

Forma Scientific, Inc.

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

Telephone: (740) 373-4763

Telefax: (740) 373-4189

Models: 3110 Series*

Single Chamber
Water Jacket Incubator

Models: 3210 Series*

Dual Chamber
Water Jacket Incubator

Operating and Maintenance Manual

Manual No: 7033110

Rev. 1

Read this Instruction Manual

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

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

Refer to the serial tag on the rear of this instruction manual

****Refer to listing of all models on Page iii.***



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

MANUAL NO. 7033110				
1	18891/SI-7832	2/23/00	Change cooling coil tubing from copper to vinyl, newloc fittings	ccs
0	--	1/3/00	Release 4, Class 100 updates	ccs
REV	ECR/ECN	DATE	DESCRIPTION	By

Single Chamber Models			
Model	CO₂ Sensor**	O₂	Voltage***
3110	T/C	No	115
3111	T/C	No	230
3120	IR	No	115
3121	IR	No	230
3130	T/C	Yes	115
3131	T/C	Yes	230
3140	IR	Yes	115
3141	IR	Yes	230

Dual Chamber Models			
Model*	CO₂ Sensor**	O₂	Voltage***
3210	T/C	No	115
3211	T/C	No	230
3220	IR	No	115
3221	IR	No	230
3230	T/C	Yes	115
3231	T/C	Yes	230
3240	IR	Yes	115
3241	IR	Yes	230

*The models listed as dual chamber models are stacked or dual chamber units. These units arrive as single stand-alone units. They may be stacked one on top of the other or configured side by side with opposite door swing.

**T/C is a thermal conductivity sensor. IR is an infrared sensor.

***All units are 50/60 Hz.

General Safety Notes used in this Manual

	<p>Important operating and/or maintenance instructions. Read the accompanying text carefully.</p> <p>Ce symbole attire l'attention de l'utilisateur sur des instructions importantes de fonctionnement et/ou d'entretien. Il peut être utilisé seul ou avec d'autres symboles de sécurité. Lire attentivement le texte d'accompagnement.</p> <p>Wichtige Betriebs- und/oder Wartungshinweise. Lesen Sie den nachfolgenden Text sorgfältig.</p> <p>Importante instrucciones de operación y/o mantenimiento. Lea el texto acompañante cuidadosamente.</p>
	<p>Potential electrical hazards. Only qualified persons should perform procedures associated with this symbol.</p> <p>Ce symbole attire l'attention de l'utilisateur sur des risques électriques potentiels. Seules des personnes qualifiées doivent appliquer les instructions et les procédures associées à ce symbole.</p> <p>Gefahr von Stromschlägen. Nur qualifizierte Personen sollten die Tätigkeiten ausführen, die mit diesem Symbol bezeichnet sind.</p> <p>Potencial de riesgos electricos. Solo personas das capacitadas deben ejecutar los procedimientos asociadas con este simbolo.</p>
	<p>Equipment being maintained or serviced must be turned off and locked off to prevent possible injury.</p> <p>Risques potentiels liés à l'énergie. L'équipement en entretien ou en maintenance doit être éteint et mis sous clé pour éviter des blessures possibles.</p> <p>Geräte, an denen Wartungs- oder Servicearbeiten durchgeführt werden, müssen abgeschaltet und abgeschlossen werden, um Verletzungen zu vermeiden.</p> <p>El equipo recibiendo servicio o mantenimiento debe ser apagado y asegurado para prevenir danos.</p>



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

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

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

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

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

Do You Need Information or Assistance on Forma Scientific Products?

If you do, please contact us 8:00 a.m. to 7:00 p.m. (Eastern Time) at:

1-740-373-4763	Direct
1-888-213-1790	Toll Free, U.S. and Canada
1-740-373-4189	FAX
http://www.forma.com	Internet Worldwide Web Home Page
fservice@forma.com	Service E-Mail Address

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

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

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

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

Regardless of your needs, our professional telephone technicians are available to assist you Monday through Friday from 8:00 a.m. to 7:00 p.m. Eastern Time. Please contact us by telephone or fax. If you wish to write, our mailing address is:

Forma Scientific, Inc.
Millcreek Road, PO Box 649
Marietta, OH 45750

International customers please contact your local Forma Scientific distributor.

Table of Contents

Section 1 - Installation and Start-Up

1.1 Name and Description of Parts	1-1
1.2 Control Panel Keys, Displays and Indicators.....	1-2
1.3 Operating the Control Panel.....	1-3
1.4 Displays.....	1-5
1.5 Installing the Unit.....	1-6
a. Choosing the Location.....	1-6
b. Stacking the Incubators	1-6
c. Preliminary Cleaning and Disinfecting.....	1-8
d. Installing the Shelves.....	1-9
e. Installing the Access Port Stopper	1-10
f. Installing the HEPA Filter.....	1-10
g. Leveling the Unit.....	1-11
h. Connecting the Unit to Electrical Power	1-12
i. Filling the Water Jacket.....	1-12
j. Filling the Humidity Pan	1-13
k. Connecting the CO ₂ Gas Supply	1-14
l. Connecting the N ₂ Gas Supply.....	1-15
1.6 Incubator Start-Up.....	1-16
a. Setting the Operating Temperature.....	1-16
b. Setting the Overtmp Setpoint	1-16
c. Setting the CO ₂ Setpoint.....	1-17
d. Setting the O ₂ Setpoint	1-18
Chart 1-1, Set Mode.....	1-19

Section 2 - Calibration

2.1 Calibration Mode2-1

- a. Calibrating the Temperature2-1
 - Temperature Stabilization Periods2-1
- b. Calibrating the Thermal Conductivity CO₂ System2-2
- c. Calibrating the T/C %CO₂ Zero to Ambient2-3
 - T/C CO₂ Sensor Stabilization Periods2-3
- d. Calibrating the T/C %CO₂ Span.....2-3
- e. Calibrating the Infra-Red CO₂ System2-4
 - IR CO₂ Sensor Stabilization Times.....2-4
- f. Calibrating the O₂ System2-5
 - O₂ Sensor Stabilization Times.....2-5
- g. Calibrating the Relative Humidity.....2-5
 - Relative Humidity Stabilization Times.....2-5

Chart 2-1, Calibrate Mode2-7

Section 3 - Configuration

3.1 Configuration Mode3-1

- a. Turning the Audible Alarm ON/OFF3-1
- b. Setting an Access Code3-1
- c. Setting a Low Temp Alarm Limit (tracking alarm).....3-1
- d. Setting a High Temp Alarm Limit (tracking alarm).....3-2
- e. Enabling Temp Alarms to Trip Contacts.....3-2
- f. Setting a Low CO₂ Alarm Limit (tracking alarm).....3-2
- g. Setting a High CO₂ Alarm Limit (tracking alarm)3-3
- h. Enabling CO₂ Alarms to Trip Contacts3-3
- i. Setting a New Zero Number for T/C CO₂ Sensors3-3
- j. Setting a New Span Number for T/C CO₂ Sensors3-4
- k. Setting a Low RH Alarm Limit3-4
- l. Enabling RH Alarms to Trip Contacts3-4
- m. New O₂ Cell3-5
- n. Setting a Low O₂ Alarm Limit (tracking alarm).....3-5
- o. Setting a High O₂ Alarm Limit (tracking alarm)3-5
- p. Enabling O₂ Alarms to Trip Contacts3-6

- q. Enabling Temp/RH to be Displayed.....3-6
- r. Enabling CO₂/O₂ to be Displayed.....3-6
- s. Enabling O₂ Compensation for T/C CO₂ Systems.....3-7
- t. Setting a RS-485 Communications Address (1535 compatible only).....3-7
- u. Selecting a Primary Tank with the Gas Guard Option3-7
- v. Enabling the Gas Guard System.....3-7
- w. Setting the Door Heat3-8
- Chart 3-1, Configure Mode (4 pages).....3-9

Section 4 - Alarms

- 4.1 Alarms (chart of alarms)4-1
- 4.2 Sensor Fault Alarms.....4-2
- 4.3 REPL O₂ SNSR (Alarm).....4-2
- 4.4 REPL IR SNSR (Alarm).....4-2
- 4.5 CO₂ SENS ERR4-2
- 4.6 IR AUTO Z ERR4-3

Section 5 - Maintenance

- 5.1 Disinfecting the Incubator Interior.....5-1
- 5.2 Cleaning the Cabinet Exterior.....5-2
- 5.3 Cleaning the Humidity Pan.....5-2
- 5.4 Reversing the Door Swing.....5-3
 - a. Reversing the Hinges for the Exterior Door5-3
- 5.5 Replacing the Fuses5-8
- 5.6 HEPA Filter Maintenance.....5-10
- 5.7 Draining Water Jacket.....5-11

Section 6 - Factory Installed Options

- 6.1 Connections to External Equipment6-1
 - a. Connecting the Remote Alarm Contacts6-1
 - b. Connecting the RS485 Interface, P/N 190523.....6-2
 - c. Connecting the Analog Output Boards.....6-2

6.2 Gas Guard for CO₂ and N₂ (190640/190642)6-4

- a. Connecting the CO₂ Gas Supplies6-5
- b. Connecting the N₂ Gas Supplies6-5
- c. Activating the Built-in Gas Guard6-6
- d. Operation of the CO₂ or N₂ Gas Guard6-6

6.3 Humidity Readout6-7

- a. Factors Affecting the Humidity Level in the Chamber.....6-7
- b. Accuracy of the Humidity Readout6-8

6.4 Factory Installed Cooling Coil6-8

- a. Installing the Cooling Coil Incubator6-10

Section 7 - Specifications

Section 8 - Spare Parts

Section 9 - Electrical Schematics

Section 10 - Warranty Information

Appendix A

Section 1 - Installation and Start-up

1.1 Name and Description of Parts

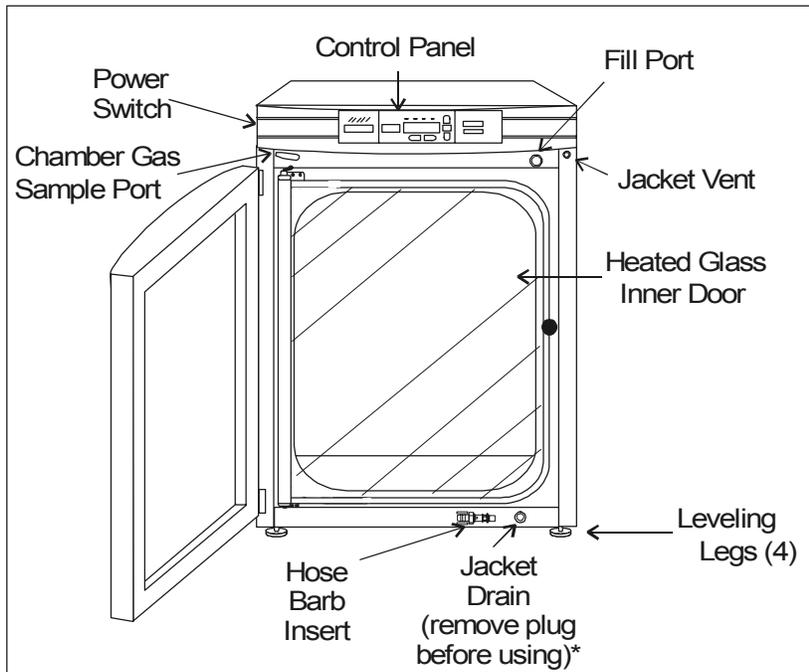


Figure 1-1
Water Jacket Incubator

- Main Power Switch
- Control Panel - Keypad, Displays and Indicators. (See 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 – Remove plug and use hose barb insert included. Plug when not in use.

- Outer Door - Reversible to opposite swing, see Section 5.4
- Heated Inner Door - Keeps chamber interior dry. Reversible to opposite swing, see Section 5.4.
- Chamber Gas Sample Port - Used for sampling chamber CO₂ content, using a Fyrite or similar instrument.

Note: The port must remain capped when not in use.

Note: The incubators are stackable. See Section 1.5.

*When setting up the incubator, install the cover plate packed inside the incubator shipping carton .

1.2 Control Panel Keys, Displays and Indicators

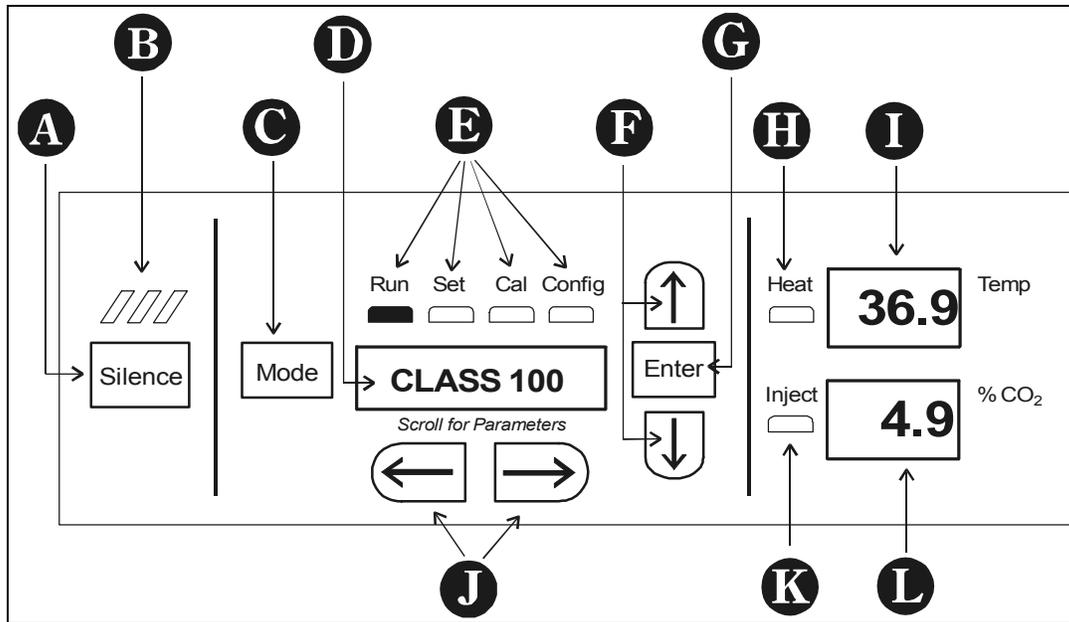


Figure 1-2 Control Panel

A SILENCE - Silences the audible alarm.

B Alarm Indicator - Pulses on/off during an alarm condition in the cabinet.

C MODE Select Switch - Used to select Run, Setpoints, Calibration and System Configuration Modes.

D Message Center - Displays system status.

E Mode Select Indicators-
Run: Run Menu
Set: Set Points Menu
Cal: Calibrate Menu
Config: Configuration Menu

F Up and Down Arrows: Increases or decreases the parameter values that are numbers, toggles the parameter values that are choices.

H Heat Indicator: Lights when power is applied to the heaters.

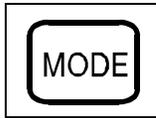
I Temp Display: Can be programmed to display temperature continuously, RH continuously (with RH option), or toggle between temperature and humidity (with RH option). See Section 3.1, Configuration.

J Scroll for Parameters Keys: Scrolls the operator through the parameters of the mode that is selected.

Ⓚ Inject Indicator: Lights when gas is being injected into the incubator. If the %CO₂/O₂ display (item L) is continuously displaying CO₂, the light indicates CO₂ injection only. If the %CO₂/O₂ display is continuously displaying O₂, light indicates N₂ injection. If the %CO₂/O₂ is toggling, either a CO₂ or N₂ injection will cause the indicator to light.

Ⓛ %CO₂/O₂ display: Can be programmed to display CO₂ continuously or O₂ continuously (on units equipped with O₂), or toggle between CO₂ and O₂ (on units equipped with O₂) See Section 3.1, Configuration.

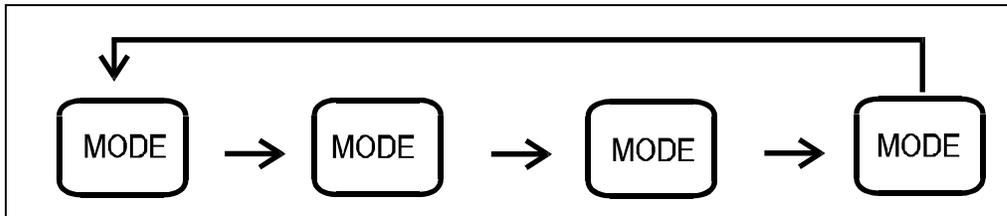
1.3 Operating the Control Panel



The Model 3110 Series water jacket incubator has four basic modes, which 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.
- **Config** is the system configuration mode that allows for custom setup of various options.

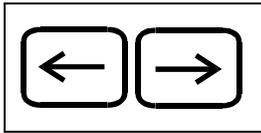
The chart on the following page shows the selections under each of the modes.



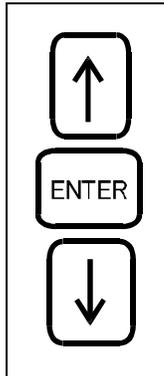
RUN	SETPOINT	CALIBRATION	CONFIGURATION
Default Mode	Temperature Overtemp CO2 <i>O2</i>	Temp Offset CO2 Zero CO2 Span <i>IR CO2 Span</i> <i>O2 Span</i> <i>RH Offset</i>	Audible Access Code Temp Lo Alarm Temp Hi Alarm Temp Relay CO2 Lo Alarm CO2 Hi Alarm CO2 Relay CO2 Z & S #'s <i>RH Lo Alarm</i> <i>RH Relay</i> <i>O2 Hi Alarm</i> <i>O2 Lo Alarm</i> <i>O2 Relay</i> <i>Display 1 Setup</i> <i>Display 2 Setup</i> <i>O2 Compensation</i> RS485 Address <i>Gas Guard Setup</i> Door Heat

****Base Unit Displays**

****Option Displays**



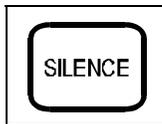
Scroll for Parameters Keys: Steps the operator through the parameters of SET, CAL and CONFIG Modes. The right arrow goes to the next parameter, the left arrow returns to the previous parameter.



Up Arrow: Increases or toggles the parameter value that has been selected in the SET, CAL, and CONFIG 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 CONFIG Modes.



Silence Key: Press to silence the audible alarm. See Section 4 for alarm ringback times.

1.4 Displays

Message Center

Displays the system status (Mode) at all times. Displays **CLASS 100** during normal operation, or alarm messages if the system detects an alarm condition. See Section 4.1, Alarms.

Upper and Lower Displays

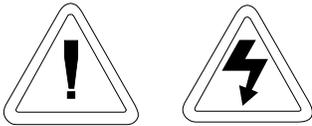
These 7-segment displays vary depending upon the options present and the configuration chosen. The upper display can display temp or RH, or toggle between them. The bottom display can display CO₂ or O₂, or toggle between them. See Section 3.1, Configuration.

1.5 Installing the Unit

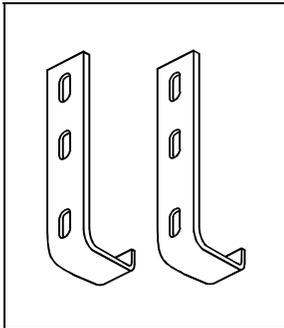
a. Choosing the Location

1. Locate the unit on a firm, level surface capable of supporting the unit's operational weight of 365 lbs. (166 kg).
2. Locate away from doors and windows and heating and air conditioning ducts.
3. Allow enough clearance behind the unit for electrical and gas hook-up.

b. Stacking the Incubators

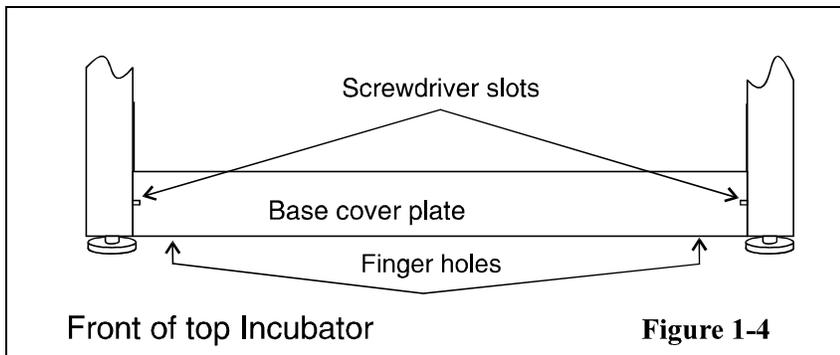


If the units have been in service, disconnect the power cord connector and drain the water jacket of the designated top unit before stacking.

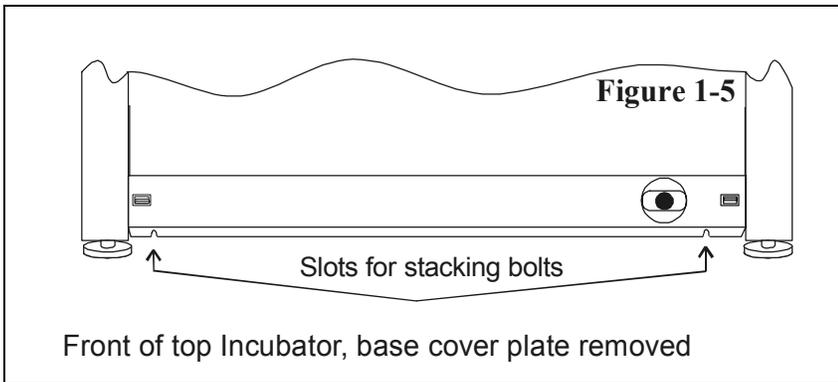


Note: Stacking brackets (shown at left) stacking bolts, washers, and bolts for stacking are included with each unit.

