



Fisher Scientific

Isotemp Plus

Chromatography and General Purpose Laboratory Refrigerator with Alarm Version Controls

Isotemp Plus Laboratory Refrigerators

<u>120V Versions</u>	<u>230V Versions</u>
13-986-120A,	13-986-120RA
13-986-122A,	13-986-122RA
13-986-124A,	13-986-124RA
13-986-126A,	13-986-126RA
13-986-128A,	13-986-128RA
13-986-128SA,	13-986-128SRA
13-986-130A,	13-986-130RA
13-986-132A,	13-986-132RA
	13-986-121A,
	13-986-123A,
	13-986-125A,
	13-986-127A,
	13-986-129A,
	13-986-129SA,
	13-986-131A,
	13-986-133A.
	13-986-121RA
	13-986-123RA
	13-986-125RA
	13-986-127RA
	13-986-129RA
	13-986-129SRA
	13-986-131RA
	13-986-133RA

Isotemp Plus Chromatography Refrigerators

Isotemp Plus Laboratory Refrigerators with Drawers

<u>120V Versions</u>	<u>230V Versions</u>
13-986-134DA,	13-986-135DA,
13-986-138DA,	13-986-139DA,
13-986-140DA,	13-986-141DA,
13-986-134DRA	13-986-135DRA
13-986-138DRA	13-986-139DRA
13-986-140DRA	13-986-141DRA

RA – Alarm controller and chart recorder

Table of Contents

Introduction	3
Safety Precautions	3
Performance Characteristics	3
Installation	4
Unpacking.....	4
Visible Loss or Damage	4
Concealed Loss or Damage.....	4
Packing List	5
Selecting a Location	5
Installing and Leveling Casters	5
Shelves.....	6
Electrical Connection	6
Controls	7
Panel Controls	7
Displays and Features	7
Key Switch.....	7
Temperature Display	8
Unit's LED's.....	8
Mode LED	8
Cool LED	8
Alarm LED's.....	8
Keypads	9
Interior Light Switch.....	9
Operation	10
Controller Set Up (Menus)	10
Temperature Setpoint	10
Units.....	11
High Alarm Set Point	11
Low Alarm Set Point	11
Alarm Silencing Menu	11
Alarm Test Menu	12
Calibration Offset.....	12
Hold-Off Time	12
Control Mode	13
Loading	13
Temperature Disregression	13
Alarms	14
Over Temp Alarm.....	14
Under Temp Alarm.....	14
Door Ajar Alarm (swinging door models only).....	14
Power Loss Alarm	15
Alarm Silencing	15
Alarm Test	15
Remote Alarm Contacts	16
Open Sensor	17
Battery Back Up	17
Calibration Using Temperature Offset	17
Convenience Outlet	18
Maintenance	19
Cabinet Cleaning	19
Cleaning the Condenser.....	19
Condensate Evaporator Pan	19
Troubleshooting	20
Replacement Parts	21
Wiring Diagrams	22
Warranty	28

Introduction



Warning

These refrigerators are not approved for the storage of blood or blood products!



DANGER: RISK OF CHILD ENTRAPMENT. BEFORE YOU THROW AWAY YOUR OLD REFRIGERATOR OR FREEZER:

- TAKE OFF DOORS
- LEAVE THE SHELVES IN THE PLACE SO THAT CHILDREN MAY NOT EASILY CLIMB INSIDE.



Warning

As a routine laboratory precaution, always wear safety glasses when working with this apparatus.



Warning

This product is NOT approved for storage of flammable or explosive materials. Also, it is NOT approved for use in hazardous locations containing explosive atmospheres.



Warning

High voltage is supplied to PC board when key switch is OFF.

This manual is intended to provide installation and operating instructions for Isotemp Plus Alarm version refrigerators. The alarm version control system provides both audible and visual indication of alarm conditions in an easy to read display format. In addition, there are two sets of contacts available for remote alarm switching.

Temperature and alarm set points are directly entered in °C or °F. More about setting up the controller and all the options are discussed in the Operating section of this manual.

Safety Precautions

Your satisfaction and safety are important to Fisher Scientific, a complete understanding of this unit is necessary to attain these objectives.

As the ultimate user of this apparatus, it is your responsibility to understand its proper function and operational characteristics. This instruction manual should be thoroughly read and all operators given adequate training before attempting to place this unit in service. Awareness of the stated cautions and warnings, and compliance with recommended operating parameters – together with maintenance requirements – are important for safe and satisfactory operation. The unit should be used for its intended application; alterations or modifications will void the Warranty.

This product is not intended, nor can it be used, as a sterile or patient connected device. It is not approved for the storage of blood or blood components. In addition, this apparatus is not designed for use in Class I, II or III locations as defined by the National Electrical Code, unless otherwise noted.

Performance Characteristics

Temperature Range: 1.0° to 12.0°C (33.8° to 53.6°F)

Temperature Stability: ±3°C

Installation

Unpacking

Save all packing material if apparatus is received damaged. This merchandise was carefully packed and thoroughly inspected before leaving our factory.

Responsibility for its safe delivery was assumed by the carrier upon acceptance of the shipment; therefore, claims for loss or damage sustained in transit must be made upon the carrier by the recipient as follows:

Visible Loss or Damage

Note any external evidence of loss or damage on the freight bill, or express receipt, and have it signed by the carrier's agent. Failure to adequately describe such external evidence of loss or damage may result in the carrier's refusing to honor your damage claim. The carrier will supply the form required to file such a claim.

Concealed Loss or Damage

Concealed loss or damage refers to loss or damage, which does not become apparent until the merchandise has been unpacked and inspected. Should either occur, make a written request to the carrier's agent within 15 days of the delivery date; then file a claim with the carrier since the damage is the carrier's responsibility.

If you follow the above instructions carefully, we will guarantee our full support of your claim to be compensated for loss from concealed damage.

