



# Series 8000WJ

## Water Jacketed CO<sub>2</sub> Incubator

Operating and Maintenance Manual 7013422 Rev. 9

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

**Thermo**  
SCIENTIFIC

Model	CO2 Sensor*	Voltage**	O <sub>2</sub>
3422	IR	115	no
3423	IR	230	no
3424	T/C	115	yes
3425	T/C	230	yes
3426	IR	115	yes
3427	IR	230	yes
3428	T/C	115	no
3429	T/C	230	no

\*T/C is a thermal conductivity sensor.

IR is an infra-red sensor.

\*\*All units are 50/60 Hz.

#### MANUAL NUMBER 7013422

9	ECNM-002047	07/15/22	Updated the Shelf count in page 7-2.	TFI
8	ECNM-001240	10/27/21	Updated Door alignment step 21 in Section Reversing the Door Swing in page 5-12.	TFI
7	ECNM-001514	06/04/21	Added EMC statements in preface page iv.	TFI
6	ECNM-000906	11/27/2020	Replaced part number 400202 with 400218 in page 8-1, updated Assembly drawing in page 8-4 and Electrical drawing in pages 9-1 to 9-6.	TFI
5	ECNM-001141	09/29/2020	Added tubing inspection details to preventive maintenance section 5-15.	TFI
4	ECNM-000651	09/11/2020	Removed Fluorinated Green Gas pare pg 7-4.	TFI
3	41345	5/02/17	Added F-gas statement	bpg
2	40087/IN-4743	7/27/16	Changed label in remote alarm artwork to note - pg 6-2	ccs
1	40639	5/20/16	Risk assessment information	ccs
0	27988/IN-4478	4/7/14	Release 7 Original	ccs
REV	ECR/ECN	DATE	DESCRIPTION	BY



**Important** Read this operating manual. Failure to read, understand and follow the instructions in this manual may result in damage to the unit, injury to operating personnel, and poor equipment performance. s

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

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

#### Intended Use:

The Thermo Scientific Water-jacketed CO<sub>2</sub> incubators listed in this manual are intended to provide an environment with controlled temperature and Carbon Dioxide (CO<sub>2</sub>) as well as elevated humidity (RH) for the growth of microbiological samples or mammalian cells for research (FDA Class 1) and IVF and other regulated cell therapy applications (FDA Class 2). Certain models also control Oxygen (O<sub>2</sub>) at suppressed levels via the injection of Nitrogen (N<sub>2</sub>).

For FDA Class 1 applications, the intended end user is a laboratory / medical professional trained in basic safety and cell culture practices. For FDA Class 2 applications, the intended user is a trained Laboratory Technician whose qualifications include a bachelor's or master's degree in chemical, physical, biological, medical technology, clinical or reproductive laboratory science from an accredited institution.

These incubators are designed to be used indoors only and operate 24 hours per day - 7 days per week, but should be shut down regularly to be disinfected to prevent contamination, maintained per the instructions for use, and calibrated to retain optimum system performance.

#### Non- intended use:

These incubators are not intended for use by the general public.

These units are not intended for use within the patient environment. These units are not to be used outdoors.

These incubators are not designed to provide O<sub>2</sub> levels greater than 20% and as such should never be connected to O<sub>2</sub> as it is highly flammable in elevated concentrations. These units are not designed to be used with Liquid CO<sub>2</sub> or Liquid N<sub>2</sub>. These units should not be used with any flammable or toxic materials.

© 2022 Thermo Fisher Scientific. All rights reserved.



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



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



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



Asphyxiation Hazard Warning. High concentrations of CO<sub>2</sub> and N<sub>2</sub> can displace oxygen and cause asphyxiation!



Lifting Hazard Warning. The incubator weighs more than 200 lbs (91 kgs). Take adequate safety measures when moving this device.



4-Person Lift. The incubator weighs more than 200 lbs (91 kgs). It is recommended that four persons lift this device together to move it.



WEEE Compliance: Thermo Fisher Scientific has contracted with companies for recycling/disposal in each EU Member State. For further information, send an email to [weee.recycle@thermofisher.com](mailto:weee.recycle@thermofisher.com).

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

## **EMC**

This equipment has been tested and found to comply with the limits for a Class A digital device. Class A covers devices for usage in all establishments other than domestic and that are not directly connected to a low voltage power supply network, which supplies domestic environment.

### **Consideration for Korea.**

EMC Registration is done on this equipment for business use only. It may cause interference when the product would be used in home.

사용자 안내문 이 기기는 업무용 환경에서 사용할 목적으로 적합성평가 를 받은 기기로서 가정용 환경에서 사용하는 경우 전파간섭의 우려가 있습니다 .

### **Consideration for EU**

This product is intended for use at locations where product is being supplied by separate power network, in most cases supplied from a high or medium voltage transformer, dedicated only for manufacturing plants or similar plants.

### **FCC Statement**

This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at his own expense.

### **Canada Statement:**

This ISM device complies with Canadian ICES-001. Cet appareil ISM est conforme à la norme NMB-001 du Canada.

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

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

1-740-373-4763

1-800-438-4851

1-877-213-8051

<http://www.thermofisher.com>

[service.led.marietta@thermofisher.com](mailto:service.led.marietta@thermofisher.com)

[www.unitylabservices.com](http://www.unitylabservices.com)

Direct

Toll Free, U.S. and Canada

FAX

Internet Worldwide Web Home Page

Tech Support Email Address

Certified Service Web Page

Our **Sales Support** staff can provide information on pricing and give you quotations. We can take your order and provide delivery information on major equipment items or make arrangements to have your local sales representative contact you. Our products are listed on the Internet and we can be contacted through our Internet home page.

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

Whatever Thermo Scientific products you need or use, we will be happy to discuss your applications. If you are experiencing technical problems, working together, we will help you locate the problem and, chances are, correct it yourself...over the telephone without a service call.

When more extensive service is necessary, we will assist you with direct factory trained technicians or a qualified service organization for on-the-spot repair. If your service need is covered by the warranty, we will arrange for the unit to be repaired at our expense and to your satisfaction.

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

Thermo Fisher Scientific (Asheville) LLC  
401 Millcreek Road, Box 649  
Marietta, OH 45750

International customers, please contact your local Thermo Scientific distributor.

## Warranty Notes

### **Information You Should Know Before Requesting Warranty Service**

- **Locate the model and serial numbers.** A serial tag is located on the unit itself.
- For equipment service or maintenance, or with technical or special application inquiries, contact your local distributor.

### **Repairs NOT Covered Under Warranty**

- **Calibration of control parameters.** Nominal calibrations are performed at the factory; typically  $\pm 1^{\circ}\text{C}$  for temperature,  $\pm 1\%$  for gases, and  $\pm 5\%$  for humidity. Our service personnel can provide precise calibrations as a billable service at your location. Calibration after a warranty repair is covered under the warranty.
- **Damage resulting from use of improper quality water, chemicals or cleaning agents detrimental to equipment materials.**
- **Service calls for improper installation or operating instructions.** Corrections to any of the following are billable services:
  - 1) electrical service connection
  - 2) tubing connections
  - 3) gas regulators
  - 4) gas tanks
  - 5) unit leveling
  - 6) room ventilation
  - 7) adverse ambient temperature fluctuations
  - 8) any repair external to the unit
- **Damage resulting from accident, alteration, misuse, abuse, fire, flood, acts of God, or improper installation.**
- **Repairs to parts or systems resulting from unauthorized unit modifications.**
- **Any labor costs other than that specified during the parts and labor warranty period, which may include additional warranty on CO<sub>2</sub> sensors, blower motors, water jackets, etc.**

# Table of Contents

<b>Section 1</b>	<b>Installation and Start-Up</b>	<b>1-1</b>
	Name and Description of Parts	1-1
	Control Panel Keys, Displays & Indicators	1-2
	Control Panel Operation	1-3
	Installing the Unit	1-4
	Stacking the Incubators	1-5
	Preliminary Cleaning	1-7
	Installing Access Port Filter	1-8
	Installing Air Sample Filter	1-8
	Installing HEPA Filter	1-8
	Installing Shelves	1-9
	Leveling the Unit	1-9
	Connecting Unit to Electrical Power	1-9
	Filling Water Jacket	1-10
	Filling Humidity Pan	1-11
	Connecting CO <sub>2</sub> or N <sub>2</sub> (O <sub>2</sub> Option) Gas Supply	1-12
	Incubator Start-Up	1-13
	Setting the Operating Temperature	1-14
	Setting the Overtemp Setpoint	1-14
	Setting CO <sub>2</sub> Setpoint	1-15
	Setting O <sub>2</sub> Setpoint	1-16
<b>Section 2</b>	<b>Calibration</b>	<b>2-1</b>
	Calibrating the Temperature	2-2
	Calibrating T/C CO <sub>2</sub> System	2-2
	Calibrating Infra-Red CO <sub>2</sub> System	2-3
	Calibrating the O <sub>2</sub> System	2-4
	O <sub>2</sub> Calibration at 20.7%	2-5
	O <sub>2</sub> Offset Calibration	2-6
	Calibrating Relative Humidity	2-6



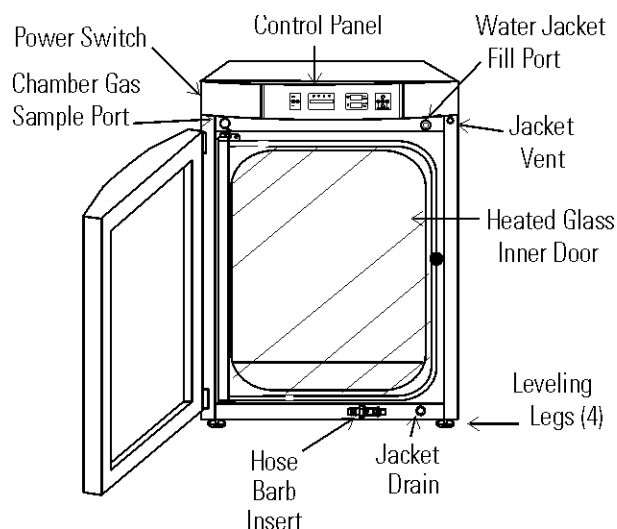
<b>Section 3</b>	<b>Configuration</b>	<b>3-1</b>
	Turning the Audible Alarm ON/OFF	3-1
	New HEPA Filter	3-1
	Setting the REPLACE HEPA Filter Reminder	3-2
	Setting an Access Code	3-2
	Setting Low Temp Alarm Limit	3-3
	Setting High Temp Alarm Limit	3-3
	Enabling Temp Alarms to Trip Contacts	3-4
	Setting Low CO <sub>2</sub> Alarm Limit	3-4
	Setting High CO <sub>2</sub> Alarm Limit	3-5
	Enabling CO <sub>2</sub> Alarms to Trip Contacts	3-5
	Setting Zero & Span #s for T/C CO <sub>2</sub> Sensors	3-6
	Setting a Low RH Alarm Limit	3-7
	Enabling RH Alarms to Trip Contacts	3-7
	Setting a Low O <sub>2</sub> Alarm Limit	3-8
	Setting a High O <sub>2</sub> Alarm Limit	3-8
	Enabling O <sub>2</sub> Alarms to Trip Contacts	3-9
	Enabling Temp/RH to be Displayed	3-9
	Enabling CO <sub>2</sub> /O <sub>2</sub> to be Displayed	3-10
<b>Section 4</b>	<b>Alarms</b>	<b>4-1</b>
	Inner Door Open Alarm	4-2
	Sensor Fault Alarms	4-2
	TSNSR1 ERR	4-2
	TSNSR2 ERR	4-2
	REPL O <sub>2</sub> SNSR	4-3
	O <sub>2</sub> SNSR ERR	4-3
	CO <sub>2</sub> SNSR ERR	4-3
	IR AUTOZ ERR	4-3
	IR AUTOZ ERR	4-4
	REPL IR SNSR	4-4
<b>Section 5</b>	<b>Routine Maintenance</b>	<b>5-1</b>
	Cleaning the Interior	5-1
	Cleaning Cabinet Exterior	5-4
	Cleaning Humidity Pan	5-4
	Replacing Fuses	5-4
	HEPA Filter Maintenance	5-6
	Replacing Air Sample Filter	5-6
	Replacing Access Port Filter	5-6
	Draining Water Jacket	5-7
	Adding /Replenishing Rust Inhibitor	5-8
	O <sub>2</sub> Sensor Fuel Cell	5-8
	Reversing the Door Swing	5-10

<b>Section 6</b>	<b>Factory Installed Options</b> .....	<b>6-1</b>
	Humidity Readout .....	6-1
	Remote Alarm Contacts .....	6-2
<b>Section 7</b>	<b>Specifications</b> .....	<b>7-1</b>
<b>Section 8</b>	<b>Spare Parts</b> .....	<b>8-1</b>
	All Models .....	8-1
	230V Units .....	8-2
	115V Units .....	8-2
	IR or O <sub>2</sub> Units .....	8-3
	Humidity Readout Units .....	8-3
	Field Installation Kits by qualified service personnel .....	8-3
	Exploded Parts Drawings .....	8-4
<b>Section 9</b>	<b>Electrical Schematics</b> .....	<b>9-1</b>
<b>Section 10</b>	<b>Warranty Information</b> .....	<b>10-1</b>



## Section 1 Installation and Start-Up

### Name and Description of Parts




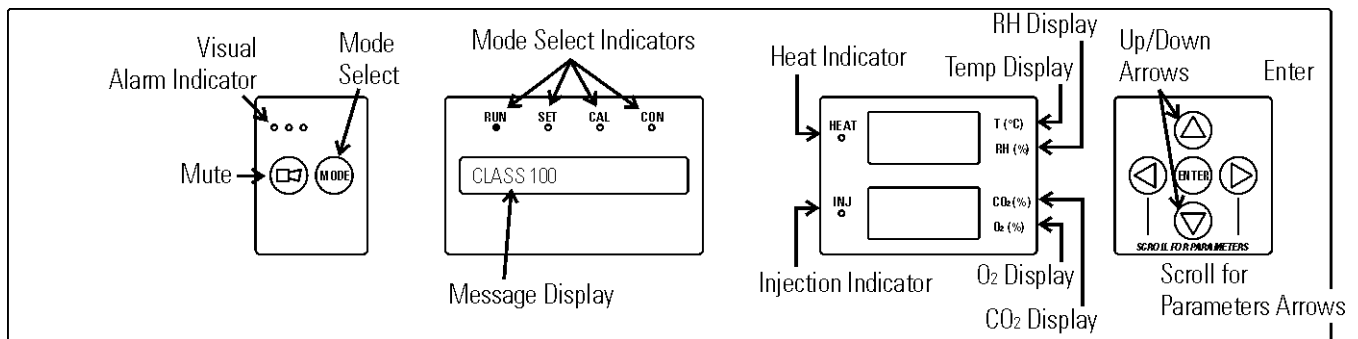
**Figure 1-1.** Water Jacket Incubator

- Heated Inner Door - Keeps chamber interior dry. Reversible to opposite swing (Section 5).
- Chamber Gas Sample Port - Used for sampling chamber CO<sub>2</sub> content, using a Fyrite or similar instrument. Cap when controlling O<sub>2</sub>.
- Main Power Switch
- Control Panel - Keypad, Displays and Indicators (Figure 1-2).
- Fill Port - Used for filling the water jacket.
- Water Jacket Vent - Do not cover! Allows air to escape from the water jacket during filling and normal expansion and contraction when the incubator heats or cools.
- Leveling Legs - Used to level the unit.
- Water Jacket Drain – Use hose barb insert included.

**Note** The incubators are stackable. Instructions follow in this section. s

## Control Panel Keys, Displays & Indicators

-  - mutes the audible alarm.
- Visual Alarm Indicator - Pulses on/off during an alarm condition in the unit.
- MODE Select Switch - Used to select Run, Setpoints, Calibration and System Configuration Modes.
- Message Display - Shows system status.
- Mode Select Indicators -
  - RUN: Run Menu
  - SET: Set Points Menu
  - CAL: Calibrate Menu
  - CON: Configuration Menu
- Up/Down Arrows: Increases or decreases number parameter values, toggles choice parameter values.
- ENTER: Accepts changes to calibration settings
- HEAT: Lights when power is applied to heaters.
- T (°C): Shows current chamber temperature in °C.
- RH (%): Shows percentage of humidity currently inside chamber.
- SCROLL FOR PARAMETERS arrows: Moves through choices in selected mode.
- INJ: Lights when gas is injected into the chamber.
- %CO<sub>2</sub>: Shows percentage of CO<sub>2</sub> in chamber.
- %O<sub>2</sub>: Shows percentage of O<sub>2</sub> in chamber, if applicable.



**Figure 1-2.** Control Panel

## Control Panel Operation

The Series 8000WJ incubator has four basic modes to allow incubator setup. The modes are as follows: Run, Setpoints, Calibration and System Configuration.

- RUN is the default mode that the incubator will normally be in during operation.
- SET is used to enter system setpoints for incubator operation.
- CAL is used to calibrate various system parameters to the customer's satisfaction.
- CON is the system configuration mode that allows for custom setup of various options.

**SCROLL FOR PARAMETERS:** Steps the operator through the parameters of SET, CAL and CON modes. The right arrow goes to the next parameter, the left arrow returns to the previous parameter.

**Up Arrow:** Increases or toggles the parameter value that has been selected in the SET, CAL and CON modes.

**ENTER:** Must press ENTER key to save to memory all changed values.

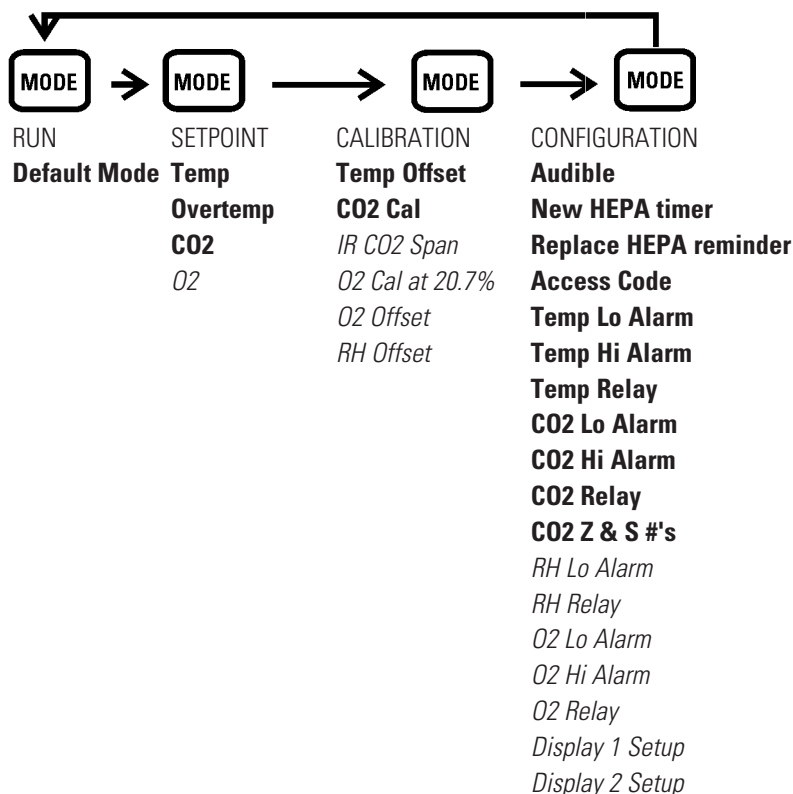
**Down Arrow:** Decreases or toggles the parameter values that have been selected in the SET, CAL and CON modes.

 **Key:** Press to mute the audible alarm. See Section 4 for alarm ringback times.

**Message Display:** Shows system status (Mode) at all times. Shows CLASS 100 or SYSTEM OK during normal operation or alarm messages if the system detects an alarm condition. See Section 4, Alarms. The message 'CLASS 100' is a timing mechanism indicating that, under normal operating conditions with the HEPA filter installed, the air inside the chamber meets the Class 100 air cleanliness standard for particulates of 0.5 micron size or larger per cubic foot of air. (For further information on the Class 100 classification of air quality, see Appendix A.)

**Upper and Lower Displays:** The upper display shows Temp and RH, depending on options installed. The lower display shows CO<sub>2</sub> and O<sub>2</sub>, depending on options installed.

## Control Panel Operation (continued)



### Base Unit Displays

*Option Displays*

## Installing the Unit

1. Locate the unit on a firm, level surface capable of supporting the unit's operational weight of 365 lbs. (166 kg) plus the estimated total sample and user equipment weight. The overall space should be well ventilated and large enough that the concentration of carbon dioxide (CO<sub>2</sub>) and/or nitrogen (N<sub>2</sub>) gases cannot accumulate, thus displacing the oxygen (O<sub>2</sub>) to dangerous levels in the room. Read Connecting CO<sub>2</sub> or N<sub>2</sub> (O<sub>2</sub> System) Gas Supply below carefully.
2. Locate the incubator away from exterior doors and windows as changes in outdoor temperatures and contact with direct sunlight can affect the anti-condensation functions of the unit.
3. Position the door opening away from forced air heating and cooling ducts as these can carry dust, dirt and other contaminants into the incubator chamber, as well as negatively affect the door opening recovery time.
4. Maintain a minimum six inch clearance behind the unit for electrical and gas hook-up. In addition, a minimum three inch ventilation clearance is needed on each side.

## Installing the Unit (continued)

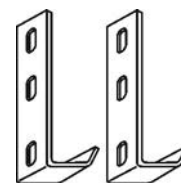
- It is recommended that the unit be placed on a manufacturer approved stand or roller base, not directly on the floor, as the air movement created each time the door is opened will pull dust, dirt and other contaminants into the incubator chamber.

**Warning** This incubator weighs 265 lbs (120 kg) before filling the water jacket. If the unit has been in service, disconnect the power cord from both the unit and the power source, turn off all gas regulators, and disconnect all tubing and any other connections from the rear of the electronics drawer before completely draining the water jacket. Do not lift the incubator by the outer door as this could damage the hinges and/or the outer door may come loose. Note that the outer door can be removed to reduce weight and make it easier to lift the unit by its corners. Have sufficient personnel available to lift the unit into place using proper lifting techniques according to local EHS policies. Lift from both sides of the unit (left and right, not front and back) holding it from the bottom. Do not push or pull the incubator along the floor as this could damage the leveling feet. s

## Stacking the Incubators

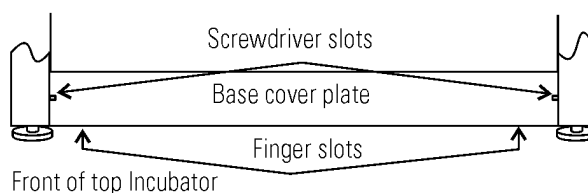
**Warning** If the units have been in service, disconnect the power cord connector and drain the water jacket of the designated top unit before stacking. s

**Note** Stacking brackets (shown at right) stacking bolts, washers, and bolts for stacking are included with each unit. s



**Figure 1-3.** Stacking brackets

- Designate one incubator to be the top unit and the other as the bottom unit. Remove the base cover plate from the top unit using the finger holes in the base or using a slotted screwdriver (Figure 1-4).



