

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

Telephone: (614) 373-4763
Telefax: (614) 373-4189

Models:
1184, 1185, 1186, 1194,
1195 and 1196

Biological Safety Cabinet
Class II, Type AB₃

Manual No. 7011186 Rev. 1

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 freezer 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 7011186				
REV	ECN	DATE	BY	DESCRIPTION
1		5-1-94	DG	Measuring blower motor voltage
0		6-1-93	DG	Minor format revision

General Safety Notes Used In This Manual

	<p>This symbol alerts the user to important operating and/or maintenance instructions. It may be used alone or with other safety symbols. Read the accompanying text carefully.</p>
	<p>Potential electrical hazards. Only qualified persons should perform the instructions and procedures associated with this symbol.</p>
	<p>Extreme temperature hazards, hot or cold. Instructions associated with this symbol should only be carried out when using special handling equipment or when wearing special, protective clothing.</p>
	<p>Potential biological hazards. Proper protective equipment and procedures must be used when following instructions associated with this symbol. Reference O.S.H.A. Regulation 1910-1030.</p>
	<p>Potentially hazardous energy. Equipment being maintained or serviced must be turned off and locked off to prevent possible injury. Reference O.S.H.A. Regulation 1910-147.</p>

Remember:

- √ 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.

TABLE OF CONTENTS

SECTION 1 - RECEIVING

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

SECTION 2 - INTRODUCTION

- 2.1 Description
- 2.2 Theory of Operation
- 2.3 Airflow Diagram (Fig. 2.0)

SECTION 3 - INSTALLATION

- 3.1 Location
- 3.2 Power Connection
- 3.3 Plumbing Connection
- 3.4 Exhaust Requirements

SECTION 4 - OPERATION

- 4.1 Control and Indicating Devices
 - a. Blower Switch
 - b. Light Switch
 - c. Ultra-Violet Light (Optional)
 - d. Alarm By-Pass Switch (sliding window models only)
 - e. Static Pressure Gauge (in. w.g.)
 - f. Blower Speed Control
 - 1. Measuring Blower Motor Voltage
 - g. Blower Motor/Lights Reset Button (15 Amp)
 - h. Receptacle Reset Button (15 Amp)
 - i. Duplex Receptacles

SECTION 4 - OPERATION (continued)

- j. Drain Valve
 - k. Service Valves
 - l. Exhaust Filter Guard
 - m. Hinged Window Assembly
 - n. Sliding Window Assembly
- 4.2 Cabinet Drawing 4.0
 - 4.3 Control Panel Drawing 4.1 (Models 1184/1186)
 - 4.4 Control Panel Drawing 4.2 (Models 1194/1196)

SECTION 5 - GENERAL CAUTIONS

- 5.1 General Caution Notes
- 5.2 Important Label Location Drawing

SECTION 6 - CABINET START-UP

- 6.1 General Recommendations
- 6.2 Use of Auxiliary Equipment in the Cabinet
- 6.3 Cabinet Check
- 6.4 Start-Up Procedure

SECTION 7 - TROUBLESHOOTING

- 7.1 Troubleshooting Guide

SECTION 8 - MAINTENANCE

- 8.1 Checking the Static Pressure Gauge "Zero"
- 8.2 Re-zeroing the Static Pressure Gauge
- 8.3 Adjusting the Damper
- 8.4 Biological Safety Cabinet Test Grids Models 1184/1186
- 8.5 Biological Safety Cabinet Test Grids Models 1186/1196

SECTION 9 - SPECIFICATIONS

- 9.1 Model 1184 (4' Cabinet with sliding window)
- 9.2 Model 1186 (6' Cabinet with sliding window)
- 9.3 Model 1194 (4' Cabinet with hinged window)
- 9.4 Model 1196 (6' Cabinet with hinged window)

SECTION 10 - ACCESSORIES

- 10.1 Accessories (Chart)

SECTION 11 - PARTS LIST

- 11.1 Model 1184
- 11.2 Model 1186
- 11.3 Model 1194
- 11.4 Model 1196

SUPPLEMENTS AND WARRANTY INFORMATION

SECTION 1 - RECEIVING

1.1 Preliminary Inspection

This safety cabinet was thoroughly inspected and carefully packed prior to shipment and all necessary precautions were taken to ensure its safe arrival. 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. Call the delivering carrier and request that their representative perform an inspection. Do not discard any of the packing material or 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.

