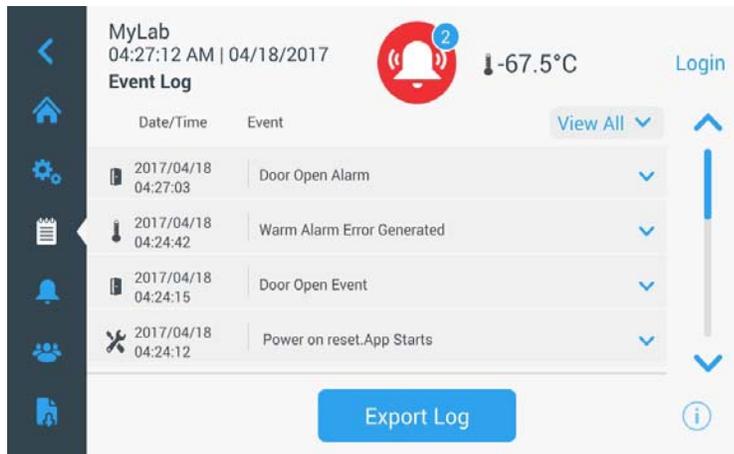
- After selecting the temperatures, tap the Save button to confirm changes.

**Note** *If the save button is not pressed, the unit will not respond to the setpoint change request.*

**Warm Alarm Test:** Tapping this button puts the system into a warm alarm test which simulates a warm alarm experience. Once this is selected, the homescreen is displayed with the current temperature readout. The temperature display will increase to the warm alarm temperature setpoint. Once the warm alarm temperature has been reached, the user is prompted to end the test.

## 9.4 Event Log

The third tab on the navigation panel is the event log that contains a record of user and system events. The Event Log screen will be displayed once the Event Log icon is pressed:



**Figure 26.** Event Log Screen

This screen displays up to two weeks of recent events, with date time stamps for each event.

The Date/Time and Event columns can be sorted in an ascending or descending order by selecting the column header.

Additional information of an individual event can be viewed by selecting the event.

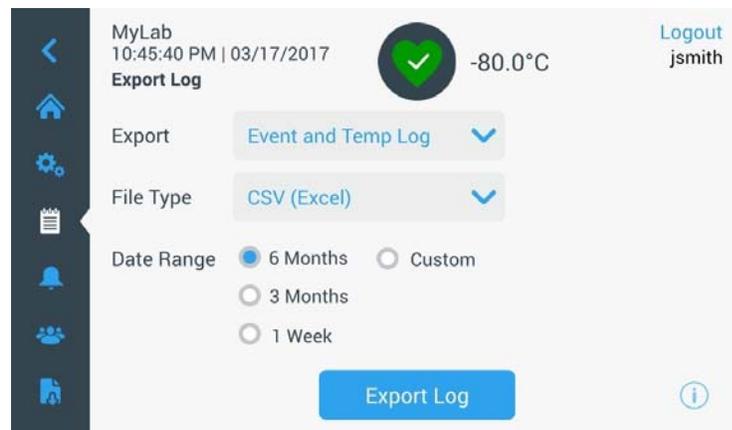
There is a drop down list based on the event types. The event types can be filtered and categorized into: Alarm, Door, User, Battery and Backup. When a filter is selected, the View All button on the right changes to Filter ON.



**Figure 27.** Event Log Screen with Filter

To export event and temperature logs,

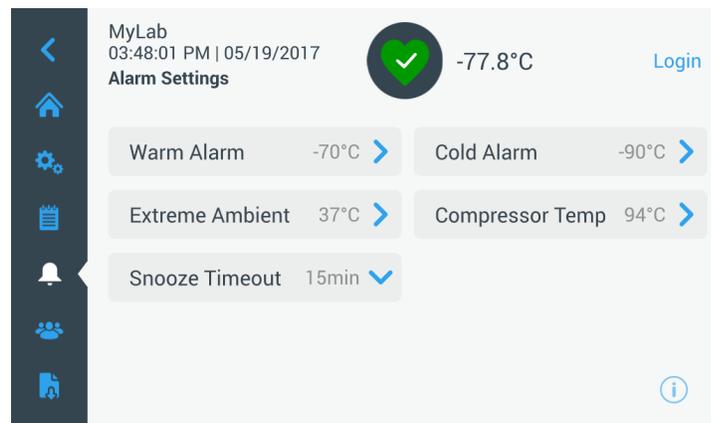
1. Choose the item to be exported from Export drop down list. The options available are: Event and Temperature Log, Event Log, Temperature Log and Reports.
2. Select the export format of the log or report.
3. A predefined or custom date range may be selected.
4. A USB drive must be inserted to store the log or report. Tap the Export Log button to download the log or report.



**Figure 28.** Export Log Screen

## 9.5 Alarms

By selecting the Alarm tab, the Alarm Settings screen is displayed. For more information on setpoints, refer to Section 9.3.1.



**Figure 29.** Alarm Settings Screen

## 9.6 Users

The User tab is used to display the Users screen. Access Mode is used to change the access mode for the system (full or secured), add a user to the system, and to import and/or export a user database. The following screen is displayed when the User icon is selected:



**Figure 30.** Users Screen

The Import button allows a user database to be imported.

**Note** *The database to be imported must be taken from another unit running the same software otherwise the system will not recognize the database.*

The user database can be exported using the Export button. A USB drive must be inserted for the data to be transferred.

### 9.6.1 Adding New Users

Select the Add User button to navigate to the Add a User screen:





**Figure 31.** Add a User Screen

Fill the information in the fields and tap Add User. An asterisk denotes a required field.

- **First Name:** Enter the user's first name.
- **Last Name:** Enter the user's last name.
- **Email:** Enter the user's email address.
- **User Name:** Enter the username as required. A default will appear based upon the email address entered.
- **Password:** Enter and confirm a user password. The password entered should be 3-16 characters long, have at least 1 uppercase, 1 number and 1 special character.
- **Phone:** Enter a user telephone number. Additional phone numbers can be added by tapping the '+' symbol.
- **User Role:** Select the access level for the user.
  - Admin: This user has access to change settings and manage profiles.
  - User Level 1: This user has access to change basic functionality such as temperature and alarm setpoints.
  - User Level 2: This user can view the temperature and alarm information but cannot change the settings.
- **Access ID:** Refer to Section 9.6.3 for more information on access cards.

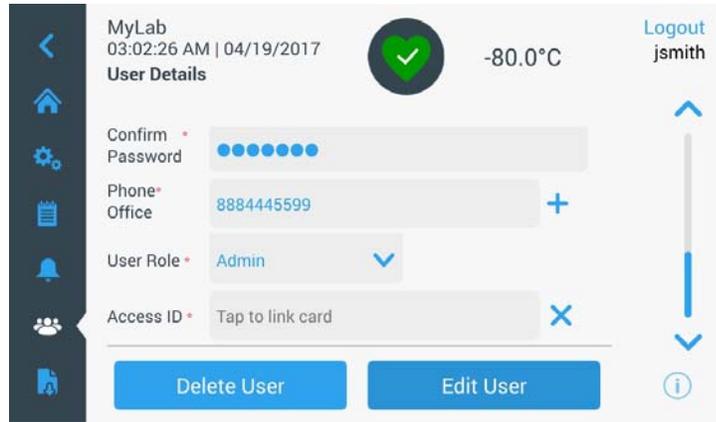
### 9.6.2 Editing and Deleting Users

To edit an existing user, tap the user entry in the Users screen and User Details screens will be displayed. Make the required edits by selecting the appropriate field and changing the information. After making the changes, tap the “Edit User” button and confirm.

