

Forma Scientific, Inc.
P.O. Box 649
Marietta, Ohio 45750

Telephone: (740) 373-4763
Telefax: (740) 373-4189

Models:

**Single Chamber
3158, 3164, 3154, 3166**

**Dual Chamber
3326, 3336, 3354**

Water Jacketed Incubator

Manual No. 7053158

Important!

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.***

NOTE:

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. 7053158			
REV	ECN	DATE	DESCRIPTION
		8/94	Corrected p. 5-10, removed reference to ringback on CO ₂ deviation.
		11/93	Manual revised to current format.

GENERAL SAFETY NOTES

The Occupational Safety and Health Administration (O.S.H.A.) has revised Section 1910-147, **The Control Of Hazardous Energy (Lockout/Tagout)**.

Hazardous energy may be: *electric, air, hydraulic, water, steam, gravity, spring & all other equally hazardous energy.*

This revised regulation, states that you will de-energize all potential sources of energy (may be more than one energy source) prior to performing service or maintenance on any equipment. It also states that a lock shall be placed on the de-energized control, along with a verified test (use of a voltmeter or other equipment) to insure no accidental starts. If you are not familiar with this regulation, review the O.S.H.A. Regulation, Section 1910-147.

In field service, full compliance with this regulation is difficult at best. Troubleshooting must often be performed with hazardous energy applied; therefore extreme caution must be followed during these troubleshooting steps. ***Only Qualified Personnel Must Perform This Work.*** This phase of the repair work must be coordinated through the customer's facilities maintenance department or designated person.

When performing service or maintenance as an outside contractor/worker, follow the **Outside Work Force's Lockout/Tagout** system. Be alert for new types of lockout/tagout devices.

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

For your safety adhere to all **DANGER** and **CAUTION** statements.

Danger: This word is used to call attention to immediate hazards of equipment or conditions which, if not avoided, could result in personal injury, loss of life or property damage.

Caution: This word is used to call attention to potential hazards of equipment or conditions which, if not avoided, could result in personal injury, loss of life or property damage.

DO YOU NEED INFORMATION OR ASSISTANCE ON FORMA SCIENTIFIC PRODUCTS?

If you do, please contact us 8:00 A.M. to 5:00 P.M. (Eastern Standard Time) at:

1-614-373-4763	Direct
1-800-848-3080	Toll Free, U.S. and Canada
1-614-373-4189	FAX
29-8205	Telex

International customers please contact your local sales group.

Forma's **Inside Sales Support Group** can provide information on pricing, give you quotations, take your order and provide delivery information on major equipment items.

The Forma **Product Service Group** can supply technical information about proper setup, operation or troubleshooting of your equipment and fill your needs for spare or replacement parts or provide you with on-site service.

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, you can count on **Life Sciences International North America (LSI/NAS)** for on-the-spot repairs by trained professional field technicians. In addition to checking the reported problem, our technicians also check basic operation, such as:

Control Calibrations	Electrical Circuits
Temperature	Recorders
CO₂ Levels	Blowers/Fans
Compressors	R.H. Levels

The following cities and their surrounding areas are covered by **LSI/NAS** personnel.

Phoenix, AZ	Chicago, IL	Cincinnati, OH
Los Angeles, CA	Boston, MA	Philadelphia, PA
San Diego, CA	Detroit, MI	Washington, PA
San Francisco, CA	St. Louis, MO	Memphis, TN
Denver, CO	Raleigh/Durham, NC	Austin, TX
Washington, DC	Newington, NH	Dallas/Ft. Worth, TX
Atlanta, GA	Newark, NJ	Seattle, WA
Des Moines, IA	New York City, NY	

Within the continental United States, service coverage is by Life Sciences International North America (LSI/NAS). They offer a Bi-Annual Performance Check which provides a complete checkout of your equipment twice a year. This keeps minor problems from becoming major ones. If you would like to know more about their Bi-Annual Performance Check, please contact LSI/NAS on their Toll-free number, 1-800-467-4627.

TABLE OF CONTENTS

SECTION 1 - RECEIVING

- 1.1 Preliminary Inspection
- 1.2 Visible Loss or Damage
- 1.3 Responsibility for Shipping Damage

SECTION 2 - UNPACKING LIST

- 2.1 Unpacking List

SECTION 3 - INTRODUCTION

- 3.1 The Water Jacket: Key Features

SECTION 4 - INSTALLATION AND START-UP

- 4.1 Location
- 4.2 Preliminary Disinfecting
- 4.3 Installing the #6 Neoprene Stopper (Access Port)
- 4.4 Installing the Shelf Brackets and Cam Latch Assembly to the Duct Sheets
 - a. Installing the Shelf Brackets
 - b. Installing the Cam Latch Assembly
- 4.5 Installing the Duct Sheets
- 4.6 Installing the Shelves
- 4.7 Leveling
- 4.8 Connecting to a Power Source

- 4.9 Preparing the Incubator for Filling
- 4.10 Filling the Water Jacket
 - a. Funnel Method
 - b. Tap Fill Method
- 4.11 Connecting the Optional Gas Guard
- 4.12 Connecting the Gas Supply without the Gas Guard
- 4.13 Filling the Humidity Reservoir or Pan
- 4.14 Setting the Chamber Temperature
- 4.15 Setting the Over Temp Safety Thermostat
- 4.16 Zeroing the CO₂ Controller
- 4.17 Setting the CO₂ Content

SECTION 5 - OPERATION

- 5.1 Operation
- 5.2 Overview of Humidification and CO₂
- 5.3 Control Panel
 - a. Fill Fitting and Vent Hole
 - b. Power Switch and Pilot Light
 - c. Circuit Breaker
 - d. Variable/37°C Switch, Temp Control and Heater Pilot Light
 - e. Gas Sample Port
 - f. Setpoint Adjustment Tool
- 5.4 (Factory-Installed Option) Gas Guard Stock #190410
 - a. Control Switch
 - b. Silence Switch
 - c. Gas Depletion Alarm
 - d. Active Tank Indicating Lights
 - e. Pressure Low Indicating Lights

5.5 CO₂ Module

- a. CO₂ Power Switch
- b. CO₂ Controller and Digital Display
- c. Audible CO₂ Alarm and Pilot Light
- d. SET/SILENCE Push Button
- e. Alarm Disable Switch
- f. CO₂ Inject Light
- g. CO₂ Zero Adjustment

5.6 Alarm/Monitor Module

- a. Over Temperature Controller and SET/SILENCE Button
- b. Over Temperature Alarm and Pilot Light
- c. Add Water Pilot Light and Audible Alarm

SECTION 6 - ROUTINE MAINTENANCE

- 6.1 Installing the Decontamination Kit
- 6.2 Disinfecting the Incubator Interior
- 6.3 Cleaning the Cabinet Exterior
- 6.4 Draining the Water Jacket
- 6.5 Changing the CO₂ Filter
- 6.6 CO₂ Test Instruments

SECTION 7 - SERVICE *(For Qualified Personnel Only)*

- 7.1 General Troubleshooting
- 7.2 CO₂ Control Calibration
 - a. Stabilizing the Incubator
 - b. Adjusting the Zero Pot
 - c. Adjusting the Span Pot
 - d. Re-checking the Zero Adjustment
 - e. Checking the CO₂ at the Desired Setpoint
- 7.3 37C° Control Calibration
- 7.4 Replacing the CO₂ Sensor

SECTION 7 - SERVICE (continued)

- 7.5 Replacing the Triac
- 7.6 Replacing the CO₂ Solenoid
- 7.7 Replacing the Temperature Control
- 7.8 Replacing the Pilot Lights
- 7.9 Replacing the Thermistor
- 7.10 Replacing the 0.3 Amp Fuse
- 7.11 Replacing the Blower Motor

SECTION 8 - SPECIFICATIONS

- 8.1 Capacity
- 8.2 Weight
- 8.3 Dimensions
- 8.4 Construction
- 8.5 Shelves
- 8.6 Alarm/Monitor Module
- 8.7 CO₂ Module
- 8.8 Temperature Control
- 8.9 Heaters
- 8.10 Blower
- 8.11 Fittings
- 8.12 Electrical Characteristics
- 8.13 Performance Data

SECTION 9 - AUXILIARY EQUIPMENT

- 9.1 Auxiliary Equipment

SECTION 10 - PARTS LIST

SECTION 11 - SCHEMATICS

SECTION 12 - WARRANTY INFORMATION

SECTION 1 - RECEIVING

1.1 PRELIMINARY INSPECTION

This incubator was thoroughly inspected and carefully packed prior to shipment and all reasonable precautions were taken to ensure safe arrival of the merchandise at its destination. Immediately upon receipt, before the unit is moved from the receiving area, carefully examine the shipment for loss or damage. Unpack the shipment and inspect both interior and exterior for any in-transit damage.

1.2 VISIBLE LOSS OR DAMAGE

If any loss or damage is discovered, note any discrepancies on the delivery receipt. Failure to adequately describe such evidence of loss or damage may result in the carrier refusing to honor a damage claim. Immediately call the delivering carrier and request that their representative perform an inspection. Do not discard any of the packing material and under no circumstances move the shipment from the receiving area.

1.3 RESPONSIBILITY FOR SHIPPING DAMAGE

For products shipped F.O.B. Marietta, Ohio, the responsibility of Forma Scientific, Inc. ends when the merchandise is loaded onto the carrier's vehicle. On F.O.B. Destination shipments, Forma Scientific's and the carrier's responsibility ends when your Receiving Department personnel sign a free and clear delivery receipt.

Whenever possible, Forma Scientific, Inc. will assist in settling claims for loss or in-transit damage.

SECTION 2 - UNPACKING LIST

2.1 UNPACKING LIST

Remove the packing box from the incubator. If the unit is to be moved by fork lift, leave the incubator on the skid until it has been moved to its designated location. A plastic bag containing the following accessories is packed inside the incubator:

Qty.	Stock #	Description	Purpose
1	190028	Decontamination Kit	Maintenance
1	180001	Polypropylene Funnel	Fill & Drain
3'	246011	Vinyl Tubing 3/8" ID	Fill & Drain
6'	246010	Vinyl Tubing 3/16" ID	CO ₂ Connection
1	285057	Extender Card	Maintenance & Calibration

Also packed with each incubator:

Qty.	Stock #	Description
24	23052	Stainless Steel Wing Nuts (#10-32)
20	55000	Stainless Steel Screws (Truss Head #10-32" x 5/8")
1	121034	Cam Latch (Nylon)
1	127019	Stainless Steel Spacer (1/4" Diameter x 9/16")
1	550017	Stainless Steel Screw (Truss Head #8-32" x 1")
1	23051	Stainless Steel Wing Nut (#8-32)
5	224140	Stainless Steel Shelves
1	600034	Snapper Hose Clamp (.375")
4	3113234	Shelf Brackets
1	3113224	Stainless Duct Channel, Top
1	3113222	Stainless Duct Sheet, Left
1	3113223	Stainless Duct Sheet, Right
1	130038	#6 Neoprene Stopper
1	190388	Telephone Line Cord, 12'
1	190392	Modular RJ-11 Jack
1	7053158	Instruction Manual

SECTION 3 - INTRODUCTION

3.1 THE WATER JACKET: KEY FEATURES

The Forma Scientific Model 3158 water jacketed incubator is designed to meet the stringent requirements of today's modern laboratory and is a combination of economy, accuracy and reliability.

The Model 3158 Incubator features:

- Ultra-flat, vibration-free shelves to provide optimum culturing conditions.
- Direct-set CO₂, temperature, and over temperature alarm set points..
- Digital readout of temperature and CO₂ levels in the chamber.
- A sealed chamber to minimize CO₂ consumption.
- A sealed water jacket to minimize water evaporation.

The 3158 water jacket incubator is designed for easy and convenient servicing by providing:

- Front access to all components so that the unit need not be moved for service, repair, or routine maintenance.
- A slide-out control panel for easy access to electrical components.
- Plug-in CO₂, and Temperature/Alarm modules to facilitate servicing or replacement.
- Self-diagnostic switches in the control modules to aid in troubleshooting the system and localizing problems.
- Minimal unit down-time to disinfect the water jacket.

- No cracks or crevices in, or around, the chamber walls to harbor hidden or hard-to-reach bacterial growth.
- Stainless steel shelves, shelf channels and duct sheets which are easily removed without tools for cleaning and disinfecting the entire interior.
- A blower wheel which is disposable and easily replaced.
- An improved water jacket drain and fill system.

Careful consideration has been given to the importance of *product protection* through the addition of:

- An audible Add Water alarm and pilot light to alert the operator when the water level in the water jacket has become too low for efficient operation.
- Audible and visual CO₂ alarms which activate when the CO₂ content in the chamber deviates 1% above or below the CO₂ control setpoint.
- An Over temperature Alarm which activates when the chamber temperature rises above the control setpoint.
- An Add Water and Over temperature silence with 30 minute ring back.
- Remote alarm contacts for Over Temperature, Add Water, Power Failure and CO₂ deviation.

SECTION 4 - INSTALLATION AND START-UP

4.1 LOCATION

Locate the incubator on a firm, level surface capable of supporting the unit.

Approximate weight with water: Single Chamber = 357 lbs (162 kg)
Dual Chamber = 714 lbs (324 kg)

The incubator should be placed in an area of the laboratory away from any centrifuges, sonicators, doors, windows and air-conditioning or heating ductwork that might produce drafts. To help prevent microbial contamination, the incubator should not be located in areas of high personnel traffic.

Adequate room must be provided behind the incubator for connections of electrical, gas and other equipment. Also, since gases are emitted from the rear of the incubator, the unit should not be installed in an unvented recess or in a poorly ventilated room.

Caution: To prevent injury to personnel and/or damage to equipment, lock inner glass door and secure outer door before tipping unit to adjust the leveling feet. Do not attempt to tilt the incubator without assistance while adjusting the leveling feet.

4.2 PRELIMINARY DISINFECTING

Before installing the duct sheets and the shelves, remove the clear plastic film from the shelf brackets and duct sheets. Forma Scientific, Inc. recommends disinfecting all interior surfaces (including both door gaskets). Rinse the surfaces with sterile distilled water (50K Ohm to 1 Meg Ohm). Also disinfect the CO₂ sensor and the blower wheel, taking care not to saturate the sensor.

The duct sheets and shelves must be washed with the same disinfectant solution and rinsed with sterile distilled water prior to their installation in the chamber. Repeat rinsing until you are satisfied that all of the disinfectant-detergent has been removed. Proceed with the installation as noted. Refer to Section 6.2 of this manual.

4.3 INSTALLING THE #6 NEOPRENE STOPPER (ACCESS PORT)

Open the incubator outer door and inner glass door. Locate the opening in the top left corner of the interior chamber. Place the beveled end of the stopper in the opening.

4.4 INSTALLING THE SHELF BRACKETS AND CAM LATCH ASSEMBLY TO THE DUCT SHEETS (Refer to Figure 4-1)

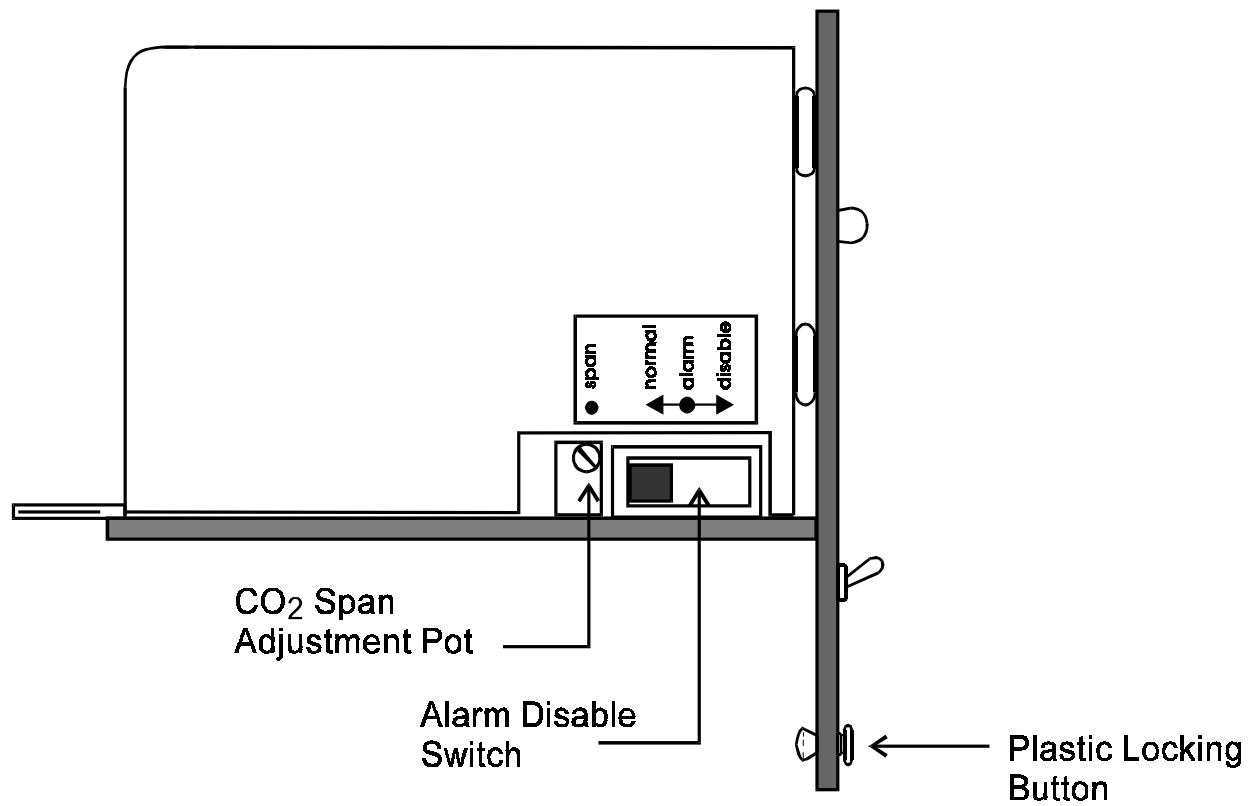
A plastic coating has been applied on the duct sheets and shelf brackets to protect the finish during shipping and handling. This plastic coating must be peeled off before the protected parts are installed.

a. Shelf Brackets

1. Locate the plastic bag containing the #10-32 x 5/8" truss head screws and wing nuts (20).
2. With the duct sheet in a vertical position, align and mount the shelf brackets to the unflanged side of the duct sheets. Wing nuts go on the flanged side.

b. Cam Latch Assembly

1. Locate the bag containing (2) #8-32 x 1" stainless steel truss head screws, (2) #8-32 wing nuts, (2) stainless steel spacers and (2) cam latches (nylon).
2. Place stainless steel spacer (sleeve) over the #8-32 x 1" screw and insert it into the opening of the cam latch.
3. Locate the five (5) small adjustment holes at the top of each duct sheet.
4. Place the cam latch on the inside (or unflanged side) of duct sheet and insert screw with spacer sleeve through the middle hole of the five hole adjustment. If the duct sheets appear to be too tight or too loose, the cam latch assembly may be re-positioned up or down one hole for proper alignment.
5. Secure screw on flanged side of duct sheet with #8-32 wing nut.
6. Repeat these procedures for the other duct sheet.



Installing the Shelf Brackets and Duct Sheets

Figure 4-1

4.5 INSTALLING THE DUCT SHEETS (Refer to Figure 4-1)

Note: The left duct sheet has large notches in both top and bottom edges. The right duct sheet has a single notch in the top edge only.