**DO NOT – FOR ANY REASON – RETURN THIS UNIT
WITHOUT FIRST OBTAINING AUTHORIZATION**

Packing List

The following items are packed in the envelope located inside the refrigerator chamber. If any of the following items are not present, report the missing item to your local Fisher representative.

1. Warranty Card
2. Instruction Manual
3. Door Lock Key
4. Power Switch Key
5. Chart Recorder Instructions (Catalog No.'s with "R")

Selecting a Location

Choose a location for the refrigerator that will provide **at least three inches of clearance** between the cabinet and any adjacent vertical surface at the sides and rear. Appropriate electrical power must be available. Locate the refrigerator within 6 feet of the power outlet so that no extension cord is required.

Installing and Leveling Casters

Isotemp Plus refrigerators and freezers come with four casters, which thread into the base of the unit, one in each corner. Use the wrench provided to thread the casters completely into the base of the unit. Back the casters in or out until the unit is level and resting on all four casters. See Figure 1.

**Note**

The refrigerator must be level in order to provide adequate condensation drainage as well as proper door alignment and operation. The refrigerator should be in its final operating location and set so that it is firmly positioned on the floor.

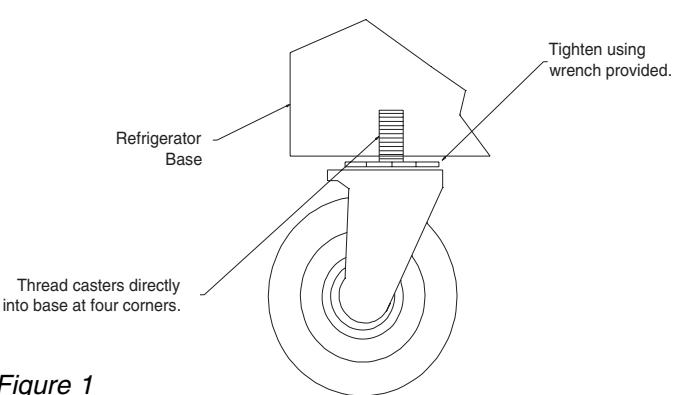


Figure 1

UNPACKING AND INSTALLATION



Caution

If the unit was tilted in excess of 30 degrees, level it and do not apply electrical power for a minimum of 12 hours.



Warning

Be sure that the power supply is the same voltage that is specified on the refrigerator's data plate.



Warning

For personal safety this unit must be properly grounded. DO NOT under any circumstance cut or remove the ground prong from the power cord. DO NOT use a plug adapter that eliminates the ground prong.

115VAC Units

NEMA 5-15P, 5-15R
13-986-120A/RA
13-986-122A/RA
13-986-124A/RA
13-986-126A/RA
13-986-128A/RA
13-986-128SA/SRA
13-986-130A/RA
13-986-132A/RA

NEMA 5-20P, 5-20R
13-986-134A/RA
13-986-136A/RA
13-986-138A/RA
13-986-140A/RA
13-986-136SA/SRA

230VAC Units

NEMA L14-20P, L14-20R
13-986-121A/RA
13-986-123A/RA
13-986-125A/RA
13-986-127A/RA
13-986-139A/RA

13-986-131A/RA
13-986-133A/RA
13-986-135A/RA
13-986-137SA/SRA
13-986-141A/RA



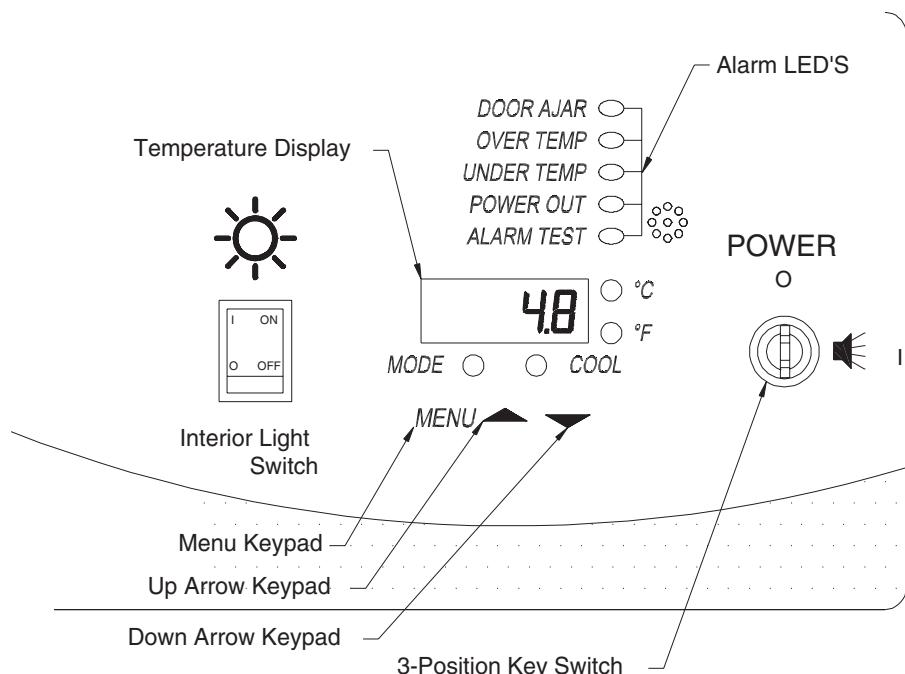
Caution

Do not use an extension cord. Use of an ungrounded cord or an overloaded circuit VOIDS the compressor warranty.

Controls

Panel Controls

Before operation, it is necessary to become familiar with the refrigerator controls. All the refrigerator controls are located on the header panel. A layout of the controller and LED's are given below.



Warning

High voltage is supplied to PC board when key switch is OFF.

Displays and Features

On the right side of the header panel is the key switch. The key switch has three positions, OFF (O), Set-up (I) and Alarm Enable (▲). With the refrigerator plugged in and the key switch in the OFF position, the displays and refrigeration system are off. The control board remains energized in order to supply a charge to the battery.