To delete a user, select the Delete User button and confirm.

**Note** *This action cannot be undone.*





**Figure 32.** User Details Screen

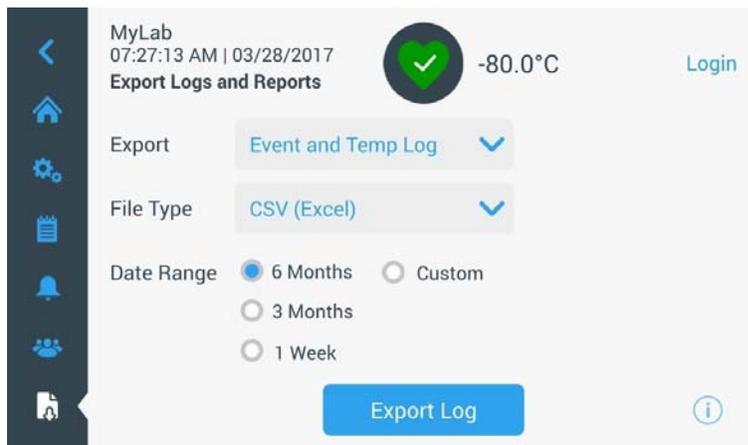
### 9.6.3 Managing Access Cards

When the Access Card system is installed, the User Details screen allows assigning cards to each user. Users may then swipe their access card to gain access to the freezer or login.

To associate an access card to a selected user, tap the Access ID field of the User Details screen (refer to Figure 32) and hold the card at the bottom center of the LCD screen. The Access ID field will automatically be populated and you will hear a beep. Only one card may be added per user.

## 9.7 Reports

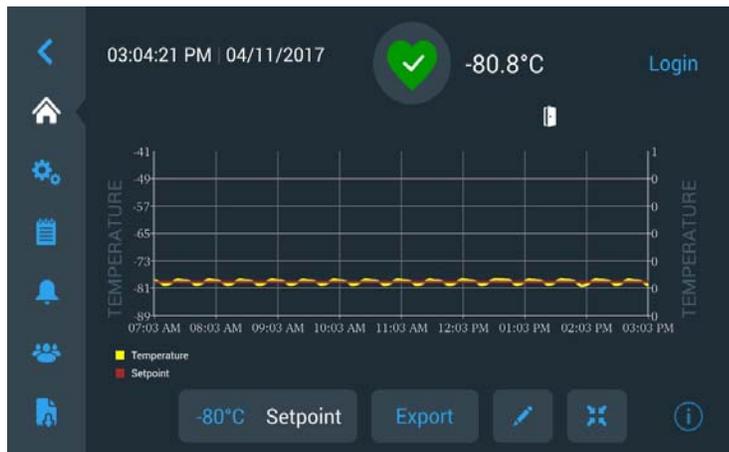
The Report tab is used to export a log or report. The Export Logs and Reports screen will be displayed when the Report icon is pressed. If Reports is selected as the Export type, the date range is limited to two options - one week (default) and one day. Refer to Section 9.4 for more information.



**Figure 33.** Export Logs and Reports Screen

## 9.8 Chart The chart displays temperature data.

There can be a maximum of two Y-axis. The left Y-axis is Primary and the right Y-axis is Secondary. Items displayed on the Primary axis are listed on the left and the ones displayed on the Secondary axis are listed on the right. Time is denoted on the X-axis.



**Figure 34.** Chart with items on Primary axis

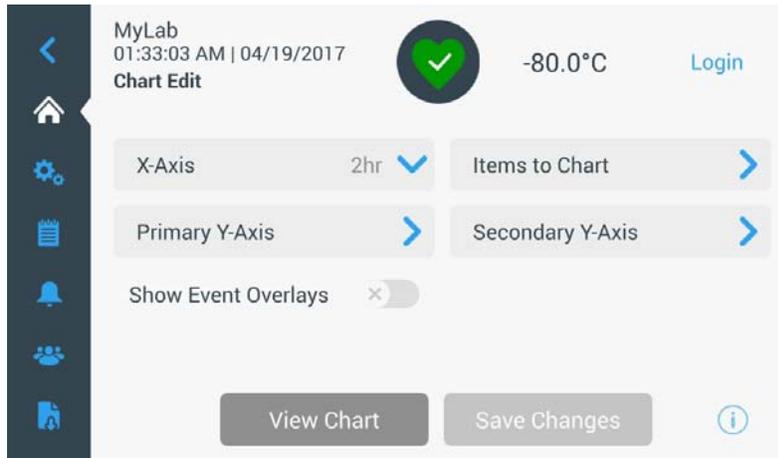


**Figure 35.** Chart with items on Primary and Secondary axes

The four buttons on the bottom of the screen are used to:

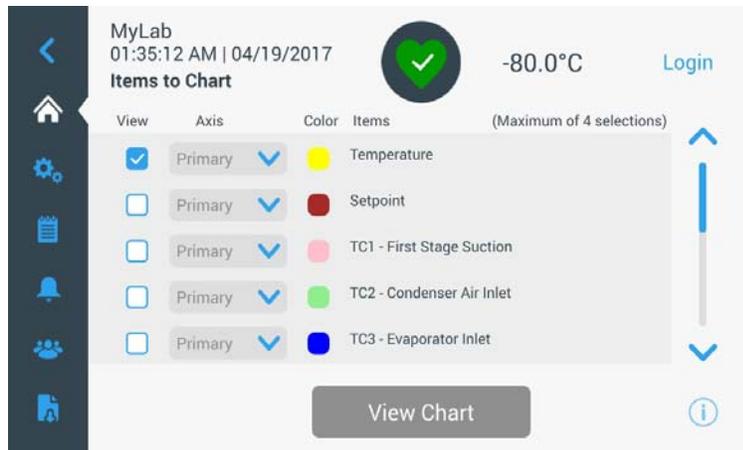
- Change the setpoint
- Export the chart
- Edit the chart, refer to Section 9.8.1 for more information
- Expand the chart to full screen or return to the default view.