1. Carefully put the right duct sheet into the incubator chamber with the flanges toward the wall.
2. Put the left duct sheet into the chamber, with the square notch at the top and the flanges toward the wall. Allow the top of the left-side duct sheet to lie diagonally across the chamber, resting upon the right-side duct sheet.
3. Hook the top channel into the top opening of the right duct sheet. The top channel must be positioned so that the brass CO₂ sample tube (mounted in the ceiling of the incubator) is aligned with the notched out area on the right rear side of the top channel. The round opening will align with the blower wheel when it is slid into place.
4. While supporting the blower channel, slide the left duct sheet up until it is vertical, making sure that the blower channel lines up into the slot on both duct sheets.
5. Turn cam latch to a vertical position (up against bottom side of top channel) to secure. If duct sheets appear to be too tight or too loose, upon installation, the cam latch assembly may be re-positioned up or down one hole for proper alignment.

4.6 INSTALLING THE SHELVES

The shelves may be placed at any level in the chamber. Slide the shelf into the shelf bracket at the desired level.

4.7 LEVELING

Check the level of the unit by placing a bubble-type gauge on one of the shelves. Turn the hex nut located on the leveling leg clockwise to lengthen the leveling leg, or raise the unit. Turning the hex nut counterclockwise will shorten the leg, or lower the unit. Be sure to level the incubator before filling the water jacket.

Caution: To prevent injury to personnel and/or damage to equipment, lock inner glass door and secure outer door before tipping unit to adjust leveling feet. Do not attempt to tilt the incubator without assistance while adjusting the leveling feet.

4.8 CONNECTING THE INCUBATOR TO A POWER SOURCE

With the incubator power switch OFF, connect the unit to an adequate power source. Refer to the serial tag located on the back of the unit or see Section 8.10 of this manual for specific power requirements.

Note: Forma Scientific, Inc. recommends that the incubator be connected to a separate electrical circuit.

4.9 PREPARING THE INCUBATOR FOR FILLING

450 ML of rust inhibitor was placed in the water jacket before the incubator was shipped. The rust inhibitor mixes with the distilled water during filling and provides a protective coating on the interior of the water jacket.

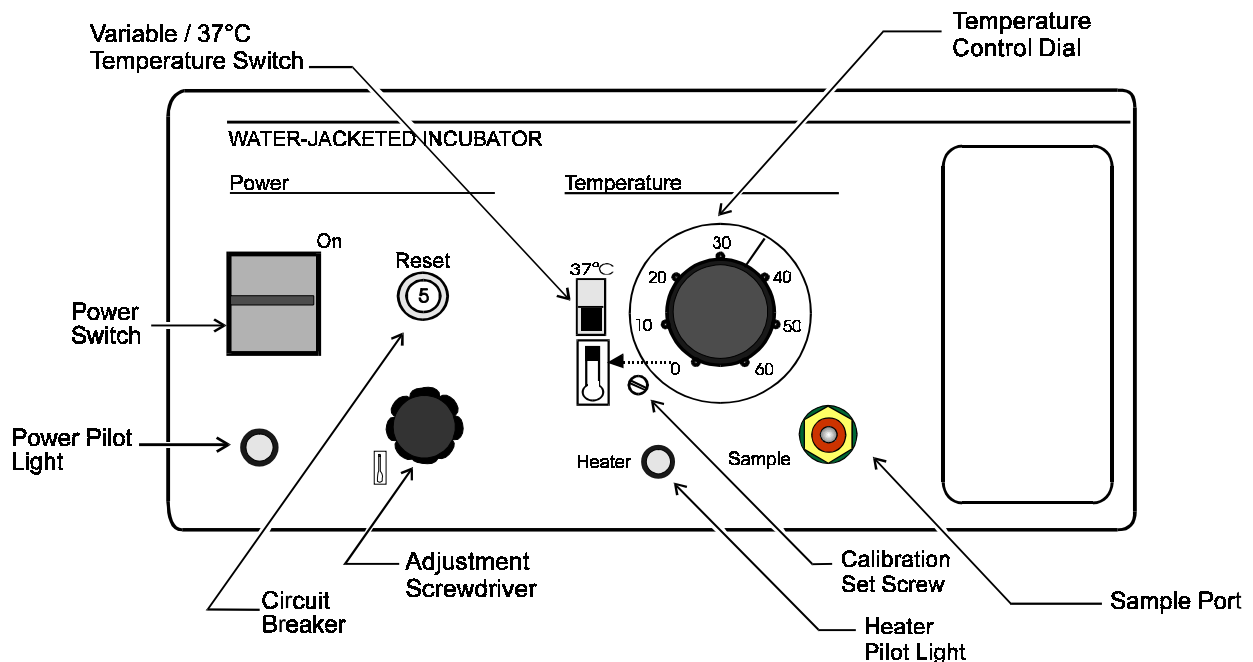
The fill fitting is used to fill the water jacket without having to move the incubator. The fill fitting is located on the center front of the unit, directly above the door. Refer to Figure 4-3.

The vent hole, located adjacent to the fill fitting allows the air displaced by water entering the jacket to escape. It also prevents distortion of the chamber by allowing air to escape as the unit expands and contracts during heating and cooling.

Caution: Do not plug this vent. A plugged vent will damage the water jacket chamber.

1. Remove the plastic protective cap from the fill fitting. Check to see that the vent hole located adjacent to the fill fitting is not covered or plugged.
2. Remove the protective metal cover from the drain fitting and valve and make sure the valve is closed (turned to the vertical position). Refer to Figure 4-4 for drain and drain valve location.
3. Set the TEMP SELECT switch to the variable position (down) and turn the temp control knob completely counterclockwise to keep the heater from coming on before the water jacket is filled. Refer to Figure 4-2 and Figure 4-3 for component locations.

4. Turn the incubator power switch ON. The ADD WATER warning light and audible alarm will come on. Pressing the SET/SILENCE button on the alarm panel will silence the audible alarm
5. Remove the adjusting screwdriver from the control panel. The black knob is the handle of the adjusting screwdriver. Refer to Figure 4-2.
6. Push the SET/SILENCE button on the Monitor Alarm Module, and adjust the over temp alarm setpoint to a setting 2° or 3° above the intended operating temperature. (Refer to Section 4.14)



Electric Power and Temperature Control Module
Figure 4-2

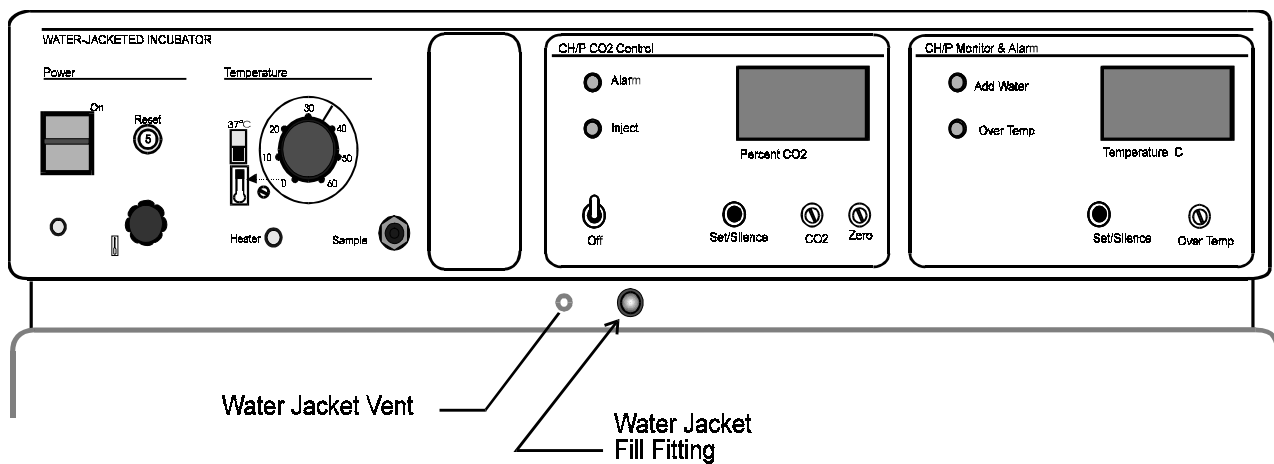
7. Press the CO₂ SET/SILENCE button and rotate the CO₂ set screw until the display reads 00.0. (Refer to Figure 4-7 for locations of the SET/SILENCE button and set screw.)

4.10 FILLING THE WATER JACKET

There are two methods of filling the water-jacket.

- a. Funnel Fill Method
- b. Tap Fill Method

Caution: *Purity of the distilled water used in the water jacket and humidifier must be within the 50K ohm to 1 Meg ohm range to protect and prolong the life of the stainless steel water jacket. The use of tap water or distilled water outside of the specified range will decrease the operating life of the unit and will void the warranty.*



Water Jacket Fill Fitting and Vent Hole
Figure 4-3

a. Funnel Fill Method (Refer to Figure 4-3)

A funnel and vinyl tubing are contained in the accessories bag included with each incubator.

1. Verify that the water jacket drain valve on the bottom of the incubator frame is closed. The valve is located behind a protective metal cover and will require two screws to be removed. The valve is closed when the valve knob is turned vertical. Refer to the illustration in Figure 4-4.

2. Attach the funnel into one end of the 3/8" I.D. vinyl tubing.
3. Remove the plastic protective cap from the fill fitting.
4. Attach the free end of the funnel tubing to the fill fitting.
5. Hold funnel above the level of the fill fitting and pour the water into the funnel until the ADD WATER alarm light goes out. The water jacket holds about 13 gallons (49 liters) of water.
6. Add one additional quart (one liter) of distilled water. The incubator is now properly filled.
7. Remove the tubing from the fill fitting and replace the plastic protective cap.

Note: Water seepage may occur from vent hole when chamber temperature increases.

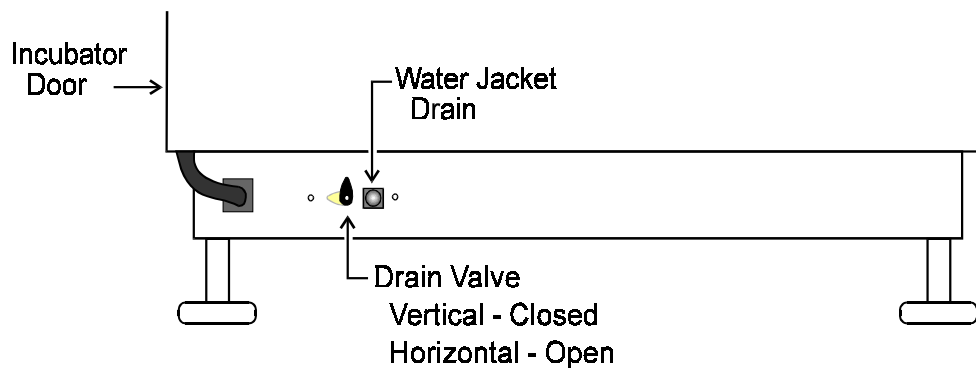


Figure 4-4
Incubator Base with Water Jacket Drain and Valve

b. Tap Fill Method (Refer to Figure 4-3)

Three feet of 3/8" ID vinyl tubing is provided in the accessories bag included with each incubator. If the distilled water outlet is more than three feet from the incubator, more tubing and a connector will be required.

1. Verify that the water jacket drain valve on the bottom of the incubator frame is closed. The valve is located behind a protective metal cover and will require two screws to be removed. The valve is closed when the valve knob is turned vertical. Refer to the illustration in Figure 4-4.

2. Remove the plastic protective cap from the fill fitting.
3. Connect the 3/8" I.D. vinyl tubing between the fitting and the distilled water tap.
4. Open the tap until water flows steadily into the water jacket.

Caution: A high rate of flow and high pressure can cause distortion to incubator chamber walls.

5. Turn the water off immediately when the ADD WATER alarm light goes out.
6. Using the funnel, add an additional quart (one liter) of distilled water.
7. Remove vinyl tubing and reinstall the plastic protective cap.

4.11 CONNECTING THE OPTIONAL GAS GUARD (Refer to Figure 4-5)

a. Connecting the Outlet of the Gas Guard to the Incubator's Gas Inlet

Before connecting the gas supply tanks, connect a piece of supply tubing between the outlet fitting and the gas inlet(s) on the back of the incubator. If both chambers are to be monitored by the gas guard both inlets need to be connected. The gas outlet is located between Tank 1 and Tank 2 inlet fittings. Check the connections for leaks.

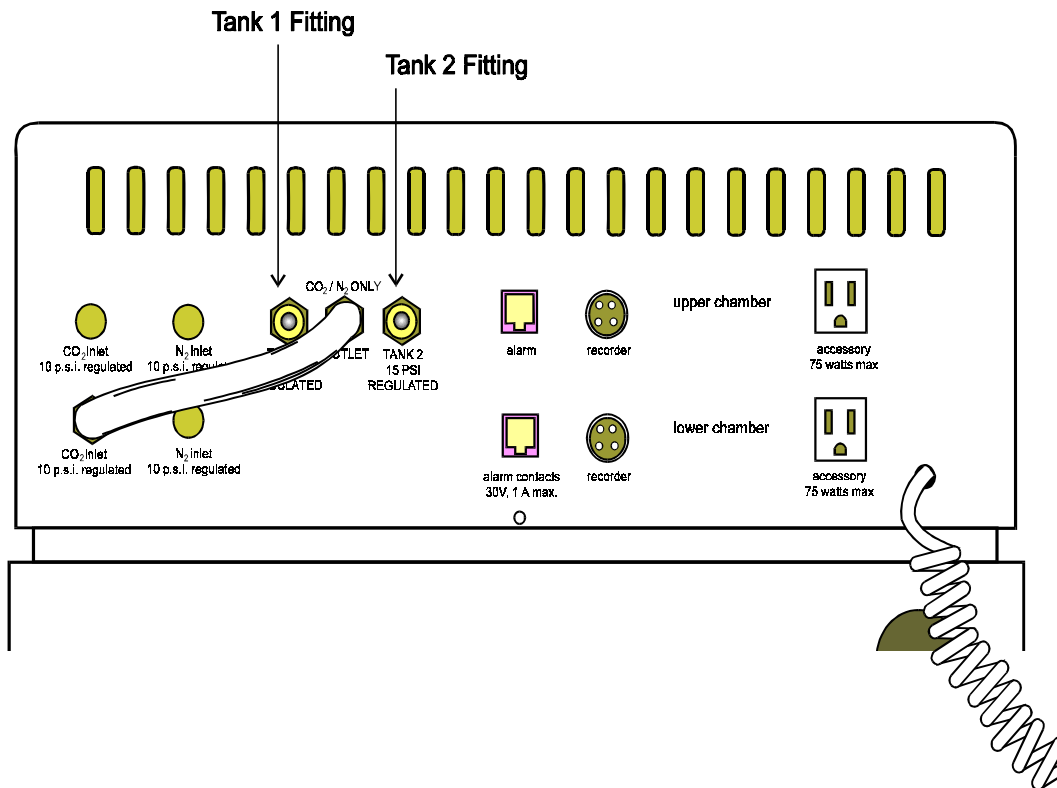
b. Connecting The CO₂ Supply Tanks

Note: Output pressure from the tanks must be regulated with a dual stage pressure gauge set at 15 PSIG. Liquid substances must be supplied in tanks *without siphon tubes* to ensure that only gas enters the incubator.

The high pressure gauge should have an indicating range of 0 to 2000 PSIG to monitor tank pressure. The low pressure gauge should have an indicating range of 0 to 30 PSIG to monitor actual input pressure to the incubator injection system. A suitable two-stage pressure regulator is available from Forma Scientific, Inc., Stock #965010.

Supply pressure lower than 13.5 PSIG will cause improper operation of the gas guard and may cause the occurrence of nuisance alarms. Pressure higher than 15 PSI will damage the CO₂ control system.

Securely attach the gas lines from the supply tank to the barbed fittings located on the back of the unit. The fittings are labeled. Check the connections for leaks.



Rear view of the control panel, stacked incubator model with Gas Guard
Figure 4-5

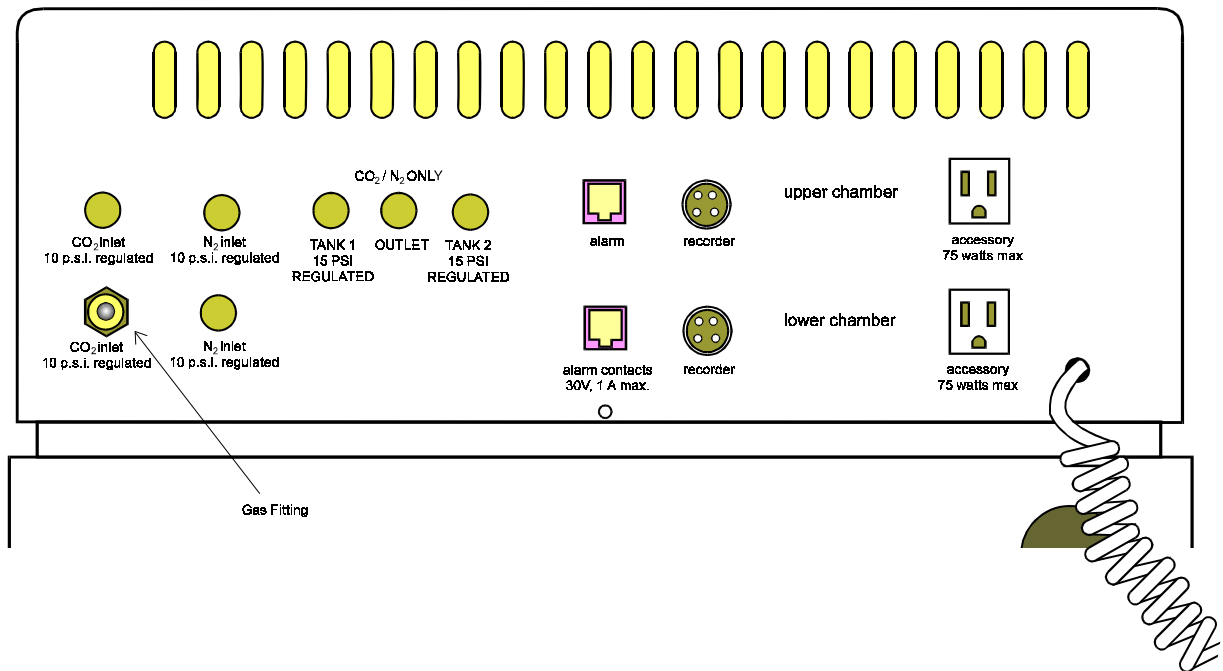
4.12 CONNECTING THE GAS SUPPLY WITHOUT GAS GUARD (Refer to Figure 4-6)

If Gas Guard is not used, the gas supply is connected directly to the CO₂ inlet.

Note: Output pressure from the tanks must be regulated with a dual stage pressure gauge set at 10 PSIG. Liquid substances must be supplied in tanks *without siphon tubes* to ensure that only gas enters the incubator.

The high pressure gauge should have an indicating range of 0 to 2000 PSIG to monitor tank pressure. The low pressure gauge should have an indicating range of 0 to 30 PSIG to monitor actual input pressure to the incubator injection system. A suitable two-stage pressure regulator is available from Forma Scientific, Inc., Stock #965010.

Securely attach the gas lines from the supply tank to the desired barbed inlet fitting(s) located on the back of the unit. The fittings are labeled. Refer to Figure 4-6. Check the connections for leaks.



Rear view of the control panel, stacked model without Gas Guard
Figure 4-6

4.13 FILLING THE HUMIDITY RESERVOIR OR PAN

Do not use plastic pans for humidification as they will have an unpredictable effect on humidity and O₂/CO₂ levels in the incubator. Use only the floor of the unit or the optional stainless steel humidity pan.

Do not use de-mineralized or de-ionized water in the humidity reservoir or pan unless it has been boiled immediately prior to use as it may be contaminated with bacteria.

Listed below are two recommended methods of providing elevated humidity in the chamber.