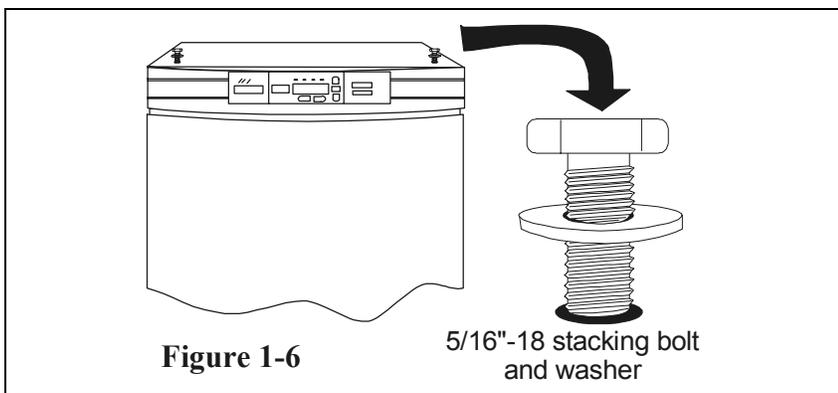
Figure 1-3 Stacking brackets



1. 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)



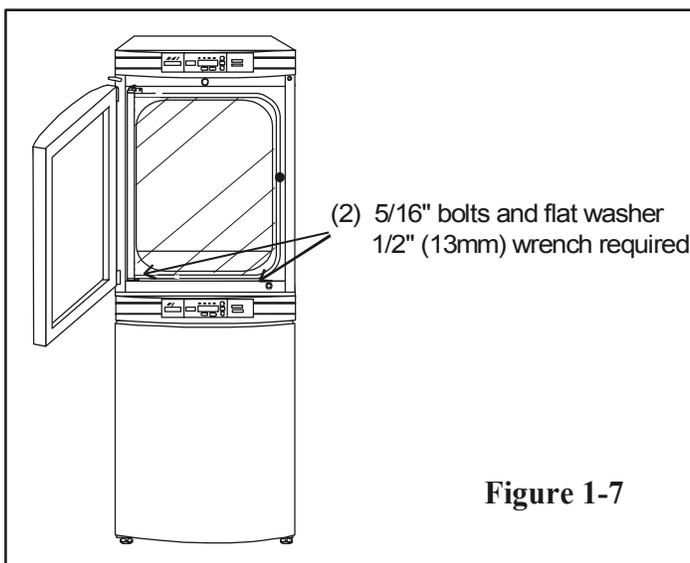
2. Note the two slots in the base of the incubator which accommodate the stacking bolts. Refer to Figure 1-5.
3. Remove the two plastic plugs from the bolt holes in the exterior top of the *bottom* unit. Install the 1/2" long 5/16-18 stacking bolts and washers into the bolt holes - do not tighten the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



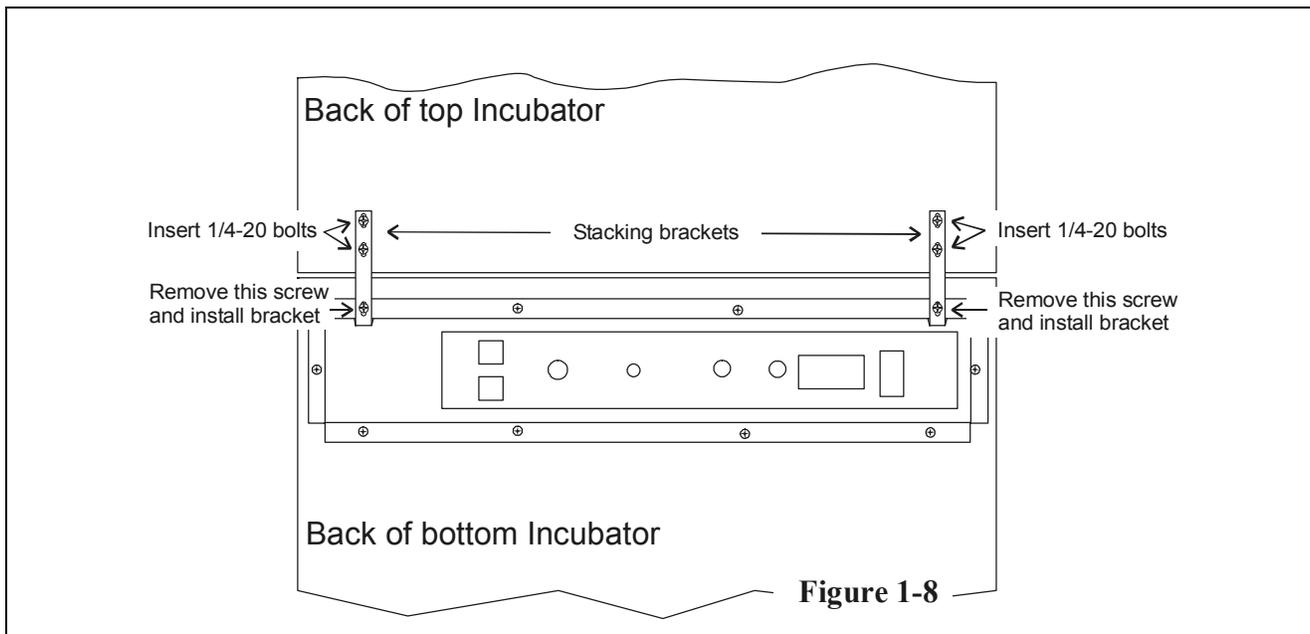
4. Unscrew and remove the leveling feet from the *top* unit and lift it onto the *bottom* unit, off-setting the base of the *top* unit approximately 2-3 inches behind the stacking bolts and washers.



This incubator weighs 265 lbs (120kg) before filling. Have sufficient personnel to lift it.



5. Align the sides of the *top* unit with the *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. Refer to Figure 1-7.
6. Remove and save the two screws from the back of the control panel on the *bottom* unit as identified in Figure 1.8.

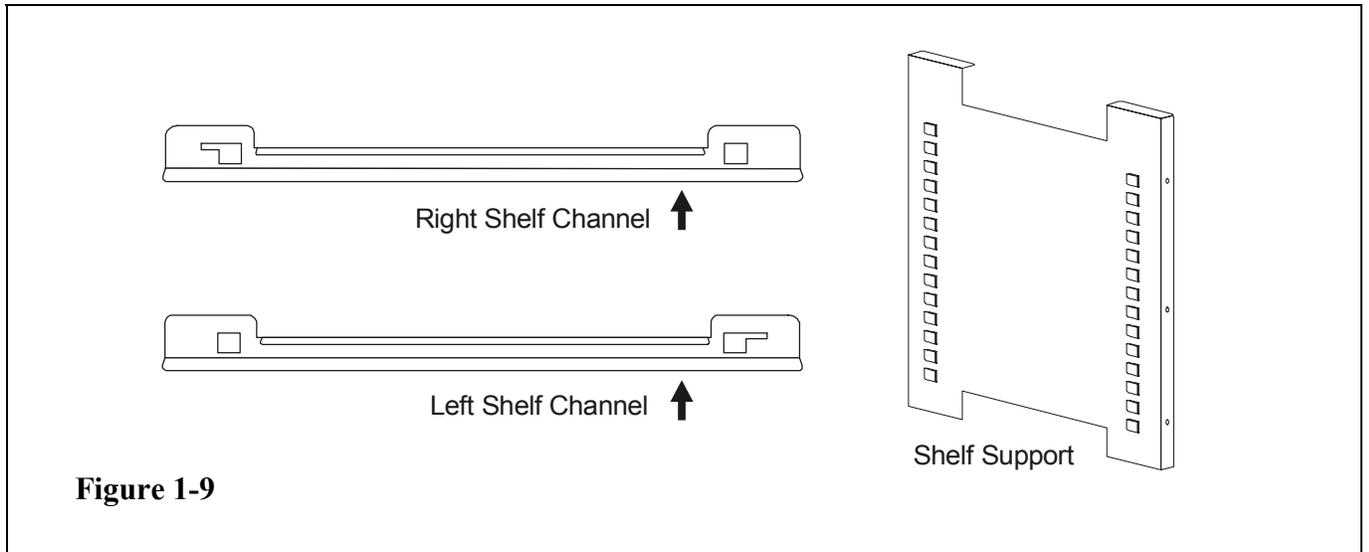


7. Insert the stacking brackets into the slots on the rear of the control panel of the *bottom* unit as shown in Figure 5-6. Align the slots in the brackets with the mounting holes on the rear of the incubators. Secure the brackets with the screws saved above and the 1/4-20 bolts provided in the stacking kit. A 7/16" wrench or socket will be 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" (13mm) wrench or suitable tool.
9. Replace the base cover on the *top* unit.
10. The stacked units are ready to be placed into service.

c. Preliminary Cleaning and Disinfecting

1. Remove the protective plastic coating on the shelf supports and air duct, if present.
2. Using a suitable laboratory disinfectant, disinfect all interior surfaces including shelves and shelf supports, door gaskets, blower wheel and CO₂ sensor. Refer to Section 5.1.

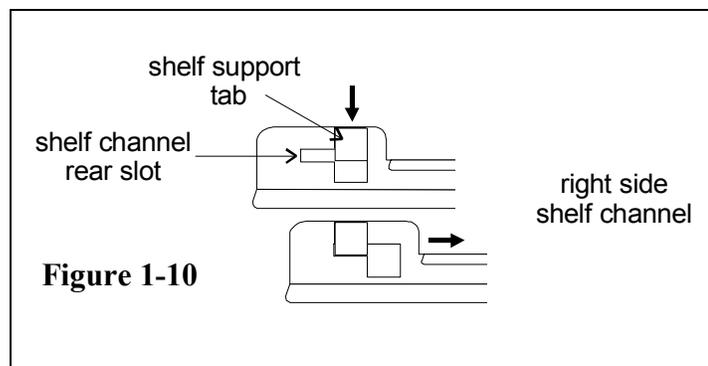
d. Installing the Shelves



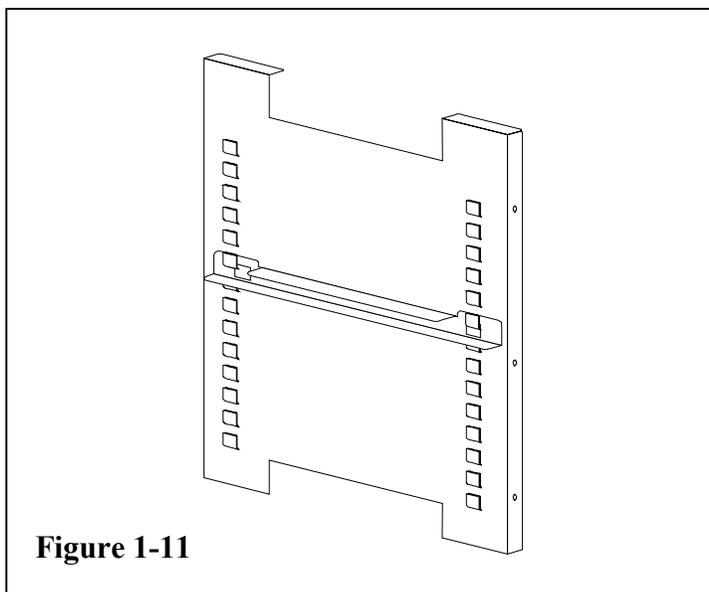
1. Install the shelf supports with the tabs facing into the center of the chamber with their slots up. There are no right side or left side supports, simply rotate one of them to fit the opposite side. Tilt the shelf supports as they are placed in the chamber so the tops fit into the top air duct, then guide them into the vertical position. Figure 1-9 shows the support as it would be oriented for the right side of the chamber.

2. Referring to Figure 1-9, note that there are left side and right side shelf channels.

3. Install the shelf channels by placing the channel's rear slot over the appropriate rear tab on the shelf support. Pull the shelf channel forward and engage the channel's front slot into the shelf support's appropriate forward tab. Refer to Figure 1-10.



- Figure 1-11 shows one of the channels installed on the right shelf support.



e. Installing the Access Port Stopper

Locate the opening in the top left corner of the interior chamber. Install the beveled end of the neoprene stopper in the opening.

f. Installing the HEPA Filter

- Remove the filter from the shipping box.
- Remove the plastic coating from the filter, using caution not to touch the filter media.
- Install the filter as shown in Figure 1-12.
- Refer to Section 5.6 for HEPA filter maintenance.



Use caution when handling the filter. The media can be damaged if it is mishandled.

To avoid damage to the incubator, do not operate the unit without the HEPA filter in place.

- Remove the filter from the shipping box.
- Remove the plastic coating from the filter using caution not to touch the filter media.
- Install the filter as shown in Figure 1-12.

Refer to Section 5.6 for HEPA filter maintenance.

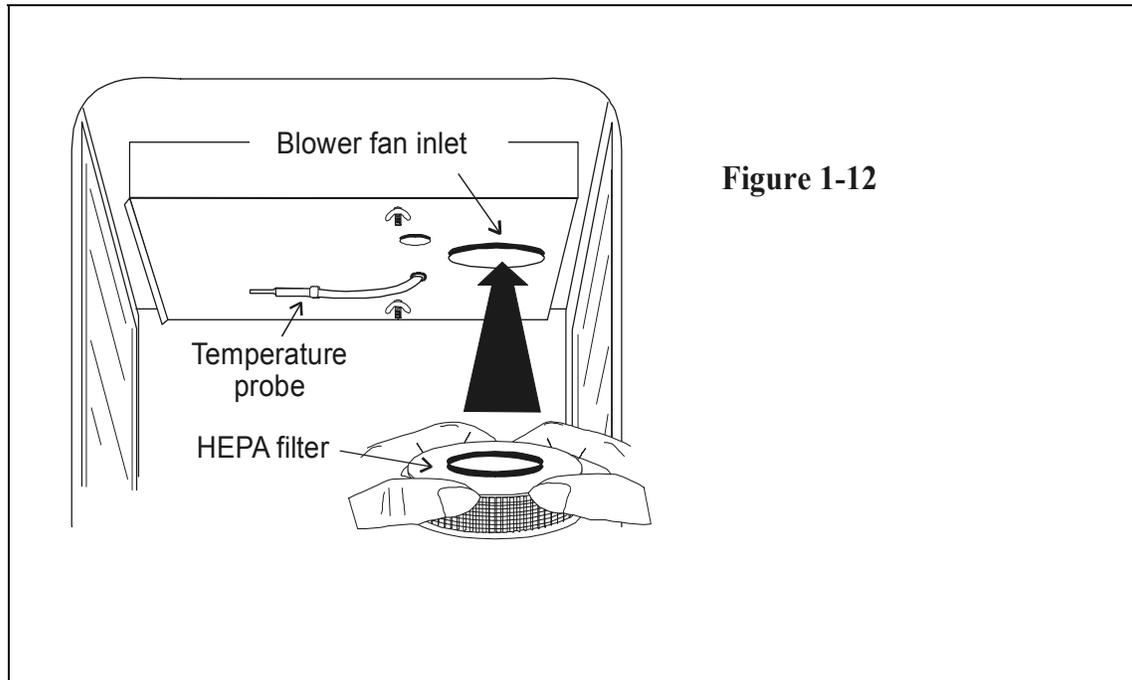


Figure 1-12

g. 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.

h. Connecting the Unit to Electrical Power

See the serial tag on the side of the unit for electrical specifications, or refer to the electrical schematics in Section 9 of this manual.



Connect the incubator to a grounded dedicated circuit only.

The power cord connector is the mains disconnect device for the incubator. Position the unit so that it can be easily disconnected.

Plug the provided power cord into the power inlet connector (See Figure 1-13) and into the grounded dedicated circuit.

Electrical Specifications

Models 3110, 3120, 3130, 3140 - 115V 50/60 Hz 3.6A 1 PH 2W

Models 3111, 3121, 3131, 3141 - 230V 50/60 Hz 2.0A 1 PH 2W

i. Filling the Water Jacket

Turn the power switch on. **ADD WATER** will appear in the message center. Press the Silence key to silence the alarm.

Note: The fill port has a plug that must be removed before filling and replaced after filling is complete.

Fill the water jacket with 11.7 gallons (43.5 liters) of water with a quality of 50K ohm to 1M ohm. Vinyl tubing and funnel for filling are included in the accessory bag shipped with the unit. Connect vinyl tubing to the fill port. 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 4.1, Table of Alarms.

j. Filling the Humidity Pan

For best operation of the incubator, sterilized distilled, demineralized or de-ionized water should be used in the humidity pan. Water purity should be in the resistance range of 50K Ohm/cm to 1M Ohm/cm, or a conductivity range of 20.0 uS/cm to 1.0 uS/cm. Refer to ASTM Standard D5391-93 or D4195-88 for measuring water purity.

Distillation systems, as well as some types of reverse osmosis water purity systems, can produce water in the quality range specified. Tap water is not recommended as it may contain chlorine, which can deteriorate the stainless steel. Tap water may also have a high mineral content, which would produce a build-up of scale in the reservoir. High purity, ultra pure or milli-q water is not recommended as it is an extremely aggressive solvent and will deteriorate the stainless steel. High purity water has a resistance of above 1M Ohm to 18M Ohm. Even high purity water can contain bacteria and organic contaminants. Water should always be sterilized or treated with a decontaminant, safe for use with stainless steel as well as safe for the product, prior to being introduced into the humidity pan.

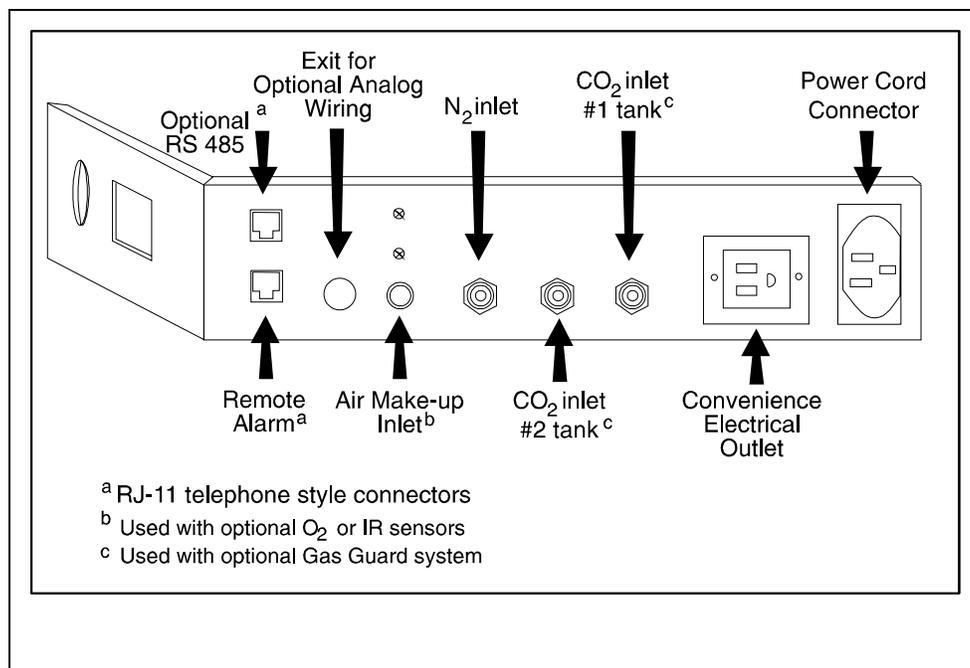


Distilled or de-ionized water used in the humidity pan must be within a water quality resistance range of 50K Ohm/cm to 1M Ohm/cm to protect and prolong the life of the stainless steel. Use of water outside the specified range will decrease the operating life of the unit and may void the warranty.

Fill the humidity pan with 0.8 gallons (3 liters) of sterile, distilled water. Place the pan directly on the incubator floor to ensure best humidity and temperature response.

Check the water level and change it frequently to avoid contamination. Do not allow the water level to fluctuate significantly. “Dry-outs” will have an adverse effect on the humidity level.

**Figure 1-13,
Rear Panel View
showing all options**



k. Connecting the CO₂ Gas Supply***Warning!******High concentrations of CO₂ gas can cause asphyxiation***

OSHA Standards specify that employee exposure to carbon dioxide in any eight-hour shift of a 40-hour work week shall not exceed the eight-hour time weighted average of 5000 PPM (0.5% CO₂). The short term exposure limit for 15 minutes or less is 30,000 PPM (3% CO₂). Carbon dioxide monitors are recommended for confined areas where concentrations of carbon dioxide gas can accumulate.

The CO₂ gas supply being connected should be industrial grade 99.5% pure and should not contain siphon tubes. Install a two-stage pressure regulator at the cylinder outlet. 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).

The incubator has serrated fittings on the back of the cabinet to connect the gas supply. Refer to Figure 1-13. The fitting is labeled CO₂ inlet #1 Tank. Make sure that the connections are secured with clamps. Check all fittings for leaks.

For units having the CO₂ Gas Guard option, refer to Section 6.2.

***Warning!***

This incubator is designed to be operated with CO₂ gas only. Connecting a flammable or toxic gas can result in a hazardous condition.

Gases other than CO₂ should not be connected to this equipment. CO₂ gas cylinders have UN1016 labeled on the cylinder and are equipped with a CGA 320 outlet valve. Check the gas cylinder for the proper identification labels. The CO₂ gas supply being connected to the incubator should be industrial grade, 99.5% pure. Do not use CO₂ gas cylinders equipped with siphon tubes. A siphon tube is used to extract liquid CO₂ from the cylinder, which can damage the pressure regulator. Consult with your gas supplier to ensure that the CO₂ cylinder does not contain a siphon tube. Gas cylinders should also be secured to a wall or other stationary object to prevent them from tipping.