**9.8.1 Editing the Chart** When the Pencil icon is selected, the Chart Edit screen is displayed:



**Figure 36.** Chart Edit Screen

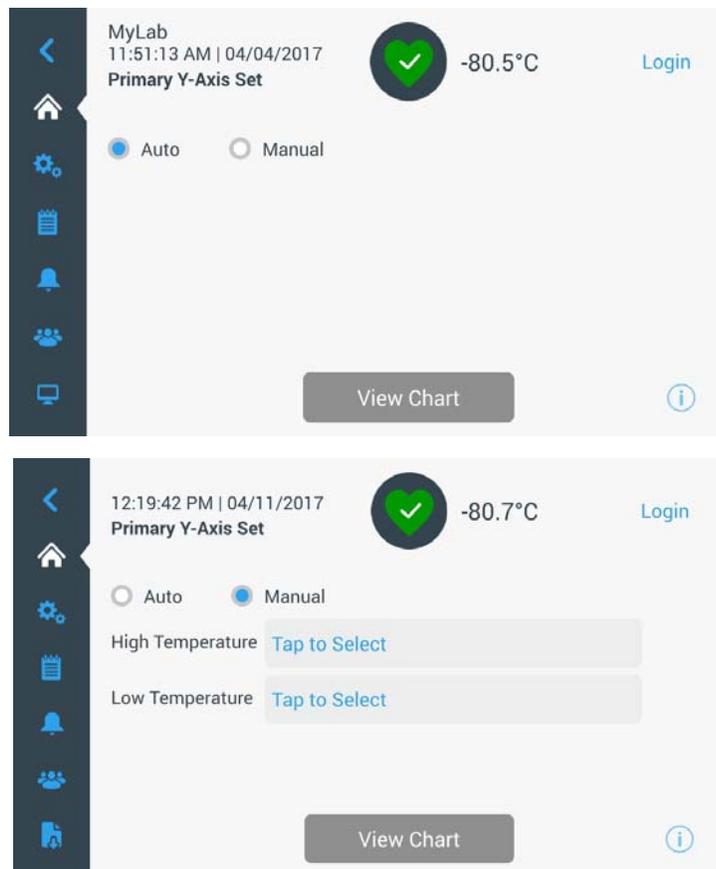
- The X-Axis button allows the user to select a time span of 2, 4, 6 or 8 hours.
- The Items to Chart button allows the user to select the items to be displayed on the chart. A maximum of four items may be selected.





**Figure 37.** Items to Chart Screen

- The Primary Y-Axis button allows the user to set the temperature range. If Auto is selected, the temperatures are automatically set. If Manual is selected, the user has to specify the high and low temperatures.



**Figure 38.** Primary Y-Axis Set screen

- The Secondary Y-Axis button allows the user to set the temperature range. If Auto is selected, the temperatures are automatically set. If Manual is selected, the user has to specify the high and low temperatures. Item associated with the Secondary Y-Axis will be displayed as a dotted line. The Axis On toggle button is used to enable the Secondary Y-axis.



**Figure 39.** Secondary Y-Axis Set Screen

- **Show Event Overlay:** Enable quick reference event icons to be displayed on the top of the chart.

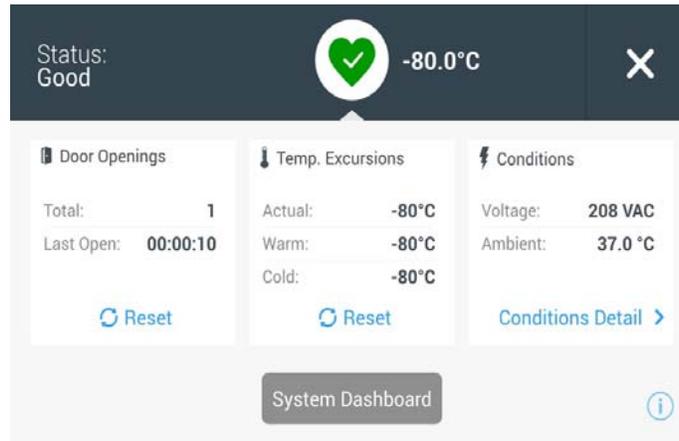
**9.8.2 Event Icons** Icons related to events will appear on the top of the chart.

Icon	Description
	Door Open
	Power Failure
	Temperature Alarm
	Service
	Multiple Events

# 10 Health Status and Alarm Management

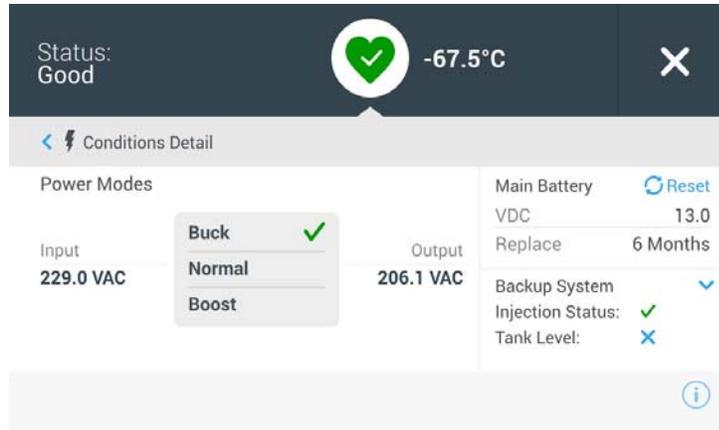
## 10.1 Health Status Overview

Selecting the Health Status (green heart) icon in the home screen displays the following screen:



**Figure 40.** Health Status Screen

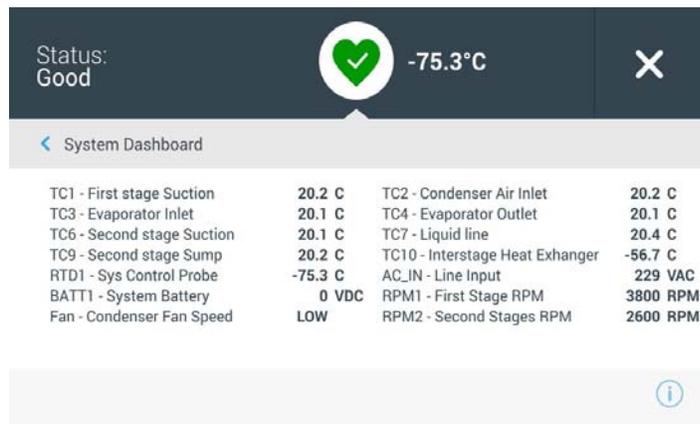
- **Door Openings:** Total displays the number of times the door has been opened since the last reset. Last Open displays the duration the door was last opened. Tap the Reset icon to change the Door Openings values to 0.
- **Temperature Excursions:** Actual displays the current cabinet temperature. Warm and Cold displays are the highest and lowest temperatures recorded since the last reset. Tap the Reset icon to set the Warm and Cold values to the current cabinet temperature.
- **Conditions:** Voltage displays the input voltage. Ambient displays the unit ambient temperature. By tapping the Conditions Detail button the following screen will be displayed:



**Figure 41.** Conditions Detail Screen

- **Power Modes:** Displays the state of the on-board voltage conditioning system.
- **Main Battery:** Displays the voltage of the main battery. There is also a 12 month countdown timer that should be reset every time the battery is replaced.
- **Backup System:** The BUS information will be displayed if a BUS is present.

Selecting the System Dashboard button in Figure 40 displays the on board sensor readings.



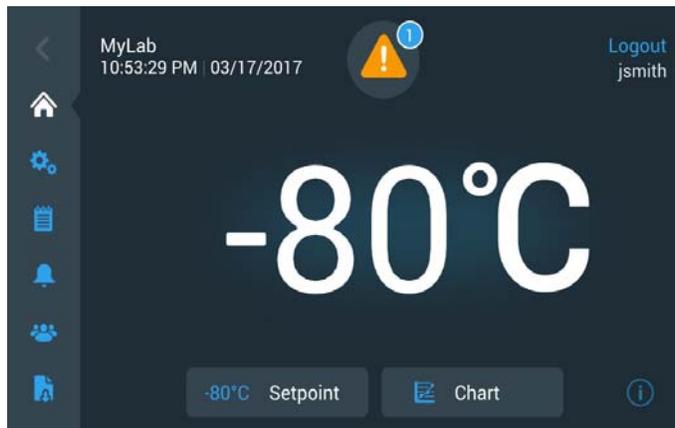
**Figure 42.** System Dashboard Screen

When a BUS is present, the BUS RTD and BUS battery values are also displayed.