1.4 Unpacking List

Included with the Installation/Operation manual are four index buttons. These buttons may be used to identify the type of service supplied to the service valves. Also included in a separate bag, is a small allen wrench (used for calibrating the Static Pressure Gauge). This allen wrench should be kept with the Manual at all times.

SECTION 2 - INTRODUCTION

2.1 Description

The 1184/1186/1194/1196 Series is a Class II, Type A/B₃ cabinet. The "Type A/B₃" designation indicates two alternative uses of the cabinet. When vented directly into the laboratory room, the unit serves a "Type A" unit. When vented to the outside atmosphere, through an in-house exhaust system, it serves as a "Type B₃" unit. Either usage of the cabinet offers both personnel and product protection.

The cabinet can be used with low-to-moderate risk hazard to the user and/or the experiment. Class 1, 2, and 3 (low-to-moderate risk) agents are described in the "Biosafety In Microbiological And Biomedical Laboratories"; CDC NIH Publication No. (NIH) 88-8395, 2nd Edition, May 1988.

The 1184/1186/1194/1196 cabinets are designed to meet the requirements of the National Sanitation Foundation Standard #49.

The cabinet is available in a sliding or hinged window version. The cabinet's window permits the user to place auxiliary equipment and research implements in the work area. The work opening must be held to 10" during all work procedures. If the window is raised higher than the designated 10", the air barrier at the front of the cabinet will be weakened and containment will be seriously impaired.

2.2 Theory of Operation

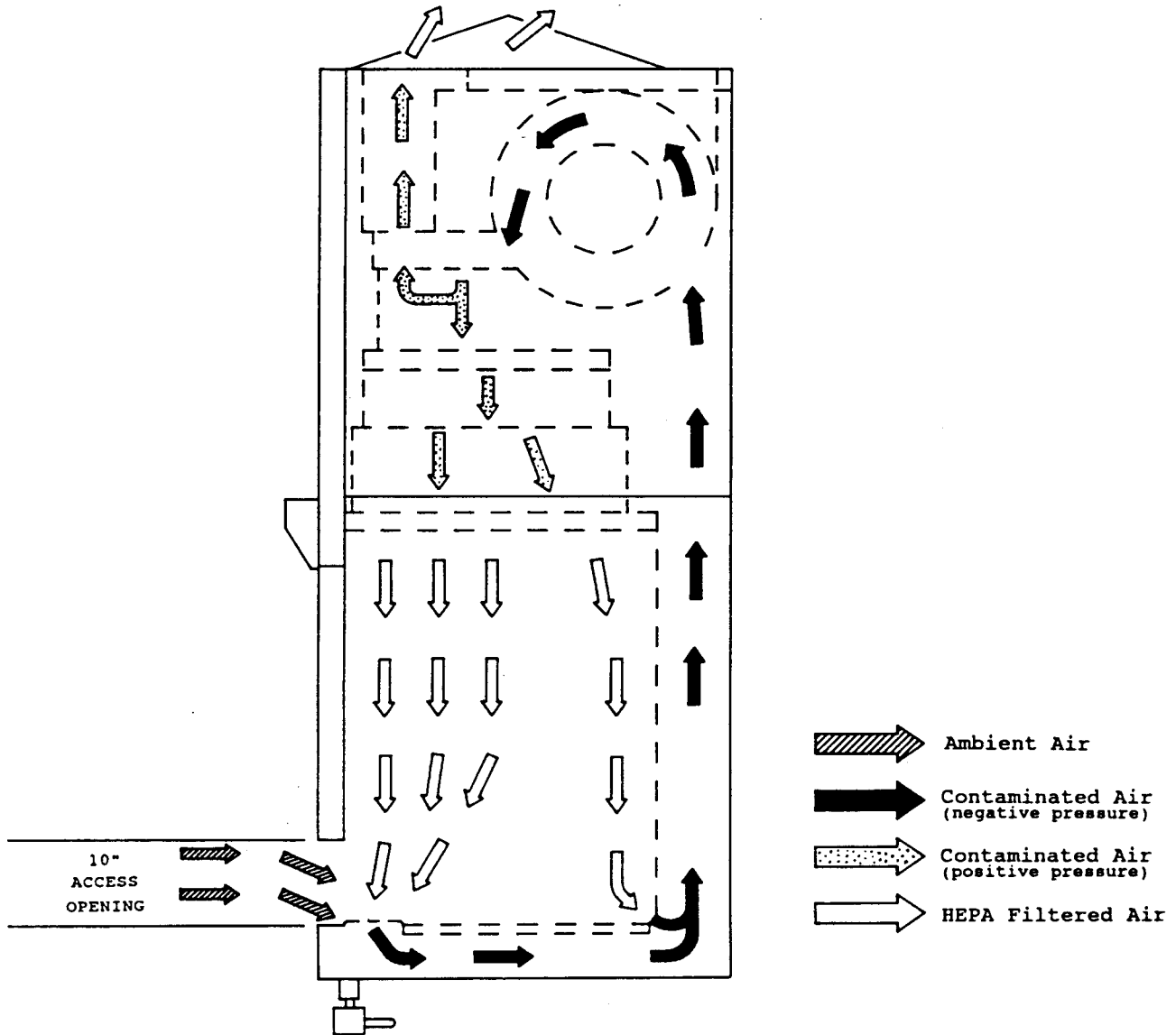
Figure 2.1 shows the general airflow pattern of the cabinet.