**Figure 1-4.** Remove Base Cover Plate



## Stacking the Incubators (continued)

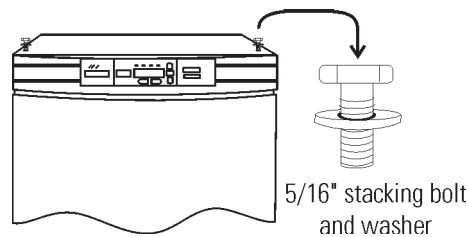
- Note the two slots in the base of the incubator which accommodate the stacking bolts. Refer to Figure 1-5.



Front of top incubator, base cover plate removed

**Figure 1-5.** Stacking Bolt Slots

- Remove the two plastic plugs from the bolt holes in the exterior top of the bottom unit. Install the 1/2" long 5/16-18 stacking bolts and washers into the bolt holes - do not tighten the bolts at this time. Refer to Figure 1-6.

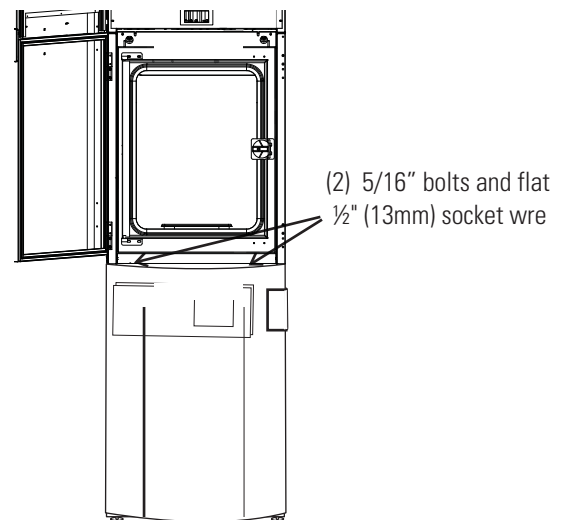


**Figure 1-6.** Bolt Holes

- Unscrew and remove the leveling feet from the top unit and lift it onto the bottom unit, off-setting the base of the top unit approximately 2-3 inches behind the stacking bolts and washers.

**Warning** This incubator weighs 265 lbs (120 kg) before filling. Have sufficient personnel available when lifting.

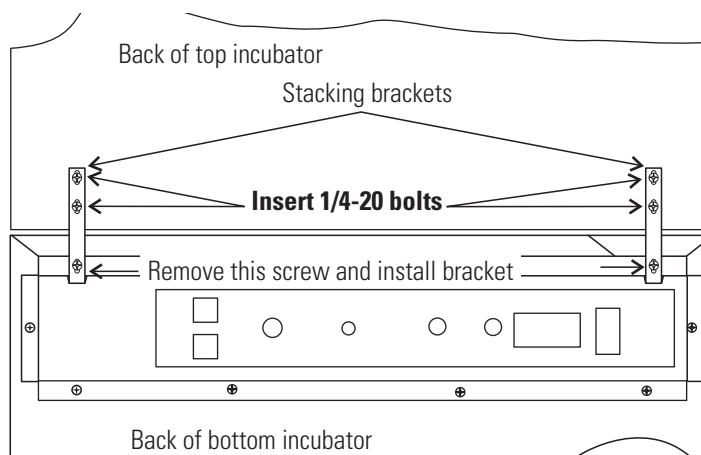
- Align the sides of the top unit and bottom unit and slide the top unit forward until the slots in the base of the top unit align with the 5/16"-18 stacking bolts in the exterior top of the bottom unit (Figure 1-7).



**Figure 1-7.** Align Slots and Bolts

## Stacking the Incubators (continued)

6. Remove and save the two screws from the back of the control panel on the bottom unit as identified in Figure 1-8.



**Figure 1-8.** Two Screws From Control Panel Back

7. Insert the stacking brackets into the slots on the rear of the control panel of the bottom unit as shown in Figure 1-5. Align the slots in the brackets with the mounting holes on the rear of the incubators. Secure the brackets with the screws saved above and the 1/4-20 bolts provided in the stacking kit. A 7/16" (11 mm) wrench or socket is required for the bolts.
8. Secure the base of the top unit to the exterior top of the bottom unit by tightening the 5/16-18 stacking bolts using a 1/2" (13 mm) wrench or suitable tool.
9. Replace the base cover on the top unit.
10. The stacked units are ready to be placed into service.

**Note** If desired, refer to center-of-gravity and attachment information in Section 7.

## Preliminary Cleaning

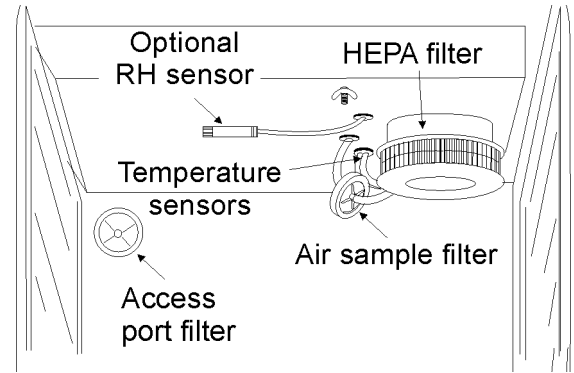
1. Remove the protective plastic coating on the shelf supports, duct sheets and air duct, if present.

**Caution** Before using any cleaning or decontamination method except those recommended by the manufacturer, users should check with the manufacturer that the proposed method will not damage the equipment. s

2. Using a suitable laboratory disinfectant, such as quaternary ammonium, thoroughly clean all interior surfaces to remove any residues which may remain from production of the incubator. Refer to Section 5.

## Installing Access Port Filter

Locate the opening in the top left corner of the interior chamber. Remove the tape from the opening on the outside of the unit. Locate the stopper with filter in the hardware bag. Install in the opening inside the chamber.



**Figure 1-9.** Filter and Sensor Locations

## Installing Air Sample Filter

1. Remove the filter from the shipping bag.
2. Separate one section of the tubing from the filter. Install this section to the fitting on the blower plate.
3. After installing the top duct, connect the filter assembly to the tubing coming through the top duct.
4. Insert the free end of the air sample filter tubing into the larger hole in the back of the blower scroll. See Figure 1-9 for completed configuration.

## Installing HEPA Filter

**Note** There are standard and Volatile Organic Chemical (VOC) HEPA filters available. Be certain to install the correct filter for the given application. See Section 8.

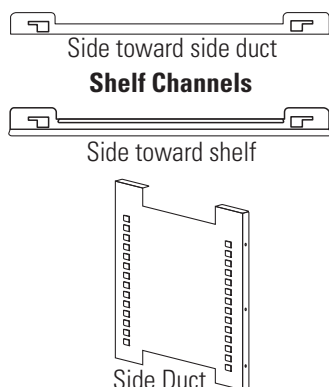
**Caution** Handle the HEPA filter very carefully as the filter media can be easily damaged. Do not touch the filter media during installation. s

1. Remove the filter from the shipping box.
2. Remove the plastic coating from the filter and inspect the filter for any visible signs of damage.
3. Install the filter as shown in Figure 1-9.
4. The default HEPA filter replacement reminder was set at the factory for 6 months. See Section 3 to change the timer value.

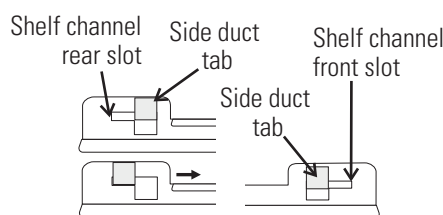
**Caution** To avoid damage to the incubator, do not operate the unit without the HEPA filter in place. If higher RH is needed and CLASS 100 air quality conditions are not required, use the restrictor plate instead of the HEPA filter in order to maintain proper air flow. s

## Installing Shelves

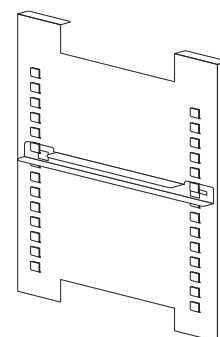
1. Install the side ducts with the tabs facing into the center of the chamber with their slots up. There are no right side or left side ducts, simply rotate one of them to fit the opposite side. Tilt the side ducts as they are placed in the chamber so the tops fit into the top air duct, then guide them into the vertical position. Figure 1-10 shows the duct as it would be oriented for the right side of the chamber. Note that there is no difference between left and right side shelf channels.



**Figure 1-10.** Shelf Channels and Side Duct



**Figure 1-11.** Slot and Tab



**Figure 1-12.** Channel Installed

2. Install the shelf channels by placing the channel's rear slot over the appropriate rear tab on the side duct. Pull the shelf channel forward and engage the channel's front slot into the side duct's appropriate forward tab. Refer to Figure 1-11.
3. Figure 1-12 shows one of the channels installed on the right side duct.

## Leveling the Unit

Check the unit for being level by placing a bubble-style level on one of the shelves. Turn the hex nut on the leveler counterclockwise to lengthen the leg, or clockwise to shorten it. Level the unit front-to-back and left-to-right.

## Connecting Unit to Electrical Power

See the serial tag on the side of the unit or the Specifications section for electrical specifications. Refer to the electrical schematics toward the end of this manual.

**Warning** Connect the incubator to a grounded dedicated circuit only. The power cord connector is the mains disconnect device for the incubator. Position the unit to allow unobstructed access to the power cord so that it can be easily disconnected in case of an emergency. s

Plug the provided power cord into the power inlet connector (Figure 1-13) on the back of the unit, and into the grounded dedicated circuit.

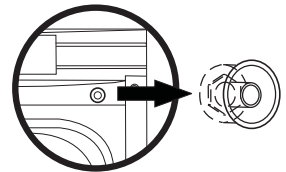
## Filling Water Jacket

**Caution** For stacked incubators, completely fill the bottom unit before filling the top unit in order to maintain the lowest possible center of gravity. s

Turn the power switch On. ADD WATER will appear in the message center. Press the Silence key to silence the alarm for 15 minutes. If the water jacket is empty, yet the message center does not display ADD WATER, then the liquid level switch could be faulty. Call Technical Services.

Remove the fill port plug and set aside.

For ease of connecting/disconnecting, use the silicone tubing provided to connect directly to the fill port. See Figure 1-1 and this detail. Do not install vinyl tubing directly onto the fill port as it may be difficult to remove. A 3/8" to 3/8" hose connector has been provided to attach lengths of vinyl tubing to the silicone fill port tubing to reach the water source, or simply attach the included funnel to manually fill using small containers.



**Detail 1**

**Caution** Chlorine is detrimental to stainless steel. Using chlorinated tap water or additives that contain chlorine will void water jacket warranty! s

For optimum incubator operation, the recommended water type is sterile distilled Type 2 water with a pH between 7 and 9, and a resistivity between 50 kΩ - 1 MΩ-cm (conductivity of 1-20 μS/cm). Reverse osmosis (RO), deionized (DI), and ultrapure Type 1 water can corrode stainless steel, copper, glass and other incubator components over time because high purity water with low ionic content and high resistivity will actively pull ions from these materials. If sterile distilled Type 2 water is not readily available, then these other water types could be used by adding the necessary amount of a sterile solution of weak sodium bicarbonate to raise the pH between 7 and 9 and lower the resistivity between 50 kΩ - 1 MΩ-cm (conductivity of 1-20 μS/cm).

**Caution** Do not block the water jacket vent! s

Carefully fill the jacket with water. Do not overfill the jacket as the excess water will pour out through the water jacket vent (see Figure 1-1).

When the jacket is full, the audible alarm will sound a continuous tone for 10 seconds and the alarm condition will be cleared. Refer to Section 4.

**Note** Model 3100 Series incubators are shipped from the factory with a rust inhibitor added to 0.2 gallons (0.8 liters) of water in the jacket. The rust inhibitor must be replenished every 2 years. See Section 5 to drain water from the jacket and add the correct proportion of rust inhibitor mixture back into the jacket.

## **Filling Water Jacket (continued)**

Disconnect the silicone tubing from the fill port, and replace the fill port plug removed above.

**Warning** Once the jacket is full, be certain to clean up any water that may have dripped down the side of the unit or spilled onto the floor during this process to prevent a possible slip hazard. s

## **Filling Humidity Pan**

**Caution** Use of chlorinated water, or decontamination products containing chlorine, will deteriorate the stainless steel and cause rust, voiding the warranty. s

Fill the humidity pan to within 1/2 inch of the top with 0.8 gallons (3 liters) of water, as described in Filling Water Jacket. If the water cannot be sterilized, a 2% solution of quaternary ammonium can be used as a disinfectant additive to control bacteria and other organic contaminants, if desired. Place the pan directly in the center of the incubator floor to ensure optimum humidity and temperature response.

For applications requiring higher humidity conditions, the pan should be placed against the left side wall of the incubator. The ductwork has been designed for this purpose. This will also accelerate humidity recovery after door openings. Even higher humidity conditions can be achieved by removing the humidity pan altogether and adding water directly to the bottom of the incubator chamber, being careful to not fill it any higher than the bottom edge of the silicone door gasket. Also the chamber gas sample port (Figure 1-1) may be capped to assist in achieving greater RH. In some ambients, this may cause condensation to form in the chamber.

**Warning** Wipe up any excessive condensation to prevent it from dripping onto the floor creating a possible slip hazard. For stacked incubators, this may also include the drip channel (just behind the face of the control panel drawer) on the bottom unit used to direct water away from the electronics drawer. s

**Warning** When the humidity pan is in place, be certain to clean up any water that may have spilled in the unit or onto the floor during this process, to prevent a possible slip hazard. For stacked incubators, this may also include the drip channel (just behind the face of the control panel drawer) on the bottom unit, used to direct water away from the electronics drawer. s

**Caution** Check the level and change the water frequently to avoid contamination. Do not allow the water level to fluctuate significantly. “Dry-outs” will have an adverse effect on the humidity level as well as CO<sub>2</sub> calibration. s

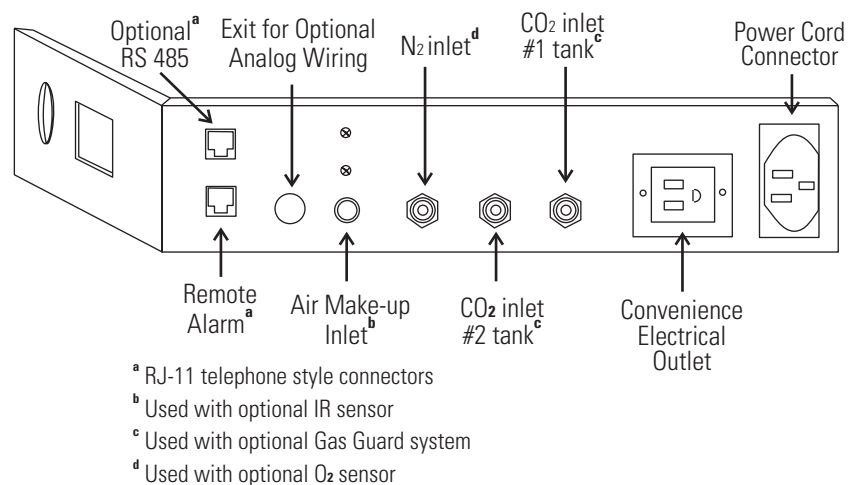
## Connecting CO<sub>2</sub> or N<sub>2</sub> (O<sub>2</sub> Option) Gas Supply

**Warning** High concentrations of CO<sub>2</sub> gas can cause asphyxiation! Refer to local EHS policies regarding employee exposure to CO<sub>2</sub>. Carbon dioxide monitors are recommended for confined areas where concentrations of carbon dioxide gas can accumulate. s

**Warning** High concentrations of N<sub>2</sub> gas can cause asphyxiation! N<sub>2</sub> and CO<sub>2</sub> can displace oxygen levels in a confined atmosphere. Refer to local EHS policies regarding employee exposure to N<sub>2</sub>. Oxygen monitors are recommended for confined areas where concentrations of N<sub>2</sub> or CO<sub>2</sub> may displace oxygen. s

**Note** For customer convenience, 12 feet of 1/4" ID vinyl tubing with two 3/8" hose clamps are included in the shipping materials for connecting the unit to the gas supply.

The incubator has 1/4" barbed fittings on the back of the cabinet to connect the gas supply. Refer to Figure 1-13. The fitting is labeled CO<sub>2</sub> Inlet #1. The N<sub>2</sub> Inlet is also labeled.



**Figure 1-13.** Back Panel and All Connections

Install 3/8" hose clamps to secure the 1/4" ID tubing to the barbed fittings on the regulator and the rear of the electronics drawer.

**Warning** Check for leaks at all connections. s

For units having the CO<sub>2</sub> or N<sub>2</sub> Gas Guard option, see Section 6.

**Warning** This incubator is designed to be operated with CO<sub>2</sub> gas, or N<sub>2</sub> in those units with an O<sub>2</sub> control system. Connecting a flammable or toxic gas can result in a hazardous condition. Gases other than CO<sub>2</sub> or N<sub>2</sub> should not be connected to this equipment. s

## Connecting CO<sub>2</sub> or N<sub>2</sub> (O<sub>2</sub> Option) Gas Supply (cont.)

The CO<sub>2</sub> gas supply being connected to the incubator should be industrial grade, 99.5% pure. Do not use CO<sub>2</sub> gas cylinders equipped with siphon tubes. A siphon tube is used to extract liquid CO<sub>2</sub> from the cylinder, which can damage the pressure regulator. Consult with your gas supplier to ensure that the CO<sub>2</sub> cylinder does not contain a siphon tube. The N<sub>2</sub> gas supply being connected to the incubator should be 99.99% pure. Do not use liquid nitrogen.

A two-stage pressure regulator is required to be installed on the outlet valve of the gas cylinder. The high pressure gauge at the tank should have 0-2000 psig range and the low pressure gauge at the incubator inlet should have a 0-30 psig range. Input pressure to the incubator must be maintained at 15 psig (103.4 kPa) maximum for proper performance of the CO<sub>2</sub> or O<sub>2</sub> control system.

If higher purity CO<sub>2</sub> is desired inside the incubator (greater than 99.5% pure), the pressure regulator should be constructed with a stainless steel diaphragm along with specifying the purity of the CO<sub>2</sub> from the gas supplier. Follow the manufacturer's instructions to ensure proper and safe installation of the pressure regulator on the gas cylinder. Gas cylinders should be secured to a wall or other stationary object to prevent tipping. Position the gas cylinders to allow unobstructed access to the pressure regulators so that they can be easily turned off in case of an emergency. Consult your facility safety officer to ensure that the equipment is installed in accordance with the codes and regulations applicable in your area.

**Note** Stored product should be protected by an activated alarm system capable of initiating a timely response 24 hours/day. These alarms provide interconnect for centralized monitoring. See Factory Installed Options for more information.

## Incubator Start-Up

Now that the incubator has been properly installed, connected to power, filled with water, humidity pan filled, and connected to gas supplies, system setpoints can be entered. The following setpoints can be entered in set mode: temperature, over temperature, CO<sub>2</sub> and O<sub>2</sub>. To enter Set Mode, press the Mode key until the Set indicator lights. Press the right and/or left arrow keys until the proper parameter appears in the message center. See Chart 1-1 for more detail.

**Caution** It is the responsibility of the user to validate proper operation of each incubator in their specific application(s) with respect to unit location, operating environment, and settings. s



## Setting the Operating Temperature

Incubator Model 3428 and 3429 have an operating temperature range of 10 to 55°C, Model 3422 and 3423 a range of 10 to 50°C and Models 3424, 3425, 3426 and 3427 a range of 10 to 45°C. The incubator is shipped from the factory with a temperature setpoint of 10°C. At this setting, all heaters and alarms are turned off.

To change the operating temperature setpoint:

1. Press the MODE key until the SET indicator lights.
2. Press the right arrow until TEMP XX.XC is displayed in the message display.
3. Press the up/down arrow until the desired temperature setpoint is displayed.
4. Press ENTER to save the setpoint.
5. Press the MODE key until the RUN Indicator lights to go to RUN mode, or right/left to go to next/previous parameter.

## Setting the Overtemp Setpoint

**Caution** In the event that the heaters are locked on as a result of a failure in the main temperature control, the independent overtemp circuit is designed as a safety to protect the incubator only. It is not intended to protect or limit the maximum temperature of the cell cultures or customer's equipment inside the incubator if an overtemp condition occurs. s

The incubator is equipped with an independent circuit that monitors the air temperature in the cabinet. Should the main temperature control fail, the overtemp circuit will disconnect power to all heaters when the chamber temperature reaches the Overtemp setpoint. When the chamber temperature falls below the Overtemp setpoint, the overtemp circuit will reconnect power to all heaters. An incubator operating in the overtemp condition will maintain the chamber temperature approximately 1°C around the Overtemp setpoint.

The factory setting for the Overtemp is 40°C. It can be set over a range of temp setpoint +0.5°C to 60°C. If the temperature setpoint is moved above the Overtemp setpoint, the Overtemp will automatically update to 1.0°C+ the temp setpoint. It is recommended that the Overtemp setpoint be 1°C over the temp setpoint.

## Setting the Overtemp Setpoint (continued)

To set the Overtemp setpoint:

1. Press the MODE key until the SET indicator lights.
2. Press the right arrow until OTEMP XX.XC is displayed in message display.
3. Press the up/down arrow until the desired Overtemp setpoint is displayed.
4. Press ENTER to save the setpoint.
5. Press the MODE key until the RUN Indicator lights to go to RUN mode or right/left to go to next/previous parameter.

## Setting CO<sub>2</sub> Setpoint

All T/C CO<sub>2</sub> cells are precalibrated at the factory at 37°C, high humidity, and 10% CO<sub>2</sub>. Therefore, if a temperature setpoint of 37°C has been entered, the humidity pan filled and the CO<sub>2</sub> control is to run between 0-10% with a T/C CO<sub>2</sub> sensor, the CO<sub>2</sub> setpoint may be entered immediately. Otherwise, it is important to allow the unit 12 hours to stabilize at the temperature setpoint before entering the CO<sub>2</sub> setpoint.

All models of the incubator have a CO<sub>2</sub> setpoint range of 0.0% to 20.0%. The incubator is shipped from the factory with a CO<sub>2</sub> setpoint of 0.0%. At this setting, all CO<sub>2</sub> control and alarms are turned off.