## 10.2 Notifications / Cautions

In the event of a notification, the green heart is replaced by a yellow triangle icon. The number in the blue circle indicates the total number of notifications.

A yellow triangle on the home screen indicates less serious caution conditions, such as extended door openings and alarms that occurred in the past.



**Figure 43.** Home Screen with Notification / Caution

Tap the yellow triangle to display additional notification details.



**Figure 44.** Notification / Caution Details

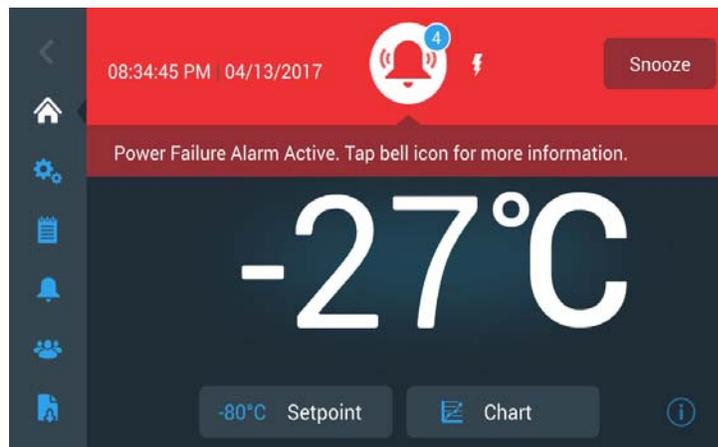
The Acknowledge button becomes highlighted in blue once a notification is selected by tapping the adjacent checkbox. The notification will be displayed until it is acknowledged. There is a link to the Event Log beside the Acknowledge button.

On the right panel, the Health Status details can also be viewed.

### 10.3 Alarms / Warning

In the event of an alarm condition, the green heart or yellow triangle is replaced by a red bell icon. A red bell alarm on the home screen indicates a serious alarm condition which must be corrected, such as a warm alarm or a power failure. The alarm can be silenced by tapping the Snooze button or the Red Bell icon which will mute the alarm for a duration set in the Snooze Timeout setting.

The number in the blue circle indicates the total number of alarms. The icon to the right of the bell helps identify the alarm type. The ticker message provides alarm details.



**Figure 45.** Home Screen with Alarm / Warning

Tap the red bell for additional information about the alarms.



**Figure 46.** Alarm / Warning Detail Screen

If an alarm and a notification occur at the same time, the alarm red bell icon is displayed and the list includes both alarms and notifications.

The Acknowledge button becomes highlighted in blue once an alarm is selected by tapping the adjacent checkbox. Active alarms cannot be cleared until the issue is addressed. There is a link to the Event Log beside the Acknowledge button.

On the right panel, the Health Status details can also be viewed.

For more information on various alarm conditions, refer to Appendix A: Alarm Summary.

## 11 Backup System (Optional)

When you purchase a built-in CO<sub>2</sub> or LN<sub>2</sub> optional backup system for the freezer, backup control is integrated into the main user interface.



**Note** Always purchase the cylinders which are equipped with siphon tubes for withdrawing liquid from the bottom of the cylinder. CO<sub>2</sub> cylinders must be kept at room temperature to function properly. LN<sub>2</sub> bottles are functional at any reasonable temperature.

### 11.1 CO<sub>2</sub> and LN<sub>2</sub> Precautions

The following are precautions for using liquid CO<sub>2</sub> and LN<sub>2</sub> backup systems.



**WARNING:** If a CO<sub>2</sub> or LN<sub>2</sub> cylinder falls and a valve is knocked off, the cylinder becomes a deadly and completely unguided missile. Transport the cylinders in a handtruck or cart with secure chain ties for the cylinder. After cylinders are connected to the equipment, securely attach them with chains to a solid, stationary object such as a building column.



**WARNING:** CO<sub>2</sub> and LN<sub>2</sub> liquids are non-poisonous but are very cold and will burn unprotected skin. Always wear protective eyewear and clothing when changing cylinders or working on the piping systems attached to an active source of liquid refrigerant.



**WARNING:** The gases produced by evaporation of CO<sub>2</sub> or LN<sub>2</sub> are non-poisonous but displace the oxygen in a confined space and can cause asphyxiation. Do not store the cylinders in subsurface or enclosed areas.



**CAUTION:** When closing the cylinder valve, make sure that the injection solenoid is energized to allow all the liquid to bleed off instead of being trapped in the supply hose. Failure to do this results in activation of the pressure relief device, which could damage the freezer and requires replacing if it is activated.



**CAUTION:** For models ordered with factory installed built-in backup systems, the flow of liquid CO<sub>2</sub> or LN<sub>2</sub> will be discontinued if the door is opened during operation of the backup system. For units operated with free-standing, field installed type backup system, the flow of liquid CO<sub>2</sub> or LN<sub>2</sub> will be discontinued upon door opening only if the switch provided with the free-standing package is installed on the freezer.

## 11.2 Installation

Field installed systems are supplied with complete installation and operating instructions. If your system is factory installed, the freezer is shipped with a coiled length of hose to connect the freezer to the bottles:

- 1/4" Flexible Hose with fittings for connection to the CO<sub>2</sub> supply.
- 1/2" Flexible Hose with fittings for connection to the LN<sub>2</sub> supply.

To install,

1. Straighten the coiled hose.
2. Connect one end to the labeled connection on the freezer.
  - Tighten the nut two flats past finger tight, approximately 120 degrees.

**Note** For CO<sub>2</sub>, remove the threaded fitting from the nut on the end of the copper tubing to access nut for connection to the freezer. Discard the threaded fitting.

3. Attach the other end to the supply bottle or building supply fitting.

- For CO<sub>2</sub>:
  - Remove Nipple from adapter (NPT Connection). Remove cable tie to release alternative nut and washer. Ensure the correct nut fitting is supplied over the nipple (US or European).
  - Add 2 wraps of Teflon tape clockwise to the 1/4" NPT fitting (on the nipple) when viewed from the threads. Tighten the NPT fittings approximately 2 turns from finger tight (approximately 720°).

**Note** The top of the nipple has a hex configuration, allowing for use of a wrench when the nut is pulled down.

- Add washer to nipple inside of nut (unless CO<sub>2</sub> supply has a built in washer).

**Note** Small raised area of washer fits into groove of nipple. The washer will feel snug when trying to shift side to side on nipple. The washers are designed for a limited number of attachments/disconnections from the supply and may wear over time. If washer appears worn and causes CO<sub>2</sub> leakage, replace washer (Part Number 45705H03).

- Wrench tighten the supply nut to the supply.
- For LN<sub>2</sub>:
  - Attach the fitting to the supply and wrench tighten.