Key Switch

Turning the key switch to the middle position energizes the circulation fans and the controller display. When the controller is in the Set Up mode, the menu items can be set to the desired values. See Controller Set Up for a description of each. Once all menu items are set, the refrigerator can then be switched to the Alarm Enable

CONTROLS



Note

The key can be pulled in the Off or the Alarm Enable position.

position. While in the Alarm Enable position, the menu items can be viewed but not changed. The key can then be pulled to secure the values entered.

Temperature Display

The temperature display indicates chamber temperature while in the control mode of operation and menu titles and values while entering menu items (see Menus).

Units LED'S

The temperature display indicates chamber temperature in both °C and °F. The current units are indicated by the corresponding LED located to the right of the temperature display.

Mode LED

The Mode LED is lit when the Menu keypad has been pressed. It indicates to the user that the controller is in the menu selection mode.

Cool LED

The Cool LED is lit when the controller calls for cooling. When this LED comes on the compressor should energize.

Alarm LED'S

Door Ajar - This LED is illuminated when a refrigerator door(s) is opened. This feature is not used in the sliding glass door models, since the circulating fans run continuously and the doors are spring loaded.

Over Temp – Indicates the current chamber temperature exceeds the high alarm temperature set point.

Under Temp – Indicates the current chamber temperature is lower than the low alarm temperature set point.

Power Out – Indicates AC power to the refrigerator is disrupted.

Alarm Test – Indicates the controller is running an alarm test.

Keypads

Menu Keypad – Pressing the Menu keypad enters the controller into the menu selection mode, designated by the illumination of the MODE LED. Repeatedly pressing the Menu keypad cycles through the list of menu items (see Menus). After the final menu, the controller returns to the control mode and the MODE LED turns off.

Up Arrow Keypad – Increases or changes menu values.

Down Arrow Keypad - Decreases or changes menu values.

Interior Light Switch

The interior lamp(s) is controlled by a rocker switch located on the header panel. This light may be operated any time the cabinet power is turned on at the keyed power switch. If the interior lamp fails, replace with the same size and wattage lamp. DO NOT USE REDUCED WATTAGE LAMPS. The reduced wattage lamps generally fail to light below 60°F (15°C).

Operation



Note

The Pout display is a normal occurrence when power is first applied to the controller. Press any key to clear it.

The refrigerator is started by inserting the power switch key, which is found in the envelope inside the unit, into the key switch. Turn the key to the middle (set up) position. The internal fans will start and the display will flash between Pout and the current chamber temperature. The Pout (power out) display indicates to the user that the controller has sensed a low voltage condition. This condition is normal during any start-up. See the Alarms section for details on the Pout message. Press any keypad to clear the Pout.

Controller Set Up (Menus)

During normal operation (control mode), the refrigerator temperature is displayed on the 4-character display. The compressor cycles on and off to control the chamber temperature between 1.0° and 12.0°C (33.8 to 53.6°F), selectable by the user. The temperature set point, high and low alarm set-points, units and display offset can be changed while the key switch is in the setup (middle) position. These values can be viewed but not changed with the key switch is in the alarm position. The menu options are explained in the following sections.

There are 8 menus to select from, each can be accessed by pressing the Menu keypad. Pressing the Menu keypad once brings up the first menu, which is the temperature set point. The displays initially flash SP followed by the last set temperature value. Pressing the Menu keypad again, accepts the value and displays the next menu, units select (CF). Each menu value is entered this way. The menus can be changed only while the key switch is in the set up (I) position. They can be viewed but not changed while the key switch is in the alarm enable (◀) position. The individual menus are described below.

Temperature Setpoint

- Temperature Set Point (SP) – Adjustable from 1.0 to 12.0°C or 33.8 to 53.6°F. Use the Up/Down Arrow keypads to change the temperature and the Menu keypad to enter the value and switch to the next menu.

Units

- Units select (C,F) – Selects the desired temperature display units in °C or °F. Use the Up Arrow keypad to change the units from C to F. Use the Down Arrow keypad to switch from F to C. Press the Menu keypad to enter the units value and continue. The appropriate units LED will indicate the selected units.

High Alarm Set Point

- High Alarm Set Point (AH) – Use the Up/Down Arrow keypad to adjust the value that triggers a high alarm condition if the chamber temperature rises above it. The high alarm temperature range is from -33.0°C (-27.4°F) to 37.0°C (98.6°F).

NOTE: The High and Low alarm setpoints can be adjusted to be within 3.5°C (6.3°F) of one another and no closer.

Low Alarm Set Point

- Low Alarm Set Point (AL) – Use the Up/Down Arrow keypad to adjust the value that triggers a low alarm condition. The low alarm condition occurs when the actual chamber temperature drops below the Low Alarm Set Point. The low alarm range is from -36.0°C (-32.8°F) to 34.0°C (93.2°F).

Alarm Silencing Menu

- Alarm Silencing (AS) – During an alarm condition, while the key switch is in the Alarm Enable Mode, this menu allows the user to silence the audible alarm for 5 minutes. After the AS stops flashing, the display will show n, press the Up Arrow keypad to change the display to y. Press the Menu Keypad and the audible alarm will turn off for 5 minutes.

CONTROLS

Alarm Test Menu

- Alarm Test (At) – Use the Up Arrow keypad to change the n to H. By subsequently pressing the Menu Keypad the controller will perform a High Alarm Test (see Alarm Test in Alarms section). Use the Down Arrow Keypad to change the display to L. Pressing the Menu Keypad will activate the Low Alarm Test. Pressing the Menu Keypad while the display shows n returns the unit to control mode (MODE LED off).

NOTE: While in any of the above menu options, the controller will wait 15 seconds for a new value to be entered. If there is no keypad activity within 15-seconds, the refrigerator will automatically shift to the Control Mode and the Mode LED will turn off.