Note: Frequent door openings will cause humidity loss from chamber and may result in desiccation of the product. Recovery time will also be affected.

a. Reservoir

The reservoir in the bottom of the incubator may be filled with at least 3/4" of sterile distilled (50K ohm to 1 Meg ohm) water.

b. Humidity Pan

The optional humidity pan (Forma Stock #237001) will hold 1.6 gallons (6 liters) of sterile distilled (50K ohm to 1 Meg ohm) water. This pan may be autoclaved. The best humidity and temperature response from the humidity pan is obtained when the pan is placed directly on the incubator floor.

The water level in the humidity reservoir should be checked frequently. If a disinfectant has been added to the water in the reservoir, it should be changed once a week to help prevent microbial contamination. If a disinfectant has not been added to the sterile distilled water, the water should be changed at least twice a week.

It is important that the water level in the reservoir or pan be kept relatively constant as fluctuations or "dry-outs" will have an adverse effect on the humidity level and CO₂ control in the chamber.

Caution: When installing the humidity pan, exercise care to avoid tearing the inner door gasket.

4.14 SETTING THE CHAMBER TEMPERATURE (Refer to Figure 4-7)

Before the initial temperature setting is made, push the SET/SILENCE button on the Alarm Monitor Module. Using the adjustment screwdriver, turn the OVER TEMP set screw until the display shows a temperature that is two to three degrees above the desired operating setpoint. The over temp safety may be reset after the chamber temperature has stabilized at setpoint.

If a chamber temperature of 37°C is desired, set the Variable/37°C switch on the Temperature Control panel to the 37°C (up) position. (Figure 4-2.)

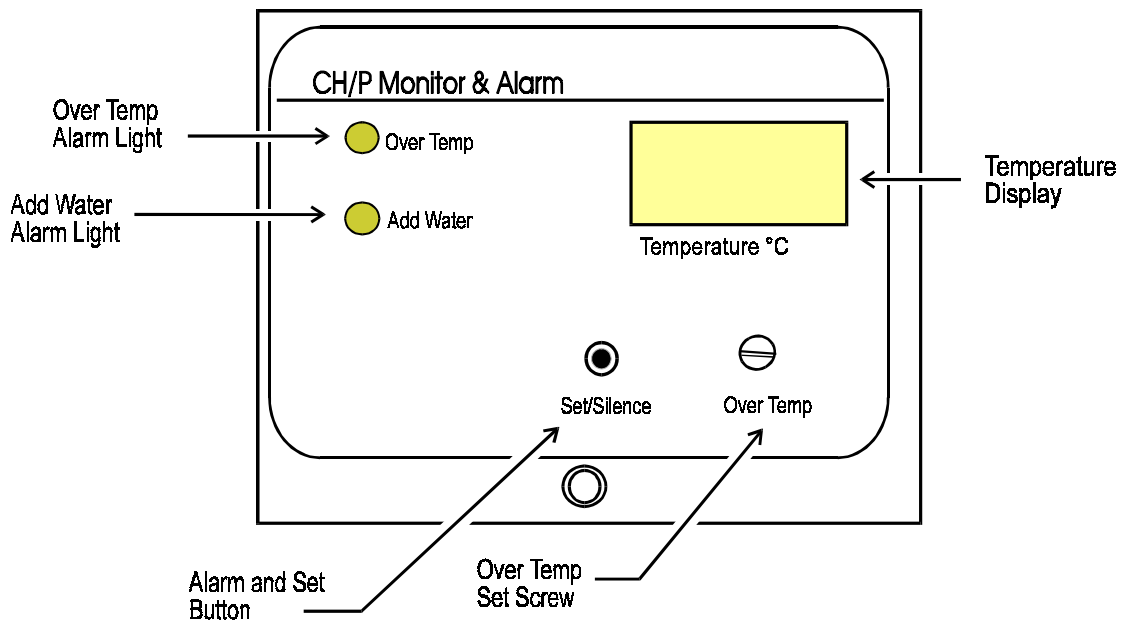
If a value other than 37°C is desired, set the switch to the Variable position (down), and set the temperature control knob to the desired setpoint. Any temperature between 5°C above ambient to 50°C may be selected.

4.15 SETTING THE OVER TEMP SAFETY THERMOSTAT (See Figure 4-7)

Once the chamber temperature has stabilized (as indicated by the digital display), the Over Temp Safety should be set as follows:

1. Push the SET/SILENCE button on the Alarm Monitor Module.
2. Using the screwdriver mounted on the control panel, turn the OVER TEMP set screw until the desired over temp alarm point is shown on the digital display. The over temp setpoint can be set within 0.1° of the operating setpoint, but it is recommended that it not be set within 0.5 of setpoint.

Note: The Over Temp Safety should be checked quarterly to insure proper operation. To check the over temp control, push the SET/SILENCE button on the Alarm Monitor Module and turn the Set screw counterclockwise until the over temp safety light is lit. Reset the Over Temp Safety after the test.



Temperature and Alarm Monitor Module, Figure 4-7

4.16 ZEROING THE CO₂ CONTROLLER (See Figure 4-8)

Important!

This adjustment is made using the CO₂ gas content of ambient air (0.03%), the most accurate standard available.

Tools Required:

Adjustment screwdriver (provided on the panel)

FYRITE CO₂ Analyzer (use only for checking) or other CO₂ measuring device.

a. Stabilizing the Incubator

1. Stabilize the incubator at the operating temperature and humidity level with no CO₂ in the interior chamber.
2. Turn off the CO₂ at the supply.
3. Fill the humidity reservoir or pan.
4. Allow the incubator temperature and humidity to stabilize. This will take a minimum of eight hours. On initial start-up, however, allow the unit three days to stabilize.

b. Adjusting the Zero Set Pot (Refer to Figure 4-8)

1. Using the small screwdriver mounted on the control panel, adjust the CO₂ control zero pot to read 00.0 on the digital display. Wait 5 minutes. Repeat if necessary until the display is stable.
2. Turn on the CO₂ at the supply.
3. Turn the CO₂ setpoint to the desired %.

c. Checking the CO₂ at the Desired Setpoint

1. Allow the incubator to reach setpoint and control (inject light will cycle) for a minimum of 30 minutes.
2. Check the CO₂ level with a FYRITE until two consecutive readings agree. If the FYRITE and display are not within plus or minus 1.0%, consult the factory.

Note: After proper zeroing, the CO₂ display will be more accurate than the FYRITE because the zero adjustment was accomplished using absolute.

4.17 SETTING THE CO₂ CONTENT

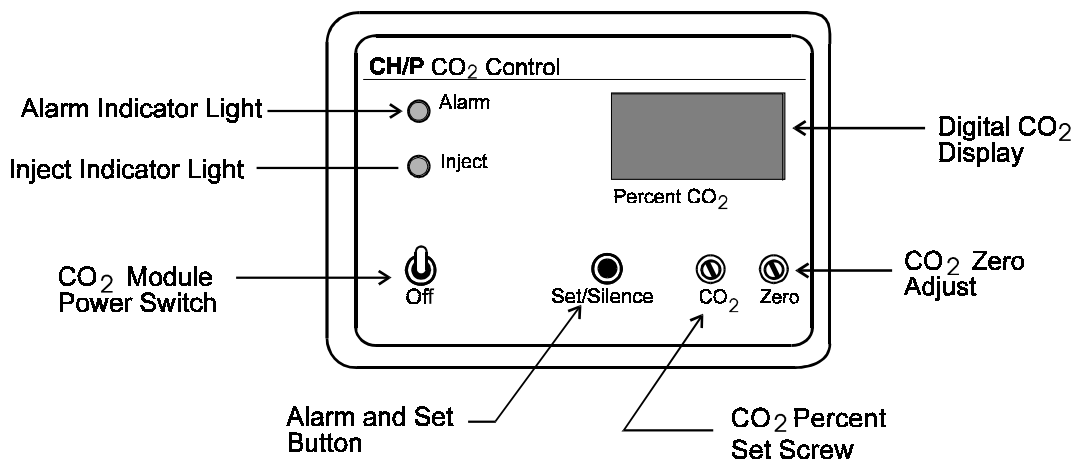
The following conditions must be satisfied before the CO₂ percentage can be set:

Allow the temperature and humidity in the chamber to stabilize. For initial settings of CO₂, it is recommended that temperature and humidity be allowed to stabilize for three days.

Check the CO₂ control zero (See Section 4.15 for detailed instructions).

To set the CO₂ percentage, press the CO₂ SET/SILENCE button and rotate the CO₂ set screw until the desired percentage is indicated on the digital display.

Note: If the unit is in Over Temp and a CO₂ injection occurs, a brief, high CO₂ percentage will appear on the digital display. This is due to the shut down of the internal fan during Over Temp. The high CO₂ percentage occurs only at the sensor. CO₂ throughout the chamber will remain normal.



CO₂ Control Module, Figure 4-8

SECTION 5 - OPERATION

5.1 OPERATION

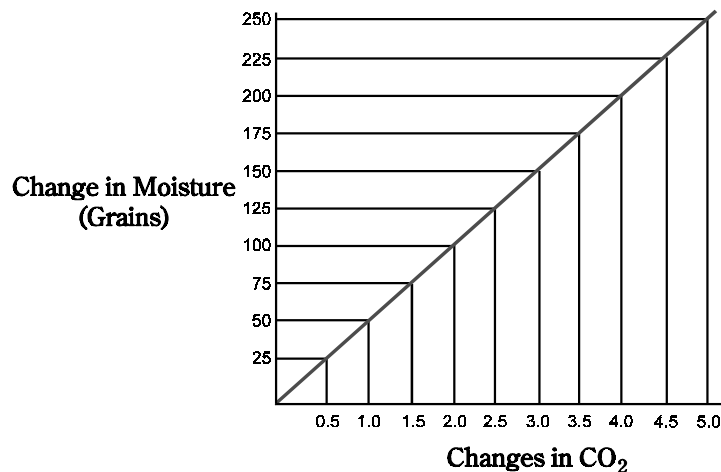
The water jacket is filled with approximately 13 gallons (49 liters) of water through the fill fitting located on the front of the unit. The water is then warmed by the chamber heater, providing very stable heating of the incubator chamber. Not only does the water stay at a constant temperature with a minimum of heater on-time, but it also acts as insulation from ambient temperature conditions.

Temperature control is maintained by a proportional, zero-switching device which provides temperature uniformity throughout the chamber. A separate and independent over temperature controller assures product safety by assuming control at the over temp setpoint if the primary controller malfunctions. Should an over temp condition develop, the monitor alarm system will alert the operator that a malfunction has occurred.

An internal blower gently circulates the air in the chamber to prevent CO₂ stratification and minimize culture desiccation.

5.2 OVERVIEW OF HUMIDIFICATION AND CO₂

Thermal conductivity of the air inside of the incubator is affected not only by the quantity of CO₂ present, but by the quantity of water vapor present in the incubator atmosphere. This effect is linearly related to the absolute humidity of the atmosphere (See figure below).



Moisture Effect on CO₂ Control Calibration
Figure 5-1

When monitoring the effects of CO₂, absolute humidity must be held constant so any change in thermal conductivity is caused only by a change in the CO₂ concentration. Under the worst circumstances, a change in absolute humidity could cause such a significant change in thermal conductivity that the controller could shift the CO₂ content by as much as 4%.

Maintaining the water level inside the incubator is a simple procedure, but an important one to keep the humidity in the incubator constant. Any water pan, used in lieu of flooding the incubator floor, must be stainless steel and at least 187 square inches of surface area. Tests indicate that smaller pans, bowls or non-metallic pans do not provide adequate humidification, which can lead to incubator humidity variations with ambient humidity shifts, resulting in CO₂ changes in the incubator.

When operating a dry incubator, as opposed to a humidified one, ambient humidity fluctuations will affect CO₂ calibrations. Since the fluctuations possible in extreme ambient changes have less effect on the total absolute humidity, the CO₂ calibration can be affected by as much as 1.5%.

When a change in humidity or temperature is needed, the CO₂ control can be easily zeroed for the new condition.

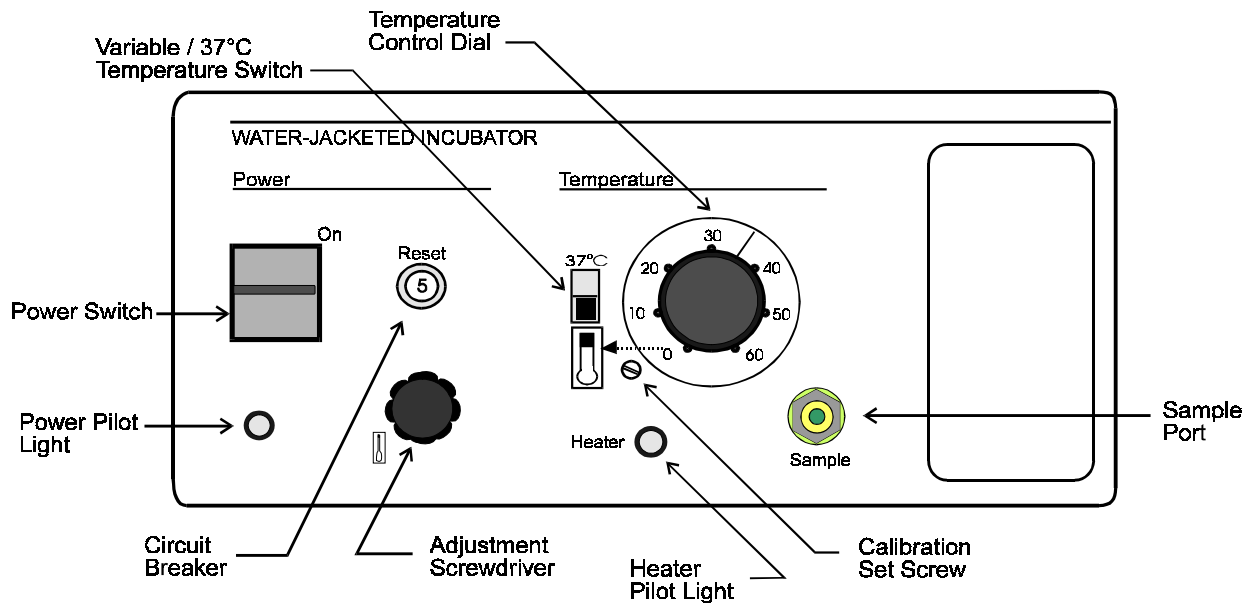
Temperature changes have little affect on CO₂ calibration, but do cause large changes in the absolute humidity which is reflected in changes in CO₂ calibration.

5.3 CONTROL PANEL (Figure 5-2)

a. Fill Fitting, Drain Fitting/Valve and Vent Hole

The fill and drain fittings are conveniently located on the front of the incubator. The water jacket fill fitting is located directly above the incubator door. The drain fitting is located below the door, behind a protective metal cover. A small valve is located adjacent to the drain fitting for convenience when draining the water jacket.

The vent hole, located adjacent to the fill fitting, allows the air displaced by water entering the jacket to escape. It also prevents distortion of the chamber by allowing air to escape as the unit expands and contracts during heating and cooling. Under no circumstances should the vent be plugged.



Incubator Control Panel

Figure 5-2

b. Power Switch and Pilot Light

The main power switch controls the ON/OFF power to the unit. The power pilot light is activated when the power switch is on, and the unit is receiving power.

c. Circuit Breaker (Reset)

The 5 amp circuit breaker for the incubator (labeled Reset) can be pushed to reset the incubator power supply within a few seconds after the breaker has tripped. If it trips a second time, the unit should be checked by a qualified electrician.

d. Variable/37°C Switch, Temp Control and Heater Pilot Light

When the Variable/37°C switch is set to the 37°C (up) position, chamber temperature will automatically be maintained at +37°C. If necessary, the 37°C control can be calibrated using the calibration screw located at the lower left side of the temperature control dial. See Section 7.3 to calibrate the 37°C setting.

When the Variable/37°C switch is set to the Variable (down) position, control is assumed by the temperature control potentiometer. The numbers (0 to 60) around the control knob indicate approximate setpoint values in degrees Centigrade. Any value between +5°C above ambient temperature and 50°C may be selected. The heater pilot light will be activated whenever the heater is energized.

e. Gas Sample Port

A sample port for checking CO₂ percentage by independent means (e.g. FYRITE or similar CO₂ measuring device).

Important!

The Sample Port should never be capped. It is the vent for the incubator chamber. Refer to Figure 5-2.

f. Setpoint Adjustment Tool

A small screwdriver, located below the circuit breaker, has been provided for setting the CO₂ and over temp setpoints. Pull out on the knob to release the screwdriver (the knob is the handle of the screwdriver).

**5.4 (FACTORY-INSTALLED OPTION) GAS GUARD STOCK #190410
(Refer to Figure 5-3)**

Forma Scientific's stock # 190410 built-in Gas Guard is a monitoring and switching device for incubator gas supplies.

The unit monitors a two cylinder gas supply. When the active supply tank is depleted, an audible alarm sounds and the gas guard automatically switches to the second tank.

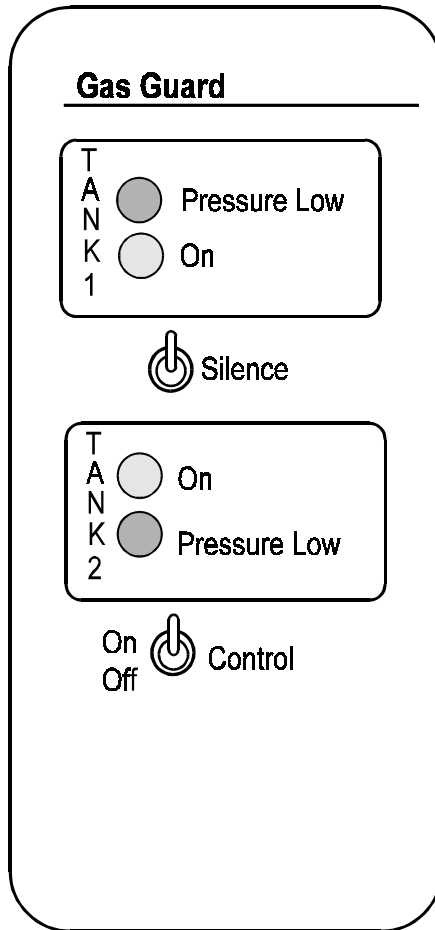


Figure 5-3
Built-in Gas Guard
(Factory Installed Option)

a. Control Switch

The CONTROL switch controls power to the gas guard control circuit board. The yellow LEDs on the front panel designate which tank is active and indicate the unit is receiving power.

b. Silence Switch

The SILENCE switch silences the alarm and selects the alternate tank. If the SILENCE switch fails to silence the alarm, it is an indication that both gas tanks are depleted or incapable of producing 15 PSIG.

Note: During initial start-up, the incubator should not be injecting gas.

c. Gas Depletion Alarm

The gas depletion alarm sounds automatically when the active supply tank becomes empty. The other tank becomes active with the sounding of the alarm.

Note: Should both supply tanks become depleted, the only way to silence the alarm is to turn the CONTROL switch to Off.

d. Active Tank Indicating Lights

The two yellow indicating lights labeled "TANK1 On" and "TANK2 On," located on the front of the control panel, indicate which of the gas tanks is active.

e. Pressure Low Indicating Lights

The two red indicating lights, labeled "Pressure Low," indicate when a supply tank is depleted. If both tanks are depleted both red indicating lights will be lit.