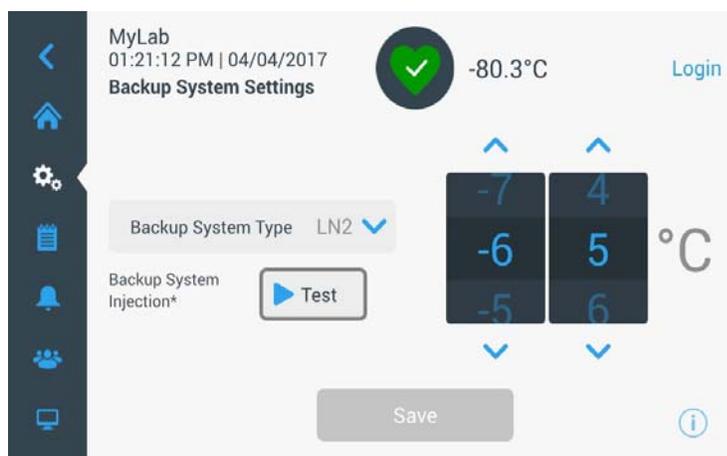


**Note** Do not twist, torque, or subject the flexible hose to sharp bends. Doing so may shorten the life of the hose.

### 11.3 Start Up

To activate the backup system:

1. Follow the instructions in Section 8 to turn on the freezer and set temperature and alarm set points.
2. Select the backup type and backup set point on the Backup Setting Screen below, which can be accessed through the Controls option in the Settings tab (refer to Figure 24).



**Figure 47.** Backup System Settings Screen

3. Tap the Save button to save the changes made.

### 11.4 Operation

When the backup system is in operation, the parameters can be viewed and configured on the settings screen.

Once the backup system has been activated, it can be tested by pressing the Test button. The system will inject as long as the button is being pressed.

The backup system can run for a minimum of 24 hours on battery power.

On average, a backup system in operation uses 8 to 10 lbs. per hour of CO<sub>2</sub> or LN<sub>2</sub> at an ambient temperature of 25°C. This rate will vary depending on setpoint, load, ambient temperature and freezer size.

## 12 Chart Recorders (Optional)

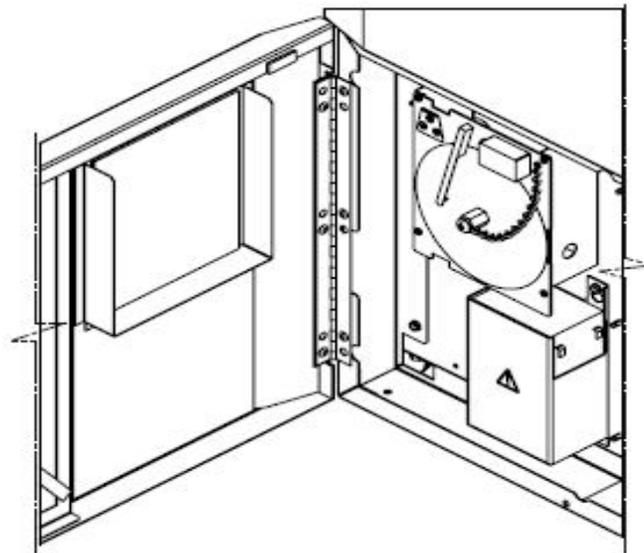
Panel-mounted six-inch seven-day recorders are available as options for all freezer models except for the smallest (300 box capacity) models.

### 12.1 Set Up and Operation

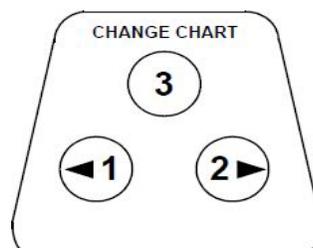
To prepare the recorder to function properly, complete the following steps:

1. Open the grille door to access the recorder.
2. Install clean chart paper (refer to Section 12.2).
3. Remove the plastic cap from the pen stylus or ink pen and close the recorder door.

Recorder operation begins when the system is powered on. The recorder may not respond until the system reaches temperatures within the recorder's range.



**Figure 48.** Chart Recorder



**Figure 49.** Chart Buttons



**CAUTION:** Do not use sharp or pointed objects to depress the chart buttons. This may cause permanent damage to the recorder.

## 12.2 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) 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 (a small groove on the left side of the recorder panel).
4. Replace the center nut and hand tighten. Press the Change Chart button again to resume temperature recording.

## 12.3 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 set point temperature. Continue steady operation for at least two hours to provide adequate time for recorder response.
2. Measure cabinet center temperature with a calibrated temperature monitor.
3. Compare the recorder temperature to the measured cabinet temperature. If necessary, adjust recorder by pressing the left (#1) and right (#2) chart buttons.

**Note** *The stylus does not begin to move until the top center button (#3) is held for five seconds.*

## 13 Maintenance and Troubleshooting



**WARNING:** Unauthorized repair of your freezer will invalidate your warranty. Contact Technical Service at 1-800-438-4851 for additional information.



**CAUTION:** Maintenance should only be performed by trained personnel.

### 13.1 Cleaning the Condenser

Clean the condenser at least every six months; more often if the laboratory area is dusty.

To clean the condenser, complete the following steps:

1. Pull the grille door open.
2. Vacuum the condenser.
3. Inspect the filter cleanliness and clean as required.
4. Close the grille door.

### 13.2 Cleaning the Condenser Filter

Clean the condenser filters every two or three months.

There are two condenser filters: a main filter and a lower filter for extra air flow into the condenser.

1. Pull the grille door open.
2. Remove the filters.
3. Shake the filters to remove loose dust, rinse the filters in clean water, shake the excess water from the filters, and replace the filters.
4. Close the grille door.

### 13.3 Gasket Maintenance

Periodically check the gaskets around the door for punctures or tears. Leaks are indicated by a streak of frost which forms at the point of gasket failure. Make sure that the cabinet is level (refer to Section 7.3 for leveling information).

Keep the door gaskets clean and frost free by wiping gently with a soft cloth.

### **13.4 Defrosting the Freezer**

Defrost the freezer once per year or whenever the ice buildup exceeds 3/8". To defrost, complete the following steps:

1. Remove all products and place in another ULT cabinet.
2. Turn off the freezer.
3. Open the outer door and all inner doors.
4. Let the freezer stand with doors open for at least 24 hours. This allows both the interior and foamed refrigerant system to warm to room temperature.
5. Dispose of the ice and wipe out any water standing in the bottom of the cabinet.
6. If there is freezer odor, wash the interior with a solution of baking soda and warm water.
7. Clean the exterior with any common household cleaner.
8. Close the doors, restart the freezer and reload. Refer to Section 5.3 to follow the instructions.

### **13.5 Battery Maintenance**

Have a certified technician replace the battery at least every 12 months. Be sure to reset the battery replacement timer when the battery is replaced.

## 14 Warranty **Be sure to register your warranty online:**

[www.thermofisher.com/labwarranty](http://www.thermofisher.com/labwarranty)

### **THERMO FISHER SCIENTIFIC USA FREEZER WARRANTY FOR Revco RLE series, Forma 89000 series, HERAFreeze HLE series and Thermo Scientific TLE series**

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

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

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

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

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

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

## 15 Warranty (International)

### **THERMO FISHER SCIENTIFIC FREEZER INTERNATIONAL WARRANTY FOR Revco RLE series, Forma 89000 series, HERAfreeze HLE series and Thermo Scientific TLE series**

The Warranty Period starts two months from the date your equipment is shipped from our facility. This allows for shipping time so the warranty will go into effect at approximately the same time your equipment is delivered. The warranty protection extends to any subsequent owner during the warranty period. Dealers who stock our equipment are allowed an additional four months for delivery and installation, providing the warranty card is completed and returned to the Technical Services Department.