To change the CO<sub>2</sub> setpoint:

1. Press the MODE key until the SET indicator lights.
2. Press the right arrow until CO<sub>2</sub> XX.X% is displayed in message display.
3. Press the up/down arrow until the desired CO<sub>2</sub> setpoint is displayed.
4. Press ENTER to save the setpoint.
5. Press the MODE key until the RUN Indicator lights to go to RUN mode or right/left to go to next/previous parameter.

## **Setting O<sub>2</sub> Setpoint**

Models 3424/3425 and 3426/3427 have a built-in O<sub>2</sub> control system. The O<sub>2</sub> setpoint range is 1.0% to 21.0%. The incubator is shipped from the factory with a O<sub>2</sub> setpoint of 21.0%. At this setting, all O<sub>2</sub> control and alarms are turned off. The gas sample port must be capped when running controlled O<sub>2</sub> levels.

To change the O<sub>2</sub> setpoint:

1. Press the MODE key until the SET indicator lights.
2. Press the right arrow until O<sub>2</sub> XX.X% is displayed in the message display.
3. Press the up/down arrow until the desired O<sub>2</sub> setpoint is displayed.
4. Press ENTER to save the setpoint.
5. Press the MODE key until the RUN Indicator lights to go to RUN mode, or right/left to go to next/previous parameter.

# Chart 1-1

## Set Mode

Press MODE to light SET

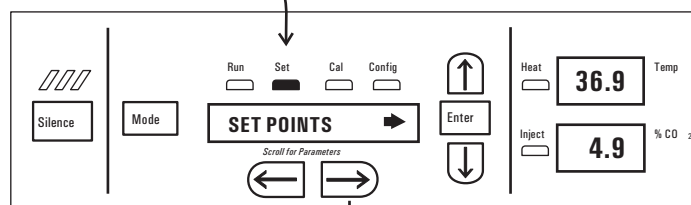
To Set:

Operating Temperature

Over Temperature

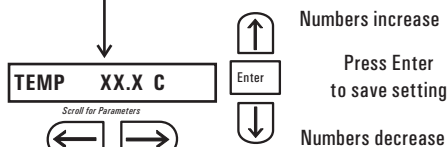
Percent CO<sub>2</sub>

Percent O<sub>2</sub>

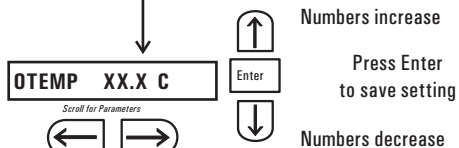


Press MODE to move to CALIBRATE mode

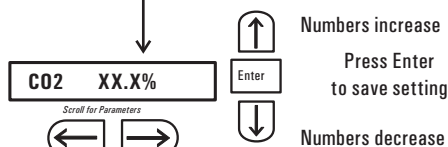
Press ← to return to previous parameter



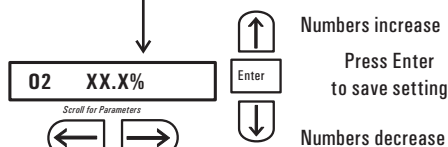
Press ← to return to previous parameter



Press ← to return to previous parameter



Press ← to return to previous parameter



**Section 1**

Installation and Start-Up

## Section 2 Calibration

After the unit has stabilized, several different systems can be calibrated. In the Calibration Mode, the air temperature, CO<sub>2</sub> reading, O<sub>2</sub> reading and RH reading can all be calibrated to reference instruments. To enter Calibration Mode, press the Mode key until the CAL indicator lights. Press the right and/or left arrow until the proper parameter appears in the message center. See Chart 2-1 for more detail.

Calibration frequency is dependent on use, ambient conditions, and accuracy required. Good laboratory practice would require at least an annual calibration check. On new installations, all parameters should be checked after the stabilization period.

Prior to calibration, the user should be aware of the following system functions. While the unit is in Calibration Mode, all system control functions will be stopped so that the unit remains stable. Readout of the system being calibrated will appear as “—” on the readout displays. If no keys are pressed for approximately five minutes while in Calibration Mode, the system will reset to Run Mode so that control functions can be reactivated.

**Caution** Before making any calibration or adjustments to the unit, it is imperative that all reference instruments be properly calibrated. It is the responsibility of the user to understand the interactive effects of temperature, CO<sub>2</sub>%, O<sub>2</sub>% and RH% on each other and the independent reference instruments, especially a Fyrite® sensor. Read and understand all reference instrument operating manuals before use. Double-check all values entered into the incubator for accuracy before completing the calibration and returning the unit to service. s

## Calibrating the Temperature

Place the calibrated instrument in the center of the chamber. The instrument should be in the airflow, not against the shelf. Before calibration, allow the cabinet temperature to stabilize.

**Start-Up** - Allow 12 hours for the temperature in the cabinet to stabilize before proceeding.

**Operating Unit** - Allow at least two hours after the display reaches setpoint for the temperature to stabilize before proceeding.

1. Press the MODE key until the CAL indicator lights.
2. Press the right arrow until TMPCAL XX.XC appears in the message display.
3. Press the up/down arrow to match the display to a calibrated instrument.
4. Press ENTER to store the calibration into memory.
5. Press the MODE key to return to RUN mode or the right/left arrow to go to the next/previous parameter.

## Calibrating T/C CO<sub>2</sub> System

Models 3424, 3425, 3428, and 3429 have a thermal conductivity (T/C) CO<sub>2</sub> sensor. Thermal conductivity of the incubator atmosphere is not only effected by the quantity of CO<sub>2</sub> present, but also by the air temperature and the water vapor present in the incubator atmosphere. In monitoring the effects of CO<sub>2</sub>, air temperature and absolute humidity must be held constant so any change in thermal conductivity is caused only by a change in CO<sub>2</sub> concentration.

Any changes to the temperature setpoint and/or humidity level control method (humidity pan location, capping ports, HEPA filter type) will necessitate a recalibration of the CO<sub>2</sub> control.

Some T/C CO<sub>2</sub> sensors go through an aging period, especially on new installations. Calibration should be checked on a weekly basis and adjusted as necessary. When stabilization occurs, checks can become less frequent.

## Calibrating T/C CO<sub>2</sub> System (continued)

**Start-up** - The CO<sub>2</sub> sensor has been calibrated at the factory for 37°.

Allow temperature, humidity and CO<sub>2</sub> levels in the chamber to stabilize at least 12 hours before checking the CO<sub>2</sub> concentration with an independent instrument.

**Presently operating** - Make sure the chamber doors are closed. Allow at least 2 hours after the temperature and CO<sub>2</sub> displays reach their setpoints for chamber atmosphere stabilization.

1. Measure the CO<sub>2</sub> concentration in the chamber through the gas sample port with an independent instrument. Take several readings to ensure accuracy.
2. Press the Mode key until the CAL indicator lights.
3. Press the right arrow until CO<sub>2</sub>CAL XX.X% is displayed in the message center.
4. Press the up/down arrow to change the display to match the independent instrument.
5. Press Enter to store calibration.
6. Press the Mode key to return to Run Mode, or the right or left arrow keys to go to the next/previous parameter.

## Calibrating Infra-Red CO<sub>2</sub> System

Models 3422, 3423, 3426, and 3427 have an infra-red CO<sub>2</sub> sensor.

Infra-red CO<sub>2</sub> sensors are not effected by chamber atmosphere temperature or humidity. However, the light detector in the sensor is effected by wide temperature changes. Therefore, changing temperature setpoints could necessitate a recalibration of the CO<sub>2</sub>. Chamber temperature should be allowed to stabilize before checking CO<sub>2</sub> concentrations with an independent instrument, especially on start-up.

All models equipped with an IR/CO<sub>2</sub> sensor have an automatic calibration that occurs every 24 hours, and lasts for 5 to 6 minutes. During automatic calibration, the CO<sub>2</sub> display is blanked out and HEPA filtered room air is pumped through the CO<sub>2</sub> sensor. A new CO<sub>2</sub> calibration value is stored in memory for use as the 0.0% CO<sub>2</sub> reference point. The keypad / control panel is locked during calibration, with the message display reading CO<sub>2</sub> AUTO CAL.



## Calibrating Infra-Red CO<sub>2</sub> System (cont.)

**Startup** - Allow the temperature and the CO<sub>2</sub> of the cabinet to stabilize at least 12 hours before proceeding.

**Operating Unit** - Allow CO<sub>2</sub> to stabilize at least 2 hours at setpoint before proceeding.

To ensure accurate calibration, the unit will not allow CO<sub>2</sub> to be spanned below 3%. If the cabinet does not contain at least 3% CO<sub>2</sub>, increase the setpoint and allow the unit to stabilize before completing this procedure.

1. Measure the CO<sub>2</sub> concentration in the chamber through the gas sample port with a Fyrite or other independent instrument. Several readings should be taken to ensure accuracy.
2. Press the MODE key until the CAL indicator lights.
3. Press the right arrow until IR CAL XX.X% appears in the message display.
4. Press the up/down arrow to adjust the display to match the independent instrument reading.
5. Press ENTER to store calibration.
6. After ENTER is pressed, the unit will go into a calibration cycle that lasts 5 to 6 minutes. Control panel is locked during this calibration cycle.
7. Press the MODE key to return to RUN mode.

## Calibrating the O<sub>2</sub> System

Models 3424, 3425, 3426 and 3427 have an O<sub>2</sub> control sensor. The sensor is a fuel cell that puts out a linear millivolt signal based on O<sub>2</sub> content of the chamber. The life expectancy for an O<sub>2</sub> sensor fuel cell is approximately 5 years (specified at an oxygen concentration of 20.7% and ambient temperature of 20°C) even if the unit is not in use. During normal incubator operation, exposure to lower concentrations of oxygen increases the life expectancy, whereas increased ambient temperatures decreases it. The cumulative effects of these conditions over time are unpredictable, but the unit is able to measure the remaining life of the O<sub>2</sub> sensor fuel cell based on its voltage output when the system is calibrated using the 20.7% method and display a sensor fault as required (Section 4); thus, it is important to calibrate the system using this method before each test experiment, or at least every 6 months.

## **Calibrating the O<sub>2</sub> System (continued)**

There are two methods available to calibrate the O<sub>2</sub> system.

- The preferred method calibrates the system to the known ambient O<sub>2</sub> value of 20.7% and checks the life of the sensor. This method should be used whenever a new sensor is installed.
- The second method available allows the system to be calibrated to an independent reference instrument by entering an offset.

### **O<sub>2</sub> Calibration at 20.7%**

1. Press the MODE key until the CAL indicator lights.
2. Press the right arrow until the display reads O<sub>2</sub> CAL@20.7%.
3. Press ENTER.
4. OPEN DOOR appears on the display. Open the outer and inner doors.
5. The display reads CALIBRATING.
6. When calibration is complete, approximately 2 minutes, an audible tone sounds and the display returns to O<sub>2</sub> CAL@20.7%.
7. The O<sub>2</sub> display will change to 20.7.
8. Press the MODE key to return to RUN mode.

A new O<sub>2</sub> span value is stored in memory for use as the 20.7% O<sub>2</sub> reference point. The keypad/control panel is "locked-up" during calibration.

## O<sub>2</sub> Offset Calibration

**Startup** - Allow the cabinet to stabilize at least 12 hours before proceeding.

**Operating Unit** - Allow O<sub>2</sub> to stabilize at least 2 hours at set point before proceeding.

Measure the O<sub>2</sub> concentration in the chamber through the gas sample port with an independent instrument. Take several readings to ensure accuracy.

1. Press the MODE key until the CAL indicator lights.
2. Press the right arrow until O<sub>2</sub> OFFS XX.X appears in message display.
3. Press the up/down arrow to adjust display to independent instrument reading.
4. Press ENTER to store calibration.
5. Press the MODE key to return to RUN mode.

## Calibrating Relative Humidity

All Series 8000WJ incubators can be equipped with an optional direct readout relative humidity sensor. This is a readout only of the chamber relative humidity. It does not provide any control of the relative humidity in the cabinet.

**Startup** - Allow 12 hours for the relative humidity and temperature in the chamber to stabilize before proceeding.

**Operating Unit** - Allow at least 2 hours after temperature display reaches setpoint for relative humidity to stabilize before proceeding.

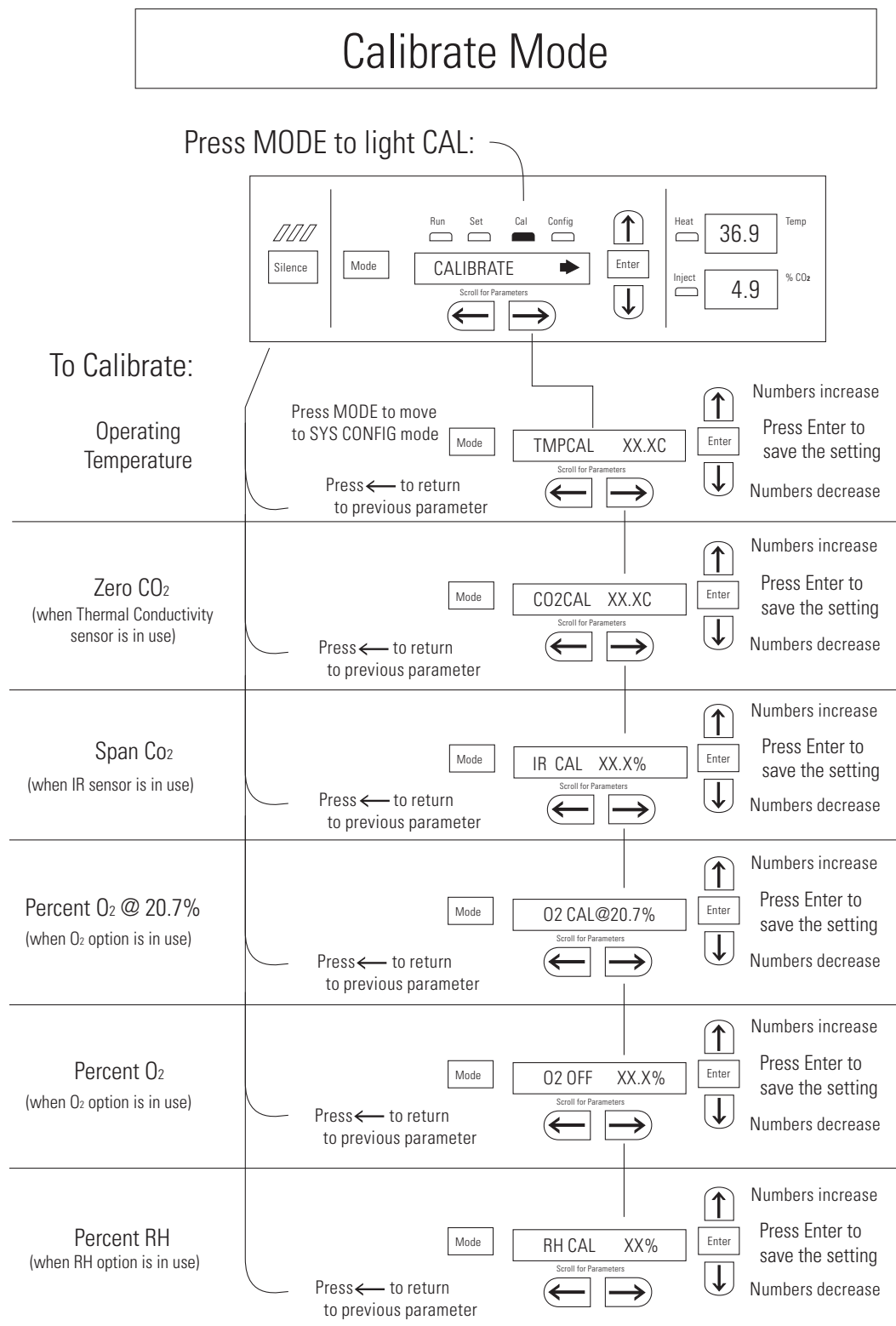
1. Press the Mode key until the CAL indicator lights.
2. Press the right arrow key until RH CAL XX% appears in message center.
3. Press the up/down arrow to match the display to the independent instrument.
4. Press Enter to store the calibration.
5. Press the Mode key to return to Run Mode.

## **Calibrating Relative Humidity (cont.)**

If a reliable RH measuring device is not available, the display may be calibrated to a typical level. This calibration method should be accurate to within 5%.

1. Follow the RH stabilization periods outlined above.
2. With a full humidity pan and stable temperature, the relative humidity in the chamber will be 95%.
3. Using Steps 3-5 of the relative humidity sensor adjustment above, adjust the display to 95%.

Chart 2-1



## Section 3 Configuration

There are many features available in Configuration mode that allow custom setup of the incubator. These features are listed and described below. All features may not be necessary in all applications, but are available if needed. To enter Configuration mode, press the MODE key until the CON indicator lights. Press the right or left arrow until the appropriate parameter appears in the message display. See Chart 3-1 for more detail.

### Turning the Audible Alarm ON/OFF

The audible alarm can be turned on or off. The factory setting is ON.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until AUDIBLE XXX is displayed in message display.
3. Press the up/down arrow to toggle AUDIBLE ON/OFF.
4. Press ENTER to save the setting.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

### New HEPA Filter

When the REPLACE HEPA reminder is displayed and the visual alarm flashes, the specified time has elapsed and the HEPA filter should be replaced. Follow the HEPA Filter Maintenance procedures in Section 5. The factory setting is 182 days (calculated from REPL HEPA XX setting of 6 months).

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until NEW HEPA is displayed in message display.

**Note** Do not press Enter if only viewing remaining days before specified filter replacement time runs out.

3. Press Enter to restart the timer and clear the REPLACE HEPA alarm. This new number is the remaining days before the specified filter replacement time runs out. For example, if 12 months was chosen in the REPL HEPA XX message screen, the NEW HEPA number would be 365 days.
4. Press the MODE key to return to RUN mode.

## Setting the REPLACE HEPA Filter Reminder

A HEPA filter replacement timer can be set for a specific amount of time, from 1 to 12 months of actual unit running time. Time will not accrue when the unit is turned off. The default time is 6 months. When the allotted time has run out, REPLACE HEPA appears in the display and the visual alarm flashes. To set the reminder, use the following procedure.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until REPL HEPA XX is displayed.
3. Press the up/down arrow to choose the number of months desired.
4. Press ENTER to save the number.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

**Note** After the reminder has been set, check the allotted time remaining by going to Configuration mode, then pressing the right arrow until NEW HEPA displays. This number is the remaining days before the filter replacement time specified runs out. For example, if 12 months was chosen in the REPL HEPA XX message screen, the NEW HEPA number would be 365 days.

## Setting an Access Code

A three-digit Access Code can be entered to avoid unauthorized personnel from changing the setpoints, calibration or configuration. A setting of 000 will bypass the access code. The factory setting is 000.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until ACC CODE XXX is displayed in the message display.
3. Press the up/down arrow to change the access code.
4. Press ENTER to save the access code.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## Setting Low Temp Alarm Limit

The low temp alarm limit (tracking alarm) is the deviation from the temperature setpoint, which will cause a low temp alarm. The low temp alarm is variable from 0.5° below setpoint to 5.0° below setpoint. The factory setting is 1.0° below setpoint. A minus sign (-) in the display indicates that the alarm setting is below the setpoint.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until TEMP LO X.XC is displayed in the message display.
3. Press the up/down arrow to change the low temp alarm limit.
4. Press ENTER to save the low temp alarm limit.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## Setting High Temp Alarm Limit

The high temp alarm limit (tracking alarm) is the deviation from the temperature setpoint that will cause a high temp alarm. It should be noted that this varies from the Overtemp setpoint, in that the Overtemp setpoint configures an independent system that monitors temperature and shuts down the system heaters if necessary. The high temp alarm limit is simply set to enable an audible and visual alarm that notifies the user of a problem. The high temp alarm is variable from 0.5° above setpoint to 5.0° above setpoint. The factory setting is 1.0° above setpoint.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until TEMP HI X.XC is displayed in the message display.
3. Press the up/down arrow to change the high temp alarm limit.
4. Press ENTER to save the high temp alarm limit.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.



## **Enabling Temp Alarms to Trip Contacts**

High and Low temperature alarms can be programmed to trip the remote alarm contacts. A setting of ON will cause this, a setting of OFF will not allow temp alarms to trip the contacts. The factory setting is ON.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until TEMP RLY XXX is displayed in the message display.
3. Press the up/down arrow to toggle the setting ON/OFF.
4. Press ENTER to save the setting.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## **Setting Low CO<sub>2</sub> Alarm Limit**

The low CO<sub>2</sub> alarm limit (tracking alarm) is the deviation from the CO<sub>2</sub> setpoint that will cause a low CO<sub>2</sub> alarm. The setpoint is variable from 0.5% CO<sub>2</sub> below setpoint to 5.0% CO<sub>2</sub> below setpoint. The factory setting is 1.0% CO<sub>2</sub> below setpoint. A minus (-) in the display indicates that the alarm setting is below the setpoint.

1. Press the Mode Key until the CON indicator lights.
2. Press the right arrow until CO<sub>2</sub> LO X.X% is displayed in the message display.
3. Press the up/down arrow to change the low CO<sub>2</sub> alarm limit.
4. Press ENTER to save the low CO<sub>2</sub> alarm limit.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## Setting High CO<sub>2</sub> Alarm Limit

The high CO<sub>2</sub> alarm limit (tracking alarm) is the deviation from the CO<sub>2</sub> setpoint that will cause a high CO<sub>2</sub> alarm. The setpoint is variable from 0.5% CO<sub>2</sub> above setpoint to 5.0% CO<sub>2</sub> above setpoint. The factory setting is 1.0% CO<sub>2</sub> above setpoint.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until CO2 HI X.X% is displayed in the message display.
3. Press the up/down arrow to change the high CO<sub>2</sub> alarm limit.
4. Press ENTER to save the high CO<sub>2</sub> alarm limit.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## Enabling CO<sub>2</sub> Alarms to Trip Contacts

High and Low CO<sub>2</sub> alarms can be programmed to trip the remote alarm contacts. A setting of ON will cause this; a setting of OFF will not allow CO<sub>2</sub> alarms to trip the contacts. The factory setting is ON.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until CO2 RLY XXX is displayed in the message display.
3. Press the up/down arrow to toggle the setting ON/OFF.
4. Press ENTER to save the setting.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## Setting Zero & Span #s for T/C CO<sub>2</sub> Sensors

If a new T/C CO<sub>2</sub> sensor is being installed, the two numbers on the factory installed sticker on the T/C cell must be entered to calibrate the CO<sub>2</sub> in the unit.