Clean, filtered air descends through the work zone with approximately 40% being discharged through the exhaust HEPA filter with the remaining air recirculating through the supply HEPA filter into the work area. Exhausted air is replaced by room air entering the system through the front access opening.

Room air entering the work zone, through the front access opening, completes the air barrier at the unit face and is responsible for the containment properties of the unit. All work must be performed beyond the intake grille, on the solid work tray.

2.3 Airflow Diagram

Figure 2-1



SECTION 3 - INSTALLATION

3.1 Location

Locate the cabinet on a firm level surface in an area of minimum temperature changes. The cabinet should be placed away from disruptive air currents caused by excessive personnel traffic, air-conditioning or heating ductwork, or laboratory windows and doors. Proper cabinet location is important, as drafts disrupt critical air flow characteristics and allow room contaminants to enter or escape the cabinet work area.

Where space permits, fourteen inches should be allowed on each side of the cabinet for maintenance. A twelve inch height should be available from the top of the cabinet to the ceiling.

3.2 Power Connection

The electrical wall outlet(s) leading to the cabinet should be accessible for electrical testing.

This cabinet is equipped with two power cords, one for the internal blower motor and lights, the other for the duplex utility outlets. The cords should be plugged into separate circuits, so if the duplex outlets overload, the cabinet itself will not shut down.

Refer to Section 9 or to the data plate on the front of the unit for electrical specifications.

3.3 Plumbing Connection

Two service valves are standard with each cabinet and are located on the right and left side of the work station. All service valves are piped within the cabinet. External connection is to 3/8" FPT coupling. Identification index buttons are supplied.

The cabinet will accommodate six service valves. The additional four service valves may be purchased from Forma Scientific, Inc.

Caution! Explosive/flammable substances should never be used in the cabinet, unless approved and monitored by a biological safety officer or other qualified individual.

3.4 Exhaust Requirements

Filtered air from the cabinet may be exhausted directly into the room or vented to the outside through an external exhaust system. Consult a biological safety officer or other qualified individual for cabinet-type exhaust requirements.

a. Direct Room Exhaust

1. Locate the exhaust filter guard on the top of the cabinet.
2. Remove and save the four acorn-type nuts and washers that secure the exhaust filter guard.
3. Remove the cardboard cover plate from the top of the cabinet.
4. Replace the exhaust filter guard with the slope toward the front and the back of the unit. Secure it with the four acorn-type nuts and washers removed in Step #2.

b. External Exhaust System (For Class II Type B₃ Only)

The intermediate connection of the cabinet to any external exhaust system (in-house) can be done by using an exhaust transition which is available from Forma Scientific, Inc. The exhaust transition requires a minimum of 16-1/2" headroom from the top of the cabinet.

Caution: The exhaust system should have safeguards against exhaust failure. It is required that a biological safety officer or industrial hygienist or other qualified individual review the agents and chemicals used inside the cabinet to determine if additional filtration treatment is necessary before venting to the atmosphere.