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

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

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

Your local Thermo Fisher Scientific 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 Thermo Fisher Scientific office or local distributor.

We're ready to answer your questions on equipment warranty, operation, maintenance, service, and special applications. Outside the USA, contact your local Thermo Fisher Scientific office or distributor for warranty information.

## Appendix A: Alarm Summary

Alarm Summary					
Alarm Type	Tone	Heart Status	Remote Alarm Event	Ringback*	Message Details
Warm Alarm Active	High	Red	Yes	Yes	<p><b>Message Header:</b> Warm Alarm</p> <p><b>Message Details:</b> The freezer temperature has exceeded the warm alarm set point of XX.X C. Prolonged door openings and warm product loading may cause warm alarms.</p> <p><b>Ticker Message:</b> Warm Alarm active. Tap bell icon for more information.</p>
Warm Alarm Inactive	None	Yellow	None	None	<p><b>Message Header:</b> Warm Alarm Recovered</p> <p><b>Message Details:</b> The freezer temperature has recovered from a warm alarm event. Select check box and tap Acknowledge to clear this notification.</p> <p><b>Ticker Message:</b> None</p>
Cold Alarm Active	High	Red	Yes	Yes	<p><b>Message Header:</b> Cold Alarm</p> <p><b>Message Details:</b> The freezer temperature has exceeded the cold alarm set point of XX.X C</p> <p><b>Ticker Message:</b> Cold Alarm active. Tap bell icon for more information.</p>
Cold Alarm Inactive	None	Yellow	None	None	<p><b>Message Header:</b> Cold Alarm Recovered</p> <p><b>Message Details:</b> The freezer temperature has recovered from a cold alarm event. Select check box and tap Acknowledge to clear this notification.</p> <p><b>Ticker Message:</b> None</p>
Door Open Ajar	High	Red	Yes	Yes	<p><b>Message Header:</b> Door Open Alarm</p> <p><b>Message Details:</b> Door open for greater than 3 minutes will cause door open alarm.</p> <p><b>Ticker Message:</b> Door Open Alarm active. Tap bell icon for more information.</p>
Door Open Recovered	None	Yellow	None	None	<p><b>Message Header:</b> Door Open Recovered</p> <p><b>Message Details:</b> Door open alarm has recovered. Select check box and tap Acknowledge to clear this notification.</p> <p><b>Ticker Message:</b> Door Open Alarm active. Tap bell icon for more information.</p>

Alarm Summary					
Alarm Type	Tone	Heart Status	Remote Alarm Event	Ringback*	Message Details
System Battery Low	None	Yellow	None	None	<p><b>Message Header:</b> System Battery Low</p> <p><b>Message Details:</b> System battery is XX.XV. Recommend replacing battery. Contact customer service for battery replacement.</p> <p><b>Ticker Message:</b> None</p>
System Battery PM Expiry	None	Yellow	None	None	<p><b>Message Header:</b> System Battery Low</p> <p><b>Message Details:</b> System battery is XX.XV. Recommend replacing battery. Contact customer service for battery replacement.</p> <p><b>Ticker Message:</b> None</p>
BUS Battery Low	None	Yellow	None	None	<p><b>Message Header:</b> BUS Battery Low</p> <p><b>Message Details:</b> BUS battery is XX.XV. Recommend replacing battery. Contact customer service for battery replacement.</p> <p><b>Ticker Message:</b> None</p>
BUS Battery PM Expiry	None	Yellow	None	None	<p><b>Message Header:</b> Battery PM Expiry</p> <p><b>Message Details:</b> Recommend to replace battery. Reset battery expiration time on XXX screen to clear this notification.</p> <p><b>Ticker Message:</b> None</p>
Extreme Ambient	None	Yellow	None	None	<p><b>Message Header:</b> Extreme Ambient Notification</p> <p><b>Message Details:</b> Ambient temperature has exceeded the Extreme Ambient set point of XX.XC. Ensure room temperature is properly controlled.</p> <p><b>Ticker Message:</b> None</p>
Control Probe Failure	High	Red	Yes	Yes	<p><b>Message Header:</b> Control Probe Failure Alarm</p> <p><b>Message Details:</b> Cannot display cabinet temperature. The freezer will continue to operate in full run mode. Contact customer service.</p> <p><b>Ticker Message:</b> Control Probe Failure Alarm active. Tap bell icon for more information.</p>
Control Probe Failure has been	None	Yellow	None	None	<p><b>Message Header:</b> Control Probe Failure Recovered</p> <p><b>Message Details:</b> Control Probe Failure Alarm has been cleared. Press the yellow triangle to clear this notification.</p> <p><b>Ticker Message:</b> None</p>

Alarm Summary					
Alarm Type	Tone	Heart Status	Remote Alarm Event	Ringback *	Message Details
Heat Exchange Probe Failure	High	Red	Yes	Yes	<p><b>Message Header:</b> Heat Exchanger Probe Failure Notification</p> <p><b>Message Details:</b> The freezer will continue to operate with current freezer set points, but cabinet temperature variation will increase. Contact customer service.</p> <p><b>Ticker Message:</b> Control Probe Failure Alarm active. Tap bell icon for more information.</p>
Heat Exchange Probe Failure has been	None	Yellow	None	None	<p><b>Message Header:</b> Heat Exchanger Probe Failure Recovered</p> <p><b>Message Details:</b> Heat Exchange Probe Failure Alarm has been cleared. Press the yellow triangle to clear this notification.</p> <p><b>Ticker Message:</b> None</p>
TC1-9 Failure	None	Yellow	None	None	<p><b>Message Header:</b> TCXX Out of Range Notification</p> <p><b>Message Details:</b> Information TC's has malfunctioned. This doesn't affect the performance of the unit. Contact service for further assistance.</p> <p><b>Ticker Message:</b> None</p>
TC1-9 has been	None	Yellow	None	None	<p><b>Message Header:</b> TCXX Out of Range Notification Recovered</p> <p><b>Message Details:</b> Information TCXX has recovered. Select check box and tap Acknowledge to clear this notification.</p> <p><b>Ticker Message:</b> None</p>
Main to UI Lost Communication	High	Red	Yes	Yes	<p><b>Message Header:</b> Lost Communication Alarm</p> <p><b>Message Details:</b> A communication error has occurred within the system. Contact customer service.</p> <p><b>Ticker Message:</b> Lost Communication Alarm Active. Tap bell icon for more information.</p>
Main to UI Lost Communication Inactive	None	Yellow	None	None	<p><b>Message Header:</b> Lost Communication Alarm Recovered</p> <p><b>Message Details:</b> A communication error has recovered. Select check box and tap Acknowledge to clear this notification.</p> <p><b>Ticker Message:</b> None</p>