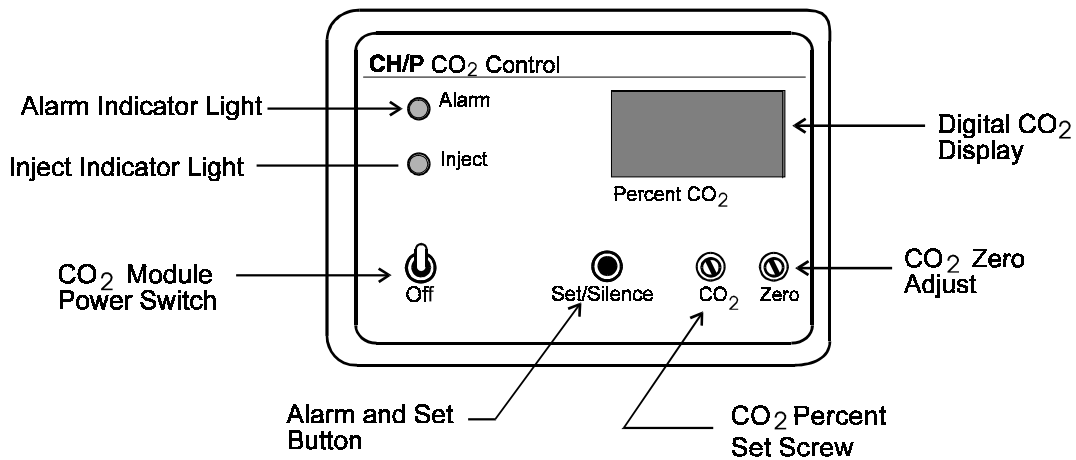
f. Start-Up Procedure

1. Turn the CONTROL switch on.
2. Select the active tank using the SILENCE switch.

5.5 CO₂ MODULE (Refer to Figure 5-4)

a. CO₂ Power Switch

The CO₂ power switch controls the electrical power to the CO₂ system, and it must be ON when the incubator is to be operated with CO₂. The switch should be turned on as soon as power is applied to the unit to allow the CO₂ system to warm up.



CO₂ Control Module, Figure 5-4

b. CO₂ Controller and Digital Display

The LCD digital readout on the CO₂ module continually displays the percent of CO₂ in the chamber. The setpoint is displayed when the CO₂ SET/SILENCE button is pushed.

The CO₂ setpoint is changed by pushing the CO₂ SET/SILENCE button and rotating the CO₂ set screw to the desired percentage.

c. Audible CO₂ Alarm and Pilot Light

The audible CO₂ alarm and pilot light are activated when the percent CO₂ deviates from setpoint by plus or minus 1% (nominal) for longer than approximately four minutes.

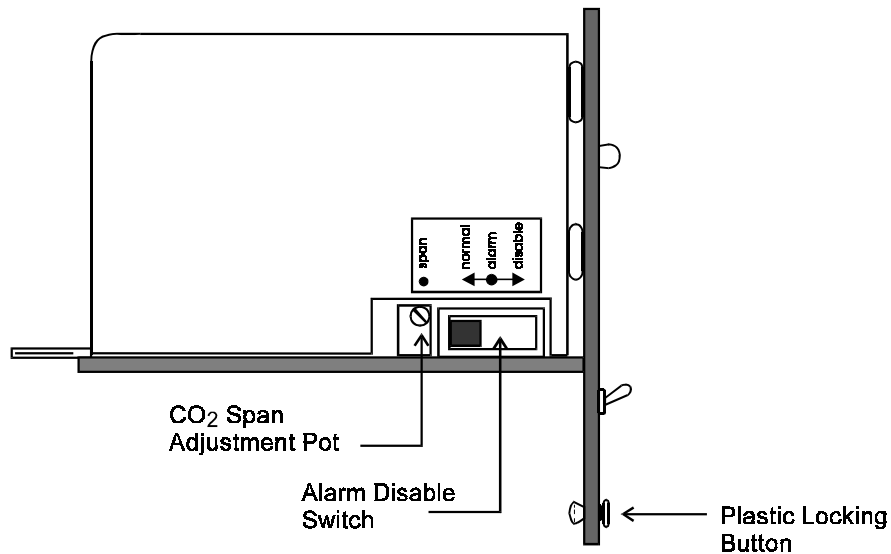
d. SET/SILENCE Push Button

When pushed, the SET/SILENCE button will silence the CO₂ alarm and extinguish the alarm light. This button must be pushed to *set* or *display* the CO₂ setpoint.

e. Alarm Disable Switch (Figure 5-5)

Note: It is necessary to pull the CO₂ module out slightly to gain access to the Alarm Disable Switch. Grasp the black plastic button on the bottom of the control module, directly beneath the Set/Silence push button. Pull on the button until the circuit board releases from its internal connector strip. The Alarm Disable Switch is located on the side of the module and is clearly marked. (See Figure 5-5.) To replace the module, push firmly until the connector strip is seated. Press the black plastic button until it snaps into place.

When the alarm switch is in the DISABLE position, the CO₂ alarm is completely disabled. When the switch is set to the NORMAL position, the alarm system is operative and can be silenced by pressing the SET/SILENCE button.



**Alarm Disable Switch on the CO₂ Control Module
Figure 5-5**

f. CO₂ Inject Light

The CO₂ inject light is activated whenever there is a demand for CO₂ to meet setpoint requirements. Since the CO₂ inject light is independent of the CO₂ alarm, it will continue to signal a need for CO₂ when the CO₂ alarm is set to either the defeat or silence position.

g. CO₂ Zero Adjustment

The CO₂ zero adjustment is used for zeroing the CO₂ controller to specific control conditions. It is the *only* user calibration adjustment on the CO₂ module. ***All internal adjustments are for qualified service personnel only.***

5.6 ALARM/MONITOR MODULE (Refer to Figure 5-6)

a. Over temperature Controller and SET/SILENCE Button

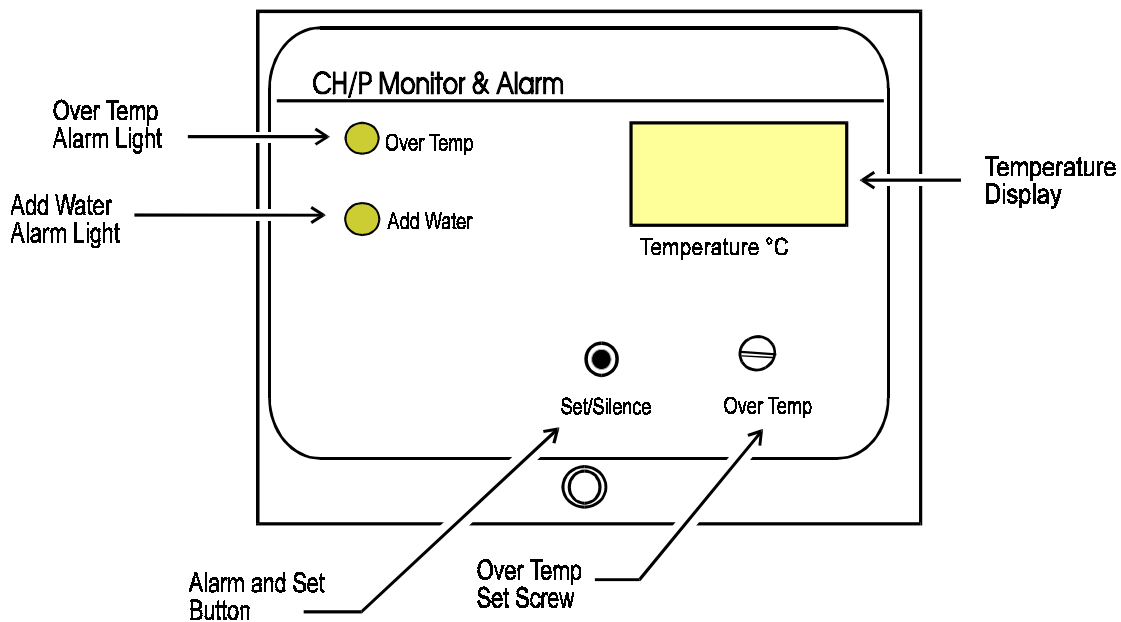
The over temperature setpoint is displayed when the SET/SILENCE button on the module is pushed. Over temp control point is adjusted by pushing the SET/SILENCE button and rotating the set screw on the module to the desired setpoint.

b. Over temp Alarm and Pilot Light

The over temperature audible alarm and pilot light will come on in the event of an over temp condition. The audible alarm can be silenced for approximately 30 minutes by pressing the SET/SILENCE button. The audible alarm will ringback every 30 minutes for as long as the over temp condition exists.

c. Add Water Pilot Light and Audible Alarm

The ADD WATER audible alarm and pilot light are activated whenever the water level in the water jacket is low. The audible alarm can be silenced for approximately 30 minutes by pressing the SET/SILENCE button. The audible alarm will ringback every 30 minutes for as long as the low water condition exists. (See Sections 4.10).



**Temperature Monitor and Alarm Control Panel
Figure 5-6**

5.7 REMOTE ALARM RELAY SYSTEM

Alarm states and alarm conditions are listed on Table 5-1. When the relay is off it is in the alarm state.

Alarms and Indicators, Table 5-1

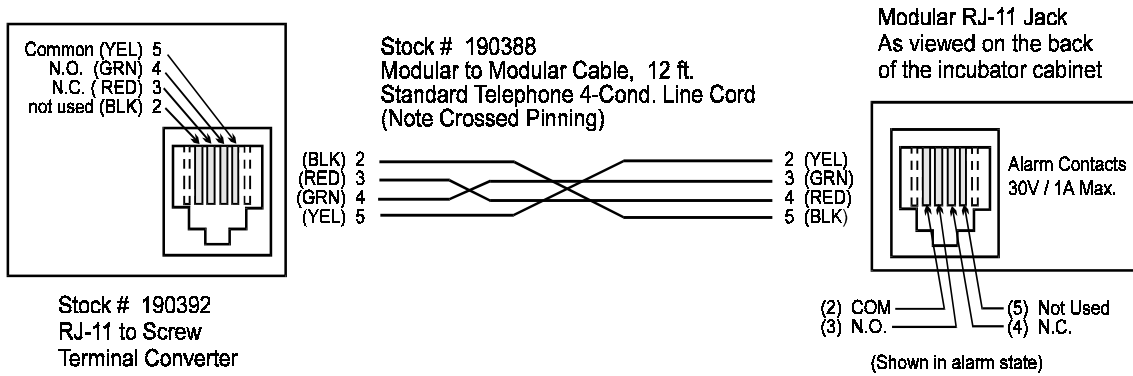
Alarm Condition	Visual Indicator	Audible Alarm	Reset Delay (Ringback)	Alarm Relay State
None	Off	Off		On
* Temp.> Over Temp Setpoint	Over Temp LED On	On	30 Minutes	Off
* Low Water in Water Jacket	Flash Add Water LED	On	30 Minutes	Off
** % CO ₂ Dev Setpoint ± 1%	CO ₂ Alarm LED On	On		Off
*** Tank1 Low	Pressure Low #1	On		Off
*** Tank2 Low	Pressure Low #2	On		Off
Power Failure	Off	Off		Off

* *Silenced audible alarm will ring back in 30 minutes if the alarm condition still exists. The Alarm LED will remain lit.*

** *Condition must exist for four minutes before alarm activates. Silenced alarm will extinguish the led alarm light, silence the audible alarm and change the alarm relay to the on state.*

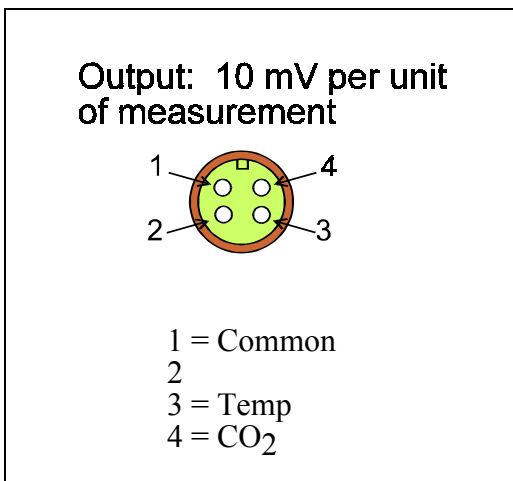
*** *When the CONTROL switch on the gas guard is turned off the indicator light will be extinguished and the audible alarm will be silenced. The built-in alarm relay on the Model 3158, however, remains active.*

A SPDT relay is provided for monitoring alarm conditions on the incubator. Connections are made by means of an RJ-11, telephone style jack, located on the rear of the cabinet. Figure 5-7 shows the connector wiring diagram. Any of the incubator alarms will activate the relay.



Remote Alarm Relay Wiring Diagram
Figure 5-7

5.8 RECORDER OUTPUT



Recorder Output Jack
Figure 5-8

Incubator temperature and CO₂ information may be directed to a recorder or other data storage device from a four-pin connector located on the back of the control panel. The connector pin numbers are identified in Figure 5-8 and are arranged as they would be seen looking at the back of the cabinet.

SECTION 6 - ROUTINE MAINTENANCE

6.1 INSTALLING THE DECONTAMINATION KIT

INSTRUCTION SHEET #7190028

INSTRUCTIONS FOR INSTALLING DECONTAMINATION KIT# 190028

1) **DISCONNECT INCUBATOR FROM POWER SUPPLY.**

2) Remove shelves, duct sheets and blower channel from incubator interior.

3) Remove blower wheel. To remove the blower wheel, grasp it as shown in Figure #1 and pull down firmly.



Figure #1

4) Match blower wheel from kit with blower wheel that was removed in Step #3. **IMPORTANT! THE BLOWER WHEEL REPLACEMENT MUST BE AN IDENTICAL MATCH.** Place new blower wheel over motor shaft and push blower wheel up against snap ring (if present). Confirm that blower wheel rotates freely.

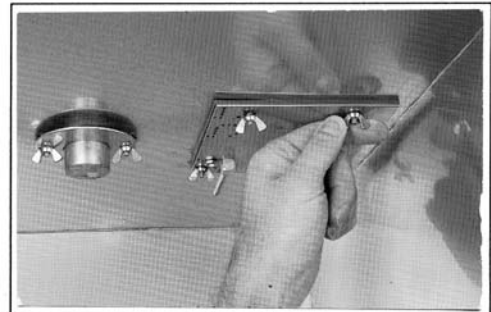


Figure #2

5) Next locate and remove the four wing nuts that secure the motor mounting plate to the incubator ceiling, shown in Figure #2.

6) Carefully pull blower motor assembly down into chamber area. **NOTE: ON SOME MODELS, THE BLOWER MOTOR ASSEMBLY MAY DROP DOWN ONLY 1-1/2 TO 2 INCHES.**

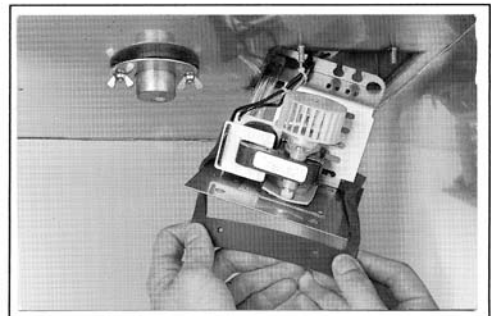


Figure #3

7) Slide the motor mounting gasket over the motor mounting plate, as shown in Figure #3. Discard gasket.

8) Install the new motor mounting gasket over the four studs, located on the incubator ceiling, as shown in Figure #4.



Figure #4

NOTE: THE MOTOR MOUNTING GASKET MUST LIE FLAT BETWEEN THE MOTOR MOUNTING PLATE AND THE INCUBATOR CEILING. TIGHTEN ALL FOUR WING NUTS FIRMLY TO ENSURE A GOOD SEAL.

THE FOLLOWING STEPS APPLY TO AUTOMATIC CO₂ AND AUTOMATIC O₂/CO₂ INCUBATORS. FOR CONSTANT FLOW INCUBATORS, SKIP TO STEP #12.

- 9) Remove the two wing nuts and large flat washer that secures the CO₂ sensor, shown in Figures #5 & #6. Allow sensor to drop down into chamber area. Disconnect sensor at electrical connector.
- 10) Peel o-ring off sensor and replace it with the new o-ring provided in kit, shown in Figure #7.
- 11) Electrically reconnect CO₂ sensor. Reinstall CO₂ sensor to original position. Tighten wing nuts firmly to ensure proper sealing.
- 12) Reinstall blower channel, duct sheets and shelves. Reconnect incubator to power supply.



Figure #5

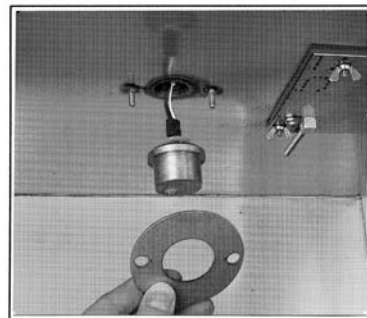


Figure #6

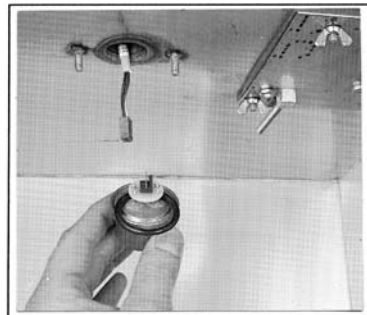


Figure #7

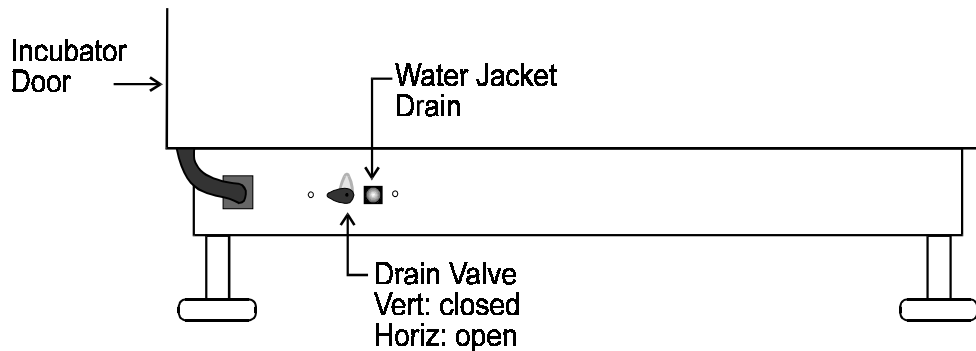
6.3 CLEANING THE CABINET EXTERIOR

The incubator exterior may be cleaned with soap and water and a general use laboratory disinfectant.

6.4 DRAINING THE WATER JACKET (Refer to Figure 6-1)

For best results, the water in the water jacket should be changed yearly.

1. Turn off the incubator and disconnect its electric plug from the power supply.
2. Remove the metal cover located below the door which protects the valve and the fill fitting. Two knurled thumb screws hold it in place.
3. Attach the 3/8" I.D. vinyl tubing to the drain fitting.
4. Route the tubing to a drain or large bucket and turn the valve to the vertical position..
5. If the water jacket is to be refilled, return the drain valve to the vertical position and review filling instructions in Section 4.10.



Incubator Base with Water Jacket Drain and Valve
Figure 6-1

6.5 CHANGING THE CO₂ FILTER (For Qualified Service Personnel Only)

Note: Under conditions of normal use, and dependent upon the purity of gas being used, the CO₂ filter should be last for about five years.

If the CO₂ filter becomes clogged, replace it with Forma stock #770001.

1. Disconnect the unit from the power supply.
2. Turn the two 1/4 turn fasteners on the control panel, grasp the CO₂ sample port, and slide the control panel out.
3. Locate the CO₂ filter at the back of the control panel. (Refer to Wiring Diagram, 140181-71-0-D for location)
4. Remove and replace the filter.

6.6 CO₂ TEST INSTRUMENTS

The best standard available for zeroing or calibrating the CO₂ controller is air, which typically contains 0.033% CO₂ by volume. Variations from this figure are insignificant.