The exhaust system must be capable of handling the volume of air passing through the cabinet.

Exhaust requirements are approximately:

Model 1186: 505-556 CFM
Model 1194: 339-369 CFM
Model 1184: 339-369 CFM
Model 1196: 510-561 CFM

SECTION 4 - OPERATION

4.1 Control and Indicating Devices (Refer To Figures 4-0, 4-1 and 4-2)

Before operating the cabinet it is recommended that the user(s) become familiar with the following items on the cabinet.

a. Blower Switch

The blower switch controls the on/off power to the internal blower.

b. Light Switch

The "dual purpose" light switch controls power to the fluorescent lamp or the optional ultra-violet lamp (both lamps are located in the work area).

Note: with the ultra-violet lamp option, the switch provides the following settings; "off"= center, fluorescent "on" top side, and ultra-violet "on" bottom side. only one lamp may be lit at one time.

c. Ultra-Violet Light (Optional)

Caution: eyes or skin should not be exposed to ultra-violet light. recommended usage is only when lab is not in use.

Cabinets may be equipped with an ultra-violet germicidal light as optional equipment. The "dual purpose" light switch labeled "Lights" controls power to the optional ultra-violet lamp.

Note: with the ultra-violet lamp option, the switch provides the following settings; "off" = center, fluorescent "on" top side, and ultra-violet "on" bottom side. only one lamp may be lit at one time.

d. Alarm By-Pass Switch (Sliding Window Models Only)

The Alarm By-Pass switch permits the operator to disable/silence the audible Window Above 10 Inch alarm for approximately five minutes. The "RED" visual indicator will remain illuminated. The alarm will "ring-back" to remind the operator that the window is still open more than 10 inches.

e. Static Pressure Gauge (In. W.G.)

The static pressure gauge, located on the control panel, measures the air pressure differential across the filters providing an indication of filter "loading". As the filters become loaded, resistance increases and the reading on the static pressure gauge increases accordingly. When the reading increases by 50% (from original measurement), cabinet airflow should be checked with a thermoanemometer, by a qualified service technician. The filters must be replaced if proper airflow cannot be obtained.

Note: The static pressure gauge should not be used as a direct measure of air flow.

f. Blower Speed Control

The blower speed control, located on the top left back side of the control panel, is used to adjust the air velocity from the internal blower motor. The blower speed is adjusted by turning the screw on the variable resistor mounted on the circuit board adjacent to the controller. (Refer to the drawing following.) A clockwise turn of the screw adjustment will increase air velocity and a counter-clockwise turn will decrease it.

The blower speed is factory set and should only be changed by a qualified technician.

Warning:

Live voltage is present on the control terminals of the switches and dials on the front of the blower panel. Use extreme care to avoid touching these controls when reaching into the drawer and making any adjustments.