Calibration Offset

The two final menu items, calibration offset and hold-off time (oS and Ho), are not accessible from the main menu. In order to view and/or change them, the Menu keypad must be pressed and held for 5 seconds. The calibration offset (oS) menu will then be displayed briefly followed by the most recent offset value.

- Display Offset (oS) – Use the Up/Down Arrow keypads to adjust the offset value. Adjustable to +/- 5.0°C (+/- 9.0°F). This value is added to the temperature display to correct for differences between actual and displayed chamber temperature. See Calibration... for details. Press the Menu keypad to enter the new value or to continue to the next menu.

Hold-Off Time

- Hold-Off (Ho) – Use the Up/Down Arrow keypads to adjust the Hold-Off time in minutes. Adjustable from 3 to 10. This value represents the delay time between compressor activations. Increasing the Hold-Off time can help prevent evaporator freeze-up by increasing the defrost cycle time. Press the Menu keypad to enter the new value and return to temperature control mode.

Control Mode

The Control Mode of operation is indicated when the Mode LED is OFF. With the desired temperature set point and alarm values entered, the refrigerator will then begin to control to the set temperature. There is a time delay between compressor activations. This delay begins the moment the controller is energized or the moment the compressor turns off during normal operation. The delay is added to provide time for the system pressure to drop before the compressor tries to start again.

Loading

Follow the guidelines below in order to provide optimal refrigerator performance:

- When loading with critical samples, verify all alarms and features are operating properly.
- Space samples uniformly in the chamber.
- Keep samples away from the back wall that may obstruct airflow.
- It is best to load the refrigerator from bottom to top. Keeping the area next to the air intake and exhaust clear will help assure proper airflow.

Temperature Digression

By pressing the Increase keypad while the refrigerator is in the Control Mode, the display will show the warmest temperature recorded since the unit was energized or the digression values were reset. Pressing the Decrease keypad will display the coldest temperature recorded since the unit was energized or the digression values were reset.

To reset the temperature digression values, simply press the Increase and Decrease keypads at the same time. The display will show Clr, indicating the controller will begin recording new warm and cold digressions once more.

Alarms

Alarm Version Refrigerators are equipped with audible and visual alarms. They also feature remote alarm contacts. The key switch must be in the Alarm Enable position for the audible alarm and remote alarm contacts to activate. Alarm messages are shown on both the temperature display and in the row of Alarm LED's located above the temperature display. The following conditions will trigger an alarm:

- Chamber temperature is greater than the high alarm set point.
- Chamber temperature is less than the low alarm set point.
- Door has been left open for 3 minutes or more.
- AC power has been interrupted to the refrigerator.
- Open sensor

NOTE: The Door Ajar Alarm is not available on the sliding glass door models.

Over Temp Alarm

If the refrigerator chamber temperature rises above the High Alarm Set Point, the display will alternately flash the refrigerator chamber temperature and "High". The Over Temp LED will be lit, the audible alarm will sound and the remote contacts will change state.

Under Temp Alarm

If the refrigerator chamber temperature falls below the Low Alarm Set Point, the display will alternately flash the refrigerator chamber temperature and "Lo". The Under Temp LED will be lit, the audible alarm will sound and the remote contacts will change state.

Door Ajar Alarm (swinging door models only)

When a swinging door is opened, the Door Ajar LED lights and the display alternately flashes between the

refrigerator temperature and "door". If a door is opened for 3 minutes, the audible alarm will sound and the remote alarm contacts change state.

Power Loss Alarm

When AC power is lost, the display will alternately flash between the current refrigerator temperature and "Pout", the Power Out LED will light, the audible alarm will sound and the remote contacts will change state. When power is restored, the display will continue flashing the "Pout" message until it is cleared by the user by pressing any of the keypads.

Alarm Silencing

The audible alarm can be silenced by deactivating it at the Alarm Silence Menu (see section earlier in this manual). The audible alarm will silence for 5 minutes. It will sound again at the end of the 5 minutes if the alarm condition persists. Alarm silencing does not affect the visual alarms.

Alarm Test

The Alarm Test feature is used to verify proper alarm operation by simulating an increase or decrease in chamber temperature. The alarm test function will work for any value of high or low alarm set points. With the key switch in the Set Up Position the remote and audible alarms are not activated during an alarm test, as is the case in an actual alarm condition.

NOTE: The key switch should be in the Alarm Enable Position in order to test all alarm features.

The procedure for testing the alarms is simple. Press the Menu Keypad repeatedly until the display shows At (alarm test). Momentarily, the display will change to n. Press the Up Arrow Keypad to change the display from n to H in order to perform a high alarm test. With the display showing H, press the Menu Keypad again and the temperature display will begin to rise until it reaches the High Alarm Set Point. Once the set point is reached, both audible and visual alarms should be activated (with the key switch in alarm enable position) and the remote alarm

CONTROLS

contacts should change state. To perform a low alarm test, simply press the Down Arrow Keypad when the display shows n. The display will then show L. Press the Menu Keypad again in order to activate the low alarm test.

Remote Alarm Contacts

In addition to the audible and visual alarms, there are two sets of remote contacts on the side of the header panel, which change state during an alarm. The upper set of contacts provides both normally open and normally closed switching. The lower set of contacts is used for the Fisher Deluxe Remote Alarm System (see Refrigerator Accessories).