**Note** For the technician's convenience, a second label containing the two numbers is supplied with the new T/C cell and should be applied to cover the original label, inside the electronics drawer. s

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until T/C ZR# XXXX is displayed in the message display.
3. Press the up/down arrow to change the zero number to match the sticker.

**Note** The '-' symbol denotes the 'minus sign' which is displayed as required by scrolling the zero number though 0.

4. Press ENTER to save the setting.

**Note** Pressing Enter will erase any zero and span offset values previously entered into the calibration menu.

2. Press the right arrow until T/C SP# XXXX is displayed in the message display.
3. Press the up/down arrow to change the span number to match the sticker.
4. Press ENTER to save the setting.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## Setting a Low RH Alarm Limit

On units that have the RH option installed, a low RH alarm limit may be entered. The low RH alarm limit is the % RH in the cabinet that will cause a low RH alarm. The setpoint is variable from setpoint 0 to 90% RH. The factory setting is 0% RH, which disables the alarm.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until RH LO XX% is displayed in the message display.
3. Press the up/down arrow to change the RH low alarm limit.
4. Press ENTER to save the RH low alarm limit.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## Enabling RH Alarms to Trip Contacts

The low RH alarm can be programmed to trip the remote alarm contacts. A setting of ON will cause this, a setting of OFF will not allow the RH alarm to trip the contacts. The factory setting is ON.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until RH RLY XXX is displayed in the message display.
3. Press the up/down arrow to toggle the setting ON/OFF.
4. Press ENTER to save the setting.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## **Setting a Low O<sub>2</sub> Alarm Limit**

On models with a O<sub>2</sub> control system, O<sub>2</sub> alarms may be configured. The low O<sub>2</sub> alarm limit (tracking alarm) is the deviation from the O<sub>2</sub> setpoint that will cause a low O<sub>2</sub> alarm. The setpoint is variable from 0.5% O<sub>2</sub> below setpoint to 5.0% O<sub>2</sub> below setpoint. The factory setting is 1.0% O<sub>2</sub> below setpoint. A minus (-) in the display indicates that the alarm setting is below setpoint.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until O2 LO -X.X% is displayed in message display.
3. Press the up/down arrow to change the low O<sub>2</sub> alarm limit.
4. Press ENTER to save the low O<sub>2</sub> alarm limit.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## **Setting a High O<sub>2</sub> Alarm Limit**

On models with a O<sub>2</sub> control system, O<sub>2</sub> alarms may be configured. The high O<sub>2</sub> alarm limit (tracking alarm) is the deviation from O<sub>2</sub> setpoint that causes a high O<sub>2</sub> alarm. The setpoint is variable from 0.5% O<sub>2</sub> above setpoint to 5.0% O<sub>2</sub> above setpoint. The factory setting is 1.0% O<sub>2</sub> above setpoint.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until O2 HI X.X% is displayed in message display.
3. Press the up/down arrow to change the high O<sub>2</sub> alarm limit.
4. Press ENTER to save the high O<sub>2</sub> alarm limit.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## Enabling O<sub>2</sub> Alarms to Trip Contacts

High and Low O<sub>2</sub> alarms can be programmed to trip the remote alarm contacts. A setting of ON will cause this, a setting of OFF will not allow the O<sub>2</sub> alarms to trip the contacts. The factory setting is ON.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until O<sub>2</sub> RLY XXX is displayed in message display.
3. Press the up/down arrow to toggle the setting ON/OFF.
4. Press ENTER to save the setting.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

## Enabling Temp/RH to be Displayed

On units that are equipped with the RH option, the upper seven-segment display on the control panel can be configured to display Temp continuously, RH continuously or toggle between Temp and RH. If the units does not have RH, the upper display will always display temperature. If temperature is set to ON and the RH is set OFF, temperature will be displayed continuously. If temperature is set to OFF and RH is set to ON, RH will be displayed continuously. If both are turned ON, the display will toggle between the two. The factory setting will default to toggle mode if the RH option is present.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until DISP TMP XXX or DISP RH XXX is displayed in the message display.
3. Press the up/down arrow to toggle the setting ON/OFF.
4. Press ENTER to save the setting.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

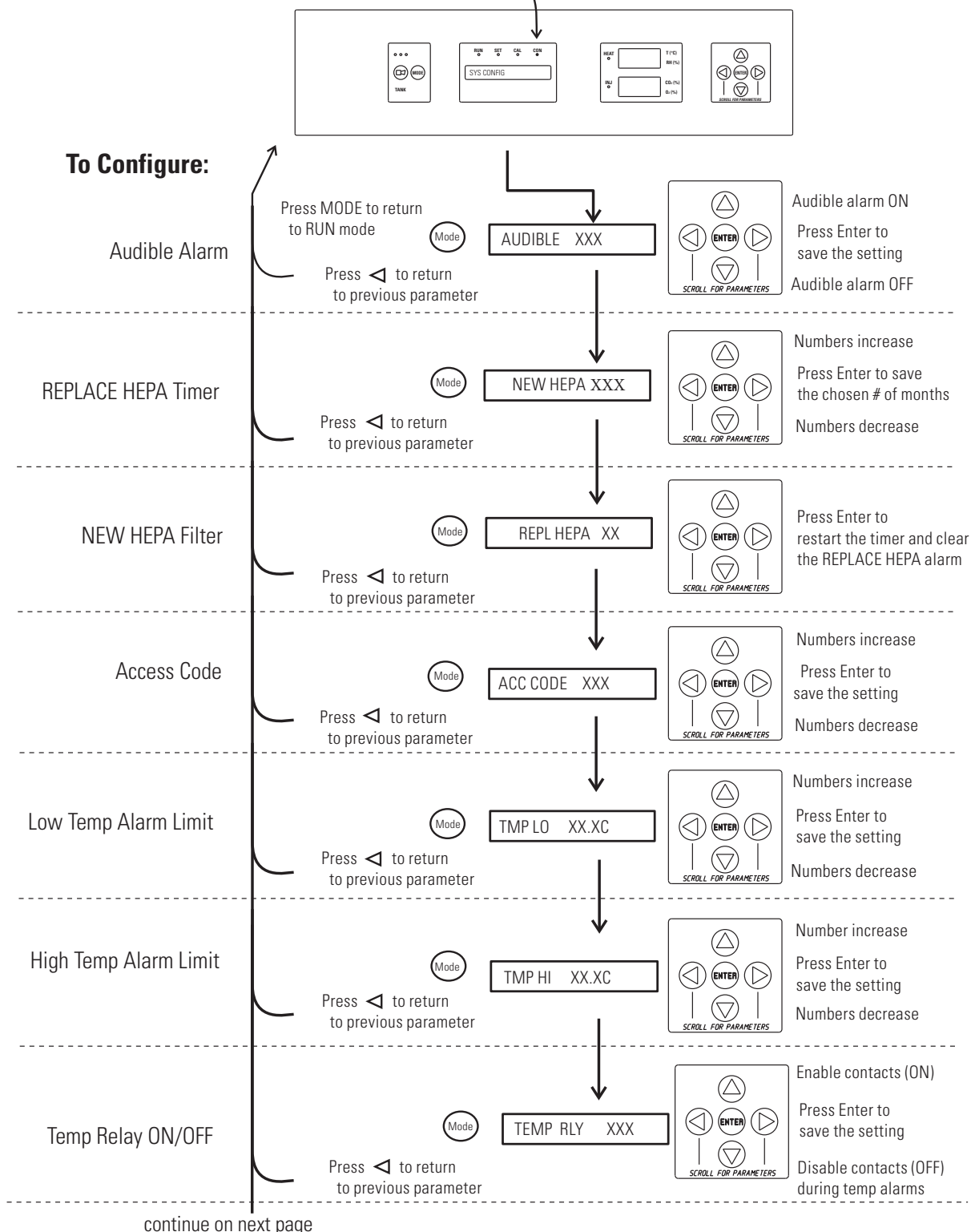
## Enabling CO<sub>2</sub>/O<sub>2</sub> to be Displayed

On models that are equipped with the O<sub>2</sub> system, the lower seven-segment display on the control panel can be configured to display CO<sub>2</sub> continuously, O<sub>2</sub> continuously or toggle between CO<sub>2</sub> and O<sub>2</sub>. If the unit does not have O<sub>2</sub>, the lower display will always display CO<sub>2</sub>. If CO<sub>2</sub> is set to ON and the O<sub>2</sub> is set OFF, CO<sub>2</sub> will be displayed continuously. If CO<sub>2</sub> is set to OFF and O<sub>2</sub> is set to ON, O<sub>2</sub> will be displayed continuously. If both are turned ON, the display will toggle between the two. The factory setting will default to toggle mode if the O<sub>2</sub> system is present.

1. Press the MODE key until the CON indicator lights.
2. Press the right arrow until DISP CO2 XXX or DISP O2 XXX is displayed in the message display.
3. Press the up/down arrow to toggle the setpoint.
4. Press ENTER to save the setpoint.
5. Press the MODE key to return to RUN mode or right/left to go to next/previous parameter.

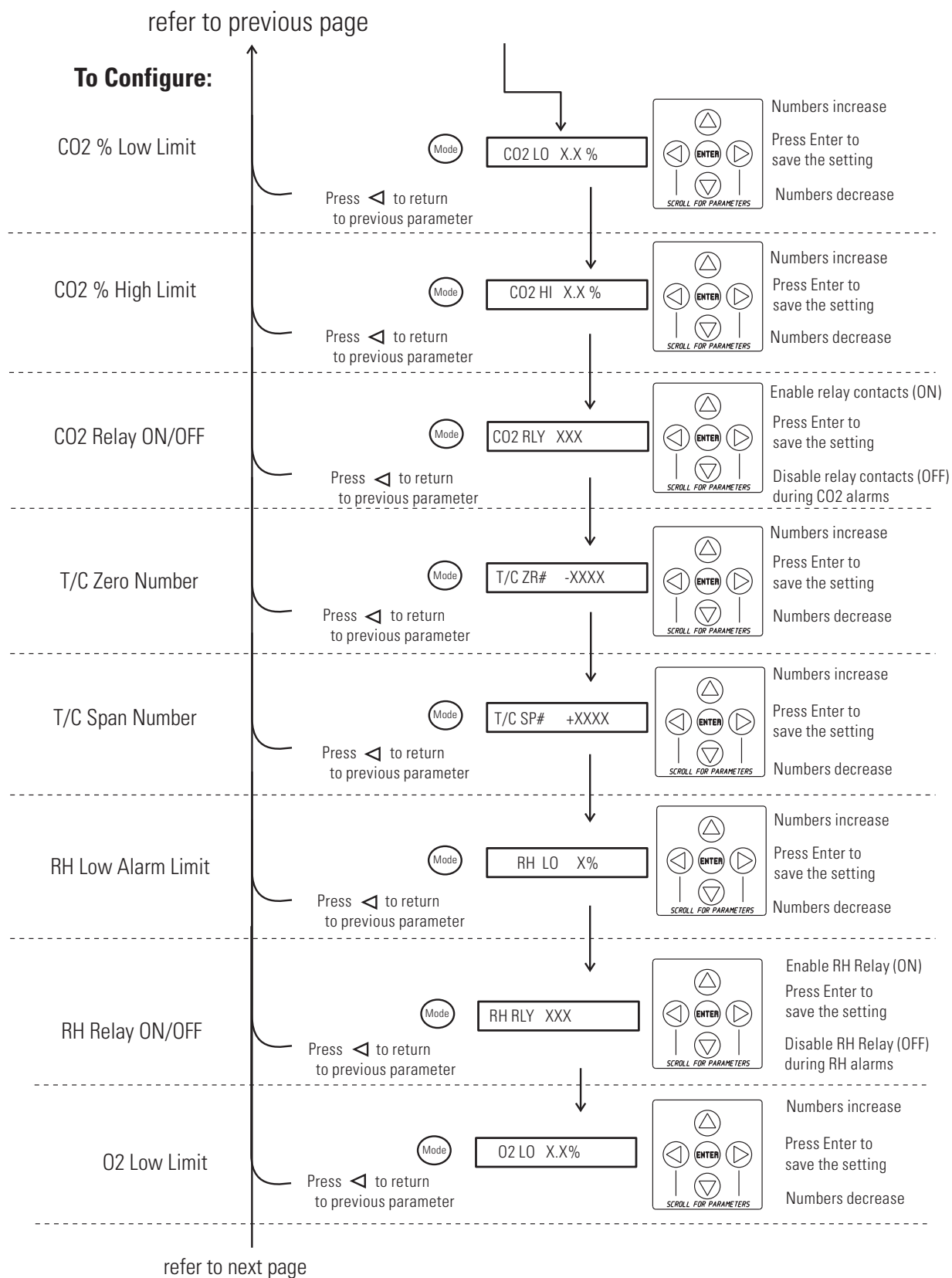
### Chart 3-1. Configuration Mode Page 1 of 3

**Press MODE to light CON:**

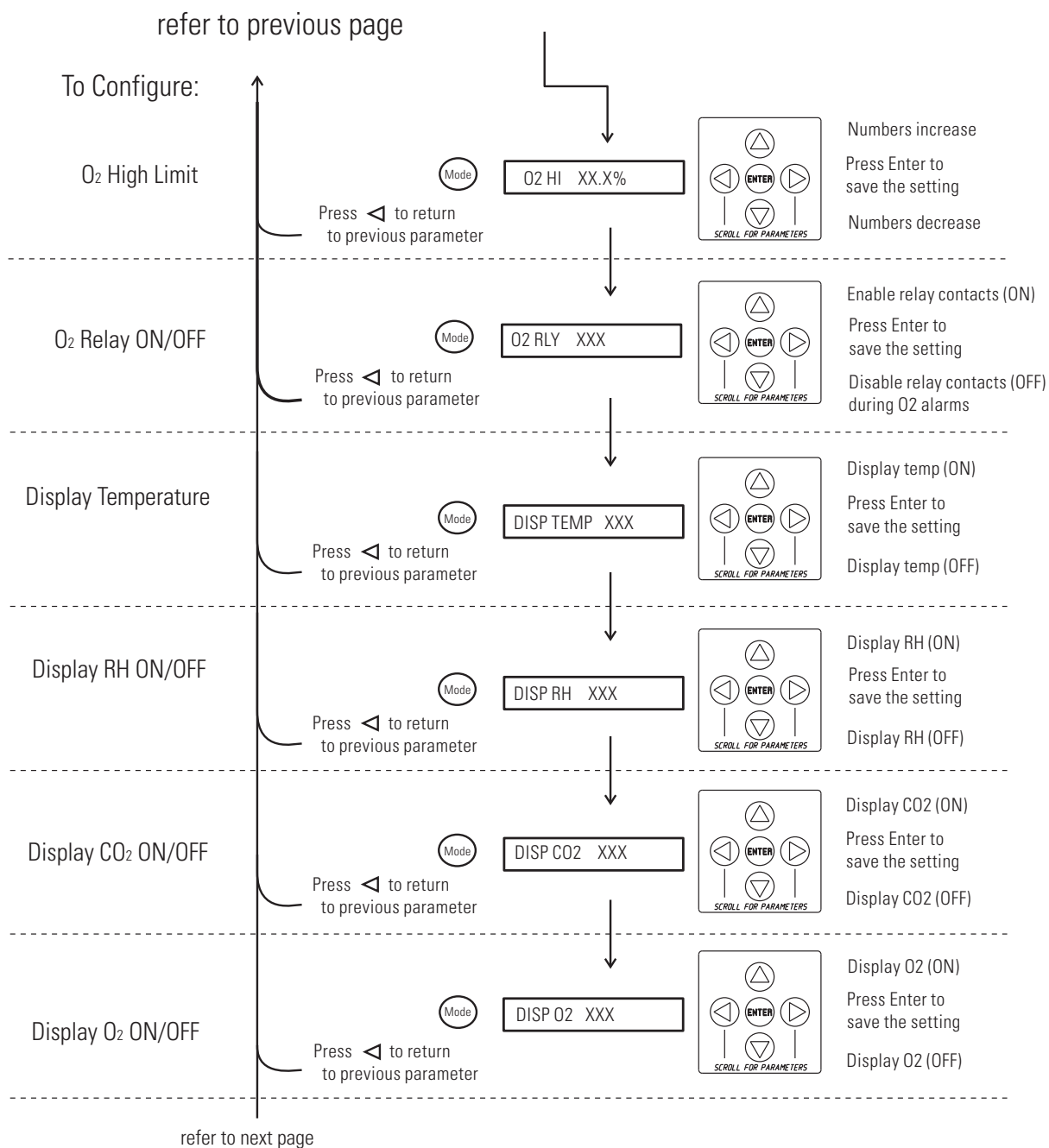




## Chart 3-1. Configuration Mode Page 2 of 3



## Chart 3-1. Configuration Page 3 of 3





## Section 4 Alarms

The Series 8000WJ incubator alarm system is shown in the table below. When an alarm is active, the message appears in the LED message display. Pressing Silence disables the audible alarm for the ringback period. However, the visual alarm continues until the incubator returns to a normal condition. The alarms are momentary alarms only. When an alarm condition occurs and then returns to normal, the incubator automatically clears the alarm condition and the display.

**Table 4-1.** Alarms

Description	Message	Delay	Ringback	Relay
No alarm condition exists	SYSTEM OK/CLASS 100	---	---	---
CO2 System Auto Calibrating	CO2 AUTO CAL	---	---	---
Temp > Otemp Setpoint	SYS IN OTEMP	0 min	15 min	Yes
Temp Control Sensor Fault (See Sect 4)	TSNSR1 ERR	0 min	15 min	No
Over Temp Sensor Fault (See Sect 4)	TSNSR2 ERR	0 min	15 min	No
CO2 Sensor Fault (See Sect 4)	CO2 SNSR ERR	0 min	15 min	No
O2 Sensor Fault (O2 option, see Sect 4)	O2 SNSR ERR	0 min	15 min	No
O2 Sensor Low (O2 option, see Sect 4)	REPL O2 SNSR	0 min	*	No
CO2 Sensor cannot be calibrated (IR option, Sect. 4)	REPL IR SNSR	0 min	15 min	No
Replace HEPA filter reminder – set time expired (See Sect. 3)	REPLACE HEPA	0 min	*	No
Water low in jacket	ADD WATER	0 min	15 min	No
Inner Door is Open	DOOR IS OPEN	15 min	15 min	No
CO2 > CO2 High Tracking Alarm	CO2 IS HIGH	15 min	15 min	Programmable
CO2 < CO2 Low Tracking Alarm	CO2 IS LOW	15 min	15 min	Programmable
TEMP > TEMP High Tracking Alarm	TEMP IS HIGH	0 min	15 min	Programmable
TEMP < TEMP Low Tracking Alarm	TEMP IS LOW	15 min	15 min	Programmable
O2 > O2 High Tracking Alarm (O2 option)	O2 IS HIGH	15 min	15 min	Programmable
O2 < O2 Low Tracking Alarm (O2 option)	O2 IS LOW	15 min	15 min	Programmable
RH < RH Low Limit Alarm (RH option)	RH IS LOW	30 min	15 min	Programmable
CO2 Auto-Zero Fault (IR option, see Sect. 4)	IR AUTOZ ERR	0 min	15 min	No
- All alarm delays and ringback times are +/- 30 seconds - *Visual alarmsonly				

When multiple alarm conditions occur, active messages are displayed in the message display one at a time, updating at 5 second intervals. Pressing Silence during multiple alarms causes all active alarms to be silenced and to ring back in 15 minutes.

The temperature alarms are disabled when Temp setpoint is 10°C. The CO<sub>2</sub> alarms are disabled when CO<sub>2</sub> setpoint is 0.0%. The O<sub>2</sub> alarms are disabled when O<sub>2</sub> setpoint is 21.0%. The RH alarm is disabled when RH setpoint is 0%.

## **Inner Door Open Alarm**

When the inner door is opened on a Series 8000WJ incubator, heat and CO<sub>2</sub> injection are disabled. The door must be securely latched for heat and CO<sub>2</sub> injection to resume after a door opening. If the door is latched yet the display still shows DOOR IS OPEN, or if the door is unlatched and the display does not show DOOR IS OPEN, the door switch could be faulty. Call Technical Services.

## **Sensor Fault Alarms**

The microprocessor in the incubator continually scans all available sensors, except the O<sub>2</sub> (see Section 4), to ensure proper operation. Should an error be detected, the incubator sounds an alarm and displays the appropriate message. If such an alarm occurs, contact your local distributor or the Technical Services department.

### **TSNSR1 ERR**

TSNSR1 ERR If the microprocessor determines that the temperature control precision thermistor is under-range or over-range, TSNSR1 ERR appears in the display, the visual alarm flashes, the audible alarm sounds, and the chamber and door heaters are disabled.

### **TSNSR2 ERR**

TSNSR2 ERR If the microprocessor determines that the over temperature precision thermistor is under-range or over-range, TSNSR2 ERR appears in the display, the visual alarm flashes, the audible alarm sounds. If the independent over temperature protection circuit determines that the thermistor is over-range, SYS IN OTEMP alternates in the display and the chamber and door heaters are disabled.

## **REPL O2 SNSR**

On units equipped with the O<sub>2</sub> system, the microprocessor checks the remaining life of the O<sub>2</sub> sensor whenever O<sub>2</sub> calibration @ 20.7% is performed. After the O<sub>2</sub> sensor declines to less than 43% (typ) of its life expectancy, REPL O<sub>2</sub> SNSR appears in the display and the visual alarm flashes. This alarm alerts the user to replace the O<sub>2</sub> sensor at the earliest convenience (Section 5). The unit will continue to function for some length of time.

## **O2 SNSR ERR**

After the O<sub>2</sub> sensor declines to the point that control cannot be accurately performed, an O<sub>2</sub> SNSR ERR alarm appears in the display, the visual alarm flashes, the audible alarm sounds, and the chamber and O<sub>2</sub> control is disabled. The O<sub>2</sub> control and alarms can be disabled by changing the setpoint to 21.0% (see Section 1) or the O<sub>2</sub> sensor fuel cell must be replaced (see Section 5).

## **CO2 SNSR ERR**

If the cables or connectors between the main microprocessor board and the CO<sub>2</sub> sensor, or between the CO<sub>2</sub> board and the sensor head on I/R CO<sub>2</sub> units become loose or disconnected, the CO<sub>2</sub> SNSR ERR alarm occurs.

## **I/R Units**

On I/R incubators, the red light on the I/R module (see Figure 5-4 for the module circuit board location) is lit continuously if communication is lost between the CO<sub>2</sub> board and the sensor head. The CO<sub>2</sub> display is locked at 00.0 without injection. Turning the incubator off and on will not clear the alarm. Only proper connections of all the components corrects the alarm.