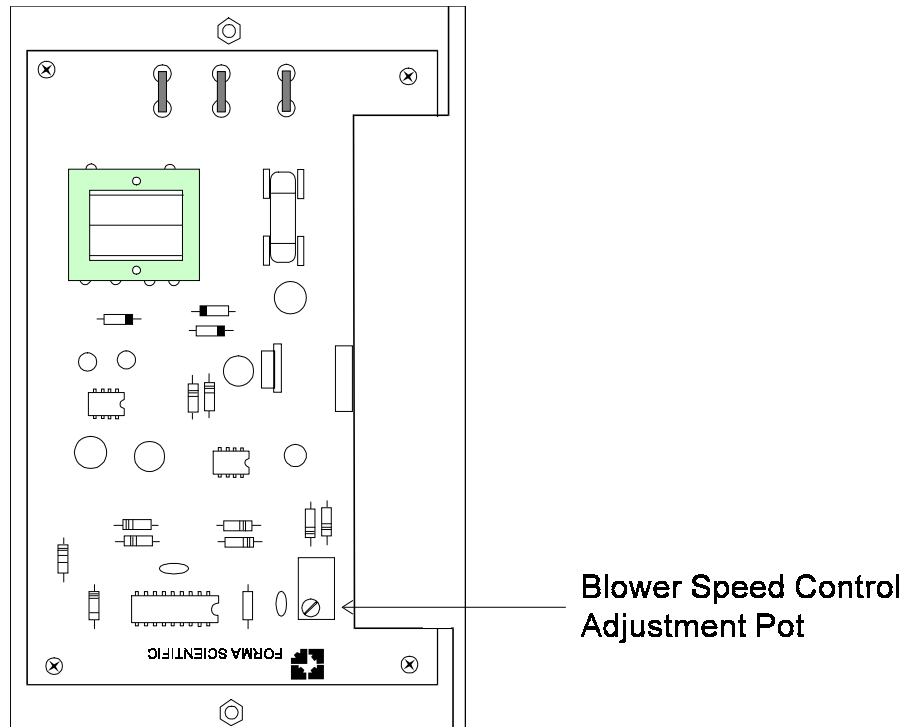


Figure 4-1
Component locations on the printed circuit control board

1. Measuring Blower Motor Voltage

Both blower motor voltage and line voltage are measured at the three terminal connectors at the top of the circuit board. Refer to Figure 4-2.

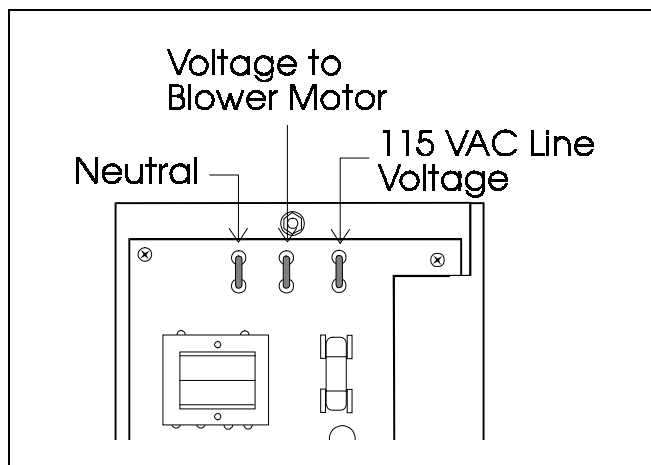


Figure 4-2
Circuit board voltage terminals

g. Blower Motor/Lights Reset Button (15 Amp)

The reset button (located on the left side of the control panel, directly below the Receptacle Reset Button) is an inline circuit breaker for the internal blower motor and lighting. If an overload condition occurs, the circuit breaker will trip, and the button will protrude from the panel. Depressing the button will reset the circuit breaker.

Note: Should an overload condition occur:

1. Turn power off to the blower.
2. Turn power off to the lighting.
3. Press the blower motor/lighting reset button.

h. Receptacle Reset Button (15 Amp)

The receptacle reset button (located directly above the Blower Motor/Lights Reset Button) is an inline circuit breaker for the receptacles only. If an overload condition occurs the circuit breaker will trip, and the button will protrude from the panel. Depressing the button will reset the circuit breaker.

Note: should an overload condition occur:

1. turn power off to appliance/apparatus.
2. unplug appliance/apparatus from receptacle.
3. turn receptacle power switch to off position.
4. depress the receptacle reset button.

i. Duplex Receptacles

A duplex receptacle (115 volts, 15 amps) is located on the left and right side wall of the work station. Power is controlled to the duplex receptacles by the receptacle switch located on the control panel.

j. Drain Valve

The drain valve, located on the right front side of the cabinet, is provided for the safe drainage of the drain pan. This valve should remain closed while work is being performed in the cabinet and should be used only in the event of a major spill.

Caution:

If an accidental spill occurs, immediately consult a biological safety officer or other qualified individual for proper procedures. To ensure the proper containment of a spill, connect a sealed hose from the drain valve to a sealed container.

k. Service Valves

Two service valves are standard with each cabinet. These valves are located on the right and left side of the work station and can be coded with the type of service that they supply. Identification index buttons are supplied.

The cabinet will support a total of six service valves. Additional service valves may be purchased from Forma Scientific Inc.

l. Exhaust Filter Guard

The exhaust filter guard, located on top of the exhaust filter, protects the exhaust air flow and also prevents the storing of objects on top the housing.

m. Hinged Window Assembly

The hinged window assembly allows the operator to raise the glass window to place items within the work area.