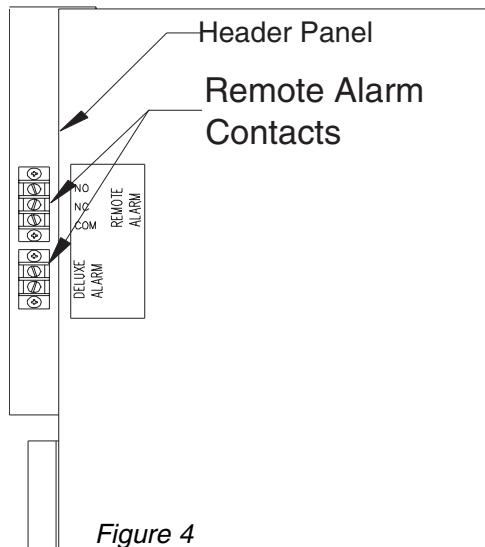


Figure 4

All contacts are capable of switching up to 24 volts AC/DC @ 2 amps. The common (C) and normally closed (NC) contacts open during an alarm. The common (C) and normally open (NO) contacts close during an alarm. The contacts for the Deluxe Remote Alarm are normally closed and open during an alarm. Figure 4 shows the location of the remote and deluxe alarm contacts.

Remote alarm contacts are active only with the key switch in the Alarm Enable position.

Open Sensor

When the controller is unable to read from the sensor, the display will show SnSr, indicating that a sensor error exists. The display will alternately flash SnSr and HigH. The audible alarm will sound, the remote alarm contacts will change state and the over temp LED will illuminate.

Battery Back Up

The Isotemp Plus Alarm Version Refrigerators are supplied with a rechargeable battery back up. The battery will continue to supply power to the controller and alarm features in the event of a power outage. A fully charged battery can continue powering the controller for up to 8 hours during a power failure.

The battery is recharged by the temperature controller during normal operation. The battery should be checked every six months and replaced when necessary (see Replacement Parts). Check the battery by unplugging the refrigerator and monitoring the controller operation. If the controller does not properly activate the alarm features for 2 hours, replace the battery. Under normal operation, the battery back up should last up to four years or 200 charge/discharge cycles.

Calibration Using Temperature Offset

In the event the refrigerator needs to be calibrated, a simple routine is available to adjust the display and control point to a referenced standard. To set a temperature offset while in the Control Mode (mode LED off), simply press and hold the Menu key for 5 seconds. The display will flash oS followed by the last temperature offset value. The factory setting is 0.0. To change the offset value, press the Increase or Decrease keypad to the value which when algebraically added to the last displayed temperature, will agree with the reference temperature. Next, press the Menu key until the Mode LED turns off.

For Example:

The display indicates 4°C but a reference thermometer in the refrigerator chamber indicates 6°C. The operator presses and holds the Menu KEY until the display shows

CONTROLS



Caution

INCREASING the offset value DECREASES the temperature at which the system controls. If the offset is inadvertently set too high, the chamber temperature can fall below zero and lead to evaporator freeze-up.



Warning

The convenience outlet is LIVE at any time the unit is connected to an electrical power source, regardless of the position of the key switch.

(oS) offset, then the value. Change the value from 0.0 to 2.0. The Menu key is then pressed repeatedly until the Mode LED turns off. Now the display indicates a chamber temperature of 6.0°C (approx.) and the controller begins to cool to the desired temperature of 4.0°C. Allow an additional 30 to 40 minutes for the refrigerator to again stabilize. If the display is still inaccurate, repeat the calibration procedure.

Convenience Outlet

(Chromatography Refrigerators Only)

Chromatography Refrigerators are equipped with a convenience outlet located in the center of the back wall. Models 140A, 140RA, 141A AND 141RA (three doors) have two convenience outlets. The power to the outlets is protected by a circuit breaker. The breaker is located just below and to the right of the outlet. Pressing the breaker button resets the breaker after a fault.

The total load to one or both convenience outlets should not exceed 5 amps.

Maintenance

**Caution**

When servicing the unit, disconnect plug from the electrical power source

**Caution**

Do not use any type of abrasive such as steel wool or fluids, such as gasoline, Naphtha or paint thinner, that could be harmful to plastic materials, door gasket and/or painted surfaces for cleaning.

Cabinet Cleaning

The cabinet interior should be cleaned frequently. Any spilled liquid should be wiped off immediately. Stains resulting from some spills can be permanent if not quickly removed. The cabinet exterior should be cleaned occasionally. A mild detergent and lukewarm water or a solution of bicarbonate of soda (1 tablespoon per gallon of water) is recommended for cleaning the interior and exterior of the cabinet. All surfaces should be rinsed and thoroughly dried.]

Cleaning the Condenser

The condenser is located in front of the compressor fan, which is on the top, left side of the swinging door model refrigerators, and behind the base grill of the sliding door model refrigerators. The condenser should be cleaned before becoming clogged with dirt/dust. Construction or other dirt causing environments may significantly increase the required frequency. **The condenser should be cleaned at least once a year.**

Condensate Evaporator Pan

The condensate evaporator pan is located behind the compressor fan, which is on the top, left side of the swinging door model refrigerators, and behind the base grill of the sliding door model refrigerators. **This pan should be cleaned at least once a year to prevent foul odors and operate efficiently.** Vacuum clean if dry, or sponge clean with soapy water.

Troubleshooting

This table is intended to assist in resolving user-correctable Refrigerator problems by relating symptoms to their likely causes. If service beyond the scope of this table is required, contact Fisher Scientific Service Division at 1-800-438-4851.

Symptom	Probable Cause	Action
Does Not Run	Unit Unplugged	Plug in Unit
	Blown fuse or tripped circuit breaker	Check fuse or circuit breaker at breaker box
Runs Continuously	Frost buildup on refrigeration coils.	Turn unit off and allow it to defrost.
	Refrigerator over-loaded.	Remove contents from top shelf of unit.
Clicking Sound	The compressor is equipped with a thermal protector. This device shuts off the compressor when it becomes too hot. A clicking sound occurring about every 30 seconds indicates this protector is working	Disconnect power and allow unit to sit for about an hour, then retry. If condition persists, call for service.
Insufficient Cooling	Set temp is too high	Reduce temperature setting, verify Cool LED turns on in 3 minutes.
	Offset (oS) value is too low.	Increase offset value.
	Condenser coil dirty	Clean condenser coil with a vacuum cleaner
	Unit frosted	Defrost unit
	Circulating fans not on.	Verify doors are completely closed.