A two-stage CO₂ pressure regulator is required to be installed on the outlet valve of the gas cylinder. Input pressure to the incubator must be maintained at 15 psig (103.4 kPa) for proper performance of the CO₂ control system. (A single stage CO₂ pressure regulator will not maintain 15 psig (103.4 kPa) to the incubator as the pressure in the CO₂ cylinder decreases; therefore, a two stage regulator is recommended.)

If higher purity CO₂ 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 CO₂ from the gas supplier. Follow the manufacturer's instructions to ensure proper and safe installation of the pressure regulator on the gas cylinder.

Consult your facility safety officer to ensure that the equipment is installed in accordance with the codes and regulations that are applicable in your area.

I. Connecting the N₂ Gas Supply

This connection applies only to those units that have an O₂ system. The N₂ gas supply being connected should be 99.99% pure. Do not use liquid nitrogen. Follow the same steps as in the section above for preparing the N₂ tank for hookup to the incubator. For units having the N₂ Gas Guard option, refer to Section 6.2. Connect the vinyl lines from the N₂ tank to the serrated fitting labeled N₂ Inlet and secure with the provided clamp. Check all fittings for leaks.

1.6 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, CO₂, and O₂. 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.

a. Setting the Operating Temperature

Incubator models 3110, 3111 have an operating temperature range of 10 to 55°C, models 3120, 3121 at 10 to 50°C, and models 3130, 3131, 3140, and 3141, at 10 to 45°C. All units require the cooling coil option to run at any temperature lower than 5°C above ambient. 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.X is displayed in the message center.
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.

b. Setting the Overtemp Setpoint



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.

The incubator is equipped with an independent circuit that monitors the air temperature in the cabinet. The independent overtemp circuit is designed as a safety for the incubator only. Should the system's temperature control fail, this circuit would cut out all heaters when the cabinet's temperature reaches the Overtemp setpoint. When an incubator is operating in an overtemp condition, the temperature control in the incubator will be $\pm 1^\circ\text{C}$ around the overtemp setpoint.

The overtemp's function is to prevent abnormally high temperatures that will occur if the heaters are locked on as a result of a failure in the main temperature control. Although the overtemp circuit will control the chamber temperature close to the overtemp setpoint, it is not intended to protect or limit the maximum temperature of the cell cultures or the equipment inside the chamber when the overtemp condition occurs.

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.

To set the Overtemp setpoint:

1. Press the Mode key until the Set indicator lights.
2. Press the right arrow until Otemp XX.X is displayed in the message center.
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.

c. Setting the CO₂ Setpoint

All T/C CO₂ cells are precalibrated at the factory at 37°C, high humidity, and 10% CO₂. Therefore, if a temperature setpoint of 37°C has been entered, the humidity pan filled, and the CO₂ control is to run between 0-10% with a T/C CO₂ sensor, the CO₂ setpoint may be entered immediately. Otherwise, it is important to allow the unit 12 hours to stabilize at the temperature setpoint before entering the CO₂ setpoint.

All models of the incubator have a CO₂ setpoint range of 0.0% to 20.0%. The incubator is shipped from the factory with a CO₂ setpoint of 0.0%. At this setting, all CO₂ control and alarms are turned off. To change the CO₂ setpoint:

1. Press the Mode key until the Set indicator lights.
2. Press the right arrow until CO₂ XX.X is displayed in the message center.
3. Press the up/down arrow until the desired CO₂ 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.

d. Setting the O₂ Setpoint

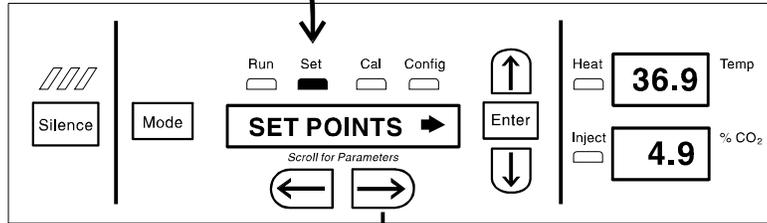
Models 3130, 3131, 3140, and 3141 of the incubator have a built-in O₂ control system. The O₂ setpoint range is 2.0% to 21.0%. The incubator is shipped from the factory with a O₂ setpoint of 21.0%. At this setting, all O₂ control and alarms are turned off. To change the O₂ setpoint:

1. Press the Mode key until the Set indicator lights.
2. Press the right arrow until O₂ XX.X is displayed in the message center.
3. Press the up/down arrow until the desired O₂ 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.

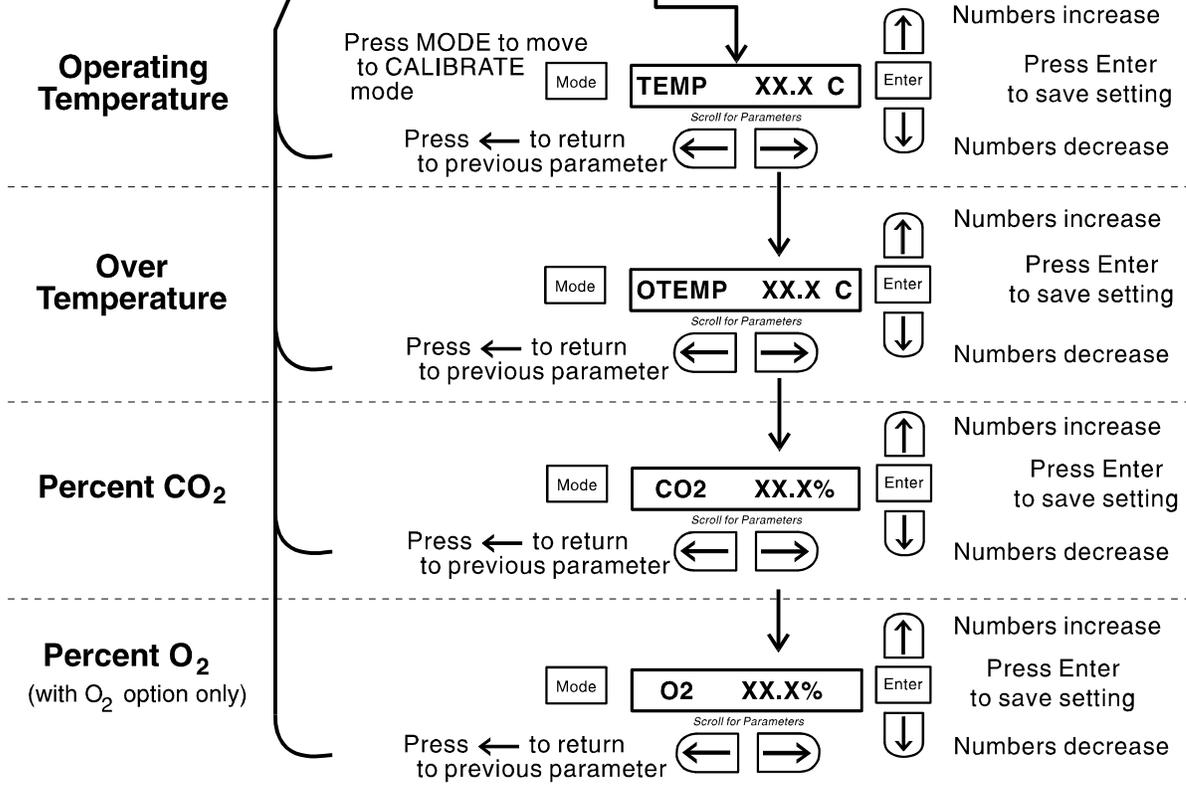
Chart 1-1

Set Mode

Press MODE to light SET



To Set:



Section 2 - Calibration

2.1 Calibration Mode

After the unit has stabilized, several different systems can be calibrated. In the Calibration Mode, the air temperature, CO₂ reading, O₂ 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 3-1 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 stabilization periods.

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.



Before making an calibration or adjustments to the unit, it is imperative that all reference instruments be properly calibrated.

a. 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.

Temperature Stabilization Periods

- **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 TEMPCAL XX.X appears in the message center.
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, or the right/left arrow to go to the next/previous parameter.

b. Calibrating the Thermal Conductivity CO₂ System

Models 3110, 3111, 3130 and 3131 have a thermal conductivity (T/C) CO₂ sensor. Thermal conductivity of the incubator atmosphere is not only effected by the quantity of CO₂ present, but also by the O₂%, the air temperature, and the quantity of water vapor present in the incubator atmosphere. In monitoring the effects of CO₂, O₂%, air temperature, and absolute humidity must be held constant so any change in thermal conductivity is caused only by a change in CO₂ concentration.

Changing temperature or O₂% set points or changing from elevated humidity levels to room ambient humidity levels would necessitate a recalibration of the CO₂ control. Models fitted with the O₂ control option, (Models 3130 and 3131) have an O₂ compensation feature. Refer to the instructions in Section 3.1.s of this manual, for enabling O₂ compensation for T/C systems. When this feature is enabled, the processor will automatically adjust the T/C CO₂ sensor for changes in the O₂ percentage.

New Installation. The CO₂ sensor has been calibrated at the factory for 37°C and elevated humidity. Allow the incubator to run at least 12 hours before checking CO₂ concentration with an independent instrument. If the incubator is operating at 37°C, a zero adjustment is all that is necessary.

1. Press the Mode key until CAL indicator lights.
2. Press the right arrow until CO2ZER0 XX.X is displayed in the message center.
3. Press the up/down arrow to change the display to match the measuring device.
4. Press Enter to store the calibration.
5. Press the Mode key to return to run or right/left arrow to go to next/previous parameter.

If the temperature setpoint is other than 37°, a zero and span adjustment is necessary. Never adjust span without first ensuring the zero is adjusted to ambient CO₂, which is approximately 0% CO₂.

c. Calibrating the T/C %CO₂ Zero to Ambient

On models with O₂ control (Models 3130 and 3131), you must disable the O₂ compensation before performing a complete CO₂ calibration. Check Section 3.1 Configuration Mode, Enabling O₂ Compensation for T/C CO₂ Systems. After completion of the CO₂ calibration, you may return to the Configuration Mode and enable the O₂ compensation feature.

T/C CO₂ Sensor Stabilization Periods

Start Up- Set CO₂ set point to 0% (which disables CO₂ valve) or shut CO₂ off at the supply tank. Allow temperature, humidity, and O₂ levels in the chamber to stabilize at least 12 hours before proceeding.

Operating Unit- Set CO₂ set point to 0% (which disables CO₂ valve) or shut off CO₂ at the supply tank. Open both doors and air-out chamber for at least 2 minutes. Close doors and allow temperature, humidity and O₂ levels to stabilize for at least 2 hours before proceeding.

1. Ensure stabilization periods as outlined above are followed.
2. Press the Mode key until the CAL indicator lights.
3. Press the right arrow until CO₂ ZERO XX.X is displayed in the message center.
4. Press the up/down arrow to adjust display to 0.0%.
5. Press Enter to store calibration.
6. Press the Mode key to return to Run Mode.

Adjust the CO₂ setpoint to the desired level. Turn on the CO₂ at the supply tank. Allow CO₂ to stabilize at least 15 minutes after CO₂ actual display reaches setpoint. Measure the chamber CO₂ content through the gas sample port with a Fyrite or other independent instrument. Several readings should be taken to ensure accuracy. If the display does not match the independent instrument reading, proceed to the CO₂ span adjustment.

d. Calibrating the T/C %CO₂ Span

To insure accurate calibration, the unit will not allow CO₂ to be spanned below 3%. If the cabinet does not contain at least 3.0% CO₂, increase the setpoint and allow the unit to stabilize before completing this procedure.

1. Ensure CO₂ zero adjustment has been properly followed before continuing.
2. Press the Mode key until the CAL indicator lights.
3. Press the right arrow until CO₂SPAN XX.X appears in the message center.
4. Press the up/down arrow to adjust the display to match the independent instrument reading.
5. Press Enter to store the calibration.
6. Press the Mode key to return to Run Mode.

e. Calibrating the Infra-Red CO₂ System.

Models 3120, 3121, 3140 and 3141 have an infra-red CO₂ sensor. Infra-red CO₂ sensors are not effected by chamber atmosphere temperature, humidity and O₂ variations. However, the light detector in the sensor is effected by wide temperature changes. *Therefore, changing temperature setpoints could necessitate a recalibration of the CO₂.* Chamber temperature should be allowed to stabilize before checking CO₂ concentrations with an independent instrument, especially on start-up.

All models equipped with an IR/CO₂ sensor have an automatic calibration that occurs every 24 hours, and lasts for 5 to 6 minutes. During automatic calibration, the CO₂ display is blanked out and HEPA filtered room air is pumped through the CO₂ sensor. A new CO₂ calibration value is stored in memory for use as the 0.0% CO₂ reference point. The keypad/ control panel is “locked up” during calibration, with the message center reading CO2 AUTO CAL.

IR CO₂ Sensor Stabilization Times

Startup- Allow the temperature and the CO₂ of the cabinet to stabilize at least 12 hours before proceeding.

Operating Unit - Allow CO₂ to stabilize at least 2 hours at set point before proceeding.

To insure accurate calibration, the unit will not allow CO₂ to be spanned below 3%. If the cabinet does not contain at least 3.0% CO₂, increase the setpoint and allow the unit to stabilize before completing this procedure.

1. Measure the CO₂ 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 center.
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 will last 5 to 6 minutes. The control panel is “locked up” during this calibration cycle.
7. Press the Mode key to return to Run Mode.

f. Calibrating the O₂ system.

If using an O₂ Fyrite, the accuracy of the instrument will be greatly affected by the concentration of CO₂ in the cabinet. Refer to the Fyrite operating manual.

Models 3130, 3131, 3140 and 3141 have an O₂ control sensor. The sensor is a fuel cell that puts out a linear millivolt signal based on O₂ content of the chamber. The fuel cell depletes over time depending on O₂ levels. The above models have an automatic calibration to maintain the O₂ calibration as the fuel cell depletes. The automatic calibration occurs every 12 hours and lasts 1 to 3 minutes. During automatic calibration, the O₂ display is locked on the existing O₂ reading and HEPA filtered room air is pumped through the O₂ sensor.

A new O₂ span value is stored in memory for use as the 20.7% O₂ reference point. The keypad/control panel is “locked-up” during automatic calibration.

O₂ Sensor Stabilization Times

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

Operating Unit - Allow O₂ to stabilize at least 2 hours at set point before proceeding.

Measure the O₂ concentration in the chamber through the gas sample port with an independent instrument. Several readings should be taken to ensure accuracy.

1. Press the Mode key until the CAL indicator lights.
2. Press the right arrow until O₂ CAL XX.X appears in the message center.
3. Press the up/downarrow to adjust display to independent instrument reading.
4. Press Enter to store calibration.
5. Press the Mode key to return to Run Mode.

g. Calibrating Relative Humidity

All 3110 Series 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.

Relative Humidity Stabilization Times

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 set point for relative humidity to stabilize before proceeding.

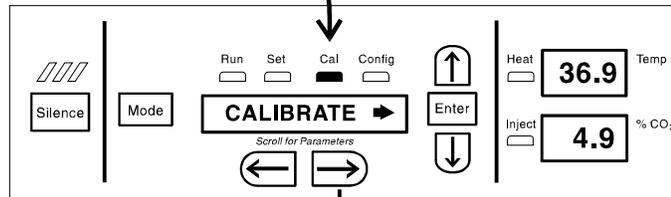
1. Place an accurate independent instrument in the center of the chamber. Allow at least 30 minutes for RH to stabilize.
2. Press the Mode key until the CAL indicator lights.
3. Press the right arrow key until RH CAL XX appears in the message center.
4. Press the up/down arrow to match the display to the independent instrument.
5. Press Enter to store the calibration.
6. Press the Mode key to return to Run Mode.

If a reliable RH measuring device is not available, you may calibrate the display to a typical level. On units with O₂ control, the nitrogen must be shut off as the incoming nitrogen lowers the optimum achievable level of RH in the chamber. Follow the RH stabilization periods outlined above. With a full humidity pan and stable temperature, the relative humidity in the chamber will be 95%. In step 5 of the relative humidity sensor adjustment, you may adjust the display to 95%. This calibration method should be accurate to within 5%.

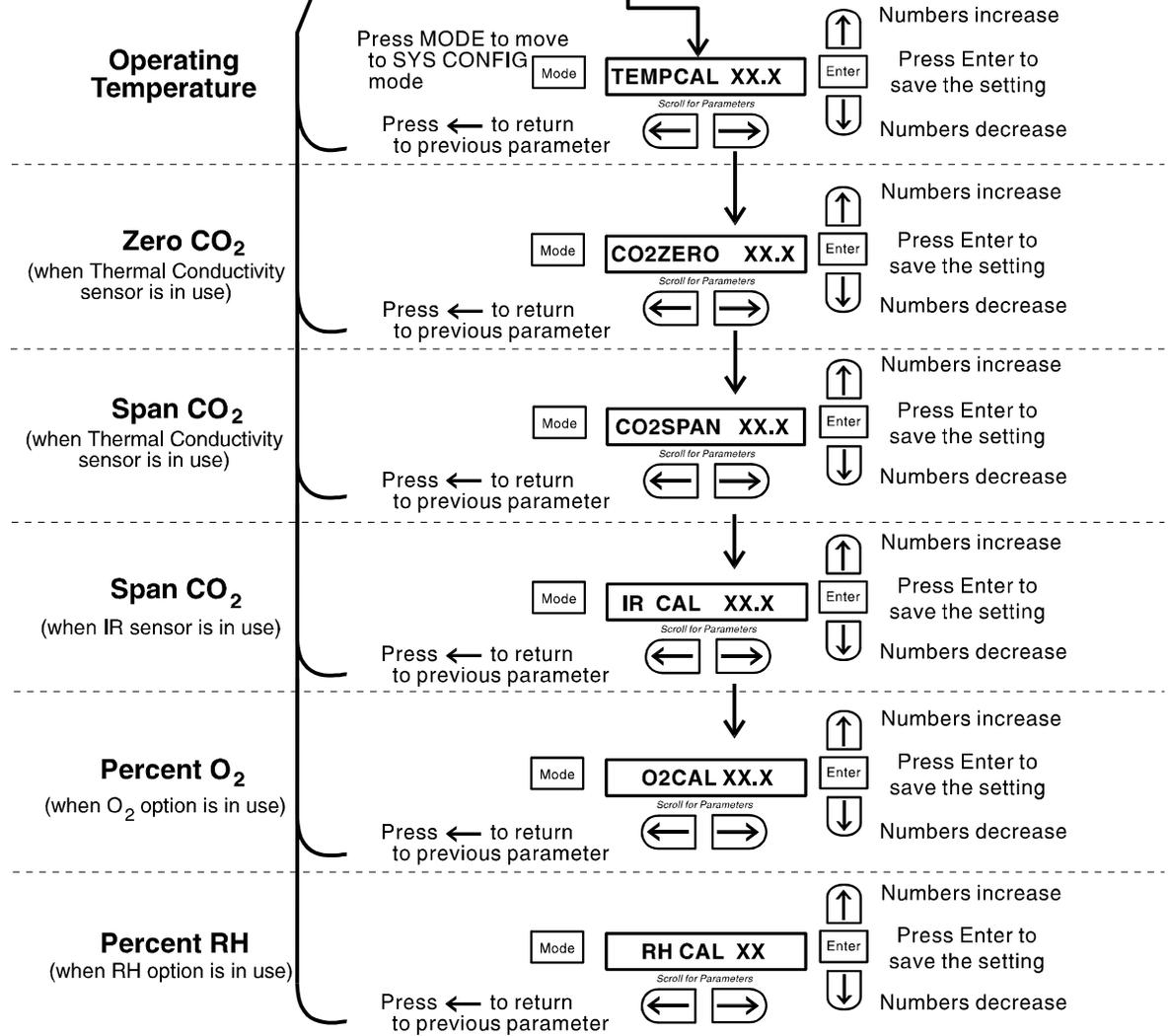
Chart 2-1

Calibrate Mode

Press MODE to light CAL



To Calibrate:



Section 3 - Configuration

3.1 Configuration Mode

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 Config indicator lights. Press the right or left arrow until the appropriate parameter appears in the message center. See Chart 3-1 for more detail.

a. 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 Config indicator lights.
2. Press the right arrow until Audible XXX is displayed in the message center.
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.

b. 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 Config indicator lights.
2. Press the right arrow until Acc Code XXX is displayed in the message center.
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.

c. Setting a Low Temp Alarm Limit (tracking alarm)

The low temp alarm limit 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 Config indicator lights.
2. Press the right arrow until Temp Lo -X.X is displayed in the message center.
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.

d. Setting a High Temp Alarm Limit (tracking alarm)

The high temp alarm limit 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 Config indicator lights.
2. Press the right arrow until Temp Hi X.X is displayed in the message center.
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.

e. Enabling Temperature 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 Config indicator lights.
2. Press the right arrow until Tmp Rly XXX is displayed in the message center.
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.

f. Setting a Low CO₂ Alarm Limit (tracking alarm)

The low CO₂ alarm limit is the deviation from the CO₂ setpoint that will cause a low CO₂ alarm. The setpoint is variable from 0.5% CO₂ below setpoint to 5.0% CO₂ below setpoint. The factory setting is 1.0% CO₂ below setpoint. A minus (-) in the display indicates that the alarm setting is below the setpoint.