## **IR AUTOZ ERR**

On incubators equipped with an IR CO<sub>2</sub> control system, calibration is performed automatically once every 24 hours using the Auto Zero feature. If a CO<sub>2</sub> correction of more than 0.45% is detected then the Auto Zero cycle is aborted and the IR AUTOZ ERR appears in the display, the visual alarm flashes, and the audible alarm sounds. The unit will continue to function using the last valid CO<sub>2</sub> correction value. The error may be cleared by performing the IR CO<sub>2</sub> control system calibration (see Section 2) or waiting 24 hours for the next Auto Zero cycle.

## **IR AUTOZ ERR**

Possible problems which will cause this alarm are:

- Auto Zero pump, orifice, filter or tubing will not allow air to the sensor.

Possibilities are:

- ~ Defective or electrically disconnected air pump
  - ~ Kinked auto zero vinyl tubing
  - ~ Disconnected tubing between air pump and sensor
  - ~ Plugged filter or orifice on auto zero assembly
  - ~ Defective auto zero circuit
- Cabinet temperature has been increased significantly from a previous setpoint. (For example, the unit was calibrated and operating at 30°C and the setpoint is increased to 50°C). In this instance, calibrating the CO<sub>2</sub> will correct the alarm.
  - There is high background CO<sub>2</sub> in the laboratory. This could be from leaks in the tank, regulator or vinyl CO<sub>2</sub> tubing.
  - High CO<sub>2</sub> sensor calibration drift occurred. This requires replacement of the sensor.

**Warning** Do not disassemble the IR CO<sub>2</sub> sensor as there are no user-serviceable adjustments. s

## **REPL IR SNSR**

If the Auto Zero feature cannot properly calibrate the sensor, then REPL IR SNSR appears in the display, the visual alarm flashes, the audible alarm sounds, and the chamber CO<sub>2</sub> control disabled. The CO<sub>2</sub> control and alarms can be disabled by changing the setpoint to 0.0% (see Section 1) or replacing the IR CO<sub>2</sub> sensor. Call Technical Services.

## Section 5 Routine Maintenance

**Caution** Before using any cleaning or decontamination method except those recommended by the manufacturer, users must check with the manufacturer that the proposed method will not damage the equipment. s

**Warning** It is the responsibility of the user to immediately clean up after all accidental spills of hazardous materials. Be certain to follow local EHS policies with regards to personal protective equipment, cleaning, and disposal. s

### Cleaning the Interior

Use an appropriate disinfectant, such as quaternary ammonium. All articles and surfaces must be thoroughly cleaned, rinsed and rough-dried.

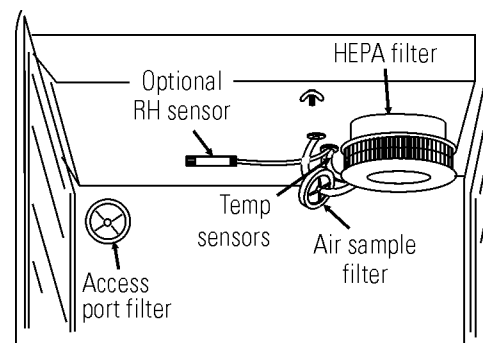
**Warning** Alcohol, even a 70% solution, is volatile and flammable. Use it only in a well ventilated area that is free from open flame. If any component is cleaned with alcohol, do not expose the component to open flame or other possible hazard. Allow the alcohol to fully dry before turning power on. s

**Warning** Do not spray the T/C sensor with flammable solutions. The internal temperature of the CO<sub>2</sub> sensor is approximately 150°C when the unit is in operation. Allow sufficient time for the sensor to cool before cleaning. s

**Caution** Do not use strong alkaline or caustic agents. Stainless steel is corrosion resistant, not corrosion proof. Do not use solutions of sodium hypochlorite (bleach), as they may also cause pitting and rusting. s

**Note** If condensation is visible on the ceiling of the chamber, yet the message center does not display ADD WATER, then the liquid level switch could be faulty. Call Technical Services.

**Warning** If the unit has been in service, disconnect the power cord from both the unit and the power source, turn off all gas regulators, and allow the unit to cool before disinfecting. s

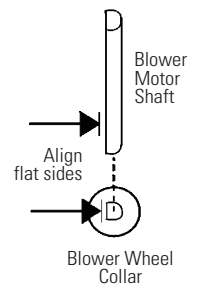


**Figure 5-1.** Sensor and Filter Locations



## Cleaning the Interior (continued)

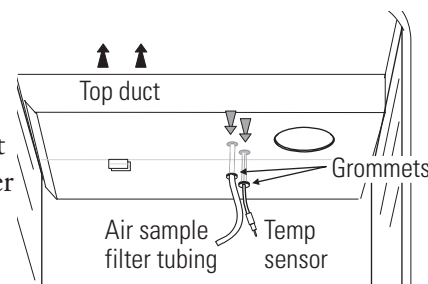
1. Remove the shelves, access port filter and side duct sheets. Remove the temperature sensors and the air sample filter tubing from the back of the blower scroll. If unit is equipped with the optional RH sensor, unfasten it from the clip on the top duct. See Figure 5-1.
2. Remove the filter from the air sample filter tubing. Carefully pull down and remove the HEPA filter.
3. Remove the wingnuts securing the top duct to the interior. Carefully slide the top duct down and off the temperature sensor, air sample filter tubing (and RH sensor, if applicable).
4. Wash the shelves, ducts, wingnuts and stopper with disinfectant and rinse with sterile water.  
**Option:** Autoclave shelves, ducts and wingnuts.
5. Remove the blower scroll by first pushing the black lever clip closest to you toward the scroll. Then turn the scroll to the right to disengage it from the blower scroll plate. Some manipulation may be required as the alignment holes are keyhole-shaped.
6. Remove the remaining wingnut, then pull down on the blower wheel. If a new wheel and scroll are to be used, discard the old ones per local EHS policies and procedures. If the old ones are to be reused, wash all parts with disinfectant and rinse with sterile water.
7. Remove the blower scroll plate by first pushing the black lever clip toward the chamber ceiling. Then turn the plate to the left to disengage it from the alignment keyholes. Clean as above, or autoclave.
8. Wash the cabinet interior with disinfectant starting at the top and working down. Wash the inner door both inside and out. The cabinet and door must be rinsed with sterile water until the disinfectant has been removed. After the cabinet has been rinsed, spray with 70% alcohol.
9. Reinstall the blower scroll plate by aligning it with the larger end of the keyholes and turning it to the left to lock it on. Pull the black lever clip downward from the ceiling.
10. Install the blower wheel onto the motor shaft, aligning the d-shaped flat sides of each. See Figure 5-2. Secure the blower wheel with the wingnut. Make sure the wheel turns freely.



**Figure 5-2.** Align

## Cleaning the Interior (continued)

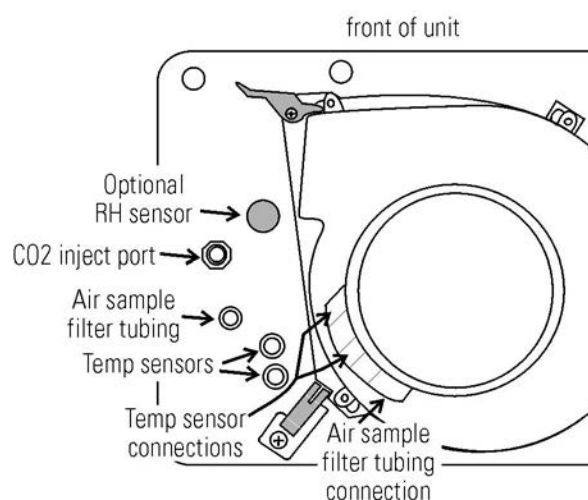
11. Locate the blower scroll over the blower wheel into the larger end of the keyholes on the scroll plate. Turn the scroll to the right to lock it into the keyholes. Pull the black lever clip closest to you toward the front of the unit.



**Figure 5-3.** Feed Through Holes

12. Install the top duct by feeding the temperature sensors, air sample tubing (and RH sensor, if applicable) through the appropriate holes in the duct as it is raised to the top of the chamber. Be careful not to pull the grommets through the duct. See Figure 5-3.
13. Locate the mounting studs and blower scroll into the appropriate holes in the top duct and install the wingnuts to secure the duct.
14. Install the air sample filter onto the top duct tubing.

15. Carefully pull the temperature sensors and air sample filter tubing down until they can be inserted approximately 1 inch into the appropriate holes in the back of the blower scroll. If applicable, place the optional RH sensor into the corresponding clip on the top duct. See Figure 5-4.



**Figure 5-4.** Underside View of Chamber Ceiling

16. Install the HEPA Filter.
17. Wipe the interior of the incubator, including the glass door and gasket, with 70% ethanol.

**Caution** Do not spray or wipe sensors.

18. Install the left and right ducts, and the access port stopper with filter, spraying each with 70% alcohol (do not saturate).
19. Install the shelves and spray with 70% alcohol.
20. Plug the incubator in and turn the power switch on. Allow the unit to run empty (heat only) for 24 hours before returning to service.

## Cleaning Cabinet Exterior

1. Use a lint-free cloth dampened in mild soapy water to clean the exterior, especially the door handle. Do not use any liquids (including spray cleaners) on the control panel; instead clean with a dry microfiber cloth.
2. Wipe the surfaces (except control panel) twice using a clean cloth dampened in clear water.
3. Wipe dry with a clean cloth.

## Cleaning Humidity Pan

Clean the humidity pan with soap and water and a general use laboratory disinfectant, such as quaternary ammonium. Rinse with sterile water and spray with 70% alcohol. The humidity pan may be autoclaved.

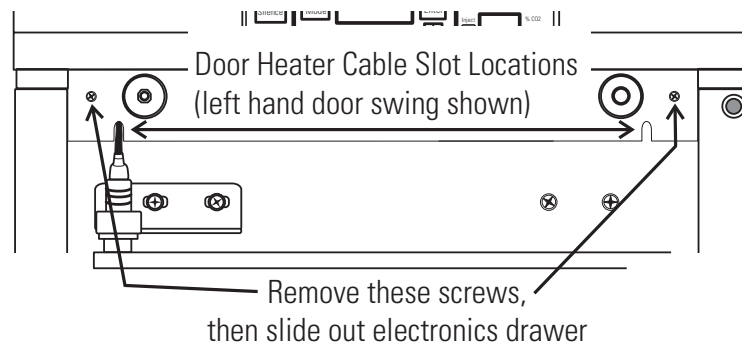
## Replacing Fuses

**Warning** The electronics drawer contains hazardous voltages. Replacing the fuses should be performed by qualified personnel only. If the unit has been in service, disconnect the power cord from both the unit and the power source, turn off all gas regulators, and disconnect all tubing and any other connections from the rear of the electronics drawer. s



There are three fuses in the incubator that can be replaced.

1. To replace a fuse, first open the exterior cabinet door.
2. Remove two screws as shown. See Figure 5-5.
3. Carefully slide out electronics drawer.



**Figure 5-5.** Screw Locations

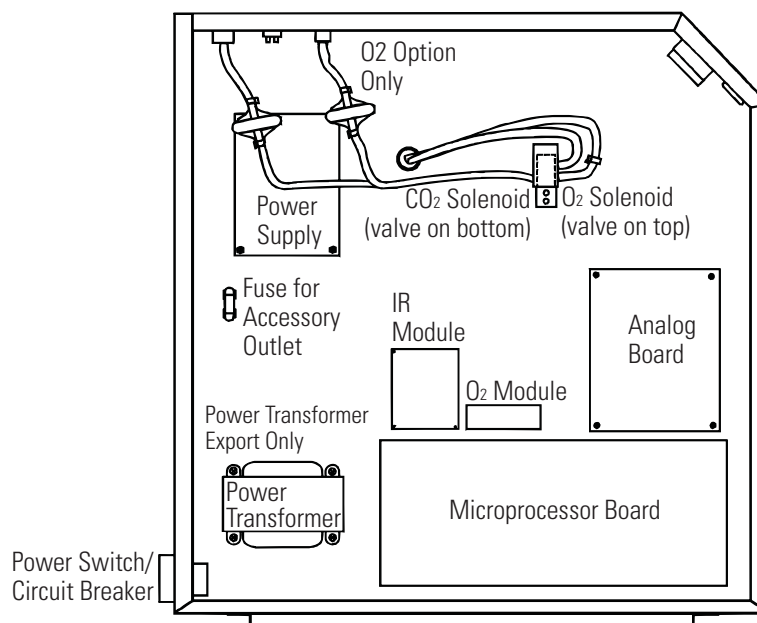
## Replacing Fuses (continued)

There are two fuses on the main microboard labeled F1 24 VAC Door Heater and F4 24 VAC Collar Heater. Refer to Figure 5-6 for the location of the main microboard. Remove the fuse and replace it with a new one of the same specification. Refer to Table 5-1.

**Table 5-1.** Fuses and Specifications

Fuse #	Manufacturer Part #	Amperage Rating	Rupture Speed	IEC Letter Code
F1	BUSS GMC-3.5A	3.5 Amp	Time-Lag	T
F4	BUSS GMC-2.5A	2.5 Amp	Time-Lag	T
115 VAC ACC	BUSS GMC-1.0A	1.0 Amp	Time-Lag	T
230 VAC ACC	BUSS BK-GMC-500mA	0.5 Amp	Time-Lag	T

The other replaceable fuse is the accessory outlet fuse mounted to the floor of the electronics drawer. To locate the fuse, refer to Figure 5-6. Remove the fuse and replace with a new one of the same specification.



**Figure 5-6.** Electronics Drawer

## **Replacing Fuses (continued)**

1. When the fuse has been replaced, slide the electronics drawer back in, being very careful to place the door heater cable back into the provided slot so that the drawer does not pinch the cable (Figure 5-5).
2. Replace the two screws removed earlier.
3. Close the exterior door.
4. Plug the power cord back into the dedicated, grounded circuit.
5. Turn on power switch. If the unit operates properly, it may now be returned to service.

There is one fuse in the incubator that is not replaceable. This fuse is intended for catastrophic failure only and is located on the power supply circuit board in the electronics drawer. See Figure 5-6. If this fuse is blown, the power supply must be replaced. Contact the factory for more information.

## **HEPA Filter Maintenance**

Replace the HEPA filter when the REPLACE HEPA reminder is displayed by following the Installing HEPA Filter procedures. Carefully pull down the HEPA filter to remove, and discard it per local EHS policies and procedures. Afterwards, clear the visible alarm and display, and reset the timer by following the NEW HEPA Filter procedures.

The REPLACE HEPA reminder can be set to alarm after a specified time from 1 to 12 months. The reminder default is the factory recommended setting of 6 months. For details, see Section 3.

## **Replacing Air Sample Filter**

The air sample filter should be replaced whenever the HEPA filter is replaced. On the inside of the chamber, inserted into the back of the blower scroll, is the air sample filter and its connecting tubing. Disengage the tubing from the back of the scroll, then remove the filter from the tubing and discard. Install the new filter. Connect it securely to the air sample filter tubing, then insert the tubing into the back of the blower scroll.

## **Replacing Access Port Filter**

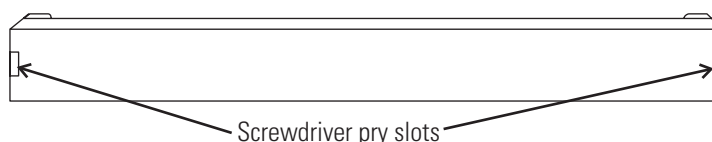
The access port filter should be replaced whenever the HEPA filter is changed. The filter is connected to the stopper in the upper left corner of the chamber back wall. Remove the filter from the connecting tube and discard it per local EHS policies and procedures. Install the new filter.

## Draining Water Jacket

**Note** Although the rust inhibitor is biodegradable, be certain to follow local EHS policies with regards to personal protective equipment, cleaning, and disposal.

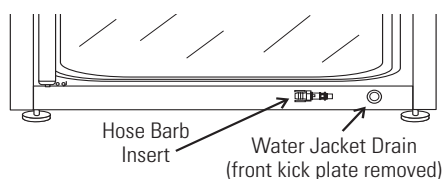
**Caution** For stacked incubators, completely drain the top unit before draining the bottom unit in order to maintain the lowest possible center of gravity.

1. Turn the unit off. Remove the plug from the power source.
2. Remove the front cover plate below the door. There are small flatblade screwdriver pry slots on each end of the plate to help remove it. See Figure 5-7.



**Figure 5-7.** Front Cover Plate Below the Door

3. If present, remove the drain plug.
4. Connect the hose barb insert to the drain on lower front of the water jacket and to the drain hose.
5. After water jacket has finished draining, remove the hose barb insert and secure on the front of the unit. See Figures 5-8 and 5-9.



**Figure 5-8.** Water Jacket Drain and Hose Barb Insert



**Figure 5-9.** Hose Barb Insert

**Warning** Once the jacket is empty, be certain to clean up any water that may have spilled onto the floor during this process to prevent a possible slip hazard.

6. Install cover plate.
7. To fill the water jacket, see Section 1. Be sure to add the rust inhibitor to the water when filling. Instructions follow in this section.

## Adding /Replenishing Rust Inhibitor

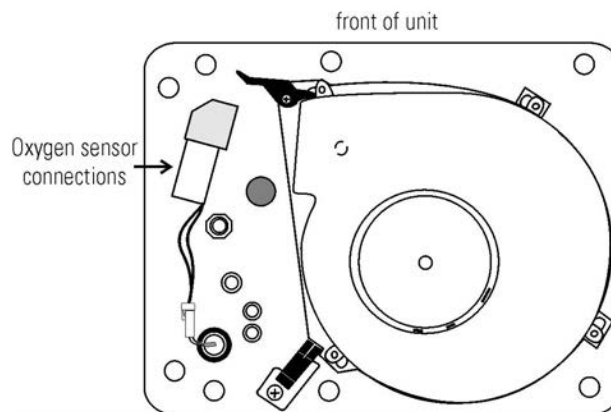
Series 8000WJ incubators are shipped from the factory with a rust inhibitor added to 0.2 gallons (0.8 liters) of the water in the jacket. The rust inhibitor must be replenished every 2 years. Mix 1 bag/bottle of the rust inhibitor with a gallon (3.8 liters) of water as described in Filling Water Jacket, Section 1. Drain 1.5 gallons (5.7 liters) of water from the jacket per Draining Water Jacket instructions above. If the message center does not display ADD WATER, then the liquid level switch could be faulty. Call Technical Services. Add the rust inhibitor mixture and continue to fill the jacket with water per Filling Water Jacket, Section 1.

Rust Inhibitor	0.5 lb. (225g) bag	1900100
----------------	--------------------	---------

## O<sub>2</sub> Sensor Fuel Cell - Models 3424, 3425, 3426, 3427

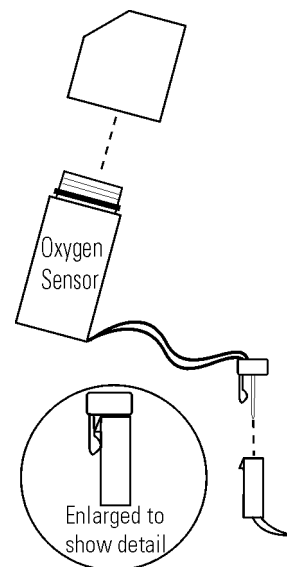
**Warning** Before removing the old O<sub>2</sub> sensor fuel cell, inspect it for any visual signs of damage. If the liquid electrolyte is leaking, follow local EHS policies with regards to personal protective equipment, cleaning and disposal. s

The O<sub>2</sub> sensor is located on the blower scroll plate in the chamber of the unit. To replace it, refer to Figure 5-10 and 5-12 and follow steps below.



**Figure 5-10.** Underside View of Top of Chamber

**Warning** If the unit has been in service, disconnect the power cord from both the unit and the power source, turn off all gas regulators and allow the unit to cool before replacing the O<sub>2</sub> sensor fuel cell. s



**Figure 5-11.** Detail

1. Remove the top duct by removing 2 wingnuts.
2. Locate the sensor on the scroll plate.

## O<sub>2</sub> Sensor Fuel Cell (continued)

**Warning** Handle the O<sub>2</sub> sensor fuel cell very carefully as it can be easily damaged. Do not install the O<sub>2</sub> Before installing the new O<sub>2</sub> sensor fuel cell, inspect it for any visual signs of damage. If the liquid electrolyte is leaking, handle per local EHS policies with regards to personal protective equipment, cleaning, and disposal. s

**Warning** Handle the O<sub>2</sub> sensor fuel cell if its shipping box is damaged or it has fallen on the floor as mechanical shock or vibration may cause irreversible internal structural damage to the O<sub>2</sub> sensor fuel cell. Do not disassemble the O<sub>2</sub> sensor fuel cell as there are no user-serviceable adjustments. If the liquid electrolyte leaks due to sensor breakage, care should be taken in handling the sensor as the solution contains lead acetate which is harmful to humans. s

**Warning** Contact with this liquid should be avoided. In case the liquid electrolyte contacts the skin or clothing, wash with soapy water and rinse generously with plain water. If the liquid electrolyte contacts the eye, flush with water for at least 15 minutes and obtain immediate medical assistance. In case of breathing in of the electrolyte, flush the nasal cavity thoroughly with water and seek immediate medical assistance. If the electrolyte is swallowed, rinse the mouth thoroughly with water and seek immediate medical assistance. s

**Warning** Before installing the new O<sub>2</sub> sensor fuel cell, inspect it for any visual signs of damage. If the liquid electrolyte is leaking, handle per local EHS policies with regards to personal protective equipment, cleaning, and disposal. s

3. Lift up slightly on the tab securing the sensor wire terminal connection. Be careful not to break the locking tab off. See enlarged detail of Figure 5-11.
4. Disengage the connectors from each other.
5. The O<sub>2</sub> sensor fuel cell is screwed into the brass fitting. Unscrew the old sensor and discard per local EHS policies and procedures.
6. Screw the new O<sub>2</sub> sensor fuel cell in firmly. Do not cross-thread!
7. Orient the connectors as shown in Figure 5-11 and connect.
8. Install the top duct and tighten the 2 wingnuts.
9. Plug the unit in. Calibrate the O<sub>2</sub> system using the O<sub>2</sub> CAL@20.7% method as described in Section 2.
10. Allow the unit to run until the temperature stabilizes. Check the O<sub>2</sub> and CO<sub>2</sub> operation and return the unit to service.