It is very important that any Forma automatic CO₂ incubator be zeroed to air (See Section 4.15). *CO₂ test instruments should be used as a secondary check or to verify the percentage of CO₂ at operating levels.*

SECTION 7 - SERVICE

7.1 GENERAL TROUBLESHOOTING

Symptom	Possible Cause
Problem 1: Pilot lights not on. Readouts are dark. CO₂ power switch is on.	<ul style="list-style-type: none"> • Unit unplugged, or main power switch <i>not on</i>. • Circuit breaker tripped. • Outlet overload protection has tripped. • No voltage at outlet. • Defective incubator wiring.
Problem 2: Pilot lights on but heat light and alarm monitor module lights out.	<ul style="list-style-type: none"> • 0.3 Amp Slo-Blo fuse in control panel is blown.
Problem 3: CO₂ system indicator and readout are dark no control of CO₂ in chamber. Switch is on.	<ul style="list-style-type: none"> • Circuit breaker tripped. • Defective switch or wiring. • Defective CO₂ module.
Problem 4: CO₂ setpoint and display agree. FYRITE reads lower.	<ul style="list-style-type: none"> • FYRITE not zeroed. • Wool filter in FYRITE is dry. • FYRITE fluid need to be changed. • Chamber absolute humidity has decreased. • Control is incorrectly zeroed. • Defective CO₂ sensor or control.
Problem 5: Alarm sounds periodically, even with the alarm disabled. CO₂ alarm light not activated.	<ul style="list-style-type: none"> • Over temp control set too low. • Shorted heater. • Defective temp control or sensor.
Problem 6: CO₂ alarm and light are on. Alarm can be reset or disabled. If reset, alarm activates again.	<ul style="list-style-type: none"> • CO₂ level has deviated more than 1% from setpoint. • CO₂ supply has been interrupted. • CO₂ inject circuit is defective. • CO₂ solenoid is defective. • CO₂ sensor fan is defective.

Symptom	Possible Cause
<p>Problem 7: Digital CO₂ display and FYRITE read more than 1% different from setpoint. CO₂ alarm not on.</p>	<ul style="list-style-type: none"> • Alarm is disabled. • Defective alarm circuit. • Defective CO₂ sampler or control.
<p>Problem 8: CO₂ display and FYRITE read 0% CO₂. Setpoint is OK. CO₂ alarm is activated. CO₂ inject light is on.</p>	<ul style="list-style-type: none"> • Loss of CO₂ supply. • Defective CO₂ solenoid. • Clogged CO₂ in-line filter.
<p>Problem 9: CO₂ display shows some random number or decimal point.</p>	<ul style="list-style-type: none"> • Defective readout board. • Defective output to readout board from CO₂ control. • Faulty interconnecting wiring.
<p>Problem 10: CO₂ display will not go to zero no matter how long the door is left open.</p>	<ul style="list-style-type: none"> • Improper procedure. Close door and allow temperature and rh to stabilize eight hours. • Defective CO₂ sensor. • Defective CO₂ control.
<p>Problem 11: Display flickers badly or counts up and down three or four digits.</p>	<ul style="list-style-type: none"> • Defective readout board. • Excessive electrical interference near the cabinet. • Faulty grounding circuit. • Sampler blower is improperly installed or is defective.

Symptom	Possible Cause
<p>Problem 12: CO₂ setpoint cannot be changed.</p>	<ul style="list-style-type: none"> • Defective CO₂ control.
<p>Problem 13: Unit cannot be zeroed.</p>	<ul style="list-style-type: none"> • Defective CO₂ potentiometer. • Defective zero control. • Defective CO₂ Sensor.
<p>Problem 14: CO₂ overshoots setpoint badly.</p>	<ul style="list-style-type: none"> • Inoperative CO₂ sampler blower. • Defective CO₂ control. • CO₂ inlet pressure too high. • Unit is in over temp.
<p>Problem 15: Actual CO₂ is higher than setpoint and readout. Re-zeroing helps for a time, then symptoms return.</p>	<ul style="list-style-type: none"> • Defective CO₂ sensor. • Incorrect calibration of CO₂ control.
<p>Problem 16: CO₂ alarm sounds while CO₂ level is controlling at setpoint.</p>	<ul style="list-style-type: none"> • Excessive RFI or EMI near unit. • Faulty grounding circuit. • Defective CO₂ control.
<p>Problem 17: CO₂ setpoint changes by itself.</p>	<ul style="list-style-type: none"> • Faulty CO₂ potentiometer. • Defective CO₂ control.

7.2 CO₂ CONTROL MODULE CALIBRATION

These calibration procedures will be necessary only under the following conditions. DO NOT calibrate the incubator for any other reasons!

- ◆ After the CO₂ controller is replaced.
- ◆ After the CO₂ sensor is replaced.
- ◆ If there is reason to believe that the controller has previously been calibrated incorrectly.

Tools Needed:

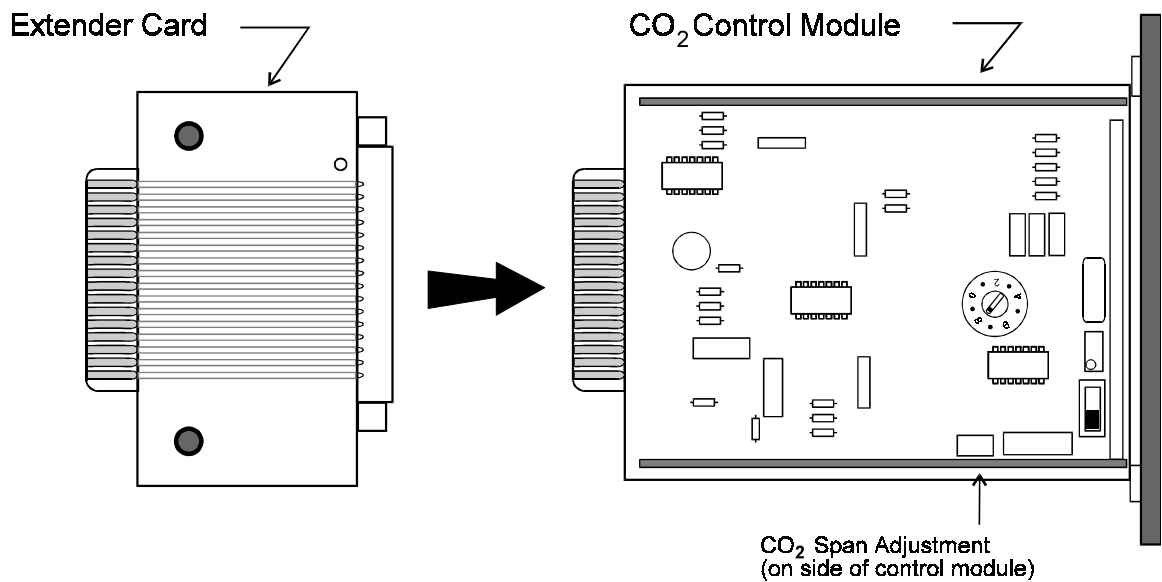
Calibration screwdriver (provided).

Circuit board extender card (provided).

FYRITE CO₂ Analyzer.

The following procedures are to be carried out in the sequence given. To prepare the incubator for calibration:

1. Turn off all electric power to the incubator.
2. Remove the CO₂ Control Module by pulling out the black plastic button locking button on the bottom of the module. (Figure 7-2.)
3. Firmly pull the button rearward until the circuit card releases from the internal connector strip.
4. Attach the extender card to the back of the control module circuit board. (Refer to Figure 7-1.)
5. Return the control module with extender card attached to the incubator control panel. Press the module firmly into the panel until the connector strip is seated. The control module will protrude a few inches from the incubator panel.



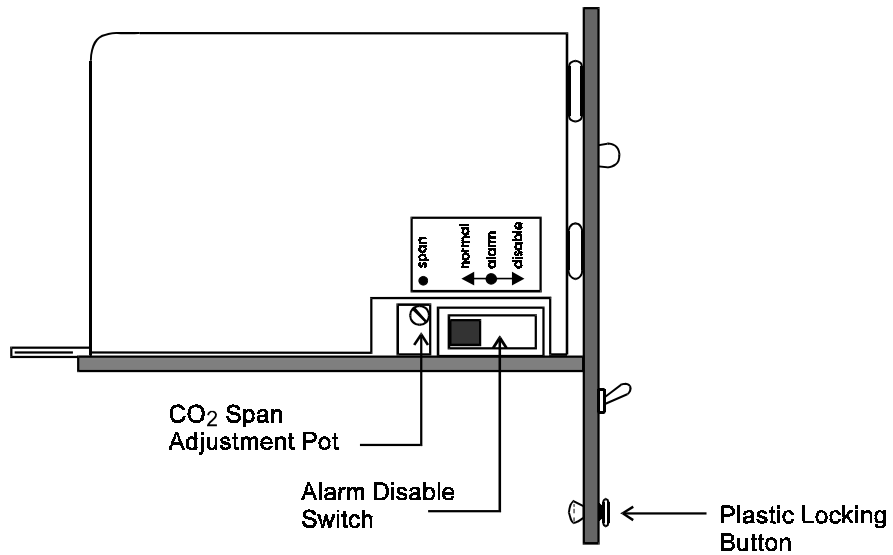
**CO₂ Control Module showing the Extender Card
and Span Adjustment Pot Location**

Figure 7-1

- a. Stabilize the Incubator at operating temperature and humidity with no CO₂ in the chamber.**
 1. Turn OFF the CO₂ supply at the source.
 2. Fill the humidity pan with sterile distilled water.
 3. Allow the chamber temperature and humidity to stabilize. This will take a minimum of eight hours if the temperature setpoint has been recently changed. Allow three days on initial start-up.
 4. Turn the span pot counterclockwise 10 turns. (Refer to Figures 7-1 and 7-2.)

b. Adjust the Zero Pot

1. Using the calibration screwdriver, adjust the CO₂ control zero pot to read 00.0 on the digital display.
2. Wait 5 minutes, and repeat as necessary until the display is stable.



**Side View of the CO₂ Control Module
Figure 7-2**

c. Adjust the Span Pot (Refer to Figure 7-2.)

1. Turn ON the CO₂ at the supply.
2. Turn the CO₂ setpoint to 10%. Allow the CO₂ to stabilize at 10% on the readout and control (inject light will cycle) for a minimum of 15 minutes.
3. Using a FYRITE or other measuring device, check the CO₂ level in the chamber until two consecutive readings agree.
4. Turn the CO₂ setpoint to 0.0% to prevent CO₂ from being injected into the chamber during the adjustment.
5. Adjust the span pot so the digital display agrees with the FYRITE reading.
6. Turn the setpoint back to 10%, and allow the CO₂ to control and stabilize for a minimum of 15 minutes.
7. Check the CO₂ in the chamber with a FYRITE or similar device until two consecutive readings agree. If the digital display is within plus or minus 1.0% of the FYRITE reading, proceed to the next step, "Recheck the Zero Adjustment." If the reading is not within plus or minus 1.0%, repeat Steps 3 through 6 of this procedure.

d. Re-check the Zero Adjustment

1. Turn OFF the CO₂ at the supply.
2. Open both doors wide for 45 seconds. Close the doors, and allow a minimum of 15 minutes for the incubator to stabilize and assure a zero CO₂ condition in the chamber.
3. If the readout is greater than 00.4, repeat the door opening for 15 seconds, and again allow the incubator to stabilize for a minimum of 15 minutes. If the display is not less than the previous reading, consult the factory. If the display now reads 00.4 or less, re-adjust the zero pot so the display reads 00.0.

e. Check the CO₂ at the desired setpoint

1. Turn ON the CO₂ at the supply.
2. Turn the CO₂ setpoint to the desired level.
3. Allow the incubator to reach setpoint and control for a minimum of 30 minutes.
4. Check the CO₂ with a FYRITE or similar device until two consecutive readings agree. If the FYRITE and display are not within 1.0%, consult the factory.

Note: After proper calibration the CO₂ display will be more accurate than the FYRITE, because the zero adjustment was made using atmospheric conditions.

Once the module is calibrated, the Extender Card must be removed and the CO₂ Control Module returned to its proper position in the incubator panel.

1. Turn off the incubator.
2. Firmly pull the Control Module rearward until it releases from the internal connector strip.
3. Remove the Extender Card.
4. Replace the Control Module into its location in the incubator panel.

7.3 37°C CONTROL CALIBRATION

If the digital display indicates a temperature other than 37°C when the chamber temperature has stabilized(temp select switch set to 37°C), it may be necessary to re calibrate the adjustment.

To calibrate:

1. Make a note of how much the display varies from 37°C.
2. Locate the 37°C calibration adjustment on the control panel. (See Figure 5-2 in Section 5.)
3. Using the calibration screwdriver, turn the calibration screw one turn clockwise for every half degree that the display is below 37°C. Turning the screw counterclockwise will lower the temperature.
4. Allow the temperature to stabilize, and check the display again. If it still varies significantly, repeat Steps 1 through 3.

Caution: Servicing should be performed by qualified service personnel only. Disconnect the incubator from the power source before starting service procedures.

7.4 REPLACING THE CO₂ SENSOR

1. Remove shelves, duct sheets, and blower channel from the chamber.
2. Locate CO₂ sensor, and remove wing nuts. The sensor will drop down.
3. There is a clip that connects sensor wiring to incubator wiring. To open the clip, pull out slightly on the clip tab to release the sensor.
4. Remove the O-ring, and place it on the new sensor.
5. Clip the new CO₂ sensor onto the incubator wiring and return the sensor to its original opening. Tighten the wing nuts securely to properly seal the O-ring.
6. After the sensor has been replaced, calibrate the CO₂ controller according to the instructions in Section 7.2.

7.5 REPLACING THE TRIAC (Refer to Wiring Diagram 140181-71-0-D)

1. Turn the incubator off and disconnect it from the power source.
2. Loosen the two 1/4-turn fasteners on the front of the control panel and pull the panel out.
3. Open the incubator outer door to gain access to the triac mounting screws from underneath the control panel.
4. Remove the two mounting screws which secure the triac.
5. Carefully disconnect the wires to the triac and note their connections.
6. Remove the triac from the unit. Note the thermal compound between the base of the triac and the floor of the control panel housing. If necessary, reapply more thermal compound before installing the new triac.
7. Install the new triac by reversing the above procedure.

7.6 REPLACING THE CO₂ SOLENOID (Refer to Wiring Diagram, 140181-71-0-D)

1. Loosen the two 1/4-turn fasteners on the front of the control panel and pull the panel out.
2. Locate the CO₂ solenoid and disconnect the tygon tubing attached to it. Note how tubing is installed so that it can be attached to new solenoid in same way.
3. Remove the screws that hold the solenoid in place.
4. Disconnect the wiring to the solenoid, making note of the connections.
5. Remove the solenoid and install the new unit by reversing the above procedure. Note the flow direction marked on solenoid.

7.7 REPLACING THE TEMPERATURE CONTROL (Refer to Wiring Diagram, 140181-71-0-D)

1. Loosen the two 1/4-turn fasteners on the front of the control panel and pull the panel out.
2. Remove the temperature control knob by loosening the two Allen head screws recessed in the knob.
3. Remove the nut on the shaft on the outside of the panel, then remove the nut on the back of the panel behind the control knob.
4. Remove the wiring to the temperature control and note the connections.
5. Remove the control and replace the new unit by reversing the above procedure.

7.8 REPLACING PILOT LIGHTS

1. Loosen the two 1/4-turn fasteners on the front of the control panel and pull the panel out.
2. Disconnect the wiring behind the defective light and break the retaining clip loose. Remove the pilot light.
3. Insert the new pilot light from the front of the panel and install the new retaining clip.
4. Reconnect the wiring.

7.9 REPLACING THE THERMISTOR

1. Loosen the two 1/4-turn fasteners on the front of the control panel and pull the panel out.
2. Locate the thermistors and remove the silastic around the thermistor cable.
3. Pull the defective thermistor out of the probe sheath.
4. Cut the wires to the defective thermistor, making note of the connections.
5. Attach the new thermistor using the electrical in-line connectors.
6. Install the new thermistor by reversing the above procedure. Take care not to damage the probe tip and make sure that the probe is fully extended in the sheath. Reseal the port with silastic or similar material.

7.10 REPLACING THE 0.3 AMP FUSE (Refer to Wiring Diagram, 140181-71-0-D)

The 3/10 amp fuse is located in the center of the control panel tray behind the CO₂ monitor module housing.

To replace the fuse:

1. Disconnect all power to the incubator.
2. Loosen the 1/4-turn fasteners on the front of the control panel and pull the panel out until the fuse holder in the center of the tray is accessible.
3. Use a fuse puller or a small screwdriver to lift one end of the fuse. Remove the fuse and discard it.
4. Insert a new 0.3 Amp fuse into the fuse holder.
5. Re-secure the control panel.
6. Connect power to incubator. Turn power switch ON and check operation of controls.

7.11 REPLACING THE BLOWER MOTOR

Replacing the Blower Motor on Single Units

1. Disconnect the unit from the power source.
2. Remove the shelves, duct sheets and blower channel from the incubator interior.
3. Remove and save the bottom blower wheel located inside the chamber by firmly pulling it down. Remove the snap ring (if present) from the motor shaft groove using needle-nose pliers.
4. Remove the four wing nuts that secure the blower mounting plate to the incubator ceiling.
5. Pull the blower motor assembly down into the chamber area.
6. Remove and save the upper blower wheel/blower blade.
7. Cut the tie wrap(s) securing the blower wires to the blower shield plate.
Note: An extra tie wrap has been provided for resealing the electrical wiring.
8. Disconnect the blower motor electrical connectors and remove the bushing securing the blower motor wiring.
9. Remove and save the two #6 nuts, washers and nylon spacers that secure the blower motor. Discard the old blower motor. *Save the #6 nuts, nylon spacers and washers.*
10. Remove the snap ring from the new blower motor shaft.
11. Install the new blower motor using the #6 nuts, nylon spacers and washers from step #9.
12. Reconnect the wires to the blower motor.
13. Install the snap bushing (located around the electrical wires from the blower motor) in the top of the blower shield plate.
14. Reinstall the new snap ring on the motor shaft groove.

15. Reinstall the lower blower wheel (inside the chamber) by pushing the blower wheel up against the snap ring.
16. Reinstall the upper blower wheel/blower blade.
17. Place the blower motor assembly into its original position.

Make sure that the vinyl tubing is not kinked and the wiring is not in contact with the top blower wheel.

Replacing the Blower Motor on the bottom Incubator of stacked units

1. Disconnect the incubator from the power source.
2. Remove the shelves, duct sheets and blower channel from the incubator interior.
3. Remove and save the bottom blower wheel (located inside the chamber) by firmly pulling it down. Remove the snap ring (if present) from the motor shaft groove using needle-nose pliers.
4. Remove the four wing nuts that secure the blower mounting plate to the incubator ceiling.
5. Disconnect the door heater plug on the top unit.
6. Remove the screws securing the base of the front cover plate from the top unit. Let cover plate hang on the left side of the cabinet.
7. Locate and disconnect the electrical connectors to the top unit's floor heater. Slide the floor heater forward and remove it.
8. Remove the duct assembly directly over the blower motor.
9. Gently pull the blower motor assembly down into the chamber area.
10. Remove and save the upper blower wheel/blower blade.
11. Cut the tie wrap(s) securing the wires from the blower motor to the blower shield plate. *An extra tie wrap has been provided for securing the electrical wiring.*

12. Disconnect the electrical connectors to the blower motor.
13. Remove the bushing securing the blower motor wiring.
14. Remove and save the two #6 nuts, washers and nylon spacers that secure the blower motor. Discard the old blower motor.
15. Remove the snap ring from new blower motor shaft.
16. Install the new blower motor using the #6 nuts, nylon spacers and washers saved in Step #14.
17. Reconnect the blower motor wires.
18. Install the snap bushing located around the wires from blower motor in the top of the blower shield plate.
19. Reinstall the new snap ring on the motor shaft groove.
20. Reinstall the lower blower wheel (inside chamber) by pushing the blower wheel up against the snap ring.
21. Reinstall the upper blower wheel/blower blade.
22. Replace the blower motor assembly, duct work and floor heater into their original positions.