1. Press the Mode Key until the Config indicator lights.
2. Press the right arrow until CO2 Lo -X.X is displayed in the message center.
3. Press the up/down arrow to change the low CO₂ alarm limit.
4. Press Enter to save the low CO₂ alarm limit.
5. Press the Mode key to return to Run Mode or right/left to go to next/previous parameter.

g. Setting a High CO₂ Alarm Limit (tracking alarm)

The high CO₂ alarm limit is the deviation from the CO₂ setpoint that will cause a high CO₂ alarm. The setpoint is variable from 0.5% CO₂ above setpoint to 5.0% CO₂ above setpoint. The factory setting is 1.0% CO₂ above setpoint.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until CO2 Hi X.X is displayed in the message center.
3. Press the up/down arrow to change the high CO₂ alarm limit.
4. Press Enter to save the high CO₂ alarm limit.
5. Press the Mode key to return to Run Mode or right/left to go to next/previous parameter.

h. Enabling CO₂ Alarms to Trip Contacts

High and Low CO₂ alarms can be programmed to trip the remote alarm contacts. A setting of ON will cause this; a setting of OFF will not allow CO₂ alarms to trip the contacts. The factory setting is ON.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until CO2 Rly XXX is displayed in the message center.
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.

i. Setting a New Zero Number for T/C CO₂ Sensors

If a new T/C CO₂ sensor is being installed, the two numbers on the factory installed sticker on the T/C cell must be entered to calibrate the CO₂ in the unit.

NOTE: For the technician's convenience, a label containing the two numbers on the T/C cell is affixed inside the electronics drawer.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until T/CZ# XXXX is displayed in the message center.
3. Press the up/down arrow to change the zero 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.

j. Setting a New Span Number for T/C CO₂ Sensors

If a new T/C CO₂ sensor is being installed, the two numbers on the factory installed sticker on the T/C cell must be entered to calibrate the CO₂ in the unit.

NOTE: For the technician's convenience, a label containing the two numbers on the T/C cell is affixed inside the electronics drawer.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until T/CS# XXXX is displayed in the message center.
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.

k. 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 will disable the alarm.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until RH Lo XX is displayed in the message center.
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.

l. 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 Config indicator lights.
2. Press the right arrow until RH Rly XXX is displayed in the message center.
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.

m. New O₂ Cell

On models equipped with the O₂ control option, the unit's microprocessor monitors the life of the O₂ sensor. When the O₂ sensor needs replaced, an alarm will sound and the message REPL O₂ SNSR will be displayed. Once a new O₂ sensor is installed, this menu will allow it to be automatically calibrated. To perform the automatic calibration:

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until NEW O₂ CELL is displayed in the message center.
3. Press Enter to start calibration.
4. Calibration will take 3-5 minutes. Once completed, press the Mode key to return to Run Mode.

n. Setting a Low O₂ Alarm Limit (tracking alarm)

On models with a O₂ control system, O₂ alarms may be configured. The low O₂ alarm limit is the deviation from the O₂ setpoint that will cause a low O₂ alarm. The setpoint is variable from 0.5% O₂ below setpoint to 5.0% O₂ below setpoint. The factory setting is 1.0% O₂ below setpoint. A minus (-) in the display indicates that the alarm setting is below setpoint.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until O₂ Lo -X.X is displayed in the message center.
3. Press the up/down arrow to change the low O₂ alarm limit.
4. Press Enter to save the low O₂ alarm limit.
5. Press the Mode key to return to Run Mode or right/left to go to next/previous parameter.

o. Setting a High O₂ Alarm Limit (tracking alarm)

On models with a O₂ control system, O₂ alarms may be configured. The high O₂ alarm limit is the deviation from the O₂ setpoint that will cause a high O₂ alarm. The setpoint is variable from 0.5% O₂ above setpoint to 5.0% O₂ above setpoint. The factory setting is 1.0% O₂ above setpoint.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until O₂ Hi X.X is displayed in the message center.
3. Press the up/down arrow to change the high O₂ alarm limit.
4. Press Enter to save the high O₂ alarm limit.
5. Press the Mode key to return to Run Mode or right/left to go to next/previous parameter.

p. Enabling O₂ Alarms to Trip Contacts

On models with an O₂ control system, O₂ alarm contacts may be configured to trip the contacts. A setting of ON will cause this, a setting of OFF will not allow O₂ alarms to trip the contacts. The factory setting is ON.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until O₂ Rly XXX is displayed in the message center.
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.

q. 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 Config indicator lights.
2. Press the right arrow until Disp Tmp XXX or Disp RH XXX is displayed in the message center.
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.

r. Enabling CO₂/O₂ to be Displayed

On models that are equipped with the O₂ system, the lower seven-segment display on the control panel can be configured to display CO₂ continuously, O₂ continuously, or toggle between CO₂ and O₂. If the units does not have O₂, the lower display will always display CO₂. If CO₂ is set to ON and the O₂ is set OFF, CO₂ will be displayed continuously. If CO₂ is set to OFF and O₂ is set to ON, O₂ 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 O₂ system is present.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until Disp CO2 XXX or Disp O2 XXX is displayed in the message center.
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.

s. Enabling O₂ Compensation for T/C CO₂ Systems

On models that have T/C CO₂ control and a O₂ control system, the T/C CO₂ reading is susceptible to O₂ levels in the chamber. The incubator is capable of compensating the CO₂ display for the varying O₂ level. A setting of ON will allow such compensation. The factory setting is OFF.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until O₂ Comp XXX is displayed in the message center.
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.

t. Setting a RS485 Communications Address (1535 compatible only)

On units that have the RS485 option, direct communication with the Forma 1535 Model alarm system can be established. Each piece of equipment connected to the 1535 must have a unique address. An address of 0-24 can be entered for the incubator. A setting of 0 is an invalid address that the 1535 will ignore. The factory setting for the RS485 address is 0.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until RS485 XX is displayed in the message center.
3. Press the up/down arrow to move the RS485 address.
4. Press Enter to save the RS485 address.
5. Press the Mode key to return to Run mode or right/left to go to next/previous parameter.

u. Selecting a Primary Tank with the Gas Guard Option

On units equipped with the Gas Guard option, a primary tank can be selected. The primary tank will be either Tank 1 or Tank 2. The factory setting is Tank 1.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until Tnk Sel X is displayed in the message center.
3. Press the up/down arrow to toggle the setting between 1 and 2 .
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.

v. Enabling the Gas Guard System

On units equipped with the Gas Guard option, the Gas Guard system may be turned ON or OFF if it is not in use. The factory setting is OFF.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until Gas Grd XX is displayed in the message center.
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.

w. Setting the Door Heat

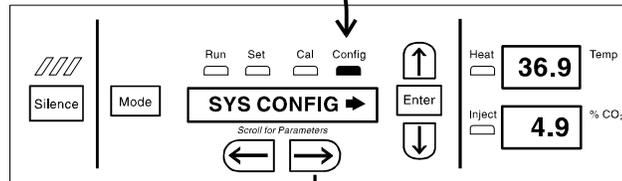
The incubator is equipped with a built-in heater on the inner glass door. The percentage of time that power is applied to this heater as compared to the main heater can be varied from 0 to 99%. The factory setting is 50%. Changing this setting can affect the unit's temperature control and uniformity. Consult the factory before changing this setting.

1. Press the Mode key until the Config indicator lights.
2. Press the right arrow until Dr Heat XX is displayed in the message center.
3. Press the up/down arrow to move the door heat setpoint.
4. Press Enter to save the door heat setpoint.
5. Press the Mode key to return to Run mode or right/left to go to next/previous parameter.

Chart 3-1

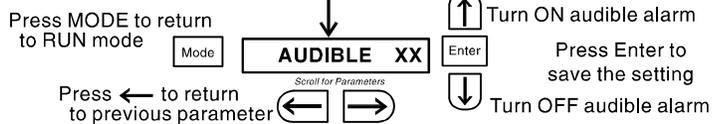
Configure Mode

Press MODE to light CONFIG

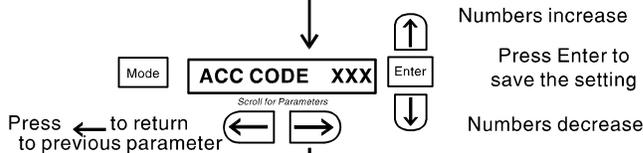


To Configure:

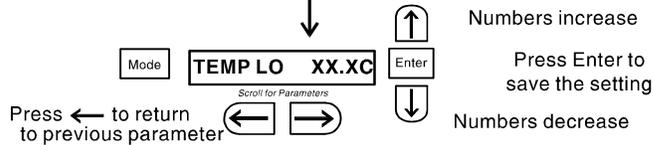
Audible ON/OFF



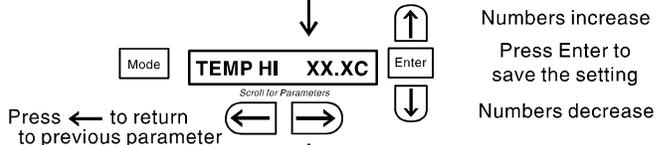
Access Code



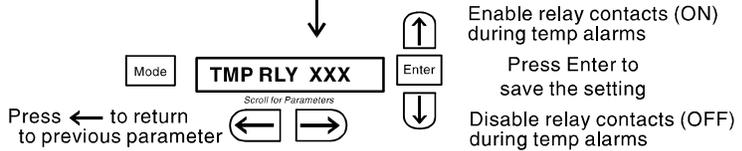
Low Temp Alarm Limit



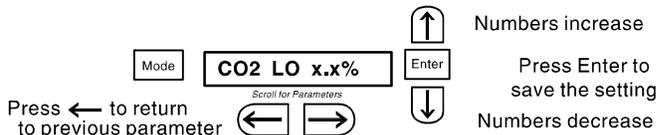
High Temp Alarm Limit



Temp Relay ON/OFF

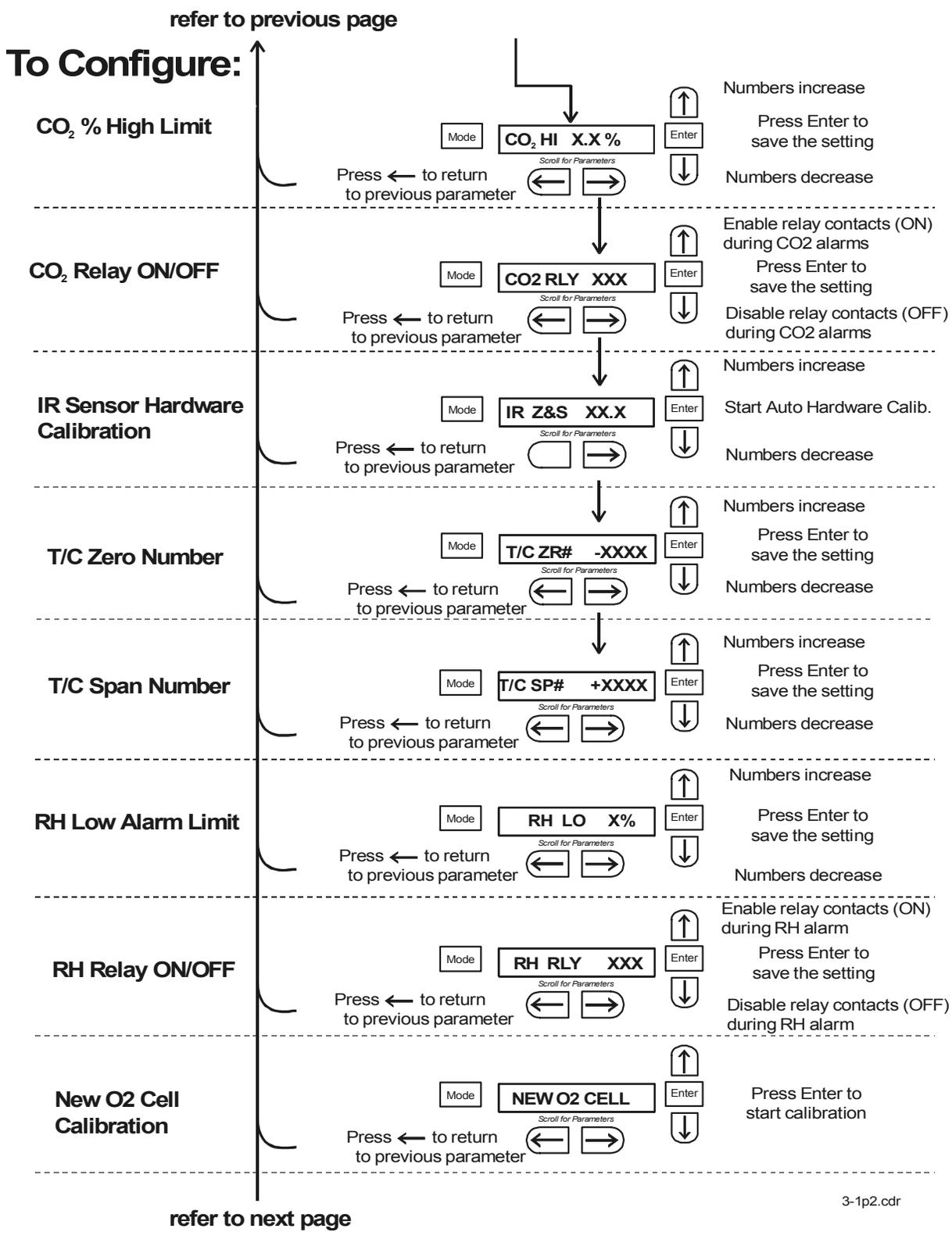


CO₂ % Low Limit

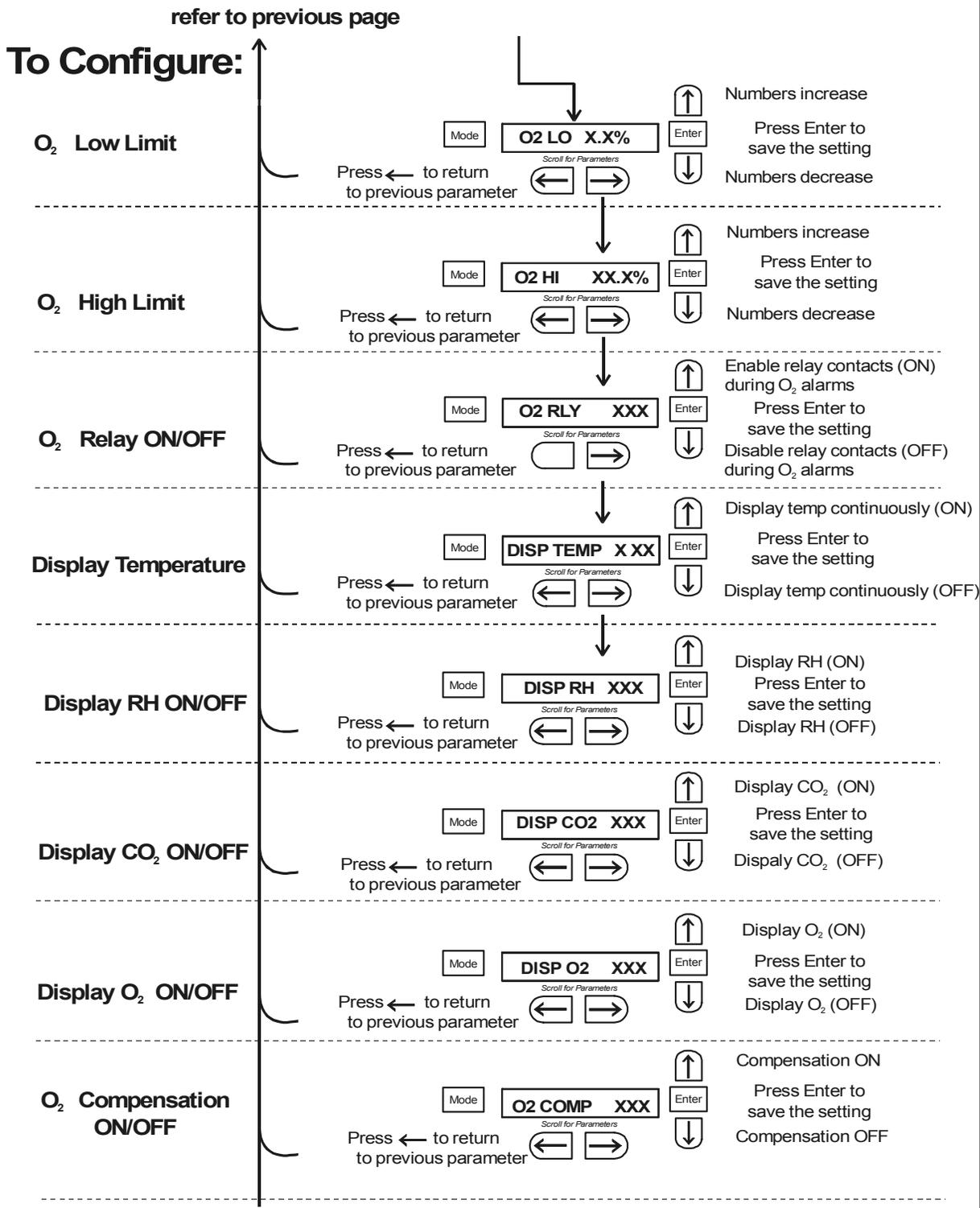


refer to next page

Configure Mode, Chart 3-1, Page 2 of 4



Configure Mode, Chart 3-1, Page 3 of 4

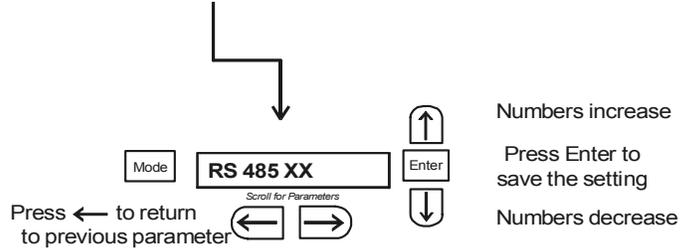


Configure Mode, Chart 3-1, Page 4 of 4

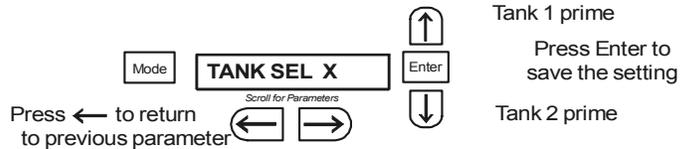
refer to previous page

To Configure:

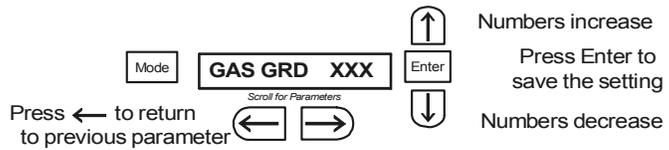
RS485



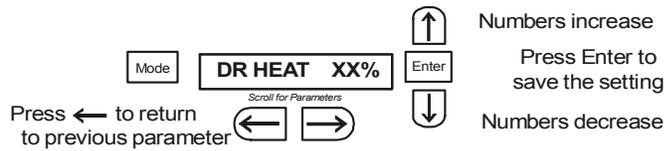
Tank Select



Gas Guard ON/OFF



Percent Door Heat



Section 4 - Alarms

4.1 Alarms

The Model 3110 Series incubator alarm system is shown in the table below. When an alarm is active, the message appears in the LED message center. 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 message center.

Description	Message	Delay	Ringback	Relay
No alarm condition exists	CLASS 100	----	----	----
CO2 System Auto Calibrating	CO2 AUTO CAL	----	----	----
Temp > Otemp Setpoint	SYS IN OTEMP	0 min.	15 min.	Yes
Water Temp Sensor Fault (See Sect 4.2)	H2O SENS ERR	0 min.	15 min.	No
Air Temp Sensor Fault (See Sect 4.2)	AIR SENS ERR	0 min.	15 min.	No
CO2 Sensor Fault (See Sect 4.5)	CO2 SENS ERR	0 min.	15 min.	No
O2 Sensor Fault (O2 option only) (See Sect 4.2)	O2 SENS ERR	0 min.	15 min.	No
O2 Sensor Low (O2 option only) (See Sect 4.3)	REPL O2 SNSR	15 min.	15 min.	No
0CO2 Sensor cannot be calibrated (IR option only) (See Sect. 4.4)	REPL IR SENSR	0 min.	15 min.	No
Water low in jacket	ADD WATER	0 min.	15 min.	No
Inner Door is Open	DOOR 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.	
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.	
O2 > O2 High Tracking Alarm (O2 option only)	O2 IS HIGH	15 min.	15 min.	Programmable ↓
O2 < O2 Low Tracking Alarm (O2 option only)	O2 IS LOW	15 min.	15 min.	
RH < RH Low Limit Alarm (RH option only)	RH IS LOW	30 min.	30 min.	Programmable
O2 Auto-Zero Fault (See Sect 4.2)	O2 AUTOZ ERR	0 min.	15 min.	No
CO2 Auto-Zero Fault (IR option only) (See Sect 4.6)	IR AUTOZ ERR	0 min.	15 min.	No
Tank 1 is low, switch to Tank 2 (Gas Guard only)	TANK1 LOW	0 min.	----	No
Tank 2 is low, switch to Tank 1 (Gas Guard only)	TANK2 LOW	0 min.	----	No
Both tanks are low (Gas Guard only)	TANK 1 and 2 LOW	0 min.	15 min.	No

- All alarm delays and ringback times are +/- 30 seconds -

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

The temperature alarms are disabled when the Temp setpoint is 10°C. The CO₂ alarms are disabled when the CO₂ setpoint is 0.0%. The O₂ alarms are disabled when the O₂ setpoint is 21.0%.