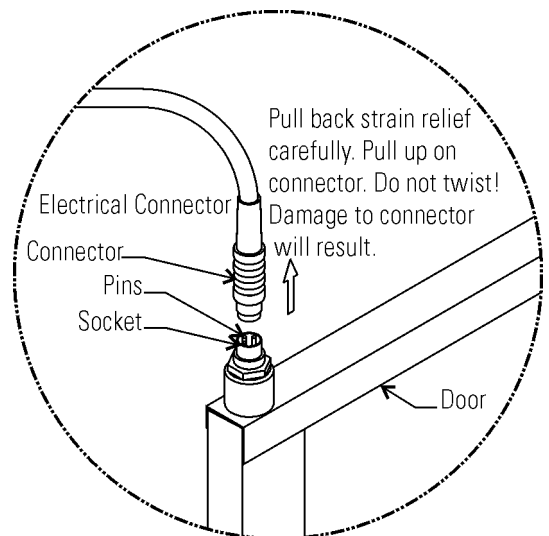


## Reversing the Door Swing

For side-by-side operation or changing lab layouts, the inner and outer doors are field-reversible. The procedure is written from the prospective of changing the door swing from a left-hand to a right-hand swing. All screw holes are pre-drilled for reversing the door. The tools required are a Phillips and a flatblade screwdriver.

**Warning** The electronics drawer contains hazardous voltages. If the unit has been in service, disconnect the power cord from both the unit and the power source, turn off all gas regulators, disconnect all tubing and any other connections from the rear of the electronics drawer, and allow the unit to cool before reversing the door swing. s

1. Referring to Illustration A following, open the outer door and remove it by lifting it off the hinge pins. Lay this door down on its face on a padded surface to prevent scratches.
2. Remove the two outer door hinges identified in the illustration.
3. Remove the four nylon screws from what will be the new locations for the inner door hinges.
4. With the Phillips screwdriver, remove and save the four nylon screws from the outer door hinge mounting holes on the right side of the door frame.
5. Remove the electrical connector on the top of the inner door hinge by carefully prying up the black strain relief.
6. Refer to Figure 5-12.  
The heater wiring connector is of yellow rubber, visible when the strain relief is moved upward as shown. See connector manufacturer's instructions in Section 11 of this manual.
7. Grasp the upper portion of the yellow connector and pull it up and out of the hinge socket.



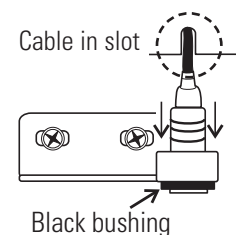
**Figure 5-12.** Connector

## Reversing the Hinges for the Exterior Door (cont.)

**Caution** The frame along the hinge side of the inner glass door has electrical connectors mounted on the frame at both the top and bottom of the door. Be careful not to damage the connectors by resting the weight of the door on them. Place the inner glass door on a flat surface where it cannot be damaged.

8. Remove the upper inner door hinge. This hinge will be inverted and become the lower inner door hinge when the hardware is reassembled. At this time, remove the 5/8" black plastic bushing from the hinge. This bushing will be inserted in the top hinge when the door is installed on the right side of the cabinet.

9. Remove the lower inner door hinge. This hinge will be inverted and become the upper inner door hinge when the hardware is reassembled. Also, remove the 1" long white plastic bushing from the center of the hinge and insert it into the hinge removed in step above.



**Figure 5-13.** Bushing

10. The shoulder of the bushing must be on top as shown in Illustration B. Install the black plastic bushing into the hinge just removed. See Figure 5-12.
11. Remove the two nylon screws opposite the door latch.
12. Remove the door strike plate and install the two nylon screws just removed (Step 11) into the vacant holes.
13. Remove the two outer Phillips screws on frame just beneath the control panel. Remove only the two outer screws and pull the electronics drawer out about one inch.

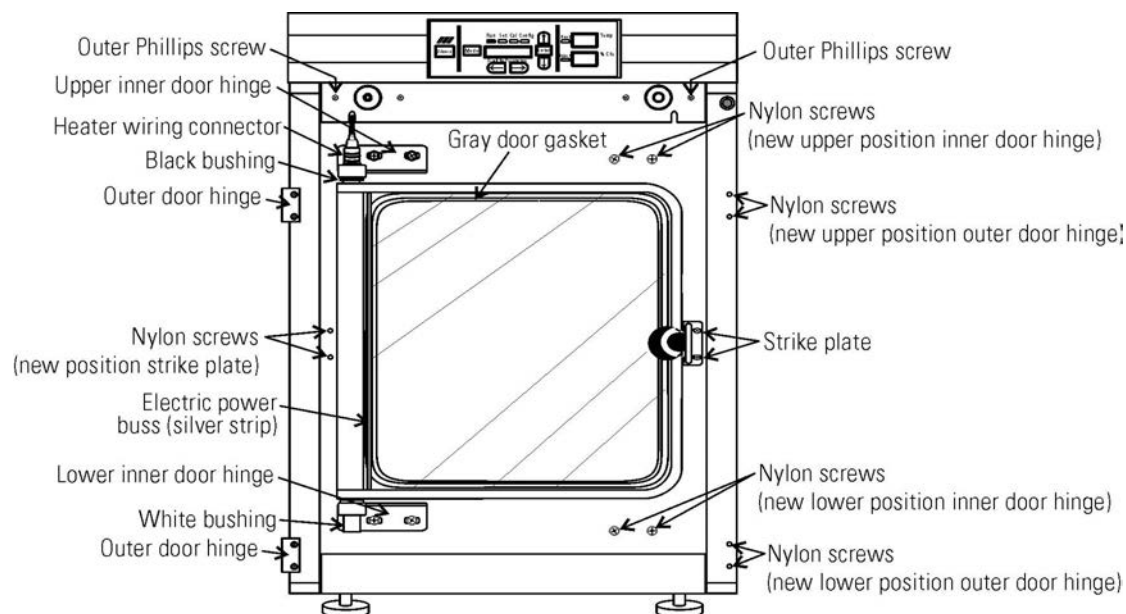
Refer to Illustration B.

14. Move the door heater cable to the slot on the right side of the electronics drawer.
15. Push any excess cable into the drawer, making sure the cable will not be damaged when the drawer is pushed back in and secured.
16. Install the lower inner door hinge in the drawing. (This was the upper inner door hinge removed in Step 8 and should have the white plastic bushing installed). Do not completely tighten the screws.
17. Insert the inner door into the lower hinge with the latch to the left. Align the door to chamber opening and place upper hinge in position (this was the lower inner door hinge removed in Step 9 with the black plastic bushing installed).
18. Insert the screws into the hinge but do not completely tighten them.

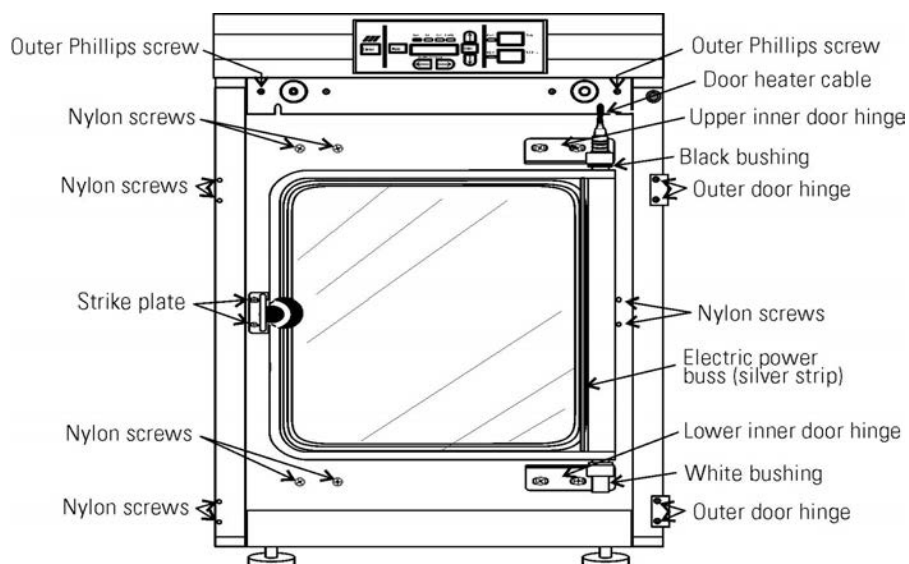
## **Reversing Hinges for the Exterior Door (continued)**

19. Match the pins on the yellow connector with the socket on the door. Press the connector completely into the socket and slide the strain relief down until no yellow is visible on the plug (Figure 5-13).
20. Slide the electronics drawer back in, being very careful to place the door heater cable back into the provided slot so that the drawer does not pinch the cable. See Figure 5-6. Secure the electronics drawer with the screws removed above.
21. Align the door by checking that the gasket is making contact with the glass door, on the latching side, top and bottom. When you close the door you can feel where the glass door is making contact with the gasket. The gasket should be making contact against the glass door and not overlapping onto the metal frame.
22. When both hinges are in place with the hinge screws still loosened, push up on the bottom hinge. This will shift both hinges and the door upward. Tighten the hinge screws.
23. Attach the strike plate, aligning it so the knob secures the door against the gasket.
24. Attach the outer door hinges.
25. Install the nylon screws in the previously upper and lower outer hinge locations.
26. Install the nylon screws in the previously upper and lower inner hinge locations. Assemble the outer door to the incubator.
27. Plug the power cord back into the dedicated, grounded circuit. Allow the unit to run until the temperature stabilizes. Check for proper unit operation, especially inner door heater and gasket seal. Clean and disinfect the unit before returning to service.

## Reversing Hinges for the Exterior Door (continued)



**Illustration A.** Disassembly



**Illustration B.** Reassembly

## **PREVENTIVE MAINTENANCE**

### **Incubators**

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

The following is a condensed list of preventive maintenance requirements. See the specified section of the operating manual for further details.

We have qualified service technicians, using NIST traceable instruments, available in many areas. For more information on Preventive Maintenance or Extended Warranties, please contact Technical Services.

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

#### **Tips for all incubators:**

- Do NOT use bleach or any disinfectant that has high chlorine content.
- Use water with characteristics as described in Filling Water Jacket.
- Avoid spraying cleaner on the CO<sub>2</sub> sensor.
- Do not use powdered gloves for tissue cultures.

### Preventive Maintenance for Water Jacket Incubators

Action	Daily	Weekly	Monthly	6 Months	Yearly	2 years
Check CO <sub>2</sub> and/or N <sub>2</sub> tank levels (tanks without optional gas guard monitor)	✓					
Inspect the inner and outer door latches and hinges for proper operation and excessive mechanical wear; inspect the line cord insulation and inner door silicone gasket for chips and cracks. Replace as required.					✓	
Check water level in humidity pan, ½" from top. See "Filling the Humidity Pan".		✓				
Check proper blower motor operation (air flow, noise, and vibration)		✓				
Verify and document CO <sub>2</sub> , humidity and temperature calibration, as applicable. See Calibration section.					✓	
Verify O <sub>2</sub> calibration, as applicable. See Calibration section				✓		
Perform a complete decontamination procedure. Wipe down interior, shelves and side ducts with disinfectant. Change or clean blower wheel and scroll. Clean top duct. Clean humidity pan. Rinse everything well with sterile water. See "Cleaning the Interior, "Cleaning the Humidity Pan".	Before first use, after every service call and between experiments to prevent cross contamination  More frequent decontamination may be required, depending on use and environmental conditions					
Change HEPA and gas filters, if applicable (or as needed). See "Installing Access Port Filter, "Installing Air Sample Filter", "Installing HEPA Filter".				✓		
Replenish rust inhibitor in water jacket. See "Adding/Replenishing Rust Inhibitor".						✓
Tubing Maintenance and Inspection: Inspect all tubing and connections & Replace any tubing that shows signs of degradation or damage.				✓		



## Section 6 Factory Installed Options

A description of the factory installed options follows.

### Humidity Readout

All Series 8000WJ incubators can be equipped with a humidity sensor to monitor the relative humidity (RH) inside the chamber. The sensor is mounted to the top air duct and provides a signal that is displayed in 1% increments on the control panel. The humidity readout can be displayed continuously or toggles with the temperature readout. In addition, a low alarm limit can be set on the humidity readout which will detect when the humidity pan runs dry. Refer to Section 3, Configuration.

Factors Affecting Humidity Level in Chamber:

- Water level in the humidity pan
- Frequency of door openings
- Humidity pan location; floor, shelf, in duct
- Flooded incubator chamber bottom
- Air leakage through the gaskets
- Gas sample port capped
- HEPA filter or air flow restrictor plate
- N<sub>2</sub> purge on incubators with O<sub>2</sub> control.
- Humidity levels in O<sub>2</sub> units will be reduced, depending on the amount of N<sub>2</sub> required to control the O<sub>2</sub> level in the chamber.

**Table 6-1.** Typical RH levels

The table at right lists some typical RH levels at different O<sub>2</sub> and CO<sub>2</sub> percentages.

O2%	CO2%	RH% (±5%)
1%	2.5%	55%
2%	5%	60%
5%	10%	75%
10%	10%	80%
21%	5%	95%

Incubators equipped with a Thermal Conductivity CO<sub>2</sub> sensor rely on a constant level of relative humidity in order to accurately measure and control the CO<sub>2</sub> concentration in the incubator.

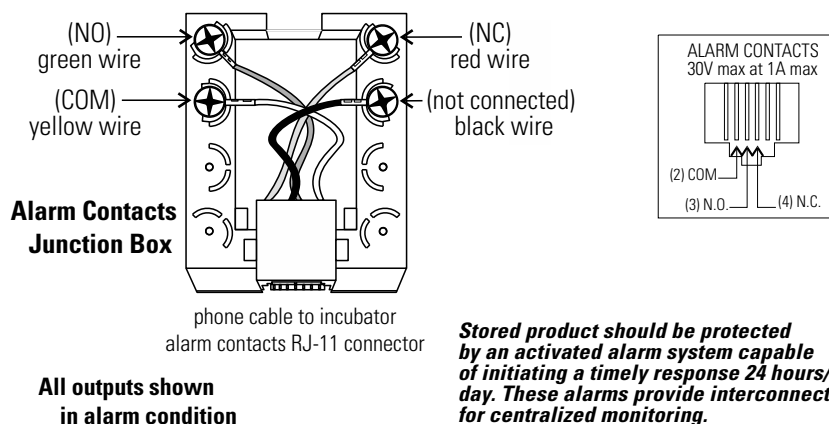
The sensor is capable of measuring relative humidity from 10% to 100% with an accuracy of ±5% above 90%. See Section 2, Calibration for details on calibrating the humidity readout.



## Remote Alarm Contacts

A set of relay contacts are provided to monitor alarms through an RJ-11 telephone style connector on the back of the cabinet. Refer to Figure 1-13 for the location of the alarm connector.

The remote alarm provides a NO (normally open) output, a NC (normally closed) output and COM (common). Refer to Figure 6-1.



**Figure 6-1.** Contacts

The contacts will trip on a power outage or an over temperature condition. The contacts may also be programmed to trip or not trip on temperature alarms, CO<sub>2</sub> alarms, O<sub>2</sub> alarms and RH alarm. See Section 3, Configuration.

## Section 7 Specifications

\*Specifications are based on a nominal voltage of 115 V in an ambient of 22°C to 25°C.

### Temperature

Control	±0.1°C
Range	+5°C above ambient to +55°C (131°F)
Uniformity	±0.2°C @ +37°C
Sensor	Precision thermistor
Tracking Alarm	User programmable high/low

### Over Temperature Protection

Sensor	Precision thermistor
Controller	Independent circuit
Setability	0.1°C
Range	Temp setpoint + 0.5C to 60C
Tracking Alarm	Not user programmable

### CO<sub>2</sub>/O<sub>2</sub>

CO <sub>2</sub> /O <sub>2</sub> Control	Better than ±0.1%
CO <sub>2</sub> Range	0-20%
O <sub>2</sub> Range	1-21%
Inlet Pressure	15 PSIG (103.4 kPa maximum)
CO <sub>2</sub> Sensor	T/C or IR
O <sub>2</sub> Sensor	Fuel Cell
Readability	0.1%
Setability	0.1%
Tracking Alarm	User programmable high/low

### Humidity

RH	Ambient to 95% @ +37°C (98.6°F)
Humidity Pan	0.8 gal. (3 liters) standard
Tracking Alarm	User programmable low only
Optional	Display in 1% increments

### **Fittings**

Fill Port	3/8" barbed
Drain Port	3/8" barbed, quick disconnect
Access Port	1-1/4" (3.18 cm) removable neoprene plug
CO <sub>2</sub> /N <sub>2</sub> Inlets	1/4" hose barbed

### **Unit Heat Load** (typical @ +37°C)

115V / 230V	344 BTUH (100 Watt)
-------------	---------------------

### **HEPA Filter**

Federal Standard 209E, Class 100 (maximum allowable number of particles 0.5 microns and larger per cubic foot of air)\*

Full recovery within 5 minutes of door opening

Replacement Reminder	User programmable
----------------------	-------------------

*\*follow manufacturer's recommendations for regular filter replacement*

### **Shelves**

Dimensions	18.5" x 18.5" (47.0 cm x 47.0 cm)
Construction	Stainless steel, perforated, electropolished
Surface area	2.4 sq. ft. (0.22 sq. m) per shelf
Max. per Chamber	38.4 sq. ft. (3.6 sq. m)
Loading	35 lbs (16kg) slide in and out, 50 lbs (23 kg) stationary
Standard	3 (shipped with unit)
Maximum	16

### **Construction**

Water Jacket Volume	11.7 gal. (44.3 liters) typ., 13.0 gal. (49.2 liters) max.
Interior Volume	6.5 cu. ft. (184.1 liters)
Interior	Type 304, mirror finish, stainless steel
Exterior	18 gauge, cold rolled steel, powder coated
Outer Door Gasket	Four-sided, molded magnetic vinyl
Inner Door Gasket	Removable, cleanable, feather-edged, silicone

## Electrical

Specifications	115 VAC, 50/60 Hz, 1 PH, 3.6 FLA
Circuit Breaker/Power Switch	6 Amp/2 Pole
Convenience Receptacle	75 Watts max. (one per chamber)
Alarm Contacts	Power interruption, overtemperature, deviation of temp., CO <sub>2</sub> , O <sub>2</sub> , and RH, customer connections through jack on back of unit. 30 V, 1 A max.

## Dimensions

Interior	21.3" W x 26.8" H x 20.0" F-B (54.1 cm x 68.1 cm x 50.8 cm)
Exterior	26.3" W x 39.5" H x 25.0" F-B (66.8 cm x 100.3 cm x 63.5 cm)

## Weight (per unit)

Net	265 lb. (120.2 kg)
Net Operational	365 lb. (165.6 kg)
Shipping	324 lb. (147.0 kg)

## Safety Specifications

Indoor Use Only	
Altitude	Up to 2000 meters
Temperature	5°C to 40°C
Humidity	Maximum 80% RH for temperatures up to 31°C, decreasing linearly to 50% RH at 40°C, non-condensing
Mains Supply Fluctuations	Not to exceed $\pm 10\%$ of the nominal voltage
Installation Category 2 <sup>1</sup>	
Pollution Degree 2 <sup>2</sup>	
Class of Equipment 1	

<sup>1</sup> Installation category (overvoltage category) defines the level of transient overvoltage which the instrument is designed to withstand safely. It depends on the nature of the electricity supply and its overvoltage protection means. For example, in CAT II which is the category used for instruments in installations supplied from a supply comparable to public mains such as hospital and research laboratories and most industrial laboratories, the expected transient overvoltage is 2500 V for a 230 V supply and 1500 V for a 120 V supply.

<sup>2</sup> Pollution Degree describes the amount of conductive pollution present in the operating environment. Pollution Degree 2 assumes that normally only non-conductive pollution such as dust occurs with the exception of occasional conductivity caused by condensation.