Caution! Make sure that the vinyl tubing is not kinked.

SECTION 8 - SPECIFICATIONS

Models 3158/3164 and 3326/3336

Specifications:

Models	3158 (3164) Single Chamber 3154 Single Chamber Right hand door swing 3326 (3336) Dual Chamber 3354 Dual Chamber Right Hand Door Swing
---------------	--

Temperature

Control	$\pm 0.1^{\circ}\text{C}$ @ $+37^{\circ}\text{C}$ (98.6°F)
Range	$+5^{\circ}\text{C}$ above ambient to $+60^{\circ}\text{C}$ (140°F)
Sensor	Thermistor
Controller	Analog electronic proportional
Setpoint	Variable analog or fixed $+37^{\circ}\text{C}$
Display	Digital LCD
Uniformity	$\pm 0.2^{\circ}\text{C}$

Temperature Alarm

Sensor	Thermistor
Controller	Independent analog electronic
Setpoint	Digital with tamper resistant - screwdriver adjust
Display	Audible/visual
Readability	0.1°C
Setability	0.1°C

CO₂

CO₂ Gas Control	Better than $\pm 0.1\%$
Range	0-20%
Inlet Pressure	10 PSIG
Filter	0.2 Micron, disposable
Sensor	Thermal conductivity
Controller	Electronic, digital
Setpoint	Tamper resistant screwdriver adjust
Display	Digital LCD

Readability 0.1%
Setability 0.1%
Alarm Differential and Delay $\pm 1.0\%$ (nominal) for approx. 4 min.

Humidity

Humidity Control Ambient or 95% RH @ +37°C (98.6°F)
Humidity Reservoir 1.6 gal. (6 liters)
Humidity Pan 1.6 gal. (6 liters) optional

Fittings

Fill Port 1/4" FPT
Drain Port 1/4" FPT
Access Port 1.25" (3.17cm) with neoprene removable plug
CO₂ Inlet 1/4" barbed

Unit Heat Load (per chamber)

115V 480 BTUH (140W)
220V 560 BTUH (160W)

Shelves

Standard 5 per chamber
Maximum 21 per chamber
Dimensions 17.75" x 17.75" (45.10cm x 45.10cm)
Construction Stainless steel, perforated, electropolished
Surface Area 2.2 sq. ft. (0.2 sq. m) per shelf
Max. Per Chamber 46.2 sq. ft. (4.2 sq. m)
Flatness $\pm .032$ " off horizontal plane

Construction

Volume 11.5 gal. (43.5 liters)
Volume Per Chamber 5.7 cu. ft. (161.4 liters)
Interior 304 2B stainless steel
Exterior Cold rolled steel
Insulation 1-1/2" (3.8cm) Fiberglass

Inner Door	1/4" (0.62cm) Fully tempered safety glass with cam action latch
Inner Door Gasket	Non-porous feather edge silicone
Outer Door Gasket	Four sided molded magnetic vinyl
Finish	Powder coated for a durable, easily maintained surface

Electrical

3154/3158	90-130 Volts 50/60HZ, 1PH, 4.2 FLA
(3164)	180-260 Volts 50/60HZ, 1PH, 2.6 FLA
3326/3354	90-130 Volts 50/60HZ, 1PH, 8.0 FLA
(3336)	180-260 Volts 50/60HZ, 1PH, 3.8 FLA
Circuit Breaker	5 Amps/chamber
Power Switch	2 Pole
Line Cord	8 ft.
Data Output	10mv/unit temperature & CO ₂ (to Recorder output connector)
Accessory Outlet	120V (75 Watts max.)
Remote Alarm Contacts	Overtemperature, add water, power failure, CO ₂ deviation

Certification

CSA	Standard C22.2 No. 151
UL	Standard No. 1262 (except 3164/3336)

Dimensions

Exterior (Single)	24.37"W x 42.75"H x 23.31"F-B (61.91 cm x 108.59 cm x 59.22 cm)
Interior	19.75"W x 26.75"H x 18.75"F-B (50.17 cm x 67.95 cm x 47.63 cm)
Exterior (Stacked)	24.37"W x 83.00"H x 23.31"F-B (61.89 cm x 210.82 cm x 59.20 cm)

Weights	Single	Dual
Net	250 lbs. (114 Kg)	500 lbs. (228 Kg)
Net Operational	357 lbs. (162 Kg)	714 lbs. (324 Kg)
Shipping		
Motor	280 lbs. (127 Kg)	600 lbs. (272 Kg)
Air	345 lbs. (156 Kg)	700 lbs. (318 Kg)
Ocean	400 lbs. (181 Kg)	792 lbs. (359 Kg)

SECTION 9 - AUXILIARY EQUIPMENT

9.1 AUXILIARY EQUIPMENT

1. Electropolished Stainless Steel Shelf

The perforated shelf is square for easy installation. 22 shelf capacity.

Catalog # 224140

2. Tissue Culture Shelves

These 11" x 16" stainless steel shelves for culture dishes are designed for stacking to increase incubator storage space. Minimum order of 6 shelves.

Catalog # 500171

3. FYRITE CO₂ Analyzer Kit

For checking chamber CO₂ (0-20%) level in the chamber. Connects to the gas sample port on the control panel. Kit comes complete with aspirator, sampling tube, and carrying case.

Catalog # 220012

4. Extra FYRITE Fluid

For replacing FYRITE tester fluid. Three bottles per carton.

Catalog # 220051

5. Digital Thermometer

Hand-held, the size of a pocket calculator, it features clear liquid crystal readout of temperature from -99.9°C to +99.9°C. Ideal for general laboratory use.

Catalog # 853227

6. Two Stage Pressure Regulator

Controls CO₂ cylinder gas pressure. First stage reduces tank pressure to pre-set intermediate level. Second stage reduces pressure to recommended incubator inlet pressure. Permits stable CO₂ flow on high or low demand through the entire cylinder supply.

Catalog # 965010

7. CO₂ Gas Guard

For use with automatic CO₂ incubators only. Protects dual tank CO₂ supply by automatically switching to another tank when one supply is exhausted. Audible alarm warns of tank depletion; reset button silences the alarm. Manual tank switch-over included.

Model # 3050

8. Replacement CO₂ Filter

Disposable 99.97% Microbiological filter to replace the inline CO₂ filter when it becomes clogged.

Catalog # 770001

9. Disinfecting Kit

Includes parts needed for improved disinfecting of the incubator. Includes disposable blower wheel, O-ring for the CO₂ sensor and motor mounting gasket.

Catalog # 190247

10. Humidity Pan

This humidity pan is made of stamped (seam-free) 304 stainless steel and will hold 6.375 quarts (6 liters) of water. Cleaning and decontamination is very easy due to the cove-shaped corners and seamless construction.

Catalog # 237001

11. Hygrometer

For measuring chamber humidity. This instrument has a 4" dial and a range of 0 to 100% rh.

Catalog # 155010

12. Glass Thermometer

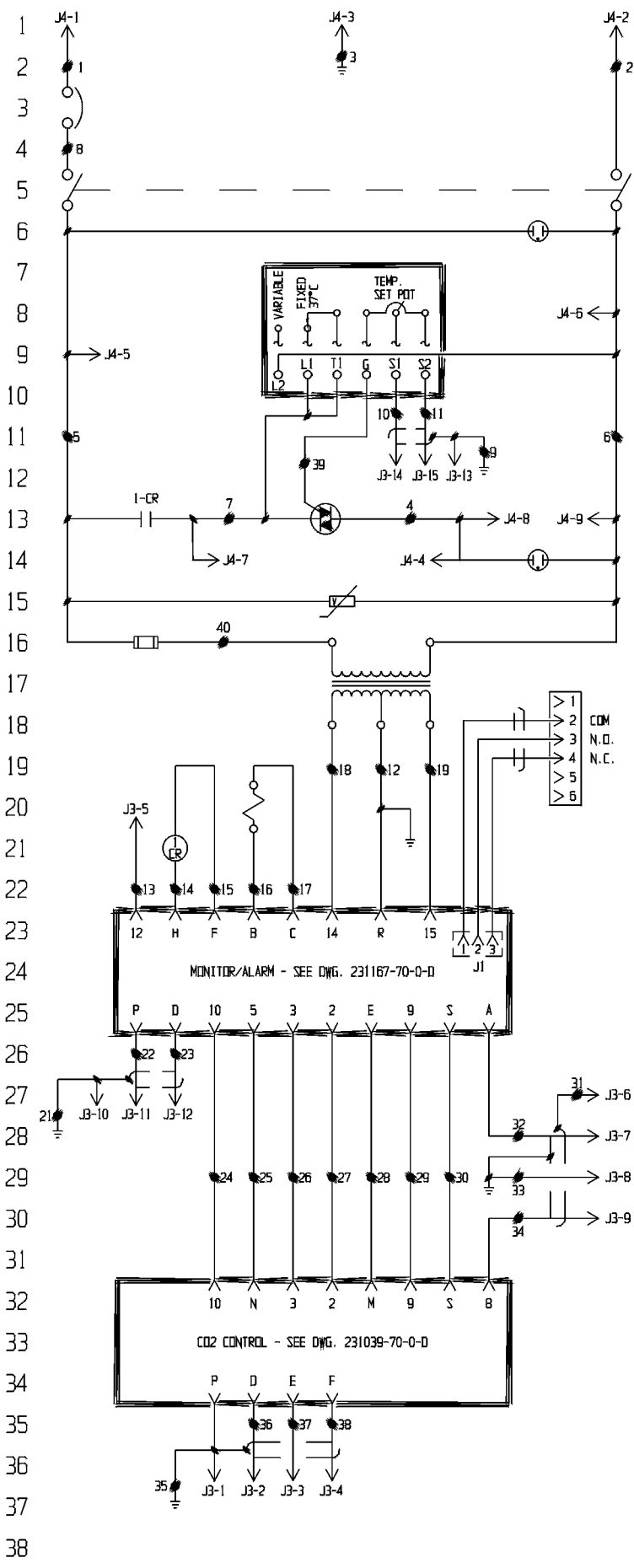
For independently measuring chamber temperature. Range from 0°C to 100°C.

Catalog # 285722

SECTION 10 - PARTS LIST

STOCK #	DESCRIPTION
770001	CO ₂ Disposable Filter
250085	CO ₂ Solenoid Valve 12VDC, 3W
300164	Over Temp Relay
431142	Four Sided Molded Magnetic Gasket
400296	Silicon Feather Gasket
230094	Glass Fuse .3A (Slow-Blow Type)
280004	#312 Green Pilot Light
280006	#312 Amber Pilot Light
111008	CLIP, Mounting, Pilot Light
285379	25A Triac
100071	Blower Wheel
227175	Blower Motor
231167	Alarm/Monitor Module
231039	CO ₂ Control Module
231047	Temperature Control
290027	1K, Thermistor Cable
285057	Extender Card Module

SECTION 11 - SCHEMATICS



CIRCUIT BREAKER, 5A
230023

POWER SWITCH
360155

POWER INDICATOR, GREEN
280004

PROPORTIONAL TEMP CONTROL
231047

TRIAC, 25A
265379

HEAT INDICATOR, AMBER
280006

METAL OXIDE VARISTOR
265823

FUSE, 0.3A TIME DELAY
230094

TRANSFORMER, 28VCT
420017

REMOTE ALARM
CONTACT OUTPUT
HARNESS ASSEMBLY
190408

CO2 SOLENOID
250085

OVERTEMP CONTROL
300164 13

CARD EDGE CONNECTOR
370065
(FOR MONITOR/ALARM MODULE)

CARD EDGE CONNECTOR
370065
(FOR CO2 CONTROL MODULE)

39 EXTERNAL CONNECTIONS

- 40 J4-1 LINE IN, 90-130VAC
- 41 J4-2 NEUTRAL IN, 90-130VAC
- 42 J4-3 GROUND
- 43 J4-4 CHAMBER HEATER, 260W; DOOR HEATER, 120W (CYCLED)
- 44 J4-5 ACCESSORY OUTLET
- 45 J4-6 NEUTRAL OUT
- 46 J4-7 FAN MOTOR; DOOR HEATER, 5W
- 47 J4-8 CHAMBER HEATER, 6W (CYCLED)
- 48 J4-9 CHAMBER HEATER, 6W (NEUTRAL)
- 49 J3-1 CO₂ SENSOR, SHIELD
- 50 J3-2 CO₂ SENSOR, REFERENCE (SEALED) CELL
- 51 J3-3 CO₂ SENSOR, COMMON
- 52 J3-4 CO₂ SENSOR, SAMPLE (OPEN) CELL
- 53 J3-5 WATER LEVEL SENSOR
- 54 J3-6 RECORDER, SHIELD
- 55 J3-7 RECORDER, TEMP OUTPUT (10mV / °C)
- 56 J3-8 RECORDER, OUTPUT COMMON
- 57 J3-9 RECORDER, CO₂ OUTPUT (10mV / %CO₂)
- 58 J3-10 OVERTEMP SENSOR, SHIELD
- 59 J3-11 OVERTEMP SENSOR
- 60 J3-12 OVERTEMP SENSOR
- 61 J3-13 CONTROL SENSOR, SHIELD
- 62 J3-14 CONTROL SENSOR
- 63 J3-15 CONTROL SENSOR
- 64
- 65
- 66
- 67
- 68
- 69
- 70
- 71
- 72
- 73
- 74
- 75
- 76

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

NUMBER	GAUGE	COLOR
1	16	BLK
2	16	WHT
3	16	GRN
4	20	RED
5	20	BLK
6	20	WHT
7	20	BLK
8	20	BLK
9	24	SHLD
10	24	BLK
11	24	WHT or CLR
12	20	GRN
13	20	BRN
14	20	BLU
15	20	BLU
16	20	ORG
17	20	ORG
18	20	YEL
19	20	YEL
20	NOT USED	
21	24	SHLD
22	24	BLK
23	24	WHT or CLR
24	20	BLK
25	20	YEL
26	20	RED
27	20	BRN
28	20	RED
29	20	ORG
30	20	GRN
31	22	SHLD
32	22	RED
33	22	BLK
34	22	WHT or CLR
35	22	SHLD
36	22	RED
37	22	BLK
38	22	WHT or CLR
39	20	YEL
40	20	BLK

} 24/2 CABLE

} 24/2 CABLE

} 22/3 CABLE

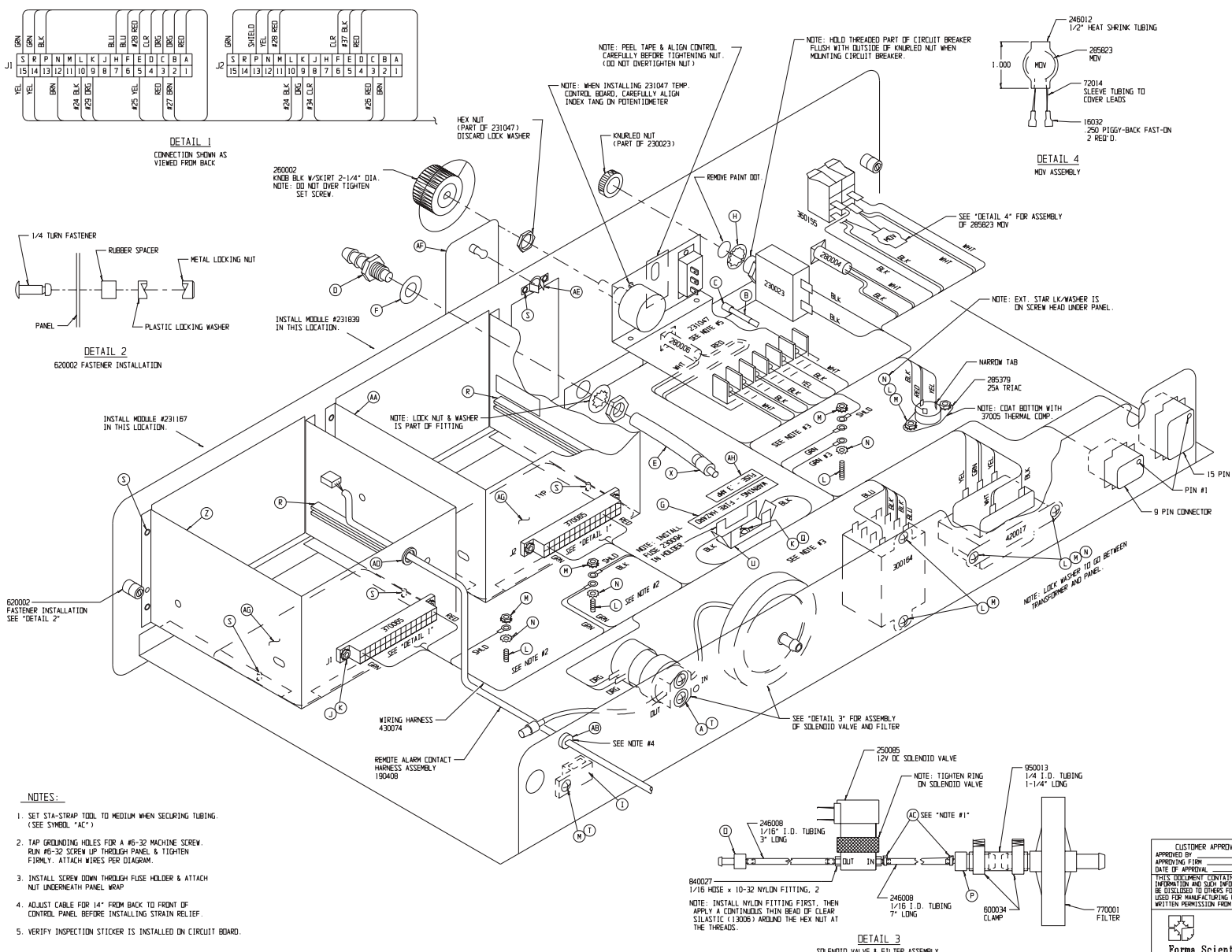
} 22/3 CABLE

NOTES:

⊗ Denotes Terminal Strip Connection	Parts List Reference Number
I-CR Last Relay Number	○ Assembly
N/A Last Terminal Number	◇ Panel
40 Last Wire Number	○ Refrigeration
	□ Wiring

CUSTOMER APPROVAL/REFERENCE						
APPROVED BY _____						
APPROVING FIRM _____						
DATE OF APPROVAL _____						
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						
0	N/A	08-19-93	GLM	KDG	RELEASED FOR PRODUCTION	
REV	ECN NO.	DATE	BY	CAD/APPD	DESCRIPTION OF REVISION	
		8-19-93	DWN	DAG	CAD	APPD LDN SCALE N/A
CUSTOMER STK NO. 140181						
JOB TITLE AUTOMATIC CO2 CONTROL PANEL						
DWG TITLE ELECTRICAL SCHEMATIC						
LOCATION			JOB NUMBER		DRAWING NUMBER	
STKND001					140181-70-0-D	





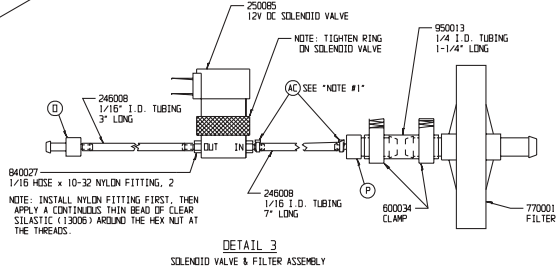
HARDWARE CHART

SYM	STOCK NO.	DESCRIPTION
A	23020	#6 FLAT WASHER
B	235013	SCREW-DRIVER
C	710002	STUD RECEIVER
D	840020	1/4" HOSE FITTING
E	950019	1/4" I. D. TUBING, 3" LONG
F	730017	3/8" BLACK FIBER WASHER
G	220437	WARNING LABEL
H	730019	7/16" INTERNAL TOOTH LOCK WASHER
I	30001	TIE WRAP ANCHOR
J	22002	#4-40 x 1/2" PHP SCREW
K	23006	#4-40 KEPS NUT
L	22049	#6-32 x 3/8" PHP SCREW
M	23001	#6-32 KEPS NUT
N	23058	#6 EXTERNAL STAR LOCK WASHER
O	840011	3/16" HS x 1/16" HS BRASS REDUCER
P	840013	1/4" HS x 1/16" HS BRASS REDUCER
Q	22130	#4-40 x 1/4" PHP SCREW
R	124001	CARD GUIDE
S	67000	COUNTERSUNK POP RIVET
T	22115	#6-32 x 1/4" PHP SCREW
U	230039	FUSE CLIP
X	840023	1/4" x 1/8" HOSE FITTING
Y		
Z	190267	TEMPERATURE MODULE BOX
AA	190268	CO ₂ MODULE BOX
AB	340036	3/8" STRAIN RELIEF
AC	30100	AUTO CABLE TIE (SEE NOTE #2)
AD	30015	SNAP BUSHING 3/4"
AE	710017	BALL STUD SPEED NUT CLIP
AF	190416	BLANK COVER PLATE
AG	190266	BOX BOTTOM
AH	220440	3 AMP LABEL

GROUND CONTINUITY CHECK LOCATIONS WITH METER, MUST MEASURE 0.1 OHMS OR LESS RESISTANCE.