4.2 Sensor Fault Alarms

The microprocessor in all 3110 series incubators continually scans all available sensors including the pump on O₂ units to ensure that they are operating properly. Should an error be detected, the incubator will sound an alarm and display the appropriate message. If such an alarm occurs, contact your local distributor or the Forma Scientific Service department at 740-373-4763 or 1-888-213-1790 (USA and Canada) or fax 740-373-4189.

4.3 REPL O2 SNSR (Alarm)

On units equipped with the O₂ system, the microprocessor continually checks the remaining life of the O₂ sensor. If the O₂ sensor functionality is declining, the REPL O2 SNSR alarm will sound. This alarm alerts the user of the need to replace the O₂ sensor at the earliest convenience. The unit will continue to function for some length of time. To order a new sensor, contact your local distributor or the Forma Scientific Service department at 740-373-4763 or 1-888-213-1790 (USA and Canada) or fax 740-373-4189.

After the O₂ sensor is replaced, see Section *New O₂ Cell*, Section 3.1.m on page 3-5 of this manual for calibrating the O₂ system.

4.4 REPL IR SNSR (Alarm)

On units equipped with an IR CO₂ control system, calibration is done automatically using an Auto Zero system. If this system cannot properly calibrate the sensor, the REPL IR SNSR alarm will sound. The unit will continue to function for some length of time. To order a new sensor, contact your local distributor or the Forma Scientific service department at 740-373-4763 or 1-888-213-1790 (USA and Canada) or fax 740-373-4189.

4.5 CO2 SENS ERR

If the cables or connectors between the main microprocessor board and the CO₂ sensor, or between the CO₂ board and the sensor head on I/R CO₂ units become loose or disconnected, the CO2 SENS ERR alarm will occur.

I/R Units

On I/R incubators, the red light on the I/R module (Refer to Figure 5-4 for the location of the module circuit board) will be lit continuously if communication is lost between the CO₂ board and the sensor head. The CO₂ display will also be locked at 00.0 without injection. Turning the incubator off and on *will not* clear the alarm. Only proper connections of all the components will correct the alarm.

4.6 IR AUTOZ ERR

On incubators equipped with I/R CO₂ control, calibration is done automatically using an Auto Zero system. Auto Zero occurs once every 24 hours. If, during the Auto Zero cycle, a CO₂ correction of more than 0.45% is detected, the IR AUTOZ ERR alarm will occur.

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 the air pump and the sensor
 - Plugged filter or orifice on the 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 CO₂ will correct the alarm.
- There is high background CO₂ in the laboratory. This could be from leaks in the tank, regulator or vinyl CO₂ tubing.
- High CO₂ sensor calibration drift occurred. This will require replacement of the sensor.

Section 5: Routine Maintenance



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.

5.1 Disinfecting the Incubator Interior



If the units have been in service, disconnect the power cord connector before disinfecting.

Use an appropriate disinfectant. All articles and surfaces to be disinfected must be thoroughly cleaned, rinsed and roughdried.



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.



Do not spray the T/C sensor with flammable solutions. The internal temperature of the CO2 sensor is approximately 150°C when the unit is in operation. Allow sufficient time for the sensor to cool before cleaning.



Do not use strong alkaline or caustic agents. Stainless steel is corrosion resistant, not corrosion proof.

Do not use solutions of sodium hypochlorite (Purex, Clorox, etc.), as they may also cause pitting and rusting.

1. Turn off the incubator power switch and unplug the power cord.
2. Remove the shelves, HEPA filter, access port stopper, left and right duct sheets, and the top duct. The top duct requires the two wing nuts to be removed.
3. Set the HEPA filter aside. Wash the shelves, ducts, wingnuts and stopper with disinfectant and rinse with sterile water. Option: Autoclave shelves, ducts and wingnuts.
4. Remove the blower scroll by first pushing the white lever clip closest to you toward the scroll. Then turn the scroll to the left to disengage it from the top plate. Some manipulation may be required as the alignment holes are keyhole-shaped.
5. Remove the blower wheel by pulling down firmly on the wheel. Remove the sample port filter. If a new blower wheel and scroll are going to be used, discard the old ones. If the old ones are being reused, wash all parts with disinfectant and rinse with sterile water.
6. 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.
7. Install the blower wheel and scroll. Make sure the wheel turns freely. Align and turn the scroll to the right to lock it into the keyholes. Pull the white lever clip toward you, away from the scroll. Install the sample filter.
8. Install the top duct, left and right ducts, and access port stopper, spraying each with 70% alcohol.
9. Install the HEPA Filter
10. Install the shelves and spray with 70% alcohol.
11. Allow the unit to run empty for 24 hours before returning to service.

5.2 Cleaning the Cabinet Exterior

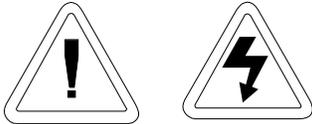
Clean the incubator exterior with a damp sponge or soft, well-wrung cloth and mild detergent dissolved in water. Dry with a soft cloth.

5.3 Cleaning the Humidity Pan

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

5.4 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 screwdriver and a flatblade screwdriver.

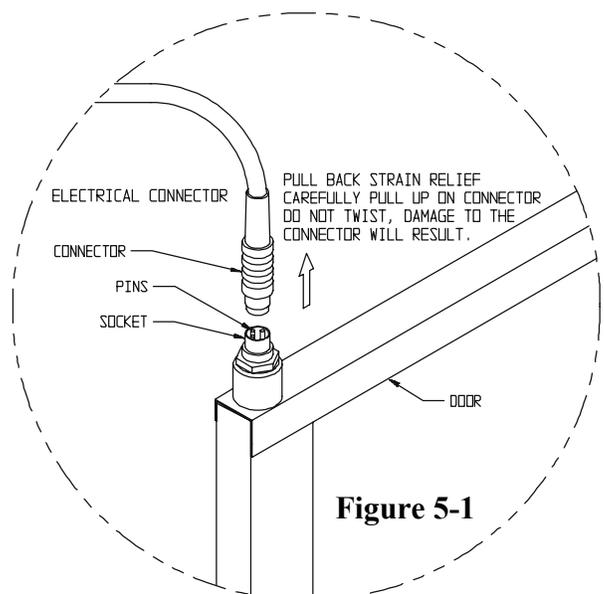


If the units have been in service, disconnect the power cord connector before reversing the door swing.

a. Reversing the Hinges for the Exterior Door

Review Illustration A on Page 5-6. The following instructions will refer to the letters within the circles.

1. Open the outer door and remove it by lifting it off its hinge pins. Lay this door down on its face on a padded surface to prevent scratches.
2. Remove the two outer door hinges identified by the “A” in the illustration.
3. Remove the four nylon screws in locations “B”, which 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, noted as locations “C” 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. Refer to item “D” in the illustration and the drawing in Figure 5-1. The heater wiring connector is of yellow rubber which should be visible when the strain relief is moved upward as shown. ***Refer to the connector manufacturer’s instructions in Appendix A of this manual.***
6. Grasp the upper portion of the yellow connector and pull it up and out of the hinge socket.



7. Remove the inner door upper hinge, shown as “E”. This hinge will be inverted and become the lower inner door hinge when the hardware is reassembled. When removing the door, set it aside on a flat surface, taking care not to damage the electrical sockets on the top and bottom of the door frame. 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.



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 before moving on to the next step.

8. Remove the lower inner door hinge, “F”. Note that 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 the step above.
9. The shoulder of the bushing must be on top as shown in Illustration B. Install the black plastic bushing into the hinge just removed and which will be the top hinge when the door is reversed. Refer also to Figure 5-2 on the next page.
10. Remove the two nylon screws at location “G”.
11. Remove the door strike plate at location “H” and install the two nylon screws just removed into the vacant holes.
12. Remove the two outer Phillips screws on the frame beneath the control panel, identified as “I” in the illustration. Remove only the two outer screws and pull the electronics tray out about one-inch.
13. Refer to Illustration B on Page 5-7:
14. Verify that the nylon screws have been installed in the vacant door strike holes, identified as “J” .
15. Move the door heater cable to the slot on the right side of the tray as shown in location “K”.
16. Push any excess cable into the tray area, making sure the wire will not be damaged when the tray is pushed back in and secured.

17. Install the lower inner door hinge, identified as location “L” in the drawing. (This was the top inner door hinge removed in Step 7 and should have the white plastic bushing installed.) Do not completely tighten the screws.
18. Insert the inner door into the lower hinge with the latch to the left. Align the door to the chamber opening and place the upper hinge in position as identified as “M”.

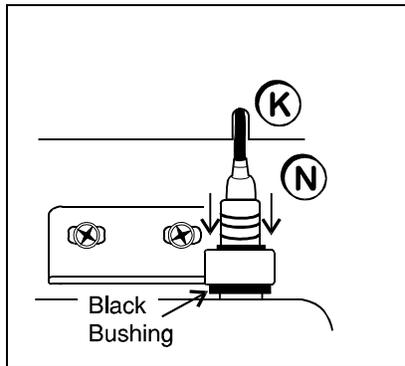


Figure 5-2

19. Insert the screws into the hinge but do not completely tighten them.
20. 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 as shown in “N” of Figure 5-2.
21. Push the power cable completely into the slot in the electronics tray. Press the tray into position and secure it with the screws at the “S” locations.
22. Align the silver power buss visible through the glass along the hinge side of the door with the gasket along the right side of the chamber opening.
23. 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.
24. Attach the strike plate at location “O”, aligning it so the knob secures the door against the gasket.
25. Attach the outer door hinges at the “P” locations.
26. Install the nylon screws at the “R” locations.
27. Install the nylon screws in the “Q” locations.
28. Assemble the outer door to the incubator and return the unit to service.

Left Hinge Door

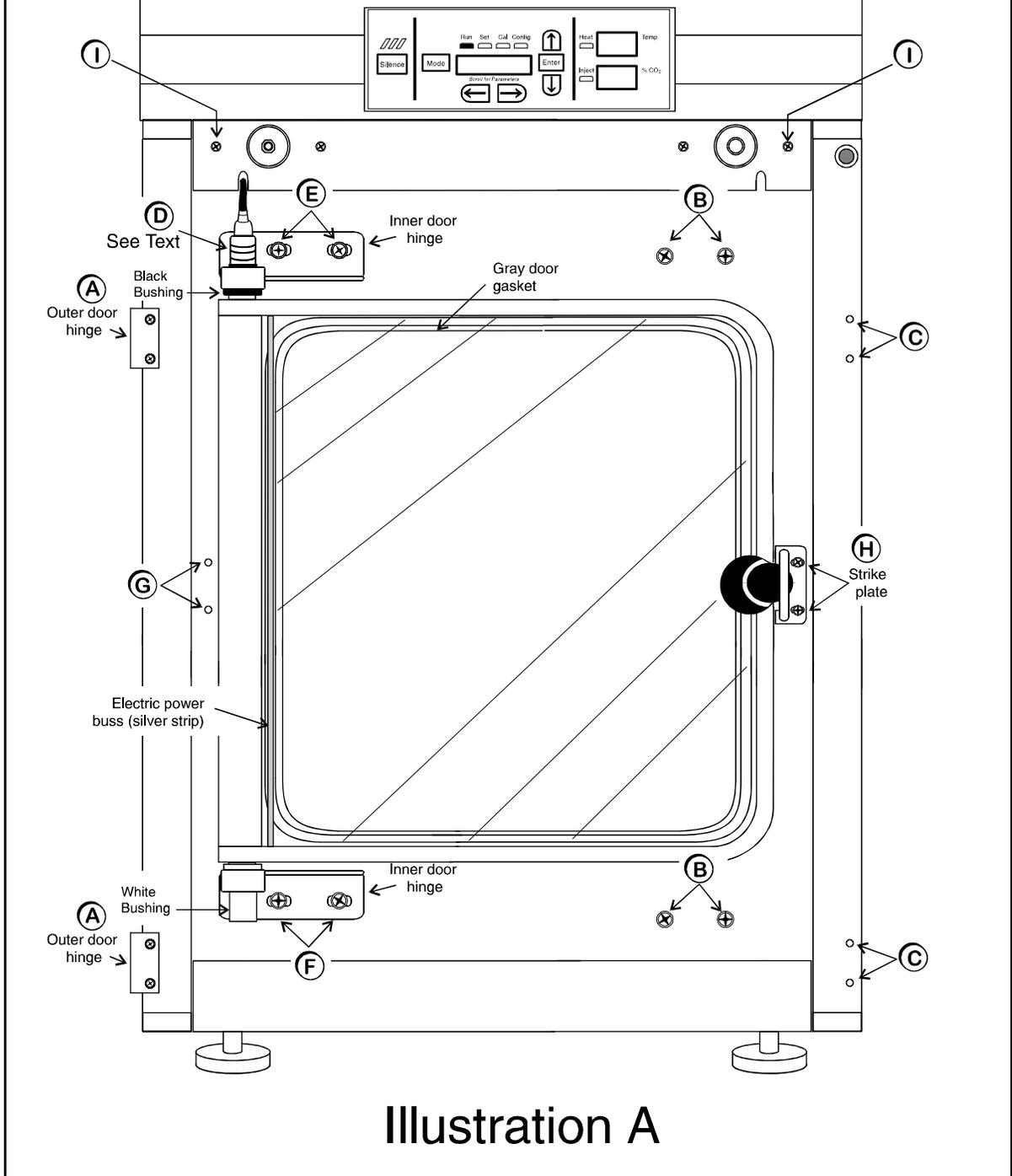


Illustration A

Right Hinge Door

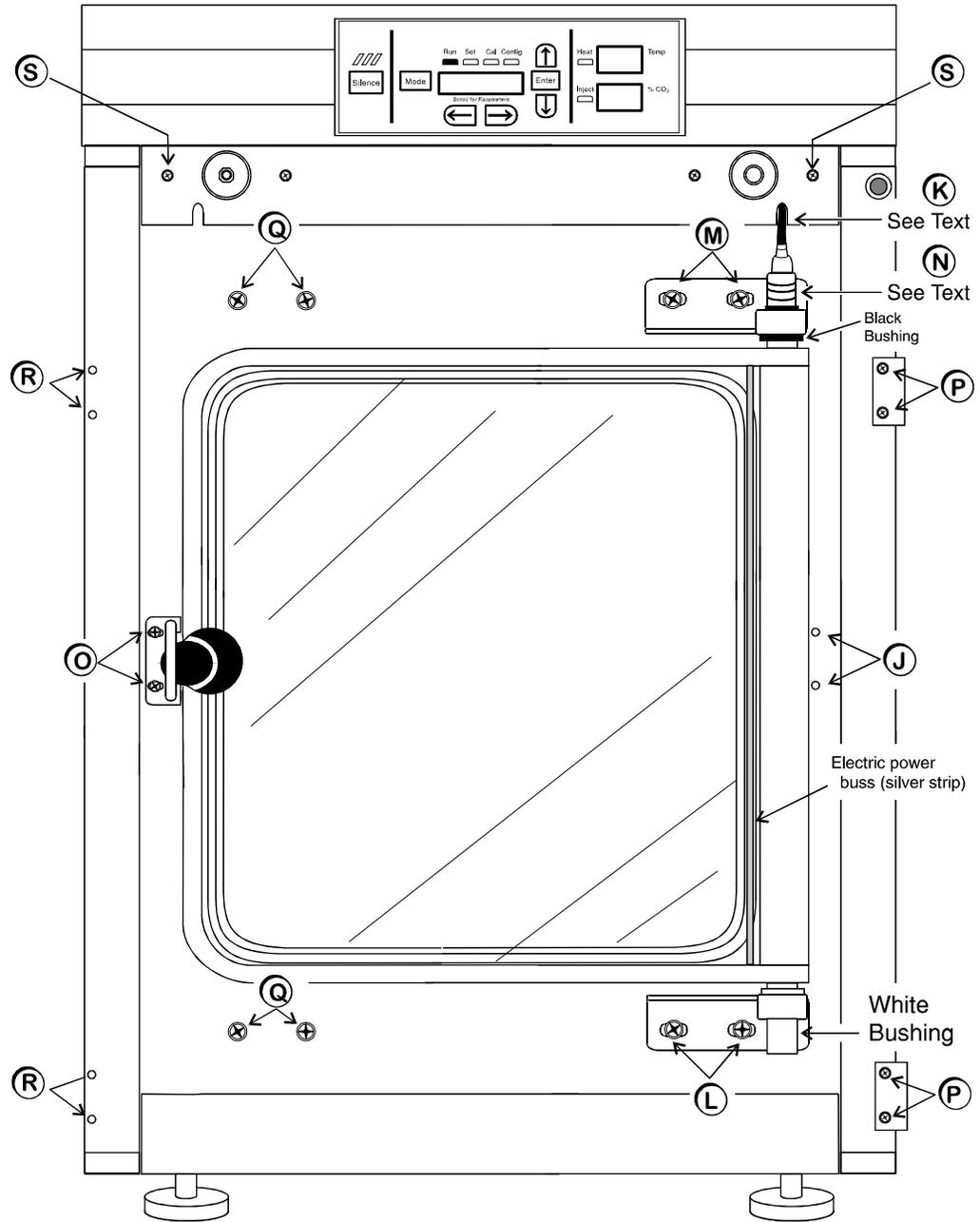
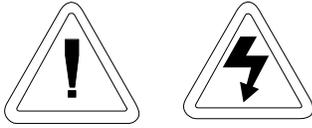


Illustration B

5.5 Replacing Fuses



***The electronics drawer contains hazardous voltages.
Replacing the fuses should be done by qualified personnel only.***

There are three fuses in the incubator that can be replaced. To replace a fuse:

1. Turn off the incubator's power switch.
2. Unplug the power cord from the wall outlet.
3. Open the exterior cabinet door.
4. Remove two screws as shown. See Figure 5-3.
5. Carefully slide out electronics drawer.

There are two fuses on the main microboard. They are labeled: F1 24 VAC Door Heater
F4 24 VAC Collar Heater

To locate the main microprocessor board, refer to Figure 5-4. Remove the appropriate fuse and replace it with a new fuse of the same specification. Refer to Table 5-1.

The other replaceable fuse is the accessory outlet fuse which is mounted to the floor of the electronics drawer. To locate the fuse, refer to Figure 5-10, and locate the accessory outlet fuse. Remove the fuse and replace with a new fuse of the same specification. Refer to Table 5-1 for the fuse specifications.

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.
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.

Table 5-1, Fuse Replacement Chart

Fuse #	Manufacturers 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

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 5-4. If this fuse is blown, the power supply must be replaced. Contact the factory for more information.

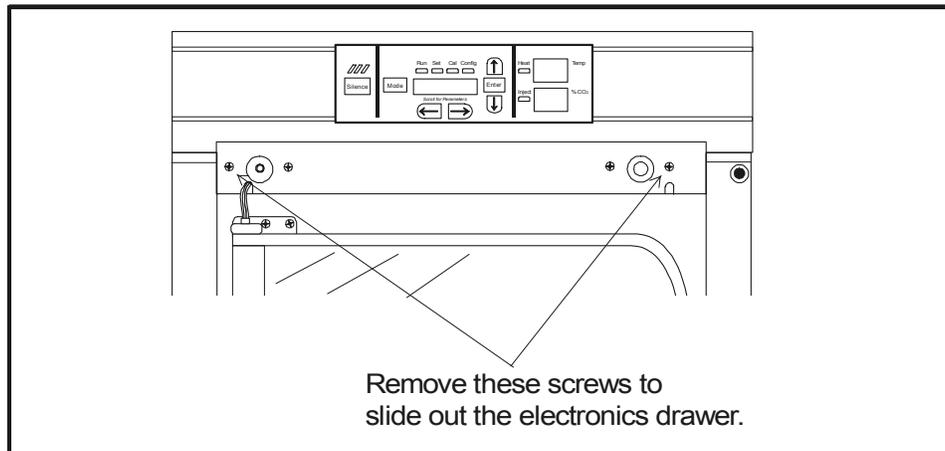


Figure 5-3

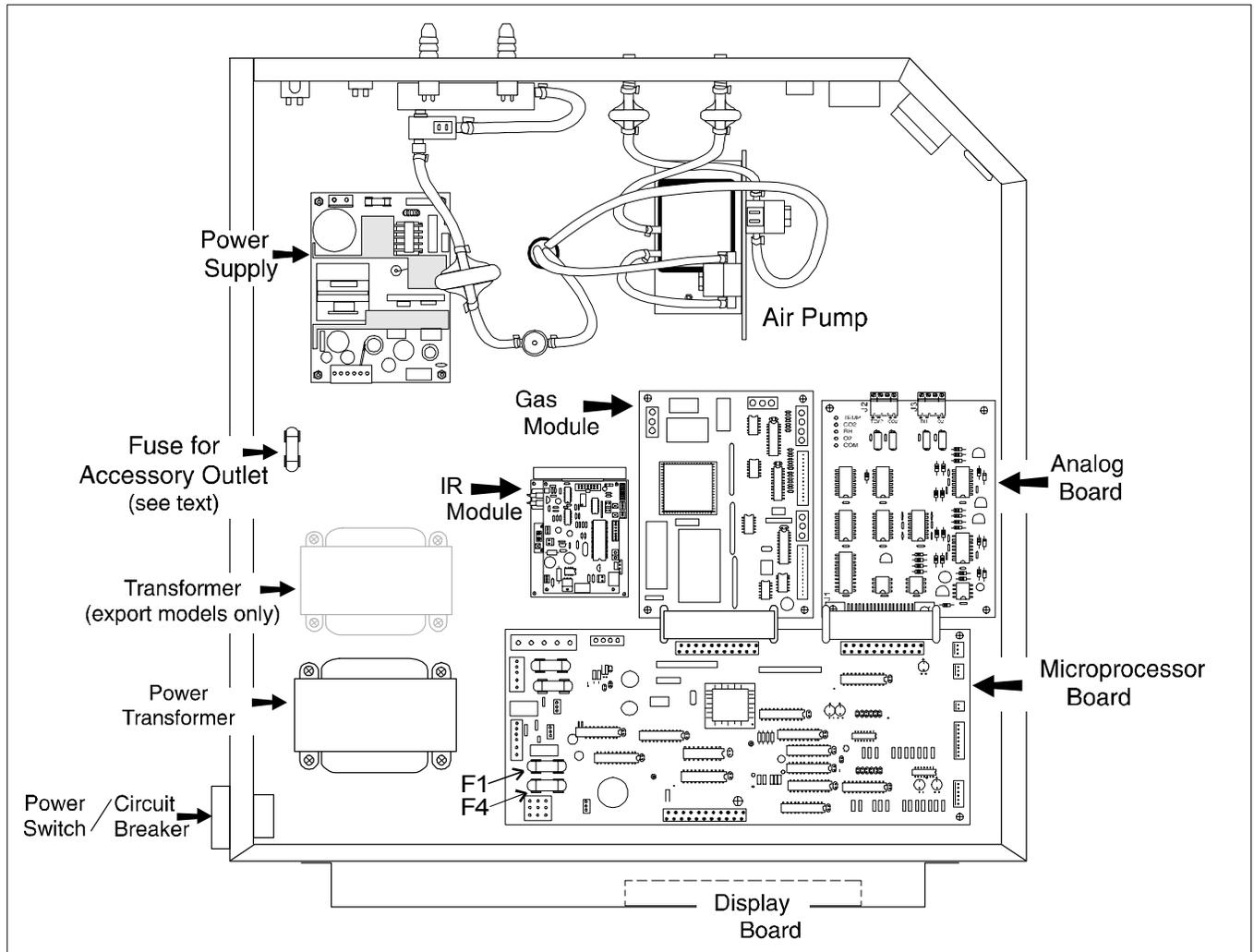


Figure 5-4
Model 3110/3210 Series Electronics Drawer

5.6 HEPA Filter Maintenance

Replace the HEPA filter routinely on an annual basis, or sooner if it changes color. The filter should also be changed each time the chamber is cleaned.