Replacement Parts

Replacements for Laboratory Refrigerator parts serviceable by the user may be ordered, by part number, from Fisher Scientific Co. at 1 800 438-4851.

Epoxy Coated Shelves

Full depth shelf for models: 13-986-120/121/126/127 13-986-130/131/132/133/134/135/138/139/140/141	13-986-210E
Half depth shelf for models: 13-986-120/121/126/127 13-986-130/131/132/133/134/135/138/139/140/141	13-986-210K
Full depth shelf for models: 13-986-122/123/124/125	13-986-210M
Half depth shelf for models: 13-986-122/123/124/124	13-986-210N
Full depth shelf for models: 13-986-128S/129S/136S/137S	13-986-210W
Half depth shelf for models: 13-986-128S/129S/136S/137S	13-986-210Y
Solid State Relay (K2)	SPN88616
Magnetic Relay (K1) SPST	SPN102259
Magnetic Relay (K1) DPDT	SPN102260
Alarm Version Temperature Controller	SPN104162
Condensate Pan Wick, 23.1 & 30 Cu. Ft.	SPN107670
Condensate Pan Wick, 46.6 & 74 Cu. Ft.	SPN107671
Condenser Pan Wick (-128S, -136S)	106961
Shelf Support	SPN103265
Chart Recorder Manual	SPN104114
Rechargeable Battery	SPN104711
Temperature Sensor	SPN104163
Drawer and Slide Assembly	SPN13999DRAW

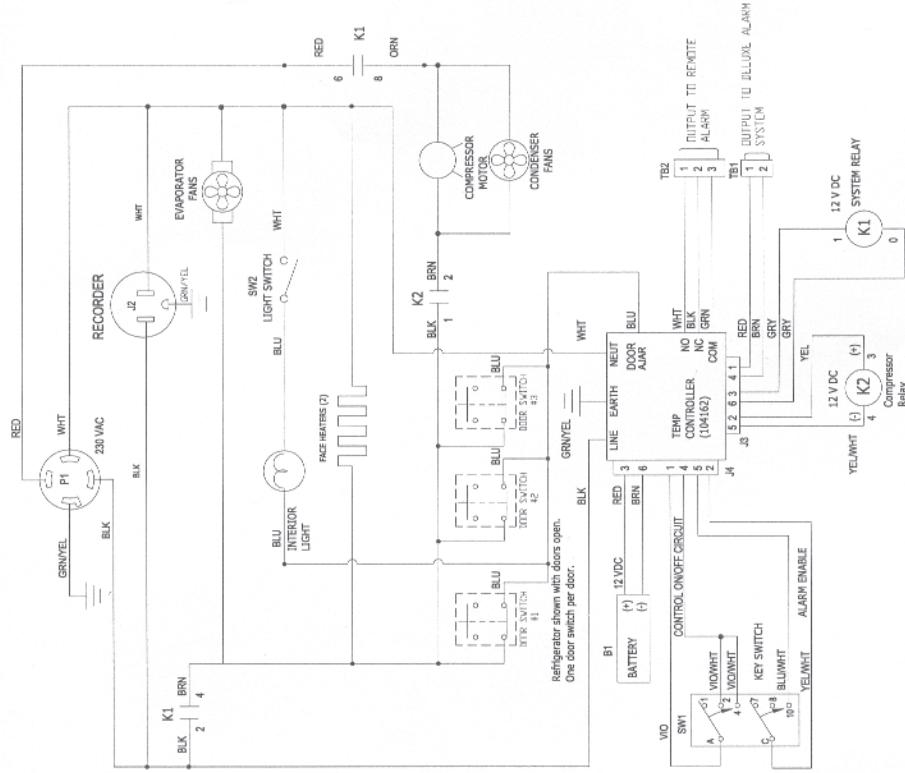
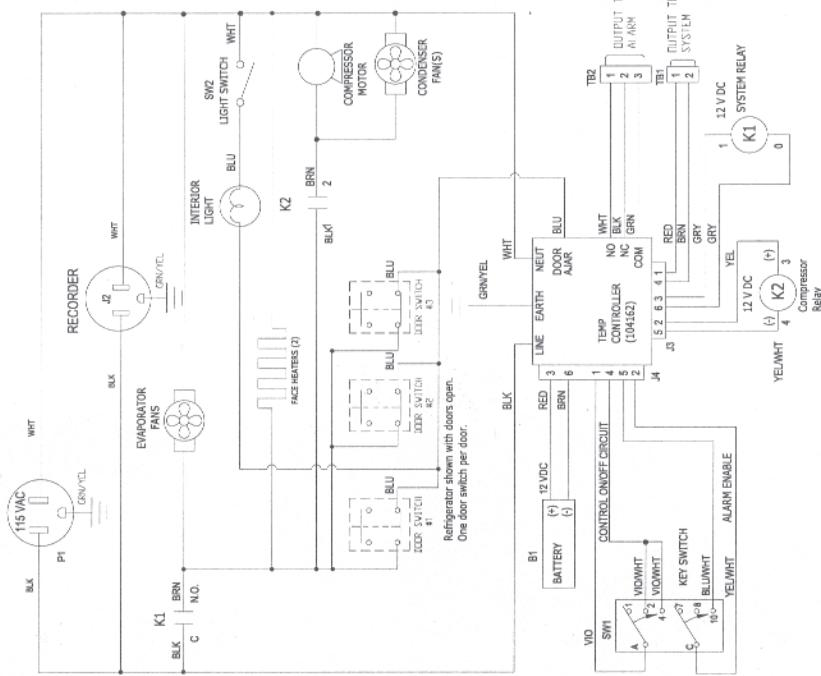
Accessories