- 1) 9 PIN CONNECTOR
PIN #3
- 2) 15 PIN CONNECTOR
PIN #1
PIN #5
PIN #10
PIN #13
- 3) CIRCUIT BREAKER (KNURED NUT)
- 4) TRIAC MOUNTING SCREWS
- 5) TRANSFORMER MOUNTING SCREWS

- NOTES:**
1. SET STA-STRAP TOOL TO MEDIUM WHEN SECURING TUBING. (SEE SYMBOL "AC")
 2. TAP GROUNDING HOLES FOR A #6-32 MACHINE SCREW. RUN #6-32 SCREW UP THROUGH PANEL & TIGHTEN FIRMLY. ATTACH WIRES PER DIAGRAM.
 3. INSTALL SCREW DOWN THROUGH FUSE HOLDER & ATTACH NUT UNDERNEATH PANEL WRAP
 4. ADJUST CABLE FOR 14" FROM BACK TO FRONT OF CONTROL PANEL BEFORE INSTALLING STRAIN RELIEF.
 5. VERIFY INSPECTION STICKER IS INSTALLED ON CIRCUIT BOARD.

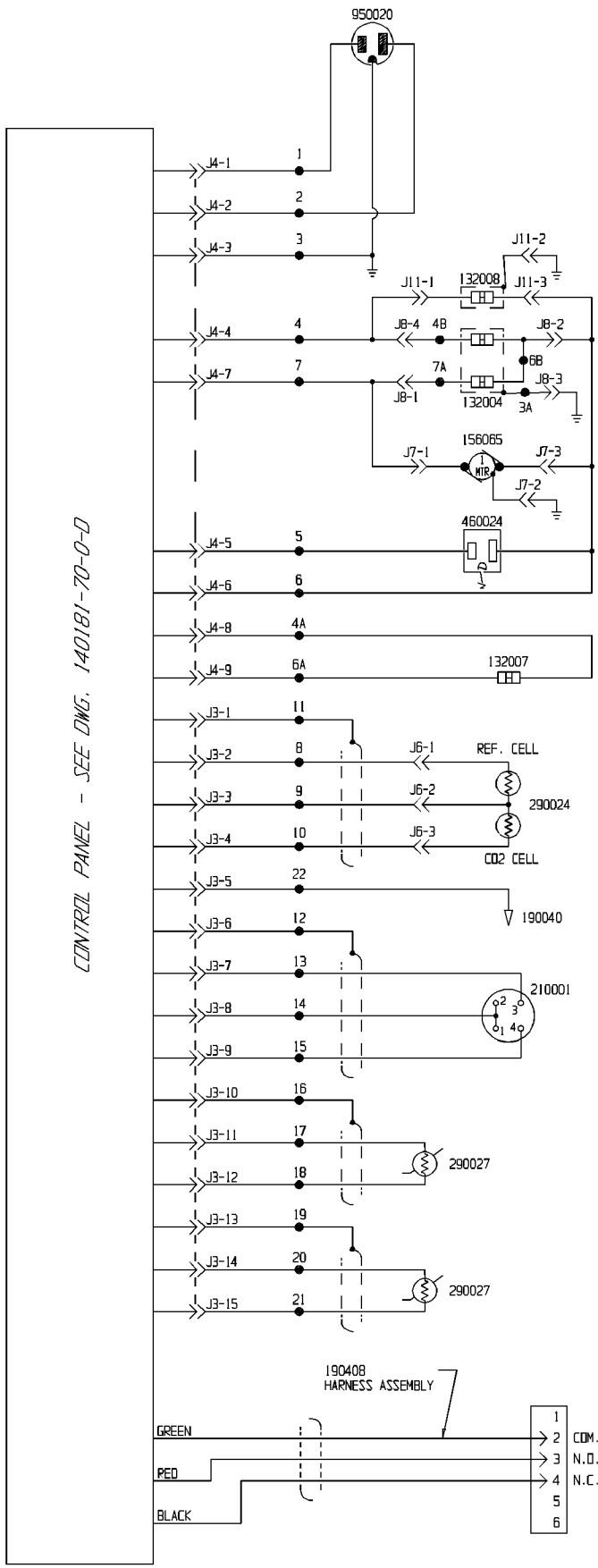


APPROVED BY	CUSTOMER APPROVAL/REFERENCE	DATE OF APPROVAL	THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM FORMA SCIENTIFIC
	6 S1-5067 03-23-95 GCF KCG		ENG. STRAIN RELIEF & REV. REF. TO NOTES
	5 S1-4893 01-04-95 BOB POK LON		REVISED LABEL, STOCK NUMBERS
	4 S1-4713 08-27-94 GCF KCG LON HOD MBR		STAR WASHER & NOTES REFILES
	3 S1-4567 09-23-94 GCF KCG LON REV.		GROUND CONTINUITY MEASUREMENT
	2 S1-4417 04-18-94 AFC KCG LON		ADD VERIFY INSPECTION STICKER NOTE
	1 S1-4300 03-18-94 GCF KCG LON		INITIAL DESIGN
			DATE 08-19-93 DWN DAG CAD POK APPD LON SCALE N.T.S.
			CUSTOMER
			JOB TITLE AUTOMATIC CO ₂ CONTROL PANEL
			DWG TITLE WIRING DIAGRAM LAYOUT
			LOCATION
			JOB NUMBER
			DRAWING NUMBER
			STKND501 140181-71-0-D

Forma Scientific
 800 469-8874, 800 469-8875, 800 469-8876
 TEL: 603 882-9400 FAX: 603 882-9400

POWER CONNECTION
 90-130VAC, 1PH, 50/60HZ, 4.2FLA

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



CONTROL PANEL - SEE DWG. 140181-70-0-D

- 16/3 LINE CORD
- CHAMBER HEATER, 260W
- ODOR HEATER, 120W
- ODOR HEATER, 5W
- BLOWER MOTOR
- ACCESSORY OUTLET
75W MAXIMUM
- CHAMBER HEATER, 6W
- REF. CELL
- CO2 SENSOR
- ADD WATER SENSOR
- RECORDER OUTPUT 10mv/UNIT
PINS 1 & 2 = OUTPUT COMMON
PIN 3 = TEMP OUTPUT
PIN 4 = CO2 OUTPUT
- OVERTEMP SENSOR
1K @ 25°C
- TEMP SENSOR
1K @ 25°C
- REMOTE ALARM
CONTACTS, 1A 30V MAX.
(SHOWN IN ALARM STATE)

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

WIRE REFERENCE CHART

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 #	GAUGE	COLOR	
1	16	BLK	
2	16	WHT	
3	16	GRN	
3A	18	GRN	
4	20	RED	
4A	18	BLK	
4B	18	RED	
5	20	BLK	
6	20	WHT	
6A	18	BLK	
6B	18	WHT	
7	20	BLK	
7A	18	BLK	
8	22	RED	} 22/3 CABLE
9	22	BLK	
10	22	WHT or CLR	
11	22	SHLD	} 22/3 CABLE
12	22	SHLD	
13	22	RED	
14	22	BLK	} 24/2 CABLE
15	22	WHT or CLR	
16	24	SHLD	
17	24	BLK	} 24/2 CABLE
18	24	WHT or CLR	
19	24	SHLD	
20	24	BLK	} 24/2 CABLE
21	24	WHT or CLR	
22	20	BRN	

CUSTOMER APPROVAL/REFERENCE		8	IN-1642	08-30-93	GLM	KDG		ADD REMOTE ALARM CONTACTS
APPROVED BY _____		7	IN-988	12-15-88	JAS	POK	LON	REVISED BLOWER MOTOR
APPROVING FIRM _____		6	IN-552	12-11-84	DWL	DWL	LON	REVISED BLOWER MOTOR
DATE OF APPROVAL _____		5	SI-390	4-18-83	DWL	DWL	LON	SEE ECR
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		4	IN-434	12-28-82	DWL	DWL	LON	ADDED VOLTAGE RANGE
		REV	ECR NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION
DATE		6-12-81	DWN	DAG	CAD	DWL	APPD	LON
CUSTOMER								
JOB TITLE		3158 AUTOMATIC CO2 INCUBATOR						
DWG TITLE		ELECTRICAL SCHEMATIC						
LOCATION		JOB NUMBER		DRAWING NUMBER				
INCUBATRO1				3158-70-0-D				

NOTES:

⊗	Denotes Terminal Strip Connection
N/A	Last Relay Number
15	Last Terminal Number
22	Last Wire Number

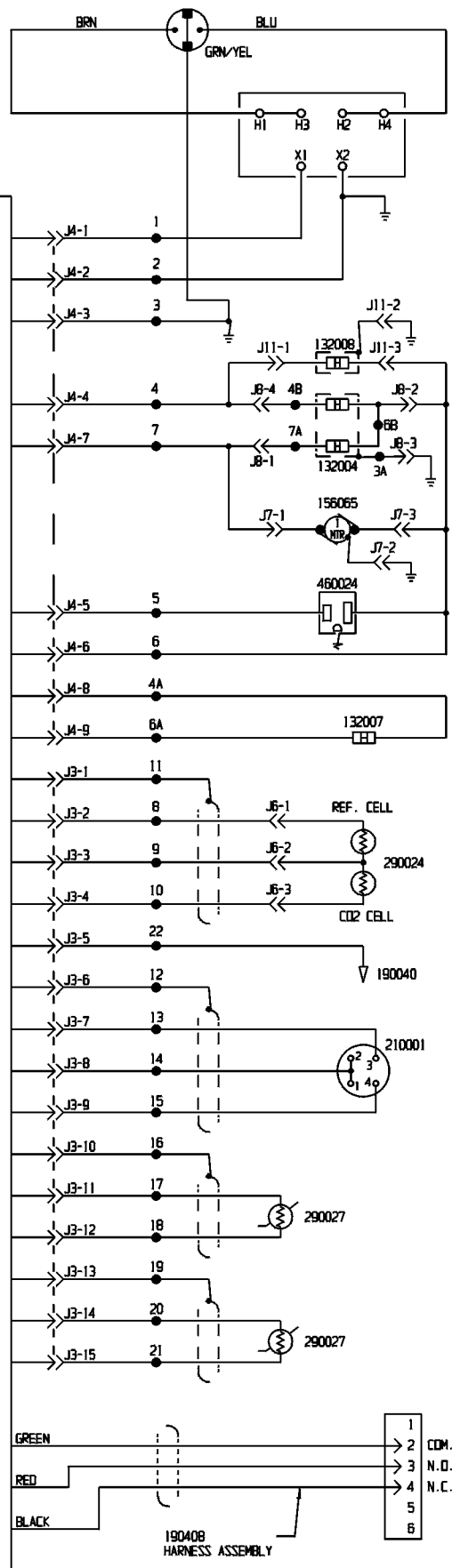


Forma Scientific
 BOX 649 HAZLETTA, OHIO 43130 TELEX 24-3394
 TOLL FREE USA 800-948-3080, OHIO 614-973-4763

POWER CONNECTION
180-260V, 1PH, 50/60HZ, 2.6FLA

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

CONTROL PANEL - SEE DWG. 140181-70-0-D



- 16/3 LINE CORD
430017
- STEPODOWN TRANSFORMER
.5 KVA
420089
- CHAMBER HEATER, 260W
- DOOR HEATER, 120W
- DOOR HEATER, 5W
- BLOWER MOTOR
- ACCESSORY OUTLET
75W MAXIMUM
- CHAMBER HEATER, 6W
- CO2 SENSOR
- ADD WATER SENSOR
- RECORDER OUTPUT 10mv/UNIT
PINS 1 & 2 = OUTPUT COMMON
PIN 3 = TEMP OUTPUT
PIN 4 = CO2 OUTPUT
- OVERTEMP SENSOR
1K @ 25°C
- TEMP SENSOR
1K @ 25°C
- REMOTE ALARM
CONTACTS, 1A 30V MAX.
(SHOWN IN ALARM STATE)

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

WIRE REFERENCE CHART

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 #	GAUGE	COLOR	
1	16	BLK	
2	16	WHT	
3	16	GRN	
3A	18	GRN	
4	20	RED	
4A	18	BLK	
4B	18	RED	
5	20	BLK	
6	20	WHT	
6A	18	BLK	
6B	18	WHT	
7	20	BLK	
7A	18	BLK	
8	22	RED	} 22/3 CABLE
9	22	BLK	
10	22	WHT or CLR	
11	22	SHLD	} 22/3 CABLE
12	22	SHLD	
13	22	RED	
14	22	BLK	} 24/2 CABLE
15	22	WHT or CLR	
16	24	SHLD	
17	24	BLK	} 24/2 CABLE
18	24	WHT or CLR	
19	24	SHLD	
20	24	BLK	} 24/2 CABLE
21	24	WHT or CLR	
22	20	BRN	

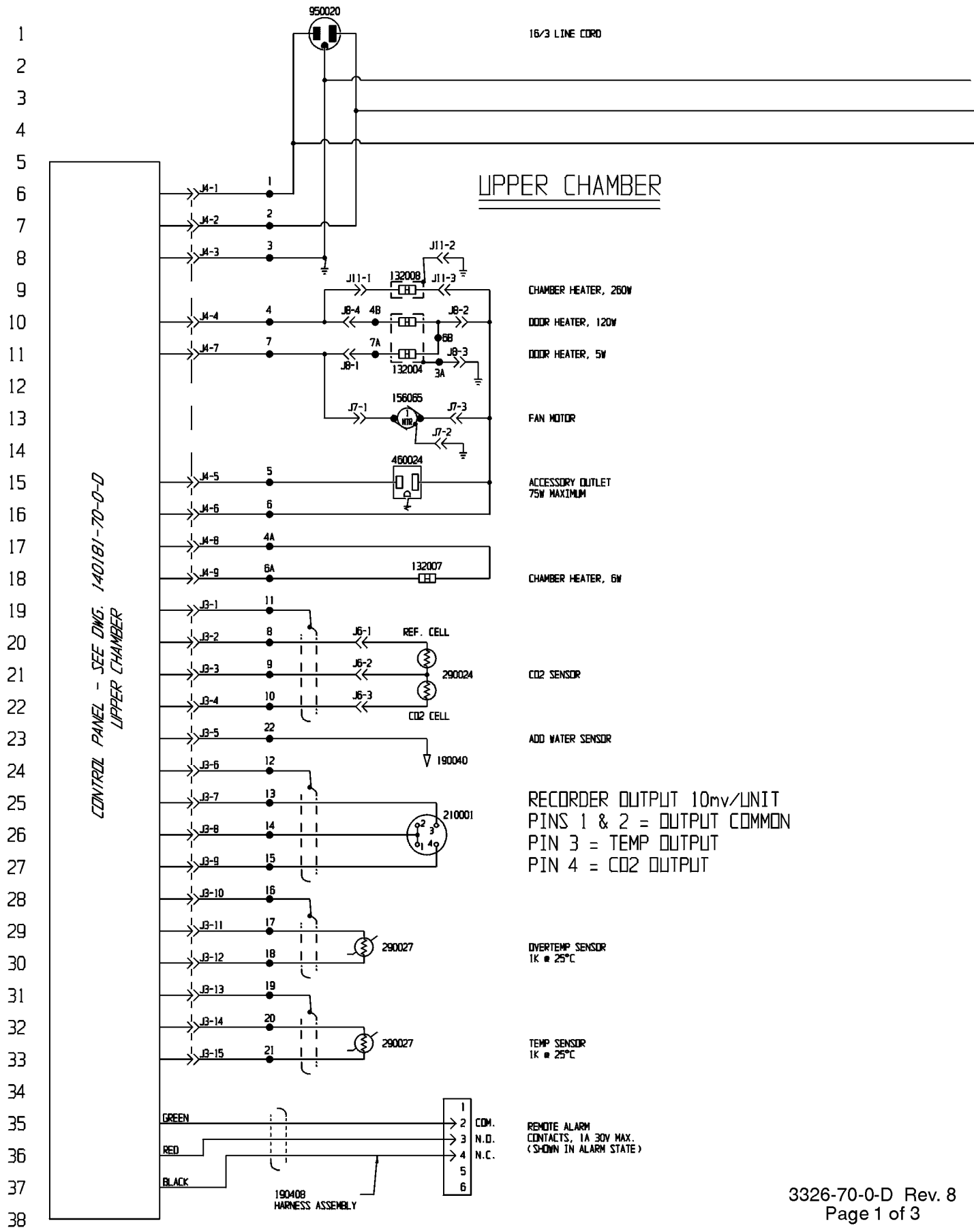
NOTES:
 ⊗ Denotes Terminal Strip Connection
 N/A Lost Relay Number
 15 Lost Terminal Number
 22 Lost Wire Number

CUSTOMER APPROVAL/REFERENCE		9	SI-4877	01-24-95	PAR	KDG		CHG. 275034 TRANSFORMER TO 420089		
APPROVED BY		8	IN-1642	08-30-93	GLM	KDG	LON	ADD REMOTE ALARM CONTACTS		
APPROVING FIRM		7	IN-988	12-15-89	JAS	POK	LON	REVISED FAN MOTOR		
DATE OF APPROVAL		6	IN-552	12-11-84	DWL	DWL	LON	REVISED FAN MOTOR		
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		5	SI-390	4-19-83	DWL	DWL	LON	REV'D. WIRING OF 6W HEATER		
REV	ECR NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION				
	DATE	6-21-81	DWN	DAG	CAD	DWL	APPD	LON	SCALE	N/A
CUSTOMER										
JOB TITLE 3164 AUTOMATIC CO2 INCUBATOR										
DWG TITLE ELECTRICAL SCHEMATIC										
LOCATION					JOB NUMBER		DRAWING NUMBER			
INCLIBATRO1							3164-70-0-D			