5.7 Draining Water Jacket

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 5-5.

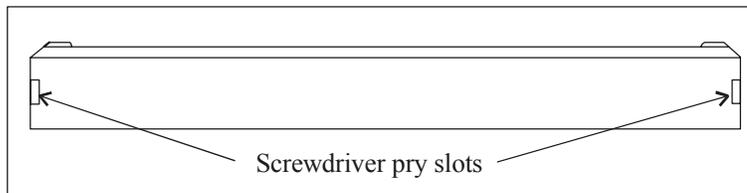


Figure 5-5

Front Cover Plate Below the Door

3. Remove the drain plug. Retain for use after draining is complete.
4. Connect the hose barb insert to the drain on lower front of the water jacket and to the drain hose.

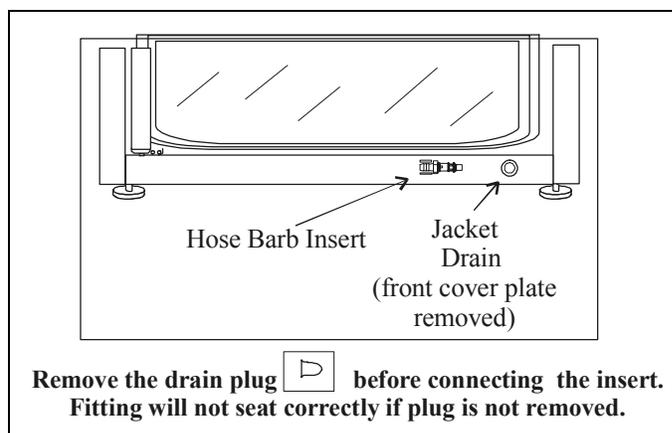


Figure 5-6

Location of Water Jacket Drain and Hose Barb Insert

5. After water jacket has finished draining, remove the hose barb insert and secure on the front of the unit. See Figure 5-7.
6. Install the plug into the drain on the incubator.
7. Install cover plate.

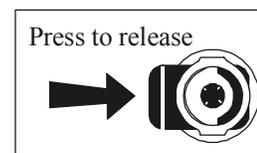


Figure 5-7

Section 6 - Factory Installed Options

6.1 Connections to External Equipment

a. Connecting the 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).

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, CO₂ alarms, O₂ alarms, and RH alarms. See Section 3.1, System Configuration.

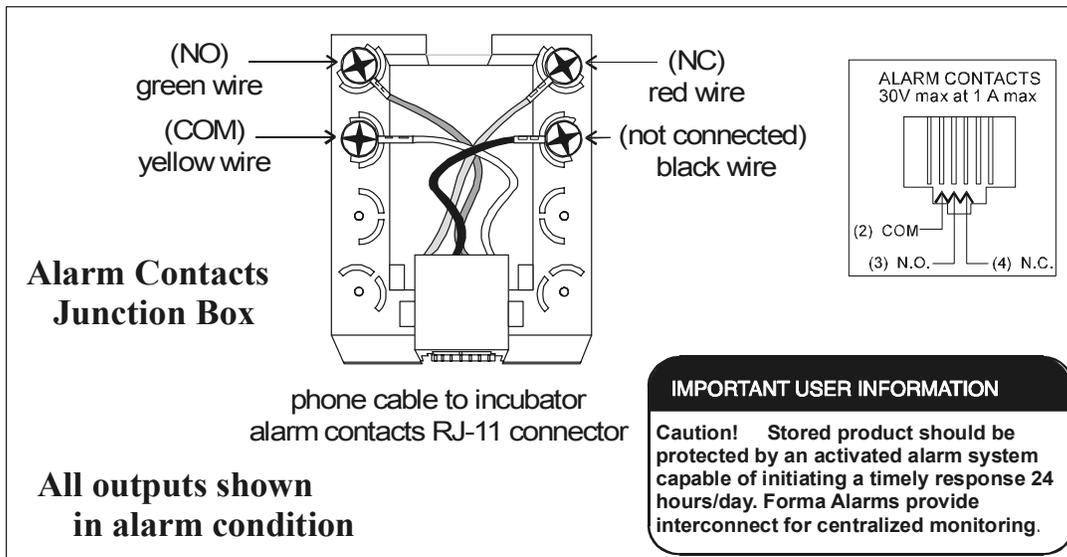
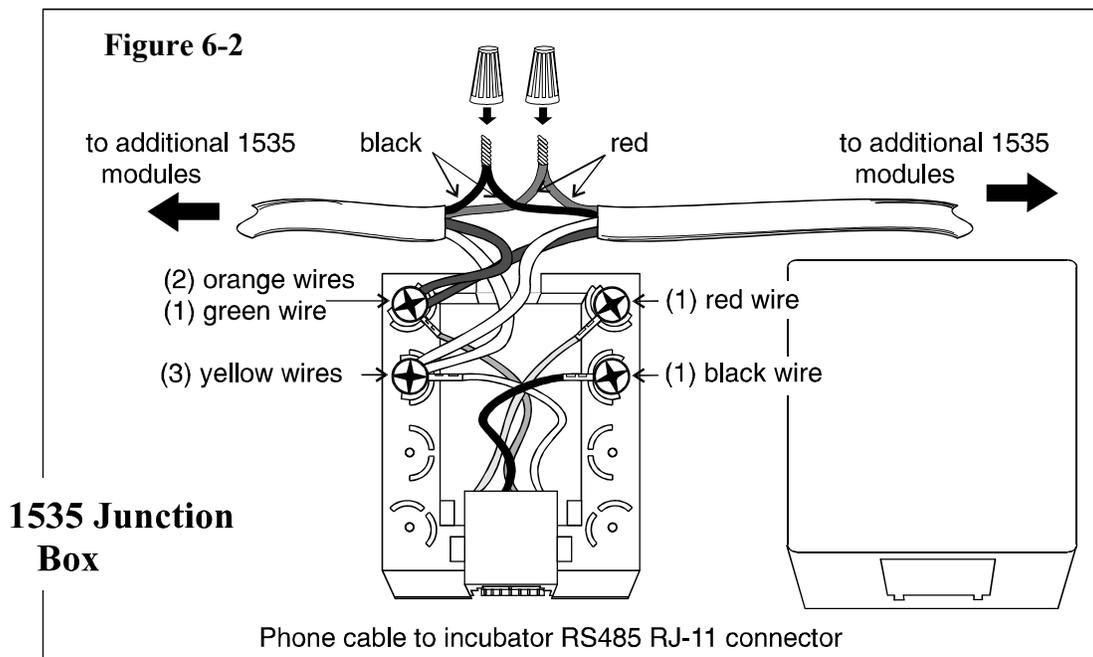


Figure 6-1

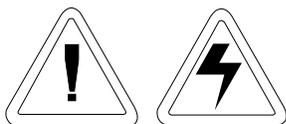
b. Connecting the RS485 Interface (190523)

All incubator models can be purchased with the RS485 communications option. This option allows the incubator to be directly connected to a Forma Model 1535 alarm system without the use of a communications module. A junction box is provided with each RS485 option. Refer to Figure 6-2 for wiring details.

To allow the incubator and the 1535 to communicate, an address must be allocated on the 1535. Refer to Section 5.8 of the 1535 operating manual. The same address number must be assigned to the incubator. Refer to Section 3.1 of this manual, System Configuration.



c. Connecting the Analog Output Boards (190512, 190543, 190544)



The electronics drawer contains hazardous voltages. Opening the drawer and/or wiring in an analog board should be done by qualified personnel only.

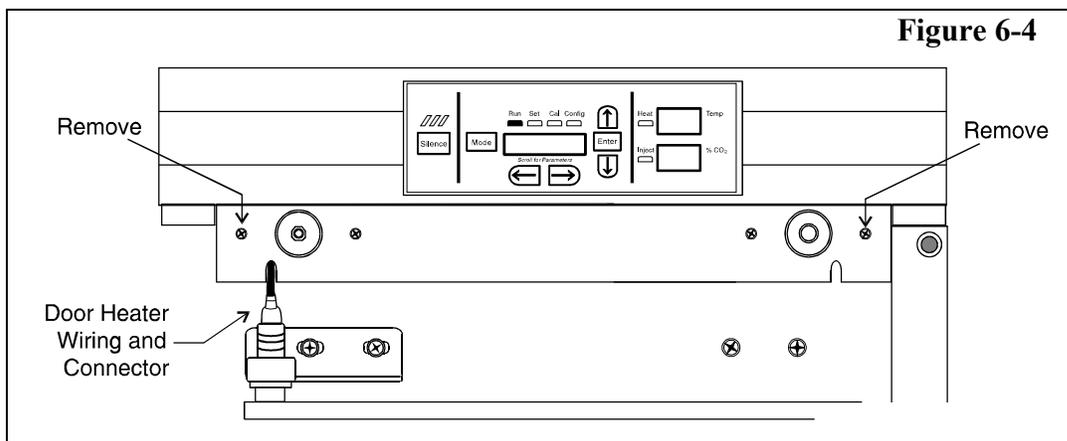
The analog output board is an option that allows the incubator to output analog signals representing the air temperature of the cabinet, the CO₂ content, the O₂ content, and the relative humidity, depending upon which systems are in the incubator. There are three different analog output board options available : 0-1V, 0-5V, or 4-20mA signals. Negative display readings will output 0V. The outputs do not have isolated grounds. Refer to Figure 6-3 for output specifications of the three boards.

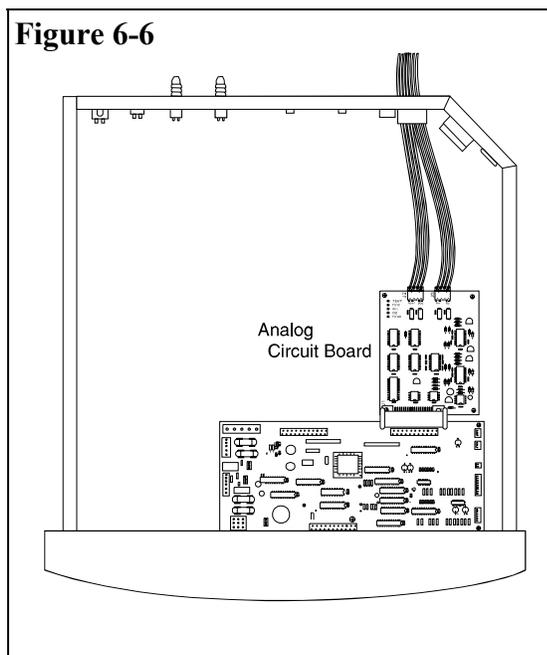
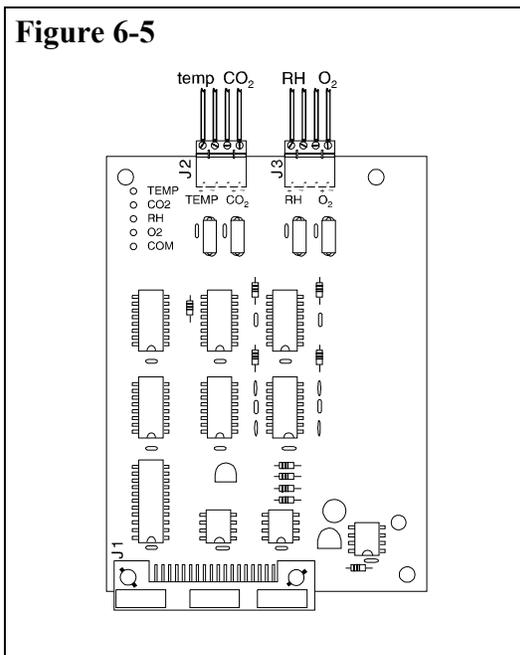
Figure 6-3 Analog Output Board Specifications

	190512 4-20mA Output Scaling 4-20 mA Equals	190544 0-1V Output Scaling 0-1 V Equals	190543 0-5V Output Scaling 0-5V Equals
Temperature	0.0-100°C	0.0-100°C	0.0-100°C
RH	0-100%RH	0-100 % RH	0-100 % RH
CO ₂	0.0-100.0% CO ₂	0-100.0 % CO ₂	0-100.0 % CO ₂
O ₂	0.0-100.0% O ₂	0-100.0 % O ₂	0-100.0 % O ₂

To wire in the analog output board, Forma Scientific recommends wire Part # 73041, which is a 22-gauge, 3-conductor wire with a shield. This is readily available from other vendors including Alpha Part #2403, and Deerborn Part # 972203.

1. Turn off the incubator’s power switch and unplug the power cord from the wall outlet.
2. Open the exterior cabinet door and remove the two screws shown in figure 6-4.
3. Carefully slide out electronics drawer and locate the Analog Output board. (Figure 6-5)
4. Each system monitored (Temp, CO₂, O₂, RH) requires two conductors. Feed the wire through the analog wiring inlet on the back of the drawer. See Figure 6-6.



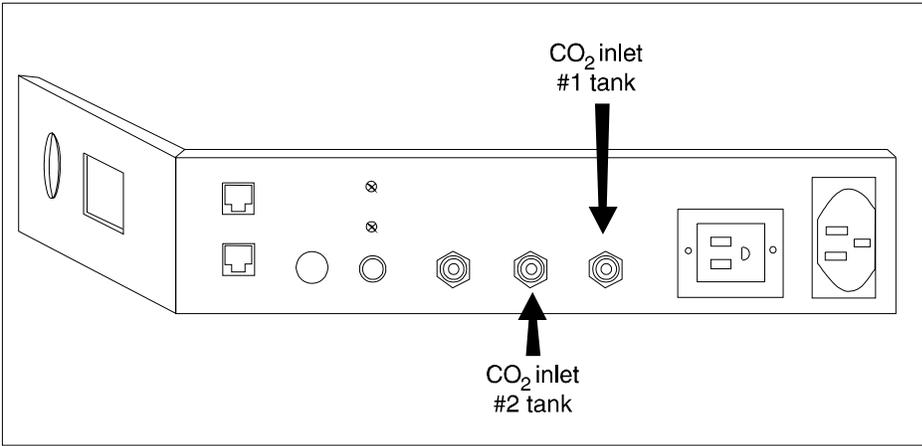


5. Strip the ends of each conductor and wire to the appropriate terminals of connectors J2 and J3 on the analog board.
6. When wiring is completed, slide the electronics drawer back in being very careful to place the door heater cable back into the provided slot so that the drawer does not pinch the cable. See Figure 6-4.
7. Replace the two screws removed earlier and return the unit to service.

6.2 Gas Guard for CO₂ or N₂ (190640/190642)

The 3110 Series incubators can be equipped with a built-in gas guard system that will operate with either a CO₂ or a N₂ gas supply. Only one gas guard can be installed on each unit. The gas guard uses two pressure switches to continuously monitor the pressures of two independent CO₂ or N₂ supplies and automatically switches from one supply to the other when the supply pressure drops below 10 psig (0.690 bar). The gas guard is not designed to be used with multiple incubators.

Both of the CO₂ or the N₂ gas supplies must be equipped with two-stage pressure regulators. The high pressure gauge at the tank should have a 0-2000 psig range and the low pressure gauge should have a 0-30 psig range. The gas supply to the incubator must be maintained at 15 psig (1.034 bar). Gas pressures below 15 psig (1.034 bar) will cause nuisance alarms to occur on incubators equipped with the built-in Gas Guard.



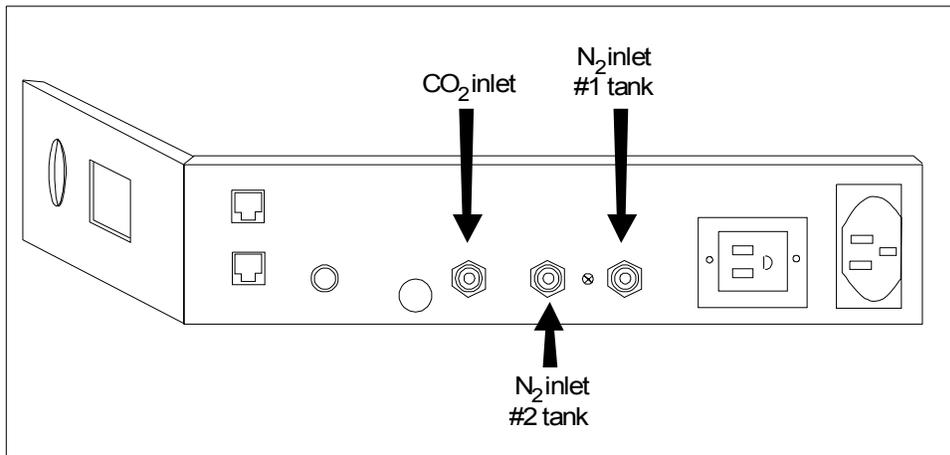
**Figure 6-7
CO₂ fittings**

a. Connecting the CO₂ Gas Supplies (refer to Figures 6-7 and 6-9):

The CO₂ inlets for a built-in gas guard are located on the rear of the electronics drawer. Using 1/4" ID tubing, connect one of the CO₂ supply tanks to the fitting labeled CO₂ Inlet #1 Tank. Connect the second CO₂ supply tank to the fitting labeled CO₂ Inlet #2 Tank. Install 3/8" hose clamps to secure the 1/4" ID tubing to the fittings on the rear of the drawer.

b. Connecting the N₂ Gas Supplies (refer to Figure 6-8):

The N₂ inlets for a built-in gas guard are located on the rear of the electronics drawer. Using 1/4" ID tubing, connect one of the N₂ supply tanks to the fitting labeled N₂ Inlet #1 Tank. Connect the second N₂ supply tank to the fitting labeled N₂ Inlet #2 Tank. Install 3/8" hose clamps to secure the 1/4" ID tubing to the fittings on the rear of the drawer.



**Figure 6-8,
N₂ fittings**

c. Activating the Built-in Gas Guard:

The built-in Gas Guard is not turned on when shipped from the factory. In addition, the Tank Select for the gas guard is specified as Tank 1 when shipped from the factory. Refer to Section 3, Configuration, to activate the Gas Guard or change the Tank Select from 1 to 2. If the operator does not want to use the Gas Guard, the incubator will function normally by supplying CO₂ (or N₂) through the CO₂ Inlet #1 Tank (or the N₂ Inlet #1 Tank).

d. Operation of the CO₂ or N₂ Gas Guard:

With the Gas Guard in operation, the incubator will use the gas supplied through CO₂ (or N₂) Inlet #1 Tank until the pressure drops below approximately 10 psig. At this time, the Gas Guard automatically switches to the gas supplied through CO₂ (or N₂) Inlet #2 Tank.

In addition, the incubator automatically changes the Tank Sel in Configure mode from 1 to 2 to indicate that the incubator is now using gas supplied through CO₂ (or N₂) Inlet # 2 Tank. If the gas supply to CO₂ (or N₂) Inlet #1 Tank is replenished, the incubator will continue to operate using the gas supplied through CO₂ (or N₂) Inlet #2 Tank unless the operator changes the Tank Select from 2 to 1 through the Configure mode. Refer to Section 3, Configuration.