Deluxe Electronic Remote Alarm System	13-986-100
Standard Remote Alarm System	13-986-32A
Freestanding Chart Recorder	13-986-202
Zinc Coated Shelves	
Full depth shelf for models: 13-986-120/121/126/127 13-986-130/131/132/133/134/135/138/139/140/141	13-986-211E
Half depth shelf for models: 13-986-120/121/126/127 13-986-130/131/132/133/134/135/138/139/140/141	13-986-211K
Full depth shelf for models: 13-986-122/123/124/125	13-986-211M
Half depth shelf for models: 13-986-122/123/124/124	13-986-211N
Full depth shelf for models: 13-986-128S/129S/136S/137S	13-986-211P
Half depth shelf for models: 13-986-128S/129S/136S/137S	13-986-211Q

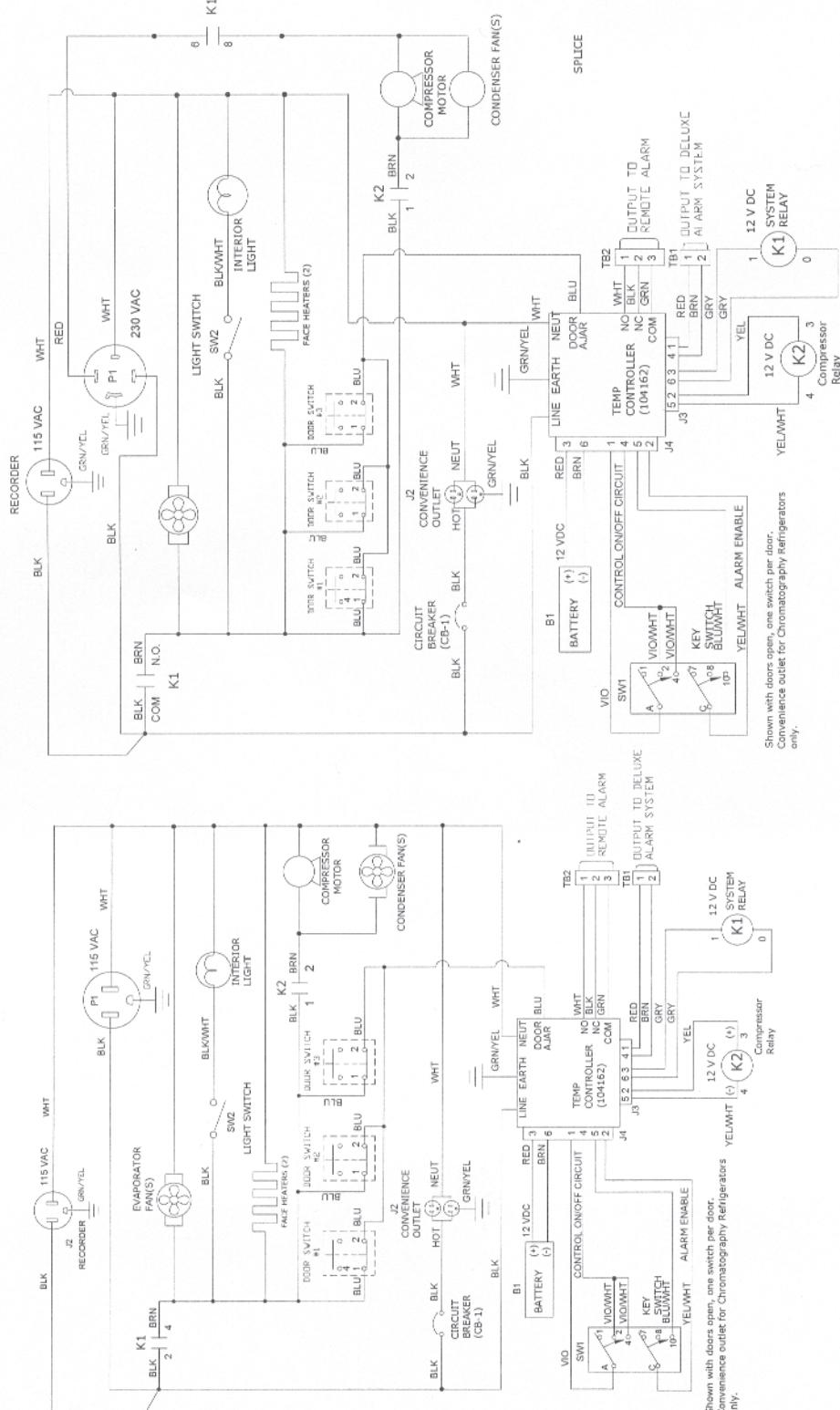
Wiring Diagrams

Swinging Door Models:
 Catalog No.'s:
 13-986-120A/RA, 122A/RA, 126A/RA, 130A/RA

13-986-121A/RA, 123A/RA, 127A/RA, 131A/RA



Sliding Glass Door Models:
 Catalog No.'s:
 13-986-128SA/SRA, 13-986-136SA/SRA, 13-986-129SA/SRA, 13-986-137SA/SRA