Caution

When work is being performed in the cabinet, the hinged window must be closed to avoid contamination to product and personnel.

n. Sliding Window Assembly

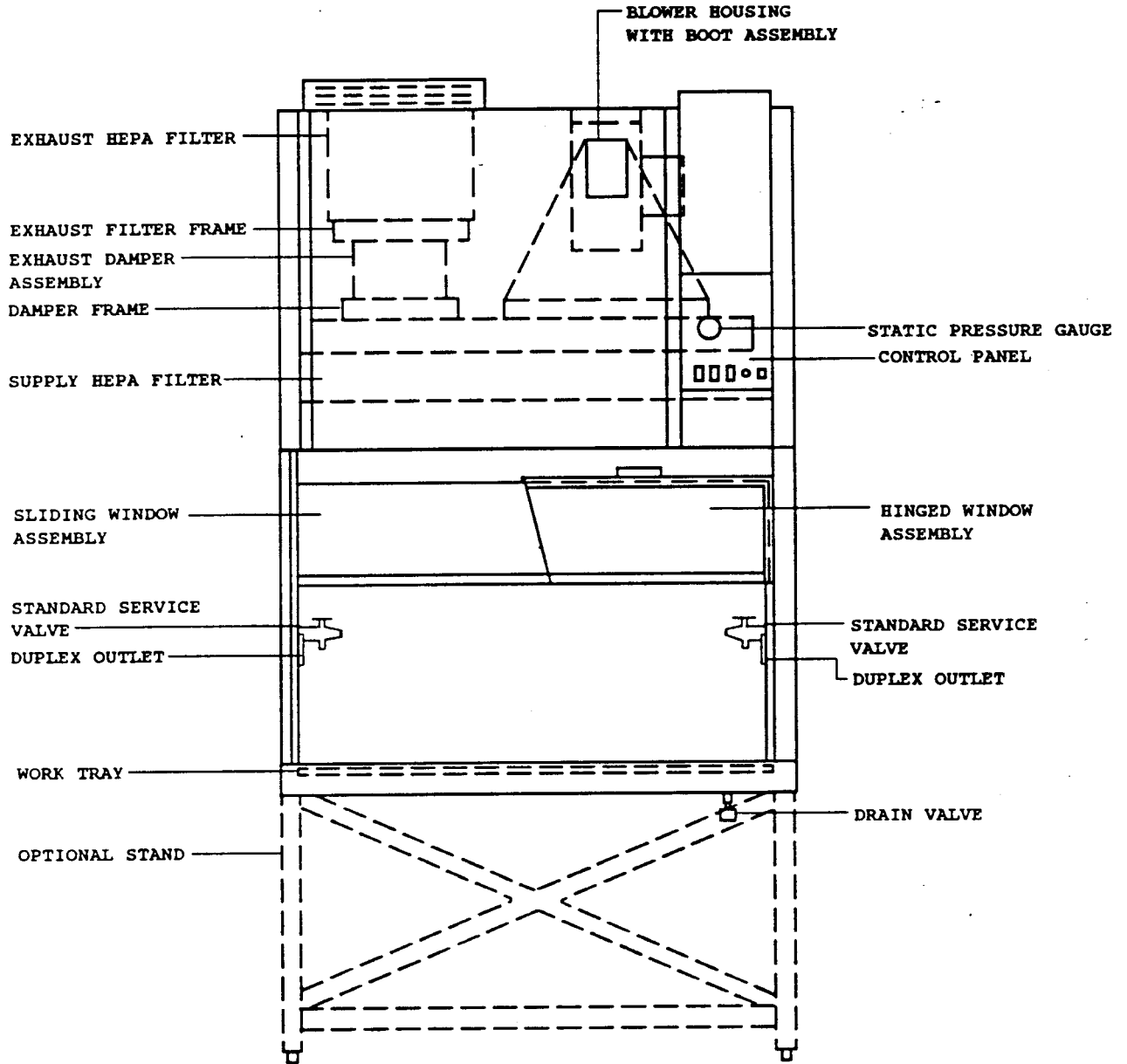
The sliding window assembly allows the operator to raise the glass window to place items within the work area.

Caution

When work is being performed in the cabinet, the sliding window must be at the 10" position to avoid contamination to product and personnel.