An audible alarm and two visual alarms occur on the control panel when the gas guard switches from one supply to the other. The audible alarm will sound until the operator presses the Silence key on the control panel. A visual alarm in the Message Center will read *Tank 1 Low* while the audible alarm is sounding, but will be removed once the operator presses the Silence key.



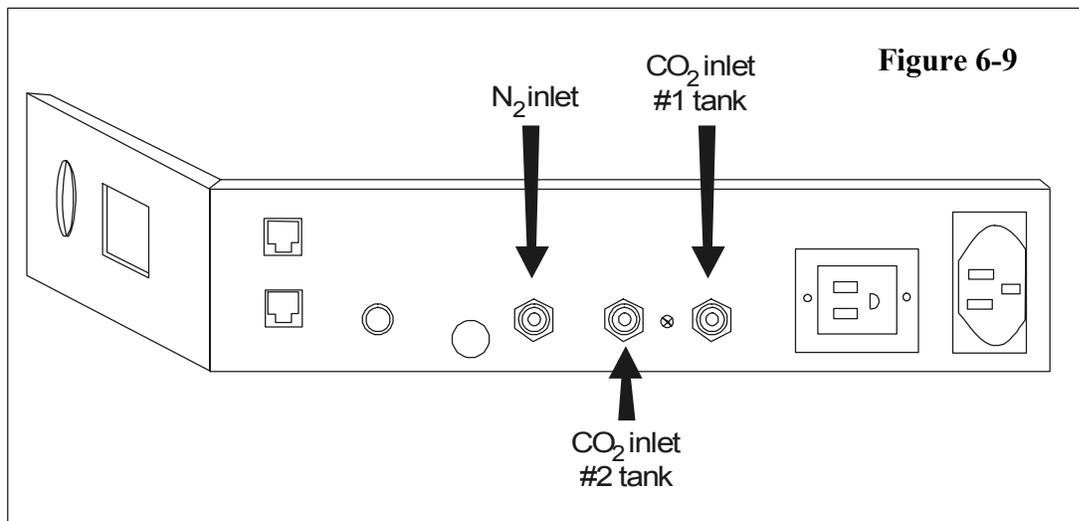
Both the audible and visual alarms described above do not ring back once the Silence key is pressed

However, there is a second visual alarm, *Tank Low*, that illuminates in red below the Silence key on the control panel when *either* of the two gas supplies fall below 10 psig (0.690 bars). The *Tank Low* remains illuminated as long as the gas guard detects a low pressure on either of the gas supply lines. This message is a reminder for the operator to replace or check for insufficient gas supply to the incubator.

If the gas guard does not detect an adequate gas supply at the CO₂ (or N₂) Inlet #1 Tank or CO₂ (or N₂) Inlet #2 Tank, a visual and audible alarm will again occur on the control panel. The visual alarm in the Message Center will read *Tank 1&2 Low*. The audible alarm will continue to ring until the Silence key is pressed. The audible alarm will ring back every 15 minutes after the alarm is silenced, if the Gas Guard continues to detect that both gas supply pressures are below 10 psig (0.690 bars).

6.3 Humidity Readout (190643)

The 3110 Series 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.



a. Factors Affecting the Humidity Level in the Chamber:

- Water level in the humidity pan
- Frequency of door openings
- Humidity pan on the bottom of the chamber versus on a shelf
- Air leakage through the gaskets or stopper in the access port
- N₂ purge on incubators with O₂ control.
- Humidity levels in O₂ units (3130, 3131, 3140, 3141) will be reduced, depending on the amount of N₂ required to control the O₂ level in the chamber.

The following table lists some typical RH levels at different O₂ and CO₂ percentages.



Incubators equipped with a Thermal Conductivity CO₂ sensor rely on a constant level of relative humidity in order to accurately measure and control the CO₂ concentration in the incubator.

Figure 6-10

%O ₂	%CO ₂	%RH (±5%)
2%	5%	60%
5%	10%	75%
10%	10%	80%
21%	5%	95%

b. Accuracy of the Humidity Readout:

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.

6.4 Factory Installed Cooling Coil

Note: For customer convenience, the following items are included in the shipping materials:
12 ft. of 3/8” I.D. vinyl tubing, 4 clamps.

The operating (setpoint) temperature range of the incubator with the cooling coil installed is from +5°C above ambient down to +15°C.

The cooling coil incubator incorporates a finned, U-shaped copper pipe installed within the water jacket. Through this pipe flows chilled water provided by a laboratory bath, typically a Forma Model 2095.

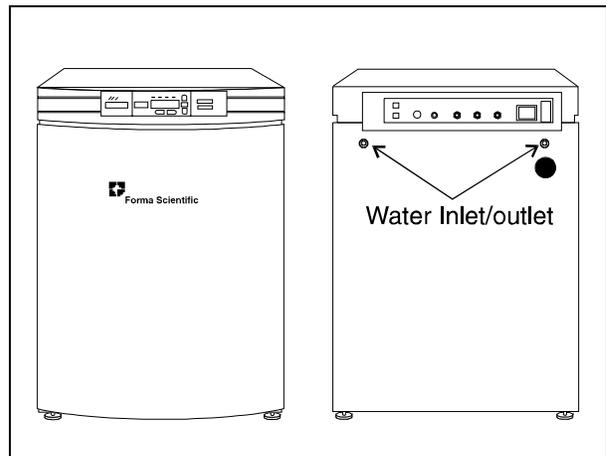


Figure 6-11

When the cooling coil is in use, several factors affect the uniformity inside the incubator chamber: the temperature difference between the operating temperature set point and the bath water temperature; the flow rate of the chilled water, and the on-time percentage of the door heaters.

As determined in carefully controlled laboratory tests, the smaller the difference between the temperature of the bath and the setpoint temperature of the incubator, the better the uniformity. However, decreasing this temperature difference does cause less control of the system because if the bath does not cool the water jacket adequately, the heaters do not cycle and the chamber temperature simply drifts with the ambient temperature of the room.

Tests have shown that as a starting point, operating the bath at 2°C to 3°C below the incubator’s operating setpoint temperature, with a cooling water flow rate of 1/2 to 1 GPM (gallons per minute), should result in good control and uniformity.

Because of the efficiency of the cooling coil design, it is possible for condensation to occur on the outside of the incubator’s water jacket when operating in certain ambient temperature and relative humidity conditions. The condensation will then saturate the fiberglass insulation between the water jacket and the incubator cabinet.

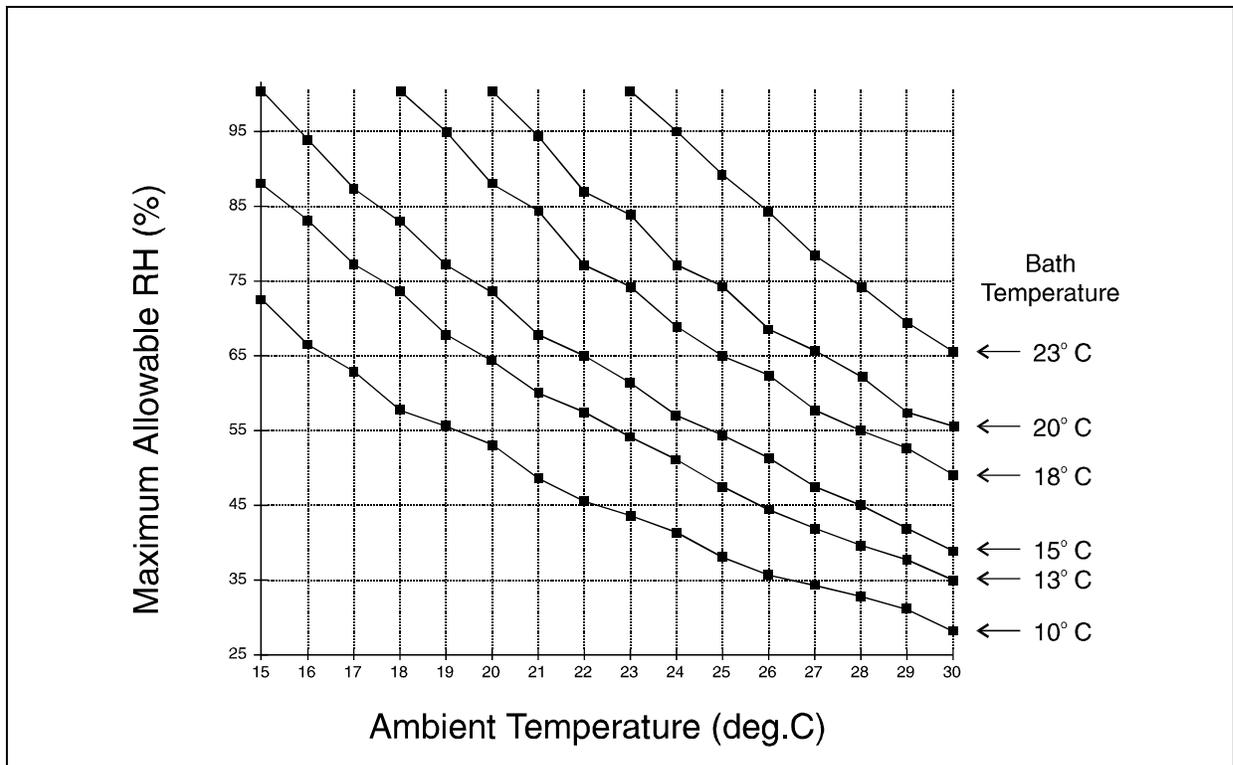


Figure 6-12
RH Limits to Avoid Water Jacket Condensation

Using psychrometric data from the Carrier Psychrometric Chart, curves of maximum allowable RH versus ambient temperature can be plotted for different incubator/bath conditions (Refer to Figure 6-12).

For example, if the bath is to be operated at 15°C and the ambient temperature is about 28°C, the RH in the room must be less than 45% to avoid condensation forming between the chamber water jacket and the outside of the cabinet.

a. Installing the Cooling Coil Incubator

Locate the (2) grey plastic hose barbs shipped inside the incubator. Insert the smooth end of the barb into the fittings on the back of the cabinet. Either can be used as the water inlet or outlet to the bath. Cutting the vinyl tubing in half, the bath can be located up to six feet from the incubator.

Section 7 - Specifications

*Specifications are based on nominal voltages of 115V or 230V in ambients 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
Tracking Alarm	±User programmable high/low

Temperature Safety	
Sensor	Precision thermistor
Controller	Independent analog electronic
Setability	0.1°C

CO₂/O₂	
CO ₂ /O ₂ Control	Better than ±0.1%
CO ₂ Range	0-20%
O ₂ Range	2-20%
Inlet Pressure	15 PSIG (1.034 bars)
CO ₂ Sensor	T/C or IR
O ₂ Sensor	Fuel Cell
Readability	0.1%
Setability	0.1%
Tracking Alarm	User programmable

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

Fittings	
Fill Port	3/8" barbed
Drain Port	1/4" barbed
Access Port	1-1/4" (3.18cm) removable neoprene plug
CO ₂ Inlet	1/4" hose barbed

Unit Heat Load	
115V/230V	344 BTUH (100 Watt)

Shelves	
Dimensions	18.5" x 18.5" (47.0cm x 47.0cm)
Construction	Stainless Steel, perforated, electropolished
Surface area	2.4 sq. ft. (0.22 sq. m)
Max. per Chamber	54.6 sp. Ft. (5.5 sq. m)
Loading	35 lb. (16kg) slide in and out 50 lb. (23kg) stationary
Standard	4
Maximum	23

Construction	
Water Jacket Volume	11.7 gal. (43.5 liters)
Interior Volume	6.5 cu. ft. (184.1 liters)
Interior	Type 304, mirror finish, stainless steel
Exterior	18 gauge, cold rolled steel
Outer Door Gasket	Four-sided, molded magnetic vinyl
Inner Door Gasket	Removable, feather-edged, silicone

Electrical	
115V Models	90-125VAC, 50/60 Hz, 1 PH, 3.6 FLA
230V Models	180-250VAC, 50/60 Hz, 1 PH, 2.0 FLA
Circuit Breaker/Power Switch	6 Amp/2 Pole
Accessory Outlet	75 Watts max. (one per chamber)
Alarm Contacts	Deviation of temp., CO ₂ , O ₂ , and RH, power fail through customer connections through jack on back of unit. 30V, 1A max.
Optional Data Outputs	RS-485,)-1V, 0-5V, 4-20mA
Installation Category	Overvoltage Category II Pollution Degree 2
Maximum Leakage Current	With ground disconnected, 0.65mA Maximum permissible leakage, 3.5mA

Dimensions	
Interior	21.4" W x 26.8" H x 20.0" F-B 54.4cm x 68.1cm x 50.8cm)
Exterior	25.9" W x 39.6" H x 24.9" F-B (63.0cm x 100.3cm x 68.6cm)

Weight (per unit)	
Net	265 lb. (120 kg)
Net Operational	365 lb. (166 kg)
Shipping	324 lb. (147 kg)

Safety Specifications	
Altitude	2000 meters
Temperature	5°C to 40°C
Humidity	80% RH at or below 31°C, decreasing linearly to 50% RH at 40°C
Mains Supply Fluctuations	Operating Voltage Range
Installation Category II ¹	
Pollution Degree 2 ²	
Class of Equipment	

¹ 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.

² 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

a. All Models

Forma Part #	Description
360171	Liquid Level Switch
103065	Feather Gasket
113002	5/16-18 Glide Foot
132046	115/230V Dual Heater
132049	Face Heater 40W, 24VAC
190630	Heated Inner Door
190618, 190619	Inner Door Hinges
190618	Right, Inner Door Hinge
700013	0.500" Flanged Nylon Bearing
990026	Door Gasket w/ Magnet
290138	Water Temp Probe, 2252 Ohm
290137	Air Temp Probe 2252 Ohm
190569	Micro Board
190509	Display Board
230153	6A Circuit Breaker/Switch
460157	Line Filter/Power Inlet
420096	130VA Transformer, Int'l 14/28V S
230135	1 AMP Fuse for 115 V Outlet
230106	1.5 Amp Fuse - Microboard
230159	3.5 Amp Fuse - Microboard
230158	2.5 Amp Fuse - Microboard
250085	Sol Valve 12V 10-32 Ports
770001	Disposable Filter 99.97
156098	Motor 2-Pole 115VAC
103075	Motor Gasket
100097	Blower Wheel 3.5x1.5 CCW
190631	Blower Scroll Assembly
760175	HEPA Filter
760142	Bacterial Air Filter
103072	Blower Plate Gasket
290090	CO ₂ Sensor Assembly
103074	CO ₂ Sensor Plate Gasket

Forma Part #	Description
130038	#6 Solid Neoprene Stopper
180001	Polypropylene Funnel
430108	Line Cord Set
110084	Drain Plug

b. Spare Parts for 230 V units (3111, 3121, 3131, 3141):

Forma Part #	Description
420097	43VA Transformer, INT. SRS
460138	Power Outlet, Snap-In Receptacle
230120	1/2 AMP Fuse for 230 V outlet

c. Spare Parts for IR or O₂ units (3120, 3121, 3130, 3131, 3140 and 3141):

Forma Part #	Description
190885	IR Sensor
190506	IR/O ₂ Circuit Board
184058	Mini Vac/Press.Pmp .75LPM
250119	AC Solenoid
770001	Filter (W/C 451)
511026	3-Hole Stopper (W/C 451)
250118	Valve, Sol, O ₂ , 12VDC, 4W
290083	O ₂ Sensor Fuel Cell
190661	IR Sensor Filter

d. Spare Parts for Gas Guards (190640/190642):

Forma Part #	Description
250121	Solenoid Valve, 3 WAY, 12VDC
360213	Pressure Switch

e. Spare Parts for Humidity Readout (190643):

Forma Part #	Description
290154	RH Sensor Assembly, 1 Ft

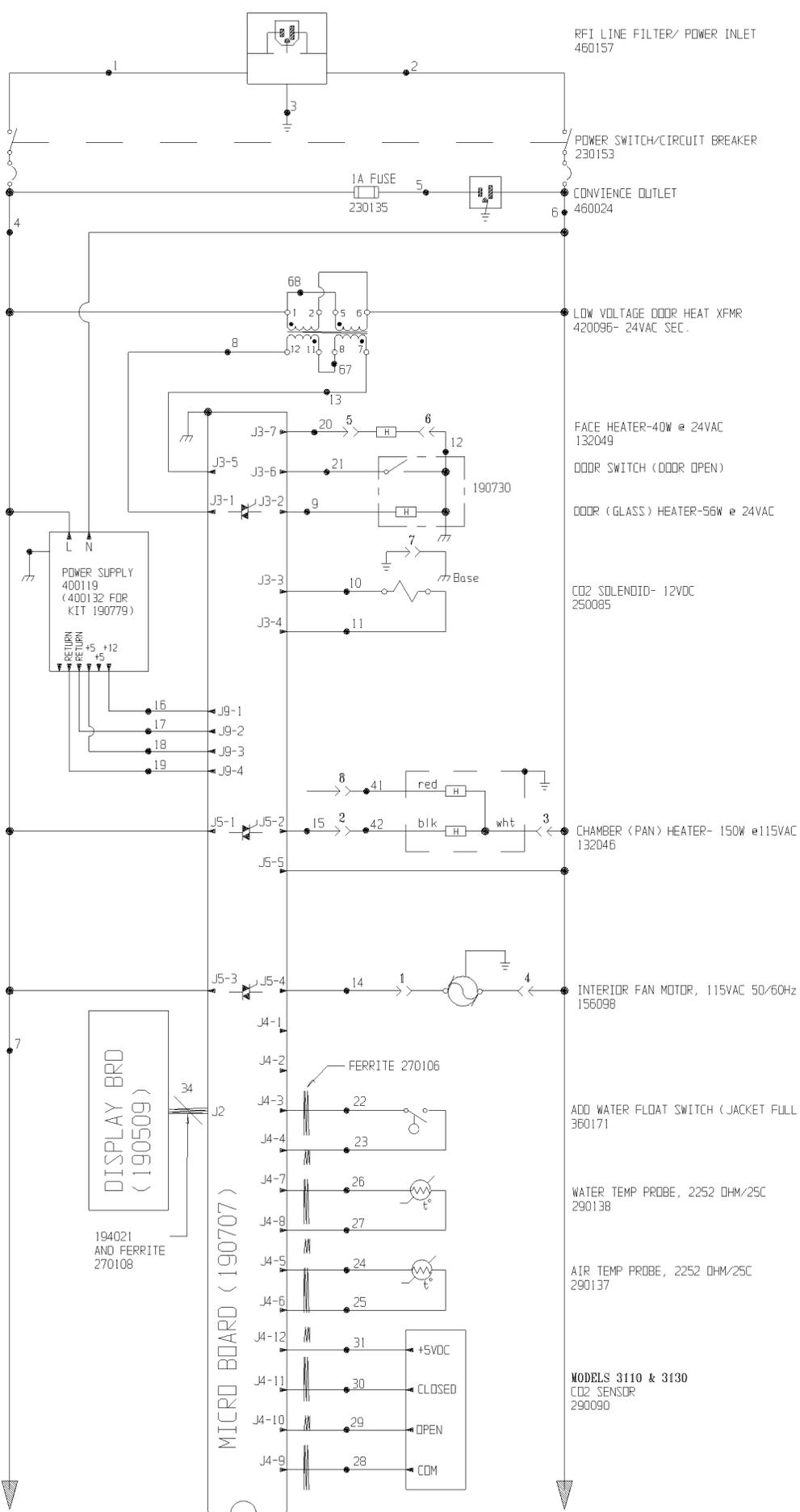
f. Spare Parts for Data Output Options:

Forma Part #	Description
190512	Analog Output Board 4-20 mA
190543	Analog Output Board 0-5V
190544	Analog Output Board 0-1V
73041	24 Gauge 3 Conductor, Analog Output Wire
190523	RS485 Output Kit