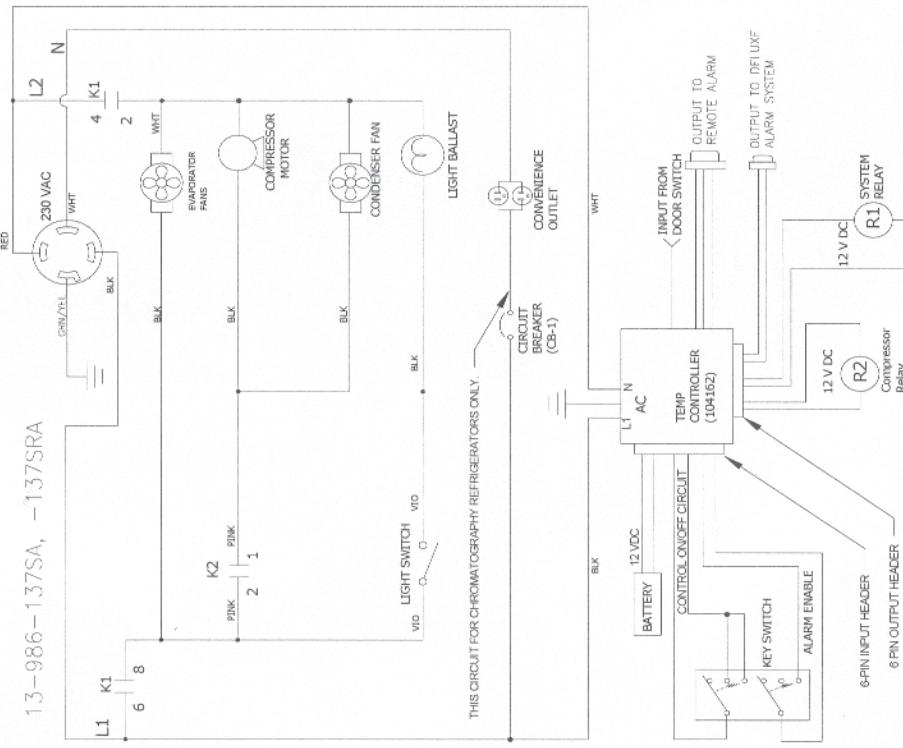
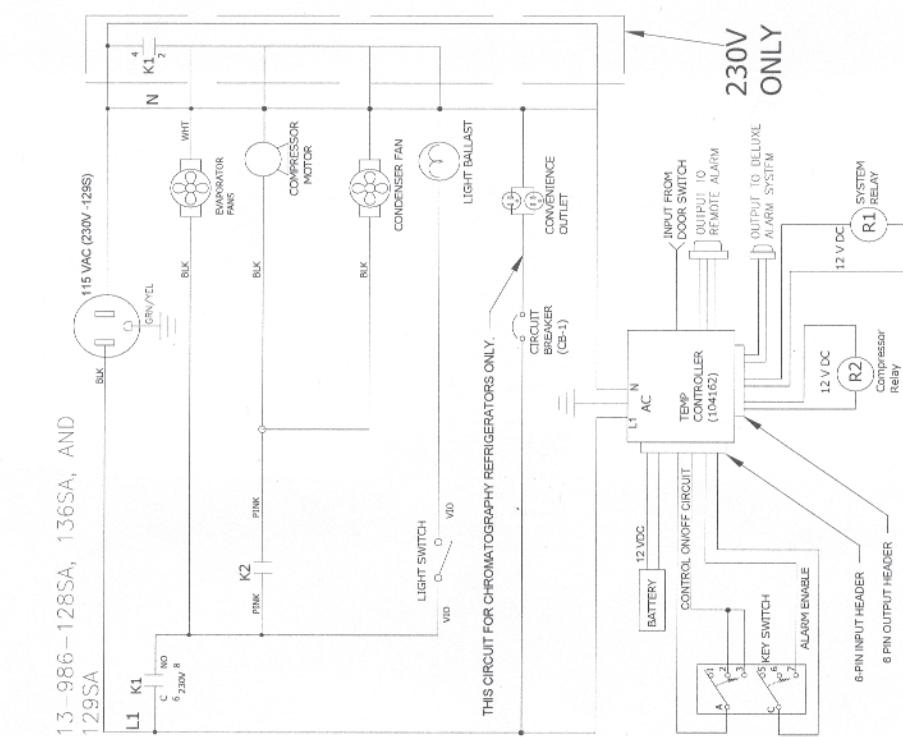


Shown with doors open, one switch per door.
 Convenience outlet for Chromatography Refrigerators
 only.

WIRING DIAGRAMS

Swinging Door Models:
Catalog No.'s:
13-986-124A/RA, 132A/RA, 134A/RA, 13-986-140A/RA

13-986-125A/RA, 133A/RA, 135A/RA, 139A/RA, 13-986-141A/RA



Warranty

Laboratory instruments and equipment manufactured by Fisher Scientific Company L.L.C. – Laboratory Equipment Division (hereinafter called “the Company”) are warranted only as stated below.

Subject to the exceptions and upon the conditions specified below, the Company agrees, at its election, to correct by repair, by replacement, or by credit to the purchaser, any defect of materials or workmanship which develops within one year (13 months for refrigerator and freezer products) from the date of purchase by the original purchaser by the Company or by an authorized dealer of the Company provided that investigation or factory inspection by the Company discloses that such defect developed under normal and proper use

The exceptions and conditions mentioned above are the following:

- a. The Company makes no warranty concerning components or accessories not manufactured by it, such as tubes, batteries, etc. However, in the event of the failure of any component or accessory not manufactured by the Company, the Company will give reasonable assistance to the purchaser in obtaining from the respective manufacturer whatever adjustment is reasonable in the light of the manufacturer's own warranty.
- b. The Company shall be released from all obligations under its warranty in the event repairs or modifications are made by persons other than its own service personnel or authorized dealer personnel unless such repairs by others are made with the written consent of the Company.
- c. **THE COMPANY MAKES NO WARRANTY OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR ANY OTHER WARRANTY, EXPRESS OR IMPLIED, EITHER IN FACT OF BY OPERATION OF LAW,...STATUTORY OR OTHERWISE.**
- d. The above warranty and the above obligations to repair, replace, or credit are complete and exclusive and the Company expressly disclaims liability for lost profits or for special, indirect, incidental, consequential, or exemplary damages of any nature whether attributable to contract, warranty, negligence, strict liability, or otherwise even if the Company has been advised of the possibility of such damages.
- e. Representations and warranties made by any person, including dealers and representatives of the Company, which are inconsistent or in conflict with the foregoing warranty shall not be binding upon the Company unless reduced to writing and signed by an officer of the Company.



Fisher Scientific