Alarm Summary					
Alarm Type	Tone	Heart Status	Remote Alarm Event	Ringback*	Message Details
BUS Lost Communication	High	Red	Yes	Yes	<p><b>Message Header:</b> Backup System Communication Alarm</p> <p><b>Message Details:</b> A communication error has occurred within the back up system. Contact customer service.</p> <p><b>Ticker Message:</b> Bus Lost Communication Alarm Active. Tap bell icon for more information.</p>
BUS Lost Communication Inactive	None	Yellow	None	None	<p><b>Message Header:</b> BUS Lost Communication Alarm Recovered</p> <p><b>Message Details:</b> A BUS communication error has recovered. Select check box and tap Acknowledge to clear this notification.</p> <p><b>Ticker Message:</b> None</p>
Failure to Reach Setpoint	None	Yellow	None	None	<p><b>Message Header:</b> Unable to Reach Setpoint Notification</p> <p><b>Message Details:</b> Door openings or product loading may cause this notification. Allow unit to stabilize. If condition persists, contact customer service</p> <p><b>Ticker Message:</b> None</p>
Compressor Temperature	None	Yellow	None	None	<p><b>Message Header:</b> Compressor Temperature Notification</p> <p><b>Message Details:</b> A refrigeration system temperature has exceeded the compressor temperature setpoint of XX.X C. Contact customer service.</p> <p><b>Ticker Message:</b> None</p>
Inefficient Buck / Boost	None	Yellow	None	None	<p><b>Message Header:</b> Inefficient buck / boost notification</p> <p><b>Message Details:</b> Freezer input voltage is out of range. Ensure proper supply voltage is applied. If unable to clear this notification, contact customer service.</p> <p><b>Ticker Message:</b> None</p>
Power Failure Alarm	High	Red	Yes	Yes	<p><b>Message Header:</b> Power Failure Alarm</p> <p><b>Message Details:</b> Unit in power failure mode. Display operating on battery power. Check unit plug, unit circuit breaker in the ON position, and supply voltage.</p> <p><b>Ticker Message:</b> Power Failure Alarm Active. Tap bell icon for more information.</p>

Alarm Summary					
Alarm Type	Tone	Heart Status	Remote Alarm Event	Ringback *	Message Details
Power Failure Inactive	None	Yellow	None	None	<p><b>Message Header:</b> Power Failure Recovered</p> <p><b>Message Details:</b> Power has been restored. Press the yellow triangle to clear this notification.</p> <p><b>Ticker Message:</b> None</p>
Clean Filter Active	None	Yellow	None	None	<p><b>Message Header:</b> Clean Filter Notification</p> <p><b>Message Details:</b> Recommend to clean the filter and condenser. If unable to clear notification, contact customer service.</p> <p><b>Ticker Message:</b> None</p>
Refrigeration System Failure	High	Red	Yes	Yes	<p><b>Message Header:</b> System Refrigeration Failure Alarm</p> <p><b>Message Details:</b> An error has occurred within the refrigeration system. Contact customer service.</p> <p><b>Ticker Message:</b> System Refrigeration Failure Alarm. Tap bell icon for more information</p>
Refrigeration System Failure Inactive	None	Red	None	None	<p><b>Message Header:</b> System Refrigeration Failure Recovered</p> <p><b>Message Details:</b> An error in refrigeration system has been recovered. Select check box and tap Acknowledge to clear this notification.</p> <p><b>Ticker Message:</b> None</p>
Wrong Power Alarm	High	Red	Yes	None, constant audible	<p><b>Message Header:</b> Wrong Power Alarm</p> <p><b>Message Details:</b> The unit has detected the wrong power connected. Please verify the proper voltage.</p> <p><b>Ticker Message:</b> Wrong Power Alarm. Tap bell icon for more information.</p>
Wrong Power Alarm Inactive	None	Yellow	None	None	<p><b>Message Header:</b> Wrong Power Recovered</p> <p><b>Message Details:</b> Wrong Power has been recovered. Select check box and tap Acknowledge to clear this notification.</p> <p><b>Ticker Message:</b> None</p>
Wrong Model Alarm	High	Red	Yes	None, constant audible	<p><b>Message Header:</b> Invalid Control Model Alarm</p> <p><b>Message Details:</b> Invalid Control Model Alarm. Contact service to ensure the correct model is selected for the system to avoid cargo loss.</p> <p><b>Ticker Message:</b> Invalid Control Model Alarm. Tap bell icon for more information.</p>

Alarm Summary					
Alarm Type	Tone	Heart Status	Remote Alarm Event	Ringback*	Message Details
Wrong Model Alarm Inactive	None	Yellow	None	None	<p><b>Message Header:</b> Invalid Control Model Recovered</p> <p><b>Message Details:</b> Invalid Control Model has been recovered. Select check box and tap Acknowledge to clear this notification.</p> <p><b>Ticker Message:</b> None</p>
Firmware Build Incompatible	High	Red	Yes	Yes	<p><b>Message Header:</b> Firmware Build Incompatible Alarm</p> <p><b>Message Details:</b> Firmware build indicates incompatibility that can result in modules to be non-coherent.</p> <p><b>Ticker Message:</b> Firmware build incompatible. Tap bell icon for more information.</p>
Firmware Build Incompatible Inactive	None	Yellow	None	None	<p><b>Message Header:</b> Firmware Build Incompatibility Recovered</p> <p><b>Message Details:</b> Firmware build Incompatibility recovered. Select check box and tap Acknowledge to clear this notification.</p> <p><b>Ticker Message:</b> None</p>

\*Ringback refers to the audible alarm activating again after the user presses the Snooze button. The time can be set in the 'Snooze Timeout' option in the Alarm Settings screen.

## Appendix B: Event Log Detail

Event Log Entry Detail	
Item	Message
System Parameter Change	<b>Header:</b> Control Settings Change <b>Role:</b> Logged in personnel or anonymous in full access mode <b>Mode:</b> Full access or secured <b>Message Detail:</b> The following parameters has been changes to: Parameter 1 from: XX.XX to XX.XX
User Interface Parameter Change	<b>Header:</b> User Interface Settings Change <b>Role:</b> Logged in personnel or anonymous in full access mode <b>Mode:</b> Full access or secured <b>Message Detail:</b> The following parameters has been changes to: Parameter 1 from: XX.XX to XX.XX
Door Open	<b>Header:</b> Door Open <b>Role:</b> Logged in personnel or blank if not HID system
Door Close	<b>Header:</b> Door Close <b>Role:</b> Logged in personnel or blank if not HID system
User Login	<b>Header:</b> User jsmith has logged in
User Logout	<b>Header:</b> User jsmith has logged out
Door History Reset	<b>Header:</b> Door History Reset <b>Role:</b> Logged in personnel or anonymous in full access mode <b>Mode:</b> Full access or secured <b>Message Detail:</b> Door Usage history has been reset
Temperature Excursion Reset	<b>Header:</b> Temperature Excursion History Reset <b>Role:</b> Logged in personnel or anonymous in full access mode <b>Message Detail:</b> Temperature Excursion history has been reset
Add User Event	<b>Header:</b> Add a User to User database <b>Role:</b> Logged in personnel or anonymous in full access mode <b>Mode:</b> Full access or secured <b>Message Detail:</b> User jsmith has been added to user database
Remove User Event	<b>Header:</b> Remove a User to User database <b>Role:</b> Logged in personnel or anonymous in full access mode <b>Mode:</b> Full access or secured <b>Message Detail:</b> User jsmith has been removed to user database
Edit User Event	<b>Header:</b> Edit a User to User database <b>Role:</b> Logged in personnel or anonymous in full access mode <b>Mode:</b> Full access or secured <b>Message Detail:</b> User jsmith details has been modified.