Section 9 - Electrical Schematics

POWER CONNECTION
115VAC, 50/60Hz, 1PH, 3.6FLA

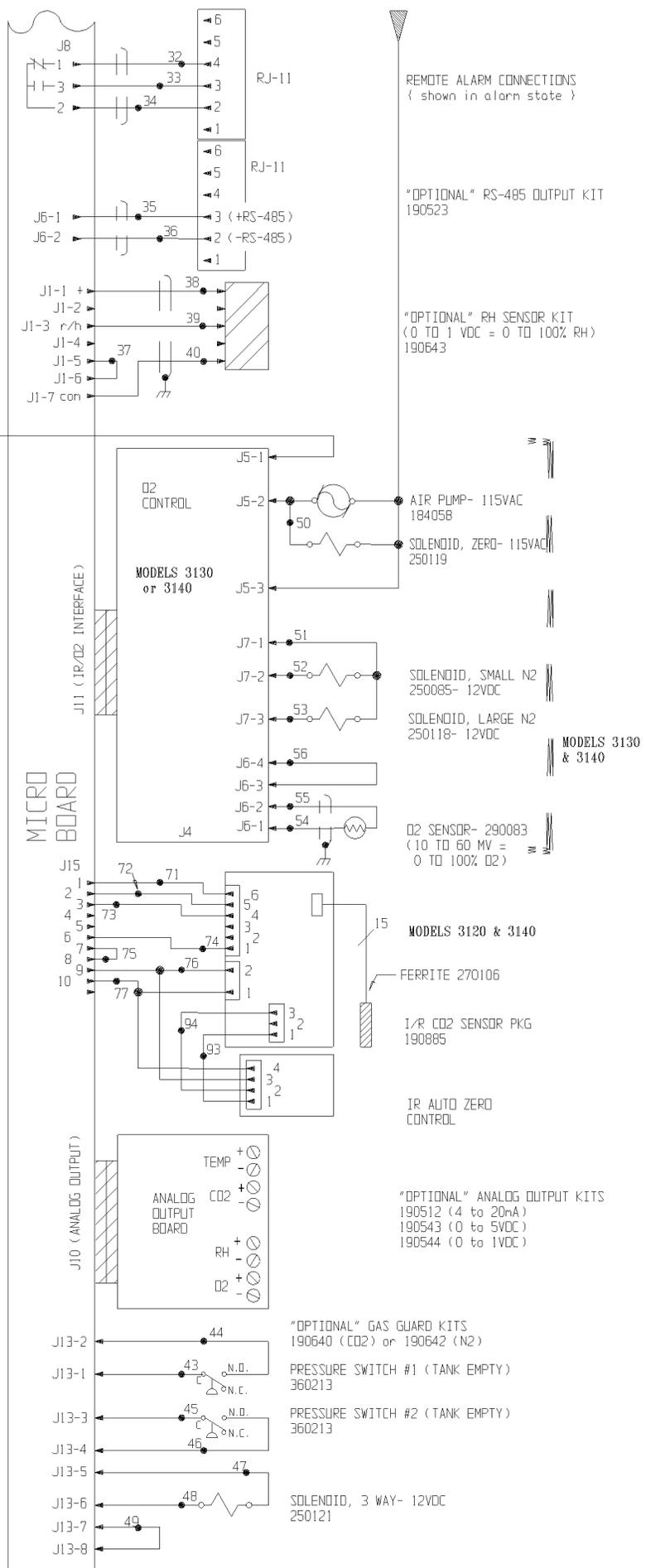
1
2
3
4
5
6
7
8
9
10
11
12
13
14
15
16
17
18
19
20
21
22
23
24
25
26
27
28
29
30
31
32
33
34
35
36
37
38



Electrical Schematic
Forma Model:
3110, 3120, 3130, 3140
Water Jacket Incubator

3110-70-0-D Rev. 6
Page 1 of 3

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



Electrical Schematic
Forma Model:
3110, 3120, 3130, 3140
Water Jacket Incubator

3110-70-0-D Rev. 6
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	NO.	GA.	COLOR
1	16	BROWN	22	22	BLUE	47	20	ORANGE	71	22	BLUE
2	16	BLUE	23	22	BLUE	48	20	ORANGE	72	22	ORANGE
3	16	GRN/YEL	24	22	RED	49	20	BLACK	73	22	YELLOW
3B	16	GREEN	25	22	RED	50	18	BROWN	74	22	BROWN
4	18	BLACK	26	22	YELLOW	51	20	YELLOW	75	22	BLACK
4A	18	BLACK	27	22	YELLOW	52	20	BROWN	76	22	RED
5	18	BLACK	28	22	GREEN	53	20	PURPLE	77	22	GREEN
6	18	WHITE	29	22	GRAY	54	22/2	RED	78	24/15	BLACK
7	18	PURPLE	30	22	PURPLE	55	22/2	BLACK	79	24/15	WHITE
8	18	BROWN	31	22	ORANGE	56	22	BLACK	80	24/15	RED
9	18	BLUE	32	22/3	BLACK	57			81	24/15	GREEN
10	18	YELLOW	33	22/3	RED	58			82	24/15	ORANGE
11	18	YELLOW	34	22/3	WHITE	59			83	24/15	BLUE
12	18	BLACK	35	22/2	RED	60			84	24/15	WHT/BLK
13	18	GREEN	36	22/2	BLACK	61			85	24/15	RED/BLK
14	18	BLUE	37	22	BLACK	62			86	24/15	GRN/BLK
15	18	ORANGE	38	22/3	BLACK	63			87	24/15	ORG/BLK
16	18	RED	39	22/3	RED	64			88	24/15	BLU/BLK
17	18	GREEN	40	22/3	GREEN	65			89	24/15	BLK/WHT
18	18	ORANGE	41	18	RED	66			90	24/15	RED/WHT
19	18	GREEN	42	18	BLACK	67	18	RED	91	24/15	GRN/WHT
20	18	ORANGE	43	20	RED	68	18	ORANGE	92	22	BLACK
21	22	BROWN	44	20	RED	69			93	22	BROWN
			45	20	BLUE	70	22	BLACK	94	22	WHITE
			46	20	BLUE						

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

NOTES:	CUSTOMER APPROVAL/REFERENCE	6 IN-2460 01-18-99 RLM/KDG LDN CHG. 184032 AIR PUMP TO 184058
⊕ Denotes Terminal Strip Connection	APPROVED BY _____	5 SI-7008 5-21-98 GJM/GJM LDN IR SENSOR REVISION
Parts List Reference Number	DATE OF APPROVAL _____	4 IN-2343 11-19-97 GLM/KDG LDN NOTED 400132 USED ON KIT
○ Assembly	THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM FORMA SCIENTIFIC	3 IN-2276 5-27-97 GJM/GJM LDN IMPLEMENT DCS IR CO2 SENSOR
∩ Panel		2 IN-2161 09-30-96 AFC/PDK LDN ADDED SENSOR REFERENCE VALUES
○ Refrigeration		REV ECN NO. DATE BY CAD APPD DESCRIPTION OF REVISION
□ Wiring		DATE 9-24-95 DWN GJM CAD GJM APPD M.H. SCALE
	CUSTOMER MODELS 3110, 3120, 3130, & 3140	JOB TITLE MODULAR WATER JACKET INCUBATOR 115 VOLT (ODM.)
	DWG TITLE ELECTRICAL SCHEMATIC	LOCATION JOB NUMBER DRAWING NUMBER
	INCUBATR	3110-70-0-D

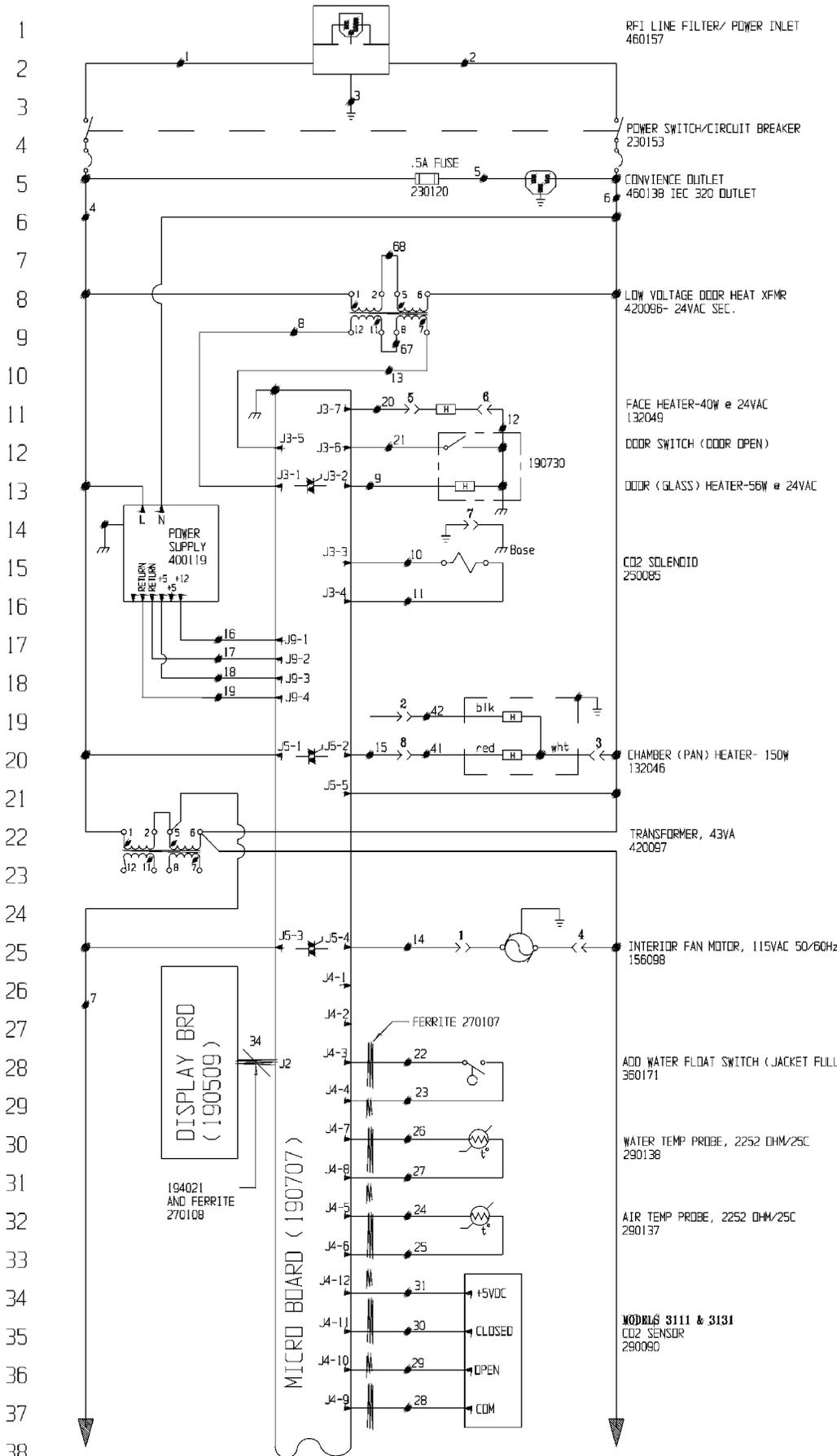
Electrical Schematic
Forma Model:
3110, 3120, 3130, 3140
Water Jacket Incubator
3110-70-0-D Rev. 6
Page 3 of 3



Forma Scientific

BOX 649 MARIETTA, OHIO 45750 TELEX 24-5394
TOLL FREE USA 800-949-3090, OHIO 614-373-4763

POWER CONNECTION
230VAC, 50/60Hz, 1PH, 2.0FLA



RFI LINE FILTER/ POWER INLET
460157

POWER SWITCH/CIRCUIT BREAKER
230153

.5A FUSE
230120

CONVIENCE OUTLET
460138 IEC 320 OUTLET

LOW VOLTAGE DOOR HEAT XFMR
420096- 24VAC SEC.

FACE HEATER-40W @ 24VAC
132049

DOOR SWITCH (DOOR OPEN)
190730

DOOR (GLASS) HEATER-56W @ 24VAC

CO2 SOLENOID
250085

CHAMBER (PAN) HEATER- 150W
132046

TRANSFORMER, 43VA
420097

INTERIOR FAN MOTOR, 115VAC 50/60Hz
156098

ADD WATER FLOAT SWITCH (JACKET FULL)
360171

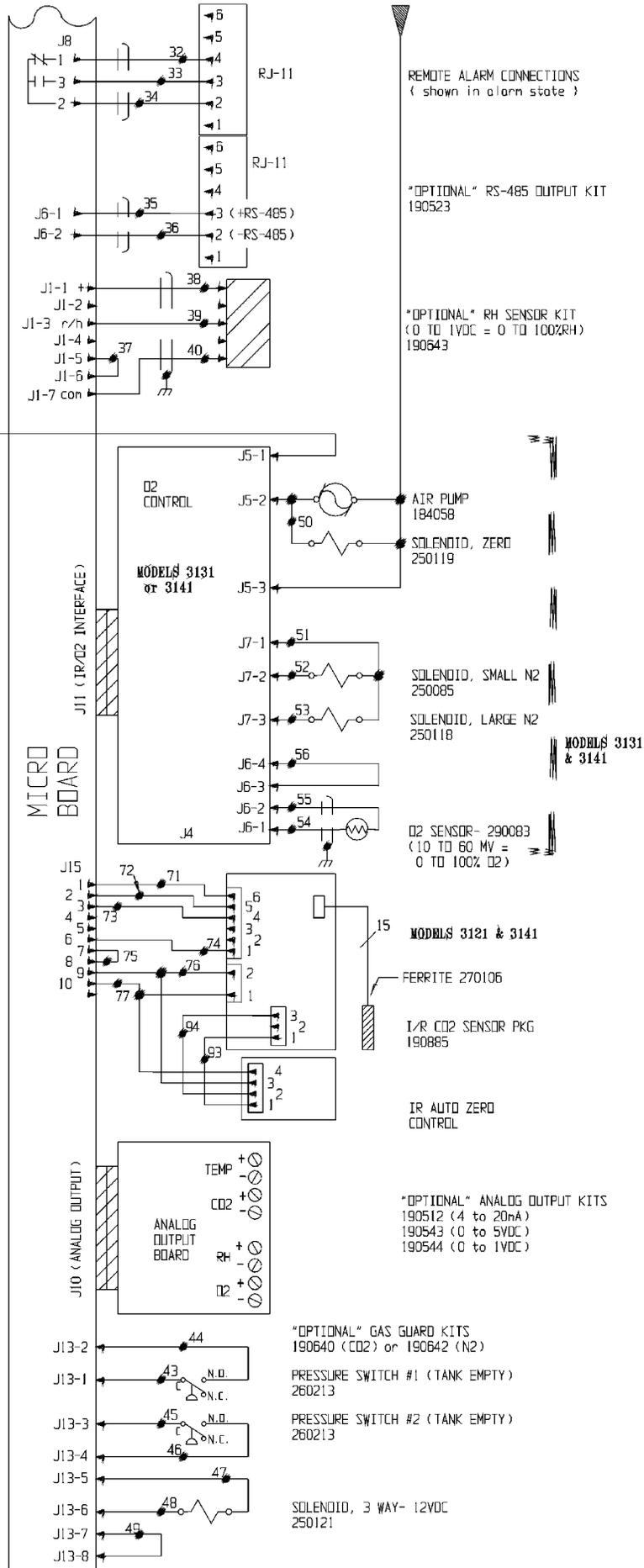
WATER TEMP PROBE, 2252 OHM/25C
290138

AIR TEMP PROBE, 2252 OHM/25C
290137

MODEL#S 3111 & 3131
CO2 SENSOR
290090

Electrical Schematic
Forma Model:
3111, 3121, 3131, 3141
Water Jacket Incubator
3111-70-0-D Rev. 5
Page 1 of 3

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



REMOTE ALARM CONNECTIONS
(shown in alarm state)

OPTIONAL RS-485 OUTPUT KIT
190523

OPTIONAL RH SENSOR KIT
(0 TO 1VDC = 0 TO 100%RH)
190643

MODELS 3131 & 3141

MODELS 3121 & 3141

OPTIONAL ANALOG OUTPUT KITS
190512 (4 to 20mA)
190543 (0 to 5VDC)
190544 (0 to 1VDC)

OPTIONAL GAS GUARD KITS
190640 (CO2) or 190642 (N2)
PRESSURE SWITCH #1 (TANK EMPTY)
260213

PRESSURE SWITCH #2 (TANK EMPTY)
260213

Electrical Schematic
Forma Model:
3111, 3121, 3131, 3141
Water Jacket Incubator

3111-70-0-D Rev. 5
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	NO.	GA.	COLOR
1	16	BROWN	22	22	BLUE	47	20	ORANGE	71	22	BLUE
2	16	BLUE	23	22	BLUE	48	20	ORANGE	72	22	ORANGE
3	16	GRN/YEL	24	22	RED	49	20	BLACK	73	22	YELLOW
3B	16	GREEN	25	22	RED	50	18	BROWN	74	22	BROWN
4	18	BLACK	26	22	YELLOW	51	20	YELLOW	75	22	BLACK
4A	18	BLACK	27	22	YELLOW	52	20	BROWN	76	22	RED
5	18	BLACK	28	22	GREEN	53	20	PURPLE	77	22	GREEN
6	18	WHITE	29	22	GRAY	54	22/2	RED	78	24/15	BLACK
7	18	PURPLE	30	22	PURPLE	55	22/2	BLACK	79	24/15	WHITE
8	18	BROWN	31	22	ORANGE	56	22	BLACK	80	24/15	RED
9	18	BLUE	32	22/3	BLACK	57			81	24/15	GREEN
10	18	YELLOW	33	22/3	RED	58			82	24/15	ORANGE
11	18	YELLOW	34	22/3	WHITE	59			83	24/15	BLUE
12	18	BLACK	35	22/2	RED	60			84	24/15	WHT/BLK
13	18	GREEN	36	22/2	BLACK	61			85	24/15	RED/BLK
14	18	BLUE	37	22	BLACK	62			86	24/15	GRN/BLK
15	18	ORANGE	38	22/3	BLACK	63			87	24/15	ORG/BLK
16	18	RED	39	22/3	RED	64			88	24/15	BLU/BLK
17	18	GREEN	40	22/3	GREEN	65			89	24/15	BLK/WHT
18	18	ORANGE	41	18	RED	66			90	24/15	RED/WHT
19	18	GREEN	42	18	BLACK	67	18	RED	91	24/15	GRN/WHT
20	18	ORANGE	43	20	RED	68	18	ORANGE	92	22	BLACK
21	22	BROWN	44	20	RED	69			93	22	BROWN
			45	20	BLUE	70	22	BLACK	94	22	WHITE
			46	20	BLUE						

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

NOTES:	CUSTOMER APPROVAL/REFERENCE	5 IN-2460 01-18-99 RLM/KDG/LDN	CHG. 184032 AIR PUMP TO 184058
<input checked="" type="checkbox"/> Denotes Terminal Strip Connection	APPROVED BY _____	4 SI-7008 5-21-98 <input checked="" type="checkbox"/> <input checked="" type="checkbox"/> LON	IR CO2 SENSOR REVISION
Last Relay Number	DATE OF APPROVAL _____	3 IN-2276 5-27-97 <input checked="" type="checkbox"/> <input checked="" type="checkbox"/> M.H.	IMPLIMENT DCS IR CO2 SENSOR
Last Terminal Number	THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM FORMA SCIENTIFIC	2 IN-2161 09-30-96 AFC/PDK/LDN	ADDED SENSOR REFERENCE VALUES
Last Wire Number		1 IN-2115 07-19-96 KDG/KDG/LDN	CORRECT I/R CO2 SENSOR
<input type="checkbox"/> Wiring		REV ECN NO. DATE BY CAD/APPD	DESCRIPTION OF REVISION
		DATE 8-24-95 DWN <input checked="" type="checkbox"/> CAD <input checked="" type="checkbox"/> APPD M.H.	SCALE
		CUSTOMER MODLS 3111, 3121, 3131 & 3141	
		JOB TITLE MODULAR WATER JACKET INCUBATOR 230 VOLT (EXPORT)	
		DWG TITLE ELECTRICAL SCHEMATIC	
		LOCATION JOB NUMBER DRAWING NUMBER	
		INCUBATR	3111-70-0-D

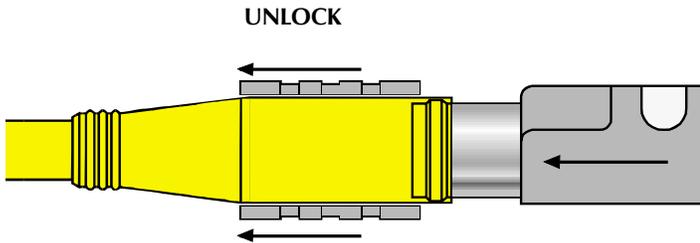
Electrical Schematic
Forma Model:
3111, 3121, 3131, 3141
Water Jacket Incubator

3111-70-0-D Rev. 5
Page 3 of 3


Forma Scientific
801 610 MARLETTA, OHIO 45750 TELEX 24-5394
TELEPHONE 614-885-3300, FAX 614-373-4763

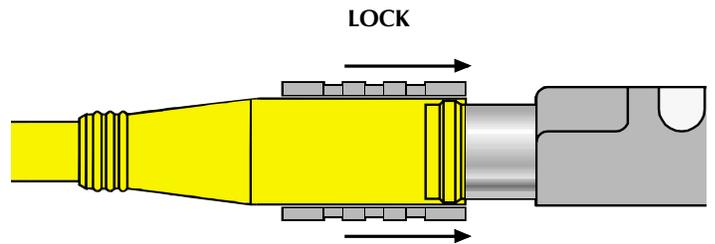
Appendix A

picofast "Snap Lock" Connector Instructions



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

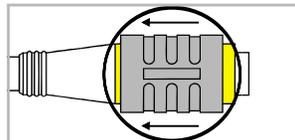
After sensor is connected, push Locking Sleeve forward to create a watertight connection. It's a snap!



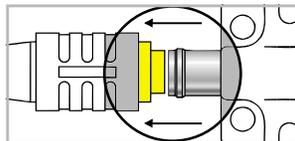
Installation Instructions

To Attach:

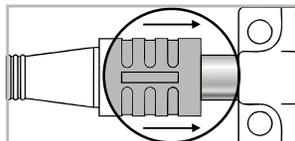
1. Be sure black locking sleeve is pulled back.



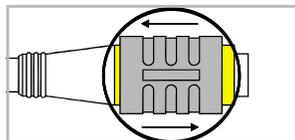
2. Line up pins and push connector onto plug. You will feel a "snap." Do not twist.



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.

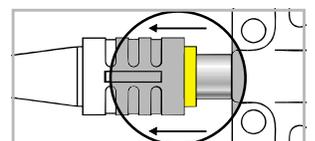


To Detach:

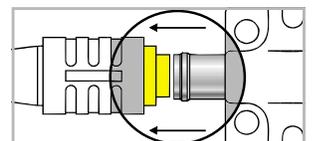
1. Do not twist.



2. Pull locking sleeve back to "unlocked" position.



3. Pull connector straight off.



4. Leave locking sleeve in unlocked position.