Forma Scientific
 BOX 640 MARICETTA, OHIO 45750 TEL# 614-24-5304
 TOLL FREE USA 800-948-3800, INTL 614-272-4763

POWER CONNECTION
90-130V, 1PH, 2W, 50/60HZ, 8FLA



CONTROL PANEL - SEE DWG. 140181-70-0-D
UPPER CHAMBER

UPPER CHAMBER

16/3 LINE CORD

CHAMBER HEATER, 260W

DOOR HEATER, 120W

DOOR HEATER, 5W

FAN MOTOR

ACCESSORY OUTLET
75W MAXIMUM

CHAMBER HEATER, 6W

CO2 SENSOR

ADD WATER SENSOR

RECORDER OUTPUT 10mv/UNIT
PINS 1 & 2 = OUTPUT COMMON
PIN 3 = TEMP OUTPUT
PIN 4 = CO2 OUTPUT

OVERTEMP SENSOR
1K @ 25°C

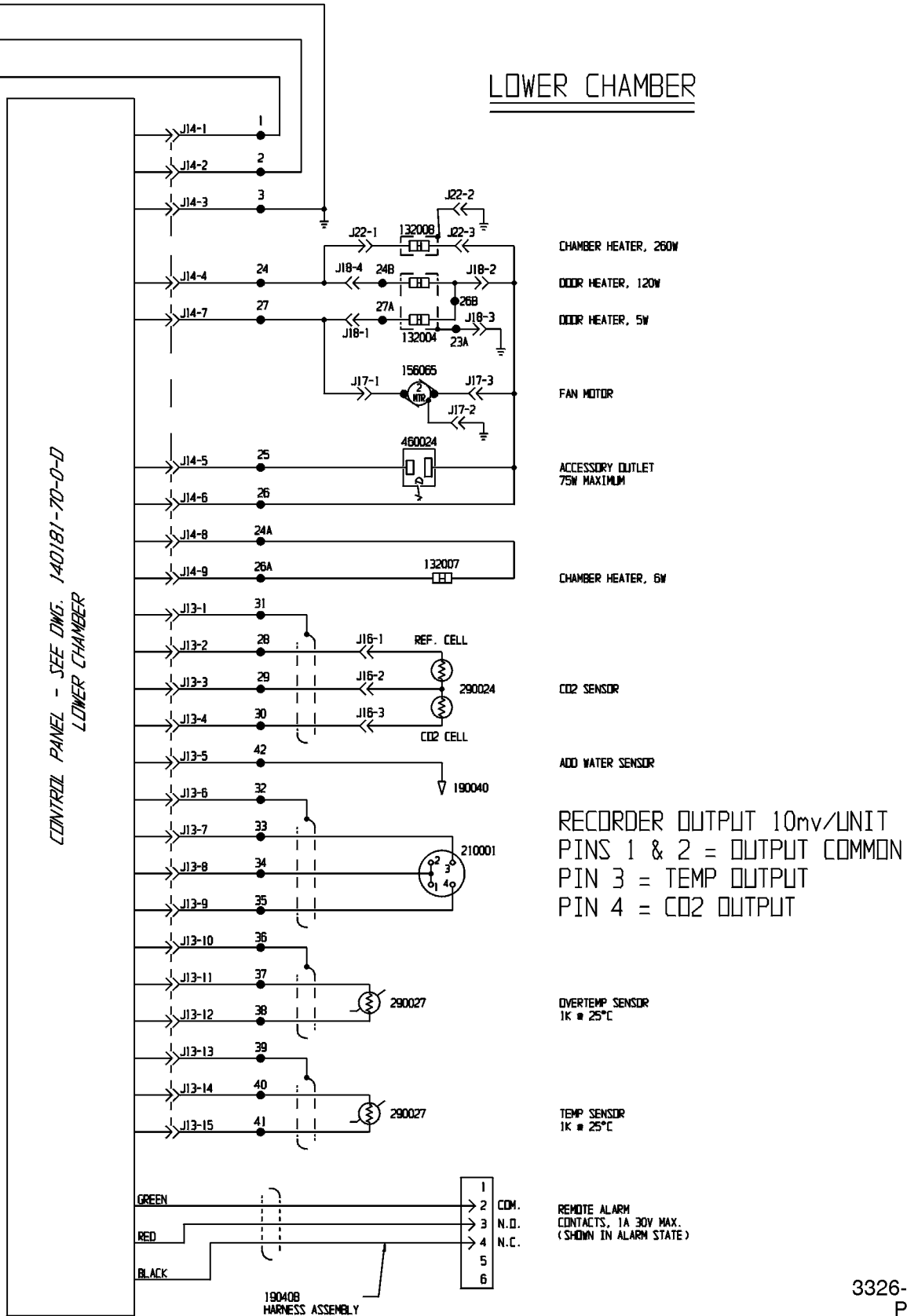
TEMP SENSOR
1K @ 25°C

REMOTE ALARM
CONTACTS, 1A 30V MAX.
(SHOWN IN ALARM STATE)

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

CONTROL PANEL - SEE DWG. 140181-70-0-D
LOWER CHAMBER

LOWER CHAMBER



- CHAMBER HEATER, 260W
- DOOR HEATER, 120W
- DOOR HEATER, 5W
- FAN MOTOR
- ACCESSORY OUTLET
75W MAXIMUM
- CHAMBER HEATER, 6W

- REF. CELL
- CO2 SENSOR
- ADD WATER SENSOR

RECORDER OUTPUT 10mv/UNIT
PINS 1 & 2 = OUTPUT COMMON
PIN 3 = TEMP OUTPUT
PIN 4 = CO2 OUTPUT

- OVERTEMP SENSOR
1K @ 25°C
- TEMP SENSOR
1K @ 25°C

REMOTE ALARM
CONTACTS, 1A 30V MAX.
(SHOWN IN ALARM STATE)

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

	U.C. WIRE #	GAUGE	COLOR	L.C. WIRE #
	1	16	BLK	1
	2	16	WHT	2
	3	16	GRN	3
	3A	18	GRN	23A
	4	20	RED	24
	4A	18	BLK	24A
	4B	18	RED	24B
	5	20	BLK	25
	6	20	WHT	26
	6A	18	BLK	26A
	6B	18	WHT	26B
	7	20	BLK	27
	7A	20	BLK	27A
	8	22	RED	28
22/3 CABLE	9	22	BLK	29
	10	22	WHT or CLR	30
	11	22	SHLD	31
	12	22	SHLD	32
22/3 CABLE	13	22	RED	33
	14	22	BLK	34
	15	22	WHT or CLR	35
	16	24	SHLD	36
24/2 CABLE	17	24	BLK	37
	18	24	WHT or CLR	38
	19	24	SHLD	39
24/2 CABLE	20	24	BLK	40
	21	24	WHT or CLR	41
	22	24	BRN	42

CUSTOMER APPROVAL/REFERENCE	8	IN-1642	08-30-93	GLM	KDG		ADD REMOTE ALARM CONTACTS
APPROVED BY	7	IN-988	12-21-89	JAS	POK	LDN	REVISED BLOWER MOTOR
APPROVING FIRM	6	IN-552	12-11-84	DWL	DWL	LDN	REVISED BLOWER MOTOR
DATE OF APPROVAL	5	SI-390	4-19-83	DWL	DWL	LDN	SEE ECR
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	4	IN-434	12-23-82	DWL	DWL	LDN	ADDED VOLTAGE RANGE
	REV	ECR NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION

NOTES:
 ⊗ Denotes Terminal Strip Connection
 N/A Last Relay Number
 15 Last Terminal Number
 42 Last Wire Number



DATE	6-15-81	DWN	DAG	CAD	DWL	APPD	LDN	SCALE	N/A
CUSTOMER									
JOB TITLE 3326 AUTOMATIC CO2 INCUBATOR									
DWG TITLE ELECTRICAL SCHEMATIC									
LOCATION					JOB NUMBER		DRAWING NUMBER		
INCUBAT01							3326-70-0-D		

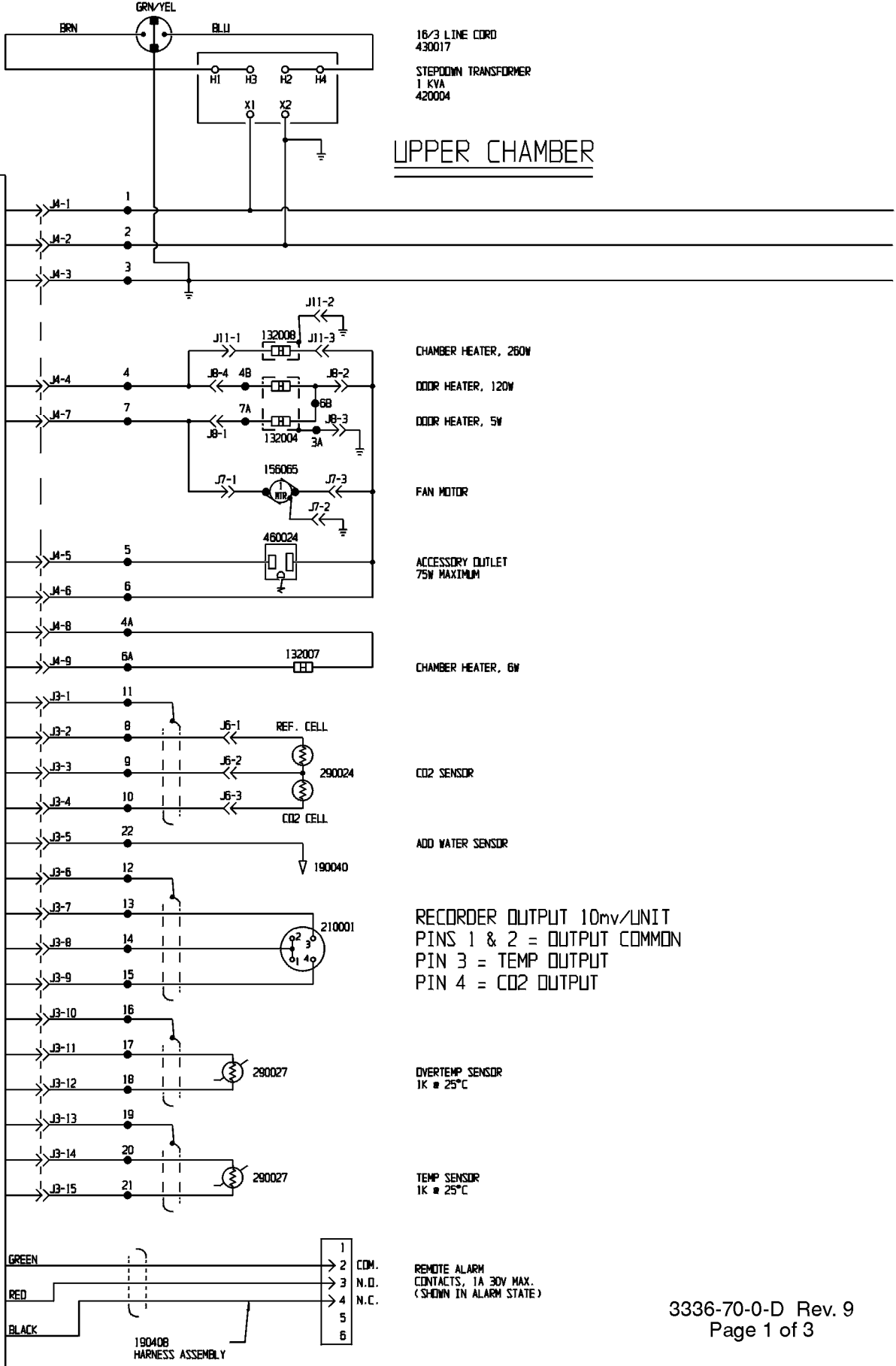
POWER CONNECTION
 180-260V, 1PH, 2W, 50/60HZ, 3.8FLA

16/3 LINE CORD
 430017
 STEPDOWN TRANSFORMER
 1 KVA
 420004

UPPER CHAMBER

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

CONTROL PANEL - SEE DWG. 140181-70-0-0
UPPER CHAMBER



CHAMBER HEATER, 260W

DOOR HEATER, 120W

DOOR HEATER, 5W

FAN MOTOR

ACCESSORY OUTLET
75W MAXIMUM

CHAMBER HEATER, 6W

CO2 SENSOR

ADD WATER SENSOR

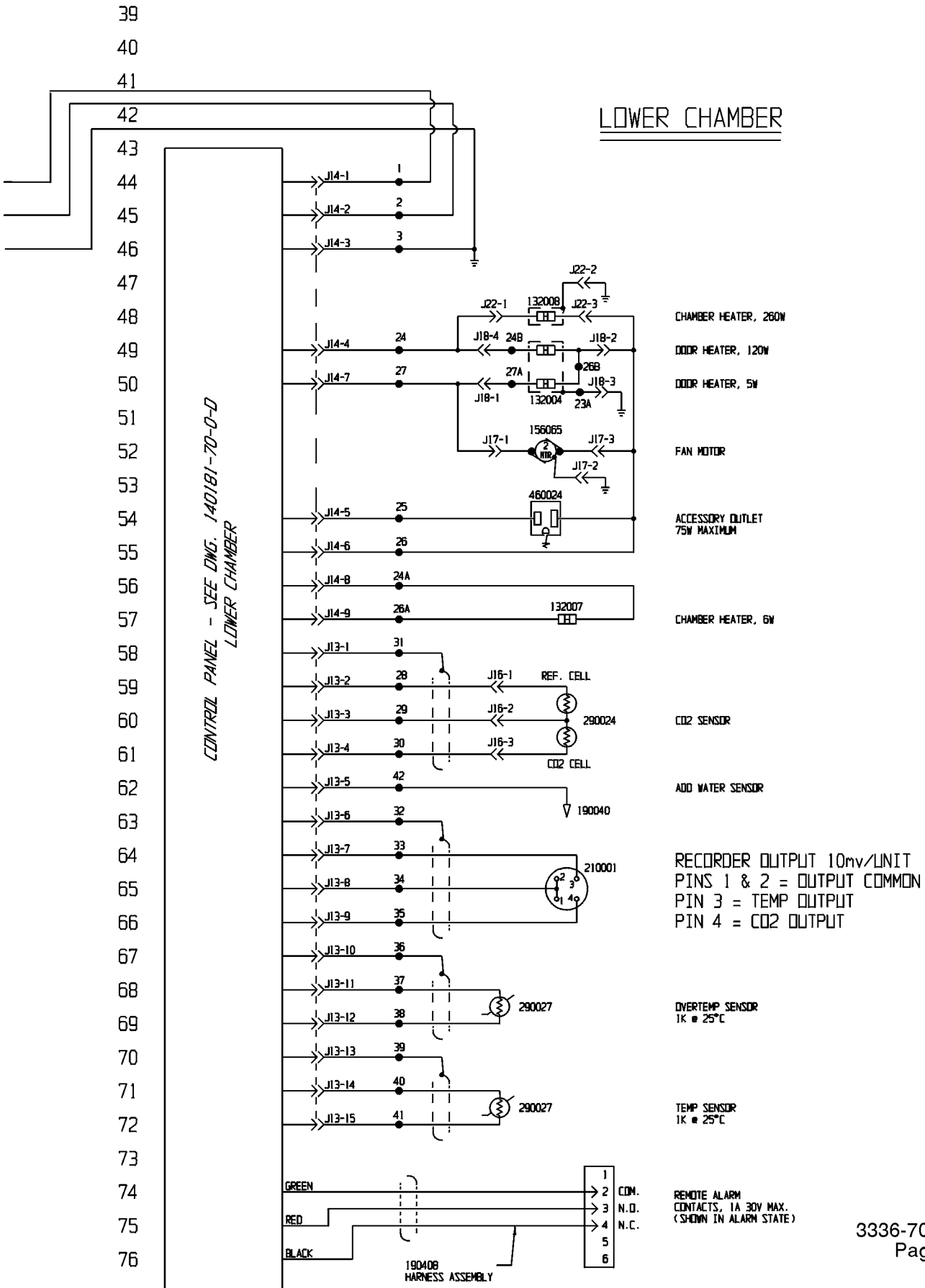
RECORDER OUTPUT 10mv/UNIT
 PINS 1 & 2 = OUTPUT COMMON
 PIN 3 = TEMP OUTPUT
 PIN 4 = CO2 OUTPUT

OVERTEMP SENSOR
 1K @ 25°C

TEMP SENSOR
 1K @ 25°C

REMOTE ALARM
 CONTACTS, 1A 30V MAX.
 (SHOWN IN ALARM STATE)

LOWER CHAMBER



CONTROL PANEL - SEE DWG. 140181-70-0-D
LOWER CHAMBER

CHAMBER HEATER, 260W

DOOR HEATER, 120W

DOOR HEATER, 5W

FAN MOTOR

ACCESSORY OUTLET
75W MAXIMUM

CHAMBER HEATER, 6W

CO2 SENSOR

ADD WATER SENSOR

RECORDER OUTPUT 10mv/UNIT
PINS 1 & 2 = OUTPUT COMMON
PIN 3 = TEMP OUTPUT
PIN 4 = CO2 OUTPUT

OVERTEMP SENSOR
1K @ 25°C


TEMP SENSOR
1K @ 25°C

REMOTE ALARM
CONTACTS, 1A 30V MAX.
(SHOWN IN ALARM STATE)

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

	U.C. WIRE #	GAUGE	COLOR	L.C. WIRE #	
	1	16	BLK	1	
	2	16	WHT	2	
	3	16	GRN	3	
	3A	18	GRN	23A	
	4	20	RED	24	
	4A	18	BLK	24A	
	4B	18	RED	24B	
	5	20	BLK	25	
	6	20	WHT	26	
	6A	18	BLK	26A	
	6B	18	WHT	26B	
	7	20	BLK	27	
	7A	20	BLK	27A	
	8	22	RED	28	}
22/3 CABLE	9	22	BLK	29	
	10	22	WHT or CLR	30	
	11	22	SHLD	31	22/3 CABLE
	12	22	SHLD	32	}
22/3 CABLE	13	22	RED	33	
	14	22	BLK	34	
	15	22	WHT or CLR	35	22/3 CABLE
	16	24	SHLD	36	}
24/2 CABLE	17	24	BLK	37	
	18	24	WHT or CLR	38	
	19	24	SHLD	39	}
24/2 CABLE	20	24	BLK	40	
	21	24	WHT or CLR	41	
	22	24	BRN	42	24/2 CABLE

CUSTOMER APPROVAL/REFERENCE	9 IN-1642 08-31-93 GLM/KDG	ADD REMOTE ALARM CONTACTS
APPROVED BY _____	8 IN-988 12-19-89 JAS/PDK LDN	REVISED FAN MOTOR
APPROVING FIRM _____	7 IN-552 12-11-84 DNL/DWL LDN	REVISED FAN MOTOR
DATE OF APPROVAL _____	6 SI-360 4-19-83 DNL/DWL LDN	SEE ECR
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	5 IN-434 12-28-82 DNL/DWL LDN	ADDED VOLTAGE RANGE
	REV ECR NO. DATE BY CAD/APPD	DESCRIPTION OF REVISION
 Forma Scientific <small>BOX 640 HOPKETTA, OHIO 45750 TELE 34-3284 TELL FREE USA 800-848-3000, OHIO 614-372-4780</small>	DATE 6-15-81 DWN DAG CAD/DWL APPD LDN SCALE N/A	
	CUSTOMER _____	
	JOB TITLE 3336 AUTOMATIC CO2 INCUBATOR	
	DWG TITLE ELECTRICAL SCHEMATIC	
	LOCATION INCUBATOR1	JOB NUMBER _____

NOTES:

Denotes Terminal Strip Connection
N/A Lost Relay Number
15 Lost Terminal Number
42 Lost Wire Number