Event Log Entry Detail	
Item	Message
Buck / Boost Change Event	<p><b>Header:</b> Buck / Boost status Change</p> <p><b>Message Detail:</b>            Buck / Boost changed from X to Y            Line Input Voltage at time of change is            Compensated Voltage at time of change is</p>
BUS Injection	<b>Header:</b> BUS Injection on or off
BUS Pressure switch	<b>Header:</b> BUS Pressure Switch active or inactive
Reset to Factory default	<p><b>Header:</b> Reset to Factory Defaults</p> <p><b>Role:</b> Logged in personnel or anonymous in full access mode</p> <p><b>Mode:</b> Full access or secured</p> <p><b>Message Detail:</b> System has been restored to factory defaults</p>
Powered Up Event	<p><b>Header:</b> Softswitch Power Up event</p> <p><b>Role:</b> Logged in personnel or anonymous in full access mode</p> <p><b>Mode:</b> Full access or secured</p> <p><b>Message Detail:</b> System has been user turned on</p>
Powered Down Event	<p><b>Header:</b> Softswitch Power Down event</p> <p><b>Role:</b> Logged in personnel or anonymous in full access mode</p> <p><b>Mode:</b> Full access or secured</p> <p><b>Message Detail:</b> System has been user turned off</p>
HID Access	<b>Header:</b> User Jsmith has accessed the unit
Invalid HID	<b>Header:</b> Invalid HID tried to access unit
Export Event	<p><b>Header:</b> (Temperature, temperature and event log or pdf report) has been exported</p> <p><b>Role:</b> Logged in personnel or anonymous in full access mode</p> <p><b>Mode:</b> Full access or secured</p> <p><b>Message Detail:</b>            (Temperature, temperature and event log or pdf report) has been exported            Date Range of export is from X to Y            File format exported is (pdf, csv or PUC)</p>
Configuration import / export	<p><b>Header:</b> Configuration import or export initiated</p> <p><b>Role:</b> Logged in personnel or anonymous in full access mode</p> <p><b>Mode:</b> Full access or secured</p> <p><b>Message Detail:</b> Following items has been imported or exported to USB</p> <ul style="list-style-type: none"> <li>- Contact info</li> <li>- User database</li> <li>- System and User Interface settings</li> <li>- Temperature and event log history</li> </ul>

# WEEE Compliance

**WEEE Compliance.** This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2012/19/EU. It is marked with the following symbol. Thermo Fisher Scientific has contracted with one or more recycling/disposal companies in each EU Member State, and this product should be disposed of or recycled through them. Further information on our compliance with these Directives, the recyclers in your country, and information on Thermo Scientific products which may assist the detection of substances subject to the RoHS Directive are available at [www.thermofisher.com](http://www.thermofisher.com) under Services & Support.

## Great Britain



**WEEE Konformität.** Dieses Produkt muss die EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2012/19/EU erfüllen. Das Produkt ist durch folgendes Symbol gekennzeichnet. Thermo Fisher Scientific hat Vereinbarungen getroffen mit Verwertungs-/Entsorgungsanlagen in allen EU-Mitgliederstaaten und dieses Produkt muss durch diese Firmen wiederverwertet oder entsorgt werden. Mehr Informationen über die Einhaltung dieser Anweisungen durch Thermo Scientific, die Verwerter und Hinweise die Ihnen nützlich sein können, die Thermo Fisher Scientific Produkte zu identifizieren, die unter diese RoHS-Anweisung fallen, finden Sie unter [www.thermofisher.com](http://www.thermofisher.com) unter Services & Support.

## Deutschland



**Conformità WEEE.** Questo prodotto deve rispondere alla direttiva dell'Unione Europea 2012/19/EU in merito ai Rifiuti degli Apparecchi Elettrici ed Elettronici (WEEE). È marcato col seguente simbolo. Thermo Fisher Scientific ha stipulato contratti con una o diverse società di riciclaggio/smaltimento in ognuno degli Stati Membri Europei. Questo prodotto verrà smaltito o riciclato tramite queste medesime. Ulteriori informazioni sulla conformità di Thermo Fisher Scientific con queste Direttive, l'elenco delle ditte di riciclaggio nel Vostro paese e informazioni sui prodotti Thermo Scientific che possono essere utili alla rilevazione di sostanze soggette alla Direttiva RoHS sono disponibili sul sito [www.thermofisher.com](http://www.thermofisher.com) in Servizi e Supporto.

## Italia



**Conformité WEEE.** Ce produit doit être conforme à la directive euro-péenne (2012/19/EU) des Déchets d'Équipements Électriques et Électroniques (DEEE). Il est marqué par le symbole suivant. Thermo Fisher Scientific s'est associé avec une ou plusieurs compagnies de recyclage dans chaque état membre de l'union européenne et ce produit devrait être collecté ou recyclé par celles-ci. Davantage d'informations sur la conformité de Thermo Fisher Scientific à ces directives, les recycleurs dans votre pays et les informations sur les produits Thermo Fisher Scientific qui peuvent aider la détection des substances sujettes à la directive RoHS sont disponibles sur [www.thermofisher.com](http://www.thermofisher.com) sous Services et Assistance.

## France



## Important

For your future reference and when contacting the factory, please have the following information readily available:

Model Number: \_\_\_\_\_

Serial Number: \_\_\_\_\_

Date Purchased: \_\_\_\_\_

The above information can be found on the dataplate attached to the equipment. If available, please provide the date purchased, the source of purchase (manufacturer or specific agent/rep organization), and purchase order number.

---

### IF YOU NEED ASSISTANCE:

Thermo Scientific products are backed by a global technical support team ready to support your applications. We also offer cold storage accessories, including remote alarms, temperature recorders and validation services. Visit [www.thermofisher.com](http://www.thermofisher.com) or call:

<b>USA/Canada</b>	+1 866 984 3766	<b>Germany international</b>	+49 6184 90 6000
<b>India toll free</b>	1800 22 8374	<b>Germany national toll free</b>	0800 1 536 376
<b>India</b>	+91 22 6716 2200	<b>Italy</b>	+32 02 95059 552
<b>China</b>	+800 810 5118 (or) +400 650 5118	<b>Netherlands</b>	+31 76 579 55 55
<b>Japan</b>	+81-120-753-670	<b>Nordic/Baltic/CIS countries</b>	+358 9 329 10200
<b>Australia</b>	+61 39757 4300	<b>Russia</b>	+7 812 703 42 15
<b>Austria</b>	+43 1 801 40 0	<b>Spain/Portugal</b>	+34 93 223 09 18
<b>Belgium</b>	+32 53 73 42 41	<b>Switzerland</b>	+41 44 454 12 22
<b>France</b>	+33 2 2803 2180	<b>UK/Ireland</b>	+44 870 609 9203
<b>New Zealand</b>	+64 9 980 6700	<b>Other Asian countries</b>	+852 2885 4613
		<b>Countries not listed</b>	+49 6184 90 6000

Thermo Fisher Scientific Inc.

275 Aiken Road  
Asheville, NC 28804  
United States  
[www.thermofisher.com](http://www.thermofisher.com)

**thermo**scientific

328398H01 Rev. C