## Section 8 Spare Parts

All Models	Part #	Description
	360171	Liquid Level Switch
	103065	Feather Gasket
	113002	5/16-18 Glide Foot
	132046	115 / 230V Dual Heater
	132049	Face Heater 40 W, 24 VAC
	1900203	Heated Inner Door
	190650	Heated Inner Glass Door Kit (8 doors)
	190619	Left Inner Door Hinge
	190646	Door Lock for Inner Glass Door
	190666	Right Hand Door Swing
	700013	0.500" Flanged Nylon Bearing
	990026	Door Gasket w/ Magnet
	290184	Temp Probe 2252 Ohm
	192057	Micro Board Assembly (for Release 7 units)
	400218	12 VDC Single Output Power Supply (for Release 7 Units)
	1900346	Display Board Assembly
	230153	6A Circuit Breaker/Switch
	460157	Line Filter/Power Inlet
	420096	130VA Transformer, Int'l, 14 / 28V S
	230159	3.5 Amp Fuse - Microboard
	230158	2.5 Amp Fuse - Microboard
	250087	Solenoid Valve Kit w/assorted fittings
	1900071	Blower Motor Replacement Kit
	100113	Blower Wheel 3.5 x 1.5 CCW
	190846	Blower Scroll Assembly
	103072	Blower Plate Gasket
	1900601	Service Kit (for P/N 290090 T/C Type CO <sub>2</sub> Sensor Assembly
	103074	CO <sub>2</sub> Sensor Plate Gasket
	965010	CO <sub>2</sub> Gas Regulator
	961027	N <sub>2</sub> Gas Regulator

## All Models (continued)

Part #	Description
950316	Wall Clamp for CO <sub>2</sub> Bottle
130097	#6 Silicone Stopper w/ 3/8" Hole
180001	Polypropylene Funnel
117044	Fill Port Cap
770001	Bacterial Air Filter (CO <sub>2</sub> line, sample and access port)
760210	10 Disposable Polypropylene Inline Filters (770001)
760175	HEPA Filter
760209	HEPA Filter Value Pack (4 pack)
1900067	Filter Replacement Kit ([2] 770001 filter and 760175 HEPA)
1900123	Air Flow Restrictor Plate
190651	Decontamination Kit
50109579	AquaTec water cleaning cell 4-pack (for water reservoirs, pans, and bottles)
3166242	Gas Shuttle Switch
3050	External Gas Guard
190884	Stainless Steel Shelf and Channels
190647	Roller Base
190648	Floor Stand
400201	40 W Switcher Kit (replace triple output power supply P/N 400119, for units prior to Release 7)
2270107	Micro Board/Power Supply Kit (to replace Micro Board P/N 191634, Release 5 & 6 units)

### 115V Units (3110, 3120, 3130, 3140)

460024	115 V Convenience Outlet
230135	1.0 A Fuse for Convenience Outlet
430108	Line Cord (North America 125V / 15A)

### 230V Units (3111, 3121, 3131, 3141)

420097	43VA Transformer, INT. SR
460138	Power Outlet, Snap-in Receptacle,
230120	0.5 A Fuse for 230 V Outlet
430109	Line Cord (Continental Europe)
430111	Line Cord (Denmark)
430112	Line Cord (Italy)
430113	Line Cord (Switzerland)
430114	Line Cord (United Kingdom)
430115	Line Cord (India)
430116	Line Cord (Israel)
430118	Line Cord (Australia)
430367	Line Cord (China)

## IR or O<sub>2</sub> Units

190885	IR Sensor
191646	O <sub>2</sub> Circuit Board
250119	AC Solenoid
770001	Bacterial Air Filter (CO <sub>2</sub> line, sample and access port)
250118	Valve, Sol, O <sub>2</sub> , 12 VDC, 4 W
290083	O <sub>2</sub> Sensor Fuel Cell
190661	IR Sensor Filter
110065	Sample Port Cap

## Humidity Readout Units

290154	RH Sensor Assembly, 1 ft.
--------	---------------------------

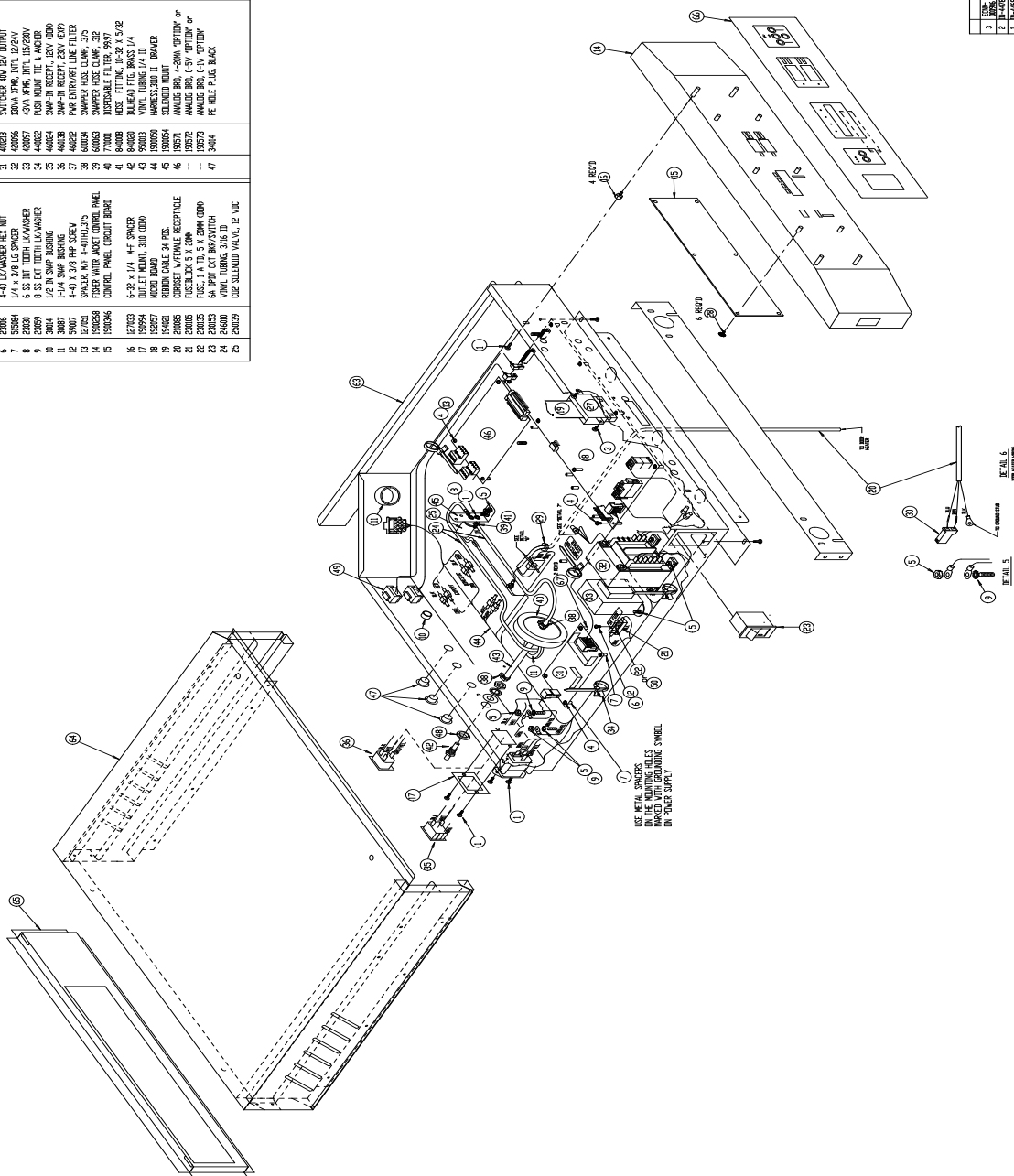
## Field Installation Kits by qualified service personnel

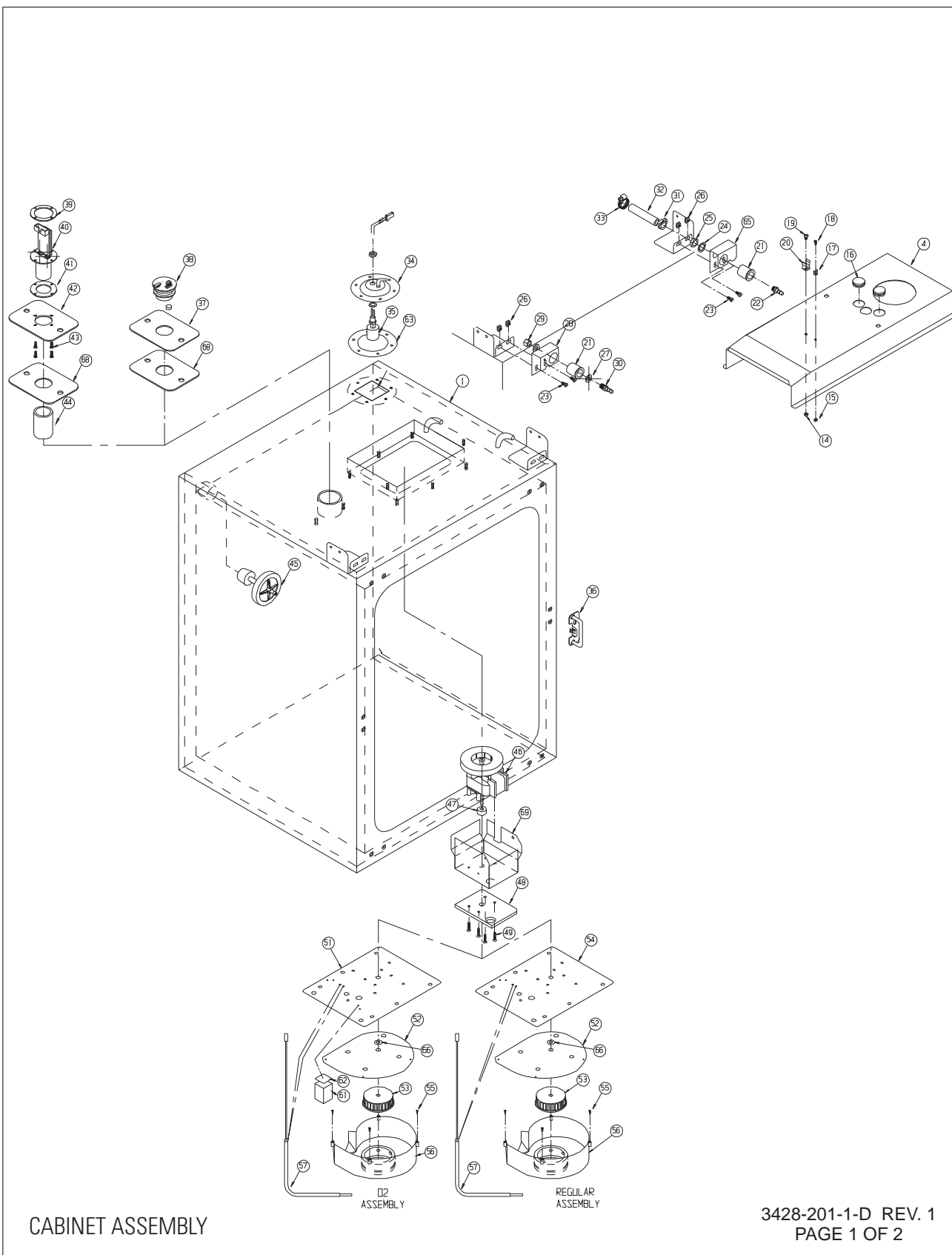
1900062	Humidity Sensor Kit
1900044	Analog Output* Kit, 4 - 20 mA
1900045	Analog Output* Kit, 0 - 5 VDC
1900046	Analog Output* Kit, 0 - 1 VDC

*\* select only one type of optional analog output*



## Spare Parts

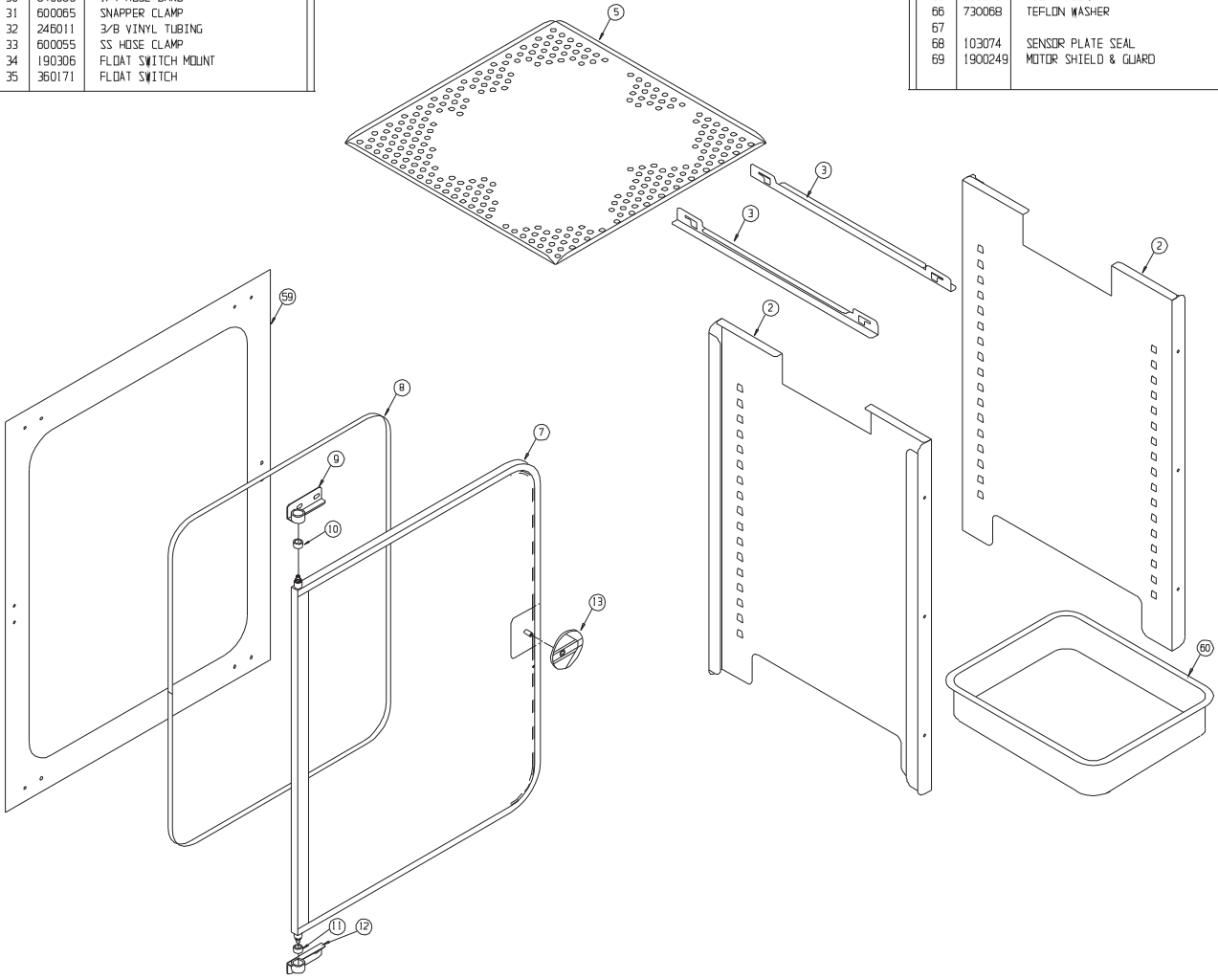
[illegible][illegible]



Section 8  
Spare Parts

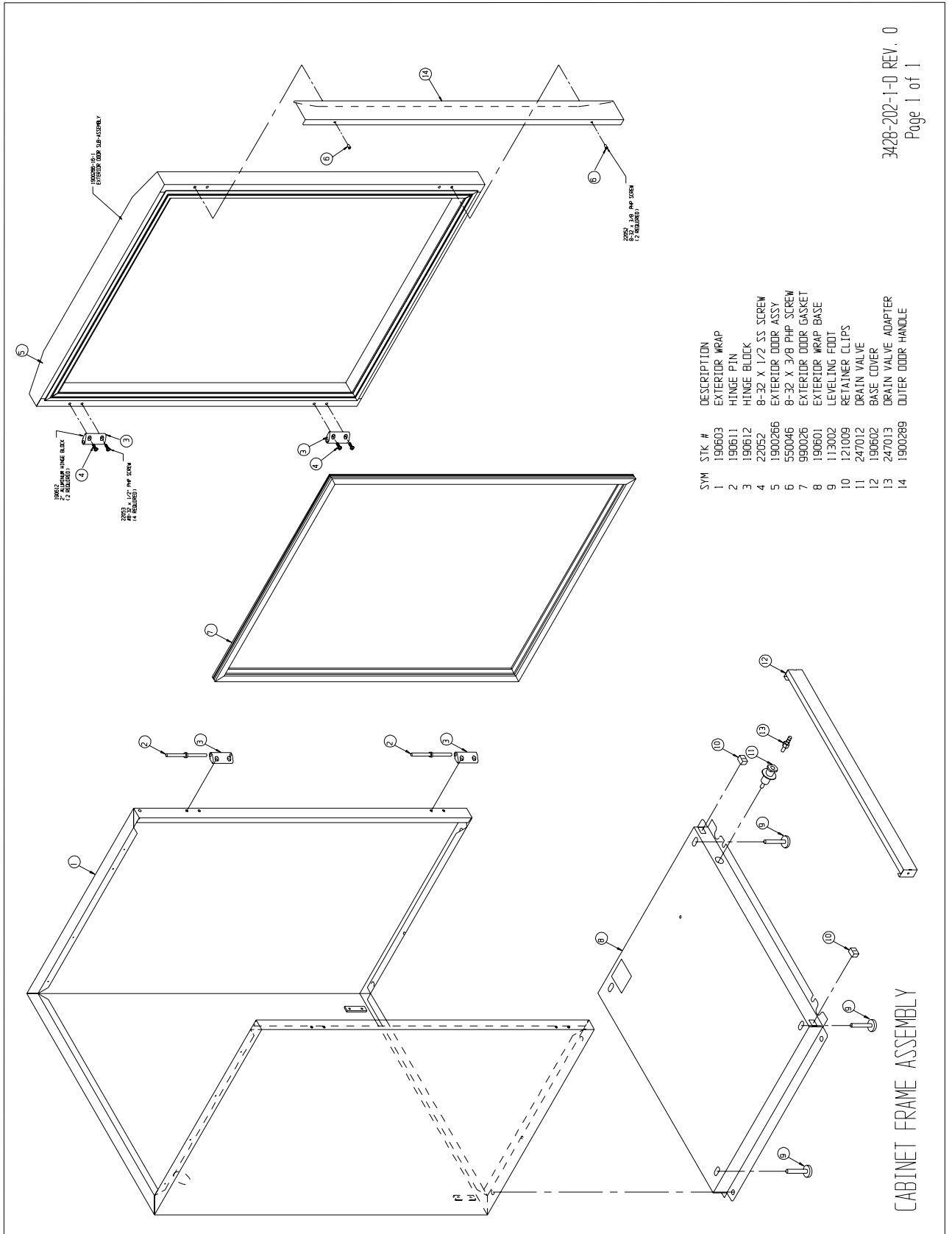
SYM	STK #	DESCRIPTION
1	1900600	WATER JACKET CHAMBER
2	190723	INTERIOR DUCT SHEETS
3	190664	SHELF CHANNELS
4	1900056	TOP DUCT(1900057 COPPER)
5	224175	SHELF
6	-----	-----
7	1900203	INNER GLASS DOOR
8	103108	INNER DOOR GASKET
9	190618	TOP INNER DOOR HINGE
10	700013	TOP INNER DOOR HINGE BUSHING
11	700046	BOTTOM INNER DOOR HINGE BUSHING
12	190619	BOTTOM INNER DOOR HINGE
13	180210	INNER DOOR KNOB
14	23010	8-32 SS NUT
15	23005	4-40 SS NUT
16	114022	15/16 BLIND GROMMET
17	111022	1/4" PROBE CLIP
18	22130	4-40 X 1/4 SS SCREW
19	22051	8-32 X 1/4 SS SCREW
20	111007	13/32 RH PROBE CLIP
21	190294	FILL PORT ACCESS
22	840043	PLASTIC HOSE BARB
23	680013	8 X 1/2 SS SCREW TYPE B
24	290128	1/2" LOCK WASHER
25	840043	NUT (PART OF #22)
26	3178097	#8 SPEED NUT
27	730060	SEAL WASHER
28	190675	LEFT MOUNT
29	840035	3/16 ADAPTER
30	840065	1/4 HOSE BARB
31	600065	SNAPPER CLAMP
32	246011	3/8 VINYL TUBING
33	600055	SS HOSE CLAMP
34	190306	FLOAT SWITCH MOUNT
35	360171	FLOAT SWITCH

SYM	STK #	DESCRIPTION
36	190634	INNER DOOR LATCH
37	190649	TC SENSOR PLATE
38	290090	TC SENSOR
39	190712	IR MOUNTING RING
40	190708	"DCS" IR SENSOR
41	103083	IR SENSOR GASKET
42	190711	IR MOUNTING PLATE
43	22130	4-40 X 1/4 SS SCREW
44	190708	IR SENSOR FILTER (PART OF ITEM 40)
45	190985	FILTERED STOPPER ASSEMBLY
46	156126	BLOWER MOTOR
47	730069	MOTOR SHAFT SEAL
48	190655	MOTOR SPACER PLATE
49	590023	8-32 X 3/4 SS SCREWS
50	-----	-----
51	1900055	O2 BLOWER PLATE ASSY
52	190893	SCROLL PLATE
53	100113	BLOWER WHEEL "D" SHAFT
54	1900053	STANDARD BLOWER PLATE
55	22134	6-32 X 1/2 "FILLISTER" HEAD SCREW
56	190846	BLOWER SCROLL
57	290184	TEMPERATURE PROBE
58	760175	HEPA FILTER
59	180174	PVC FACE FRAME
60	237016	HUMIDITY PAN
61	1900058	BRASS O2 BLOCK
62	100064	BRASS BLOCK SEAL
63	103032	FLOAT SWITCH GASKET
64	23024	3/8 FLAT WASHER
65	190675	RIGHT MOUNT
66	730068	TEFLON WASHER
67	-----	-----
68	103074	SENSOR PLATE SEAL
69	1900249	MOTOR SHIELD & GUARD

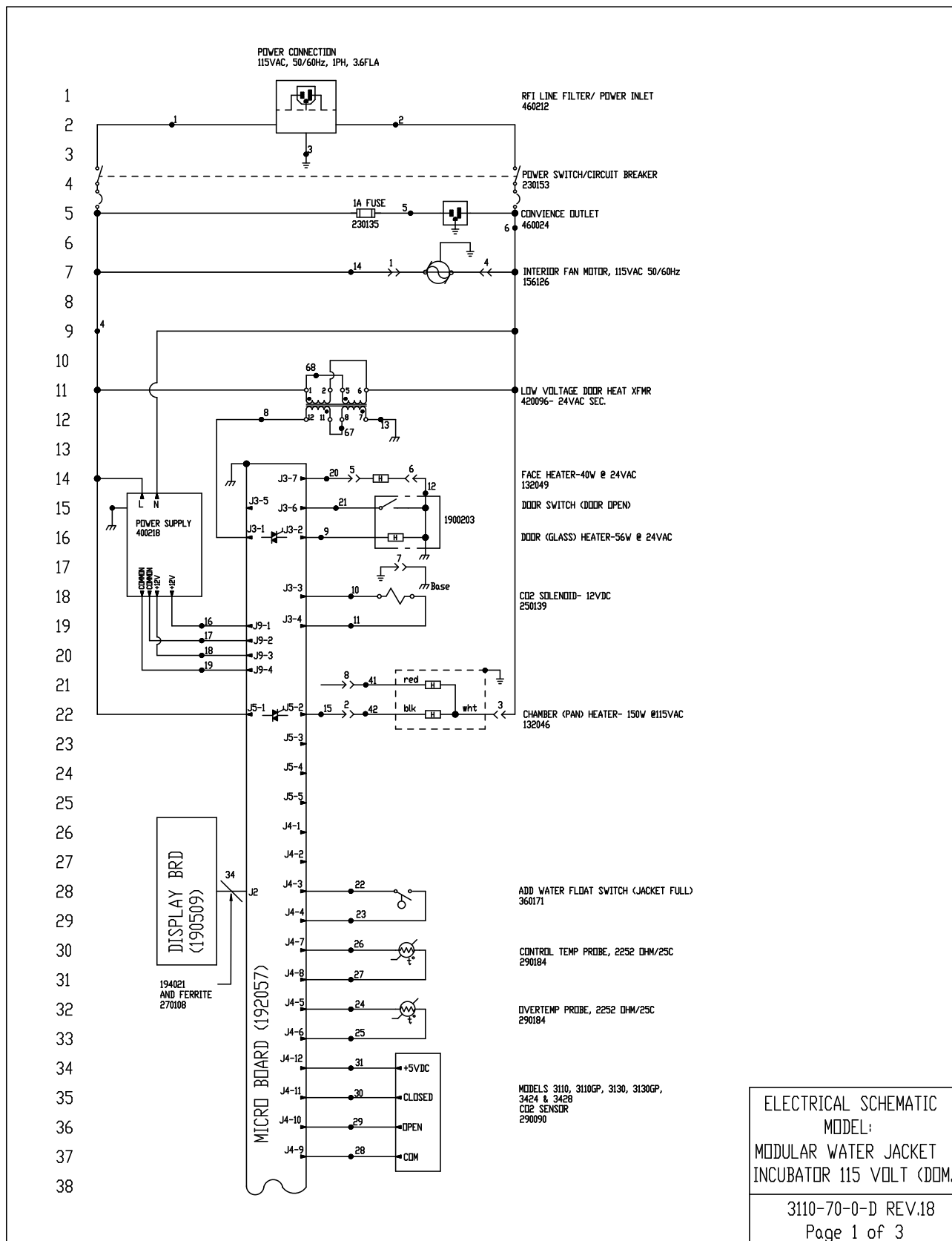


CABINET ASSEMBLY

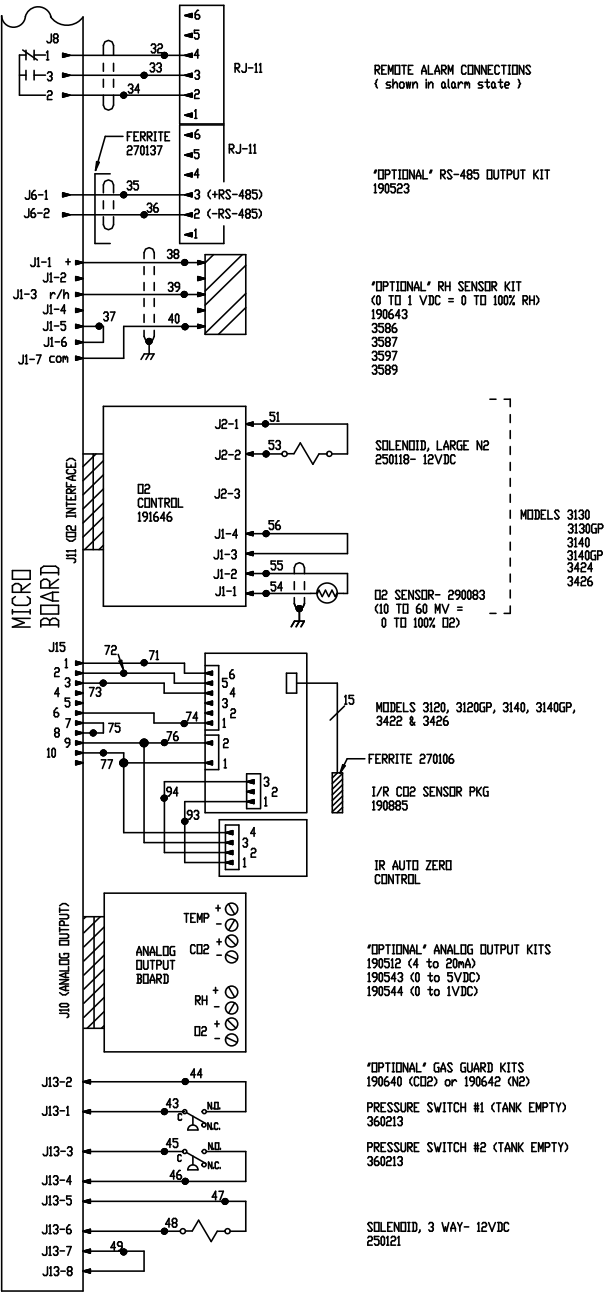
3428-201-1-D REV. 1  
PAGE 2 OF 2







39  
40  
41  
42  
43  
44  
45  
46  
47  
48  
49  
50  
51  
52  
53  
54  
55  
56  
57  
58  
59  
60  
61  
62  
63  
64  
65  
66  
67  
68  
69  
70  
71  
72  
73  
74  
75  
76



ELECTRICAL SCHEMATIC  
MODEL:  
MODULAR WATER JACKET  
INCUBATOR 115 VOLT (DOM.)  
3110-70-0-D REV.18  
Page 2 of 3

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

WIRE REFERENCE CHART								
NO.	GA.	COLOR	NO.	GA.	COLOR	NO.	GA.	COLOR
1	16	BROWN	26	22	YELLOW	52	N/U	
2	16	BLUE	27	22	YELLOW	53	20	PURPLE
3	16	GRN/YEL	28	22	GREEN	54	22/2	RED
38	16	GREEN	29	22	GRAY	55	22/2	BLACK
4	18	BLACK	30	22	PURPLE	56	22	BLACK
5	18	BLACK	31	22	ORANGE	57		
6	18	WHITE	32	22/3	BLACK		N/U	
7	N/U		33	22/3	RED	66		
8	18	BROWN	34	22/3	WHITE	67	18	RED
9	18	BLUE	35	22/2	RED	68	18	ORANGE
10	18	YELLOW	36	22/2	BLACK	69	N/U	
11	18	YELLOW	37	22	BLACK	70	N/U	
12	18	BLACK	38	22/3	BLACK	71	22	BLUE
13	18	GREEN	39	22/3	RED	72	22	ORANGE
14	18	BLUE	40	22/3	GREEN	73	22	YELLOW
15	18	ORANGE	41	18	RED	74	22	BROWN
16	18	RED	42	18	BLACK	75	22	BLACK
17	18	GREEN	43	20	RED	76	22	RED
18	18	ORANGE	44	20	RED	77	22	GREEN
19	18	GREEN	45	20	BLUE	78		
20	18	ORANGE	46	20	BLUE		24/15	I/R CABLE
21	22	BROWN	47	20	ORANGE	91		
22	22	BLUE	48	20	ORANGE	92	22	BLACK
23	22	BLUE	49	20	BLACK	93	22	BROWN
24	22	RED	50	N/U		94	22	WHITE
25	22	RED	51	20	YELLOW			

N/U= Not Used

SENSOR REFERENCE VALUES

CO2 (290090) DIFFERENCE VOLTAGE OF 3-6MV/%CO2  
J4-9 & J4-10 TO J4-9 & J4-11

RH (190643) J1-7 & J1-1 = 12VDC  
J1-7 & J1-3 = 10MV/%RH

O2 (290083) J6-1 & J6-2 = 12MV @ 21%O2

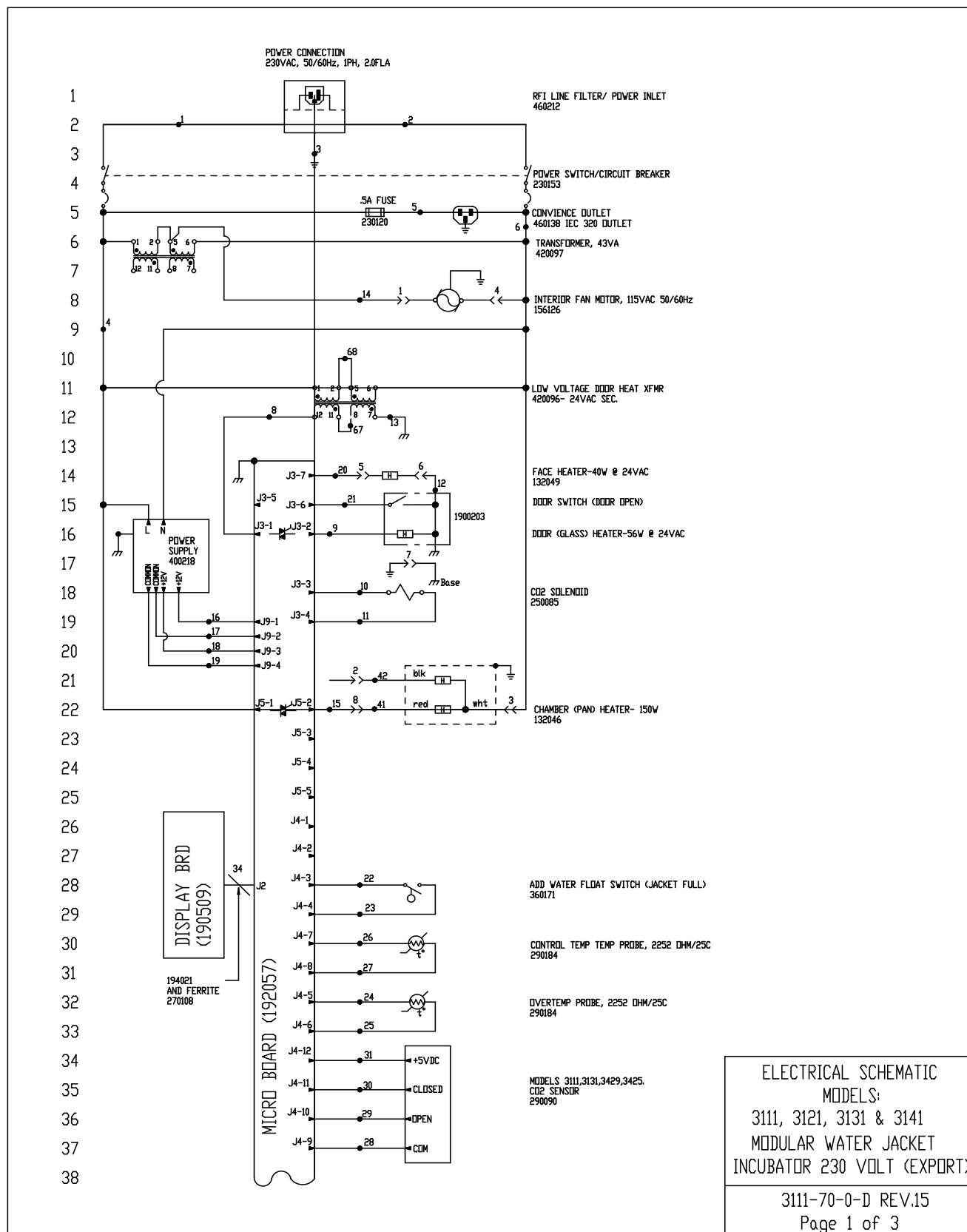
MODEL	DESCRIPTION
3110	WATER JACKET INCUBATOR - TC CO2
3120	WATER JACKET INCUBATOR - IR CO2
3130	WATER JACKET INCUBATOR - TC O2
3140	WATER JACKET INCUBATOR - IR O2
3428	WATER JACKET INCUBATOR - TC CO2
3422	WATER JACKET INCUBATOR - IR CO2
3424	WATER JACKET INCUBATOR - TC O2
3426	WATER JACKET INCUBATOR - IR O2
3110GP	WATER JACKET INCUBATOR - TC CO2
3120GP	WATER JACKET INCUBATOR - IR CO2
3130GP	WATER JACKET INCUBATOR - TC O2
3140GP	WATER JACKET INCUBATOR - IR O2

NOTES:	
⊗ Denotes Terminal Strip Connection	Parts List Reference Number
Last Relay Number	○ Assembly
Last Terminal Number	◇ Panel
Last Wire Number	○ Refrigeration
→ >— Denotes Pin & Socket Connection	□ Wiring

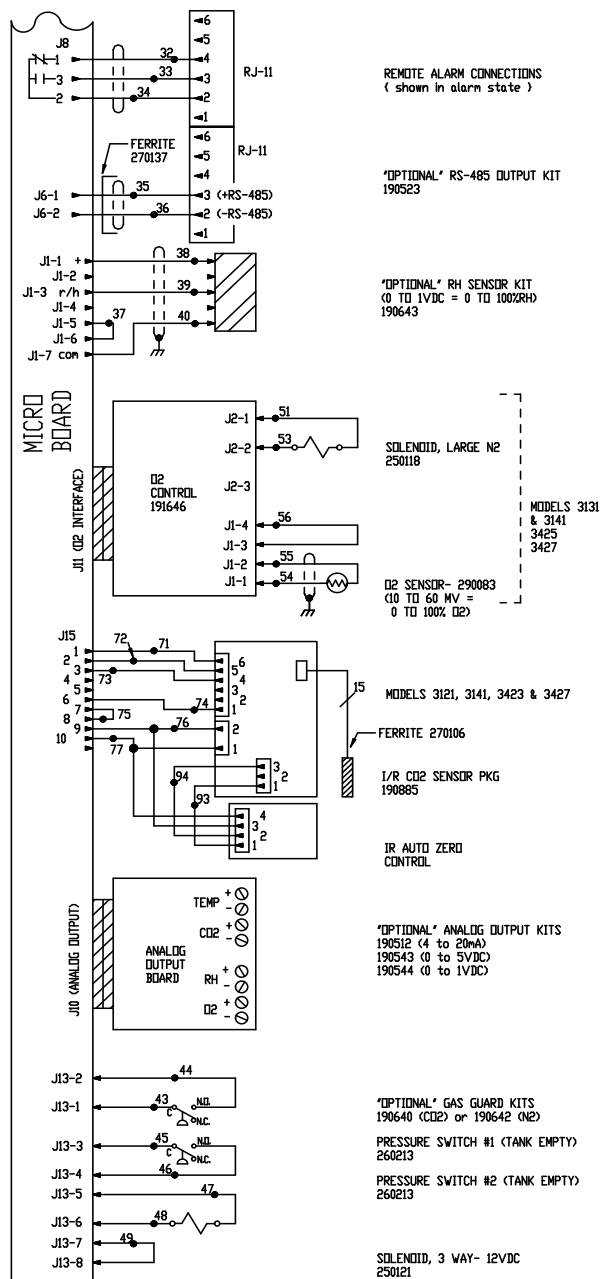
ELECTRICAL SCHEMATIC  
MODEL:  
MODULAR WATER JACKET  
INCUBATOR 115 VOLT (DOM.)  
3110-70-0-D REV.18  
Page 3 of 3



**Section 9**  
Electrical Schematics



39  
40  
41  
42  
43  
44  
45  
46  
47  
48  
49  
50  
51  
52  
53  
54  
55  
56  
57  
58  
59  
60  
61  
62  
63  
64  
65  
66  
67  
68  
69  
70  
71  
72  
73  
74  
75  
76



ELECTRICAL SCHEMATIC  
MODELS:  
3111, 3121, 3131 & 3141  
MODULAR WATER JACKET  
INCUBATOR 230 VOLT (EXPORT)

3111-70-0-D REV.15  
Page 2 of 3

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

WIRE REFERENCE CHART								
NO.	GA.	COLOR	NO.	GA.	COLOR	NO.	GA.	COLOR
1	16	BROWN	26	22	YELLOW	52	N/U	
2	16	BLUE	27	22	YELLOW	53	20	PURPLE
3	16	GRN/YEL	28	22	GREEN	54	22/2	RED
3B	16	GREEN	29	22	GRAY	55	22/2	BLACK
4	18	BLACK	30	22	PURPLE	56	22	BLACK
5	18	BLACK	31	22	ORANGE	57		
6	18	WHITE	32	22/3	BLACK		N/U	
7	N/U		33	22/3	RED	66	↑	
8	18	BROWN	34	22/3	WHITE	67	18	RED
9	18	BLUE	35	22/2	RED	68	18	ORANGE
10	18	YELLOW	36	22/2	BLACK	69		N/U
11	18	YELLOW	37	22	BLACK	70	22	BLACK
12	18	BLACK	38	22/3	BLACK	71	22	BLUE
13	18	GREEN	39	22/3	RED	72	22	ORANGE
14	18	BLUE	40	22/3	GREEN	73	22	YELLOW
15	18	ORANGE	41	18	RED	74	22	BROWN
16	18	RED	42	18	BLACK	75	22	BLACK
17	18	GREEN	43	20	RED	76	22	RED
18	18	ORANGE	44	20	RED	77	22	GREEN
19	18	GREEN	45	20	BLUE	78		I/R
20	18	ORANGE	46	20	BLUE		24/15	CABLE
21	22	BROWN	47	20	ORANGE	91	↑	
22	22	BLUE	48	20	ORANGE	92	22	BLACK
23	22	BLUE	49	20	BLACK	93	22	BROWN
24	22	RED	50	N/U		94	22	WHITE
25	22	RED	51	20	YELLOW			

N/U= Not Used

SENSOR REFERENCE VALUES

CO2 (290090) DIFFERENCE VOLTAGE OF 3-6MV/%CO2  
J4-9 & J4-10 TO J4-9 & J4-11

RH (190643) J1-7 & J1-1 = 12VDC  
J1-7 & J1-3 = 10MV/%RH

O2 (290083) J6-1 & J6-2 = 12MV @ 21%O2

MODEL	DESCRIPTION
3111	WATER JACKET INCUBATOR - TC CO2
3121	WATER JACKET INCUBATOR - IR CO2
3131	WATER JACKET INCUBATOR - TC O2
3141	WATER JACKET INCUBATOR - IR O2
3429	WATER JACKET INCUBATOR - TC CO2 (THERMO 8000)
3423	WATER JACKET INCUBATOR - IR CO2 (THERMO 8000)
3425	WATER JACKET INCUBATOR - TC O2 (THERMO 8000)
3427	WATER JACKET INCUBATOR - IR O2 (THERMO 8000)
3111GP	WATER JACKET INCUBATOR - TC CO2 GENERAL PURPOSE
3121GP	WATER JACKET INCUBATOR - IR CO2 GENERAL PURPOSE
3131GP	WATER JACKET INCUBATOR - TC O2 GENERAL PURPOSE
3141GP	WATER JACKET INCUBATOR - IR O2 GENERAL PURPOSE

NOTES:		ELECTRICAL SCHEMATIC MODELS: 3111, 3121, 3131 & 3141 MODULAR WATER JACKET INCUBATOR 230 VOLT (EXPORT)  3111-70-0-D REV.15 Page 3 of 3
⊗	Denotes Terminal Strip Connection	
	Last Relay Number	
	Last Terminal Number	
	Last Wire Number	
→↘	Denotes Pin & Socket Connection	
	Parts List Reference Number	
	○ Assembly	
	○ Panel	
	○ Refrigeration	
	□ Wiring	

## THERMO FISHER SCIENTIFIC STANDARD PRODUCT WARRANTY

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

During the first year, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo's expense, labor included. The Watlow EZ-ZONE PM controller is covered for one additional year for repair or replacement (parts only), provided the unit has not been misapplied. Installation and calibration are not covered by this warranty agreement. The Technical Services Department must be contacted for warranty determination and direction prior to performance of any repairs. Expendable items, glass, filters and gaskets are excluded from this warranty.

Replacement or repair of components parts or equipment under this warranty shall not extend the warranty to either the equipment or to the component part beyond the original warranty period. The Technical Services Department must give prior approval for return of any components or equipment. At Thermo's option, all non-conforming parts must be returned to Thermo Fisher Scientific postage paid and replacement parts are shipped FOB destination.

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

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

If equipment service is required, please call your Technical Services Department at 1-800-438-4851 (USA and Canada) or 1-740-373-4763. We're ready to answer your questions on equipment warranty, operation, maintenance, service and special application. Outside the USA, contact your local distributor for warranty information.



Rev. 5 8/13

## THERMO FISHER SCIENTIFIC INTERNATIONAL DEALER WARRANTY

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

During the first year, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo's expense, labor excluded. The Watlow EZ-ZONE PM controller is covered for one additional year for repair or replacement (parts only), provided the unit has not been misapplied. Installation and calibration are not covered by this warranty agreement. The Technical Services Department must be contacted for warranty determination and direction prior to performance of any repairs. Expendable items, glass, filters, reagents, tubing, and gaskets are excluded from this warranty.

Replacement or repair of components parts or equipment under this warranty shall not extend the warranty to either the equipment or to the component part beyond the original warranty period. The Technical Services Department must give prior approval for return of any components or equipment. At Thermo's option, all non-conforming parts must be returned to Thermo postage paid and replacement parts are shipped FOB destination.

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

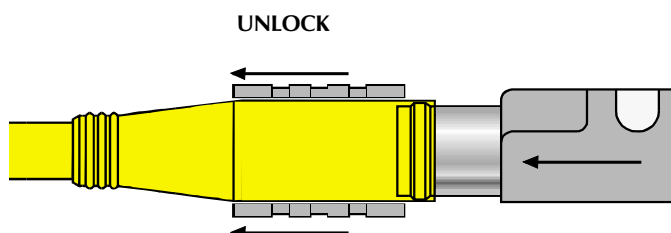
Your local Thermo Sales Office is ready to help with comprehensive site preparation information before your equipment arrives. Printed instruction manuals carefully detail equipment installation, operation and preventive maintenance.

Contact your local distributor for warranty information. We're ready to answer your questions on equipment warranty, operation, maintenance, service and special application.



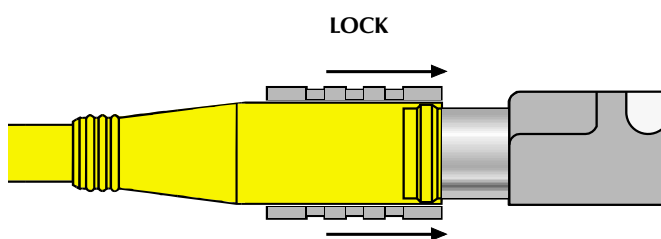
Rev. 5 8/13

## *picofast* “Snap Lock” Connector Instructions



With TURCK’s patented Locking Sleeve pulled back, any *picofast* sensor slides on without any difficulty.

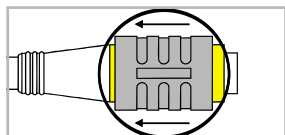
After sensor is connected, push Locking Sleeve forward to create a watertight connection. It’s a snap!



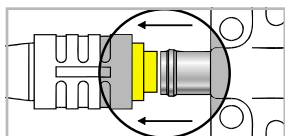
## Installation Instructions

### To Attach:

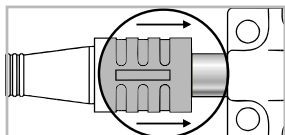
1. Be sure black locking sleeve is pulled back.



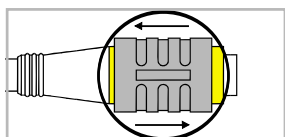
2. Line up pins and push connector onto plug. You will feel a “snap.” Do not twist.



3. Push locking sleeve forward on connector until it is flush with the front of the connector.

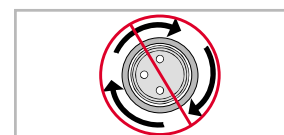


4. If sleeve is difficult to slide on a new connector, “exercise” it a few times. Do not use tools.

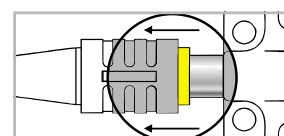


### To Detach:

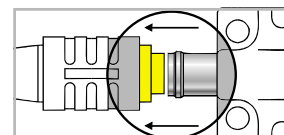
1. Do not twist.



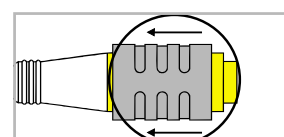
2. Pull locking sleeve back to “unlocked” position.



3. Pull connector straight off.



4. Leave locking sleeve in unlocked position.



**thermoscientific.com**

© 2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

---

**Thermo Fisher Scientific (Asheville) LLC**

401 Millcreek Road  
Marietta, Ohio 45750  
United States

**Thermo**  
S C I E N T I F I C  
A Thermo Fisher Scientific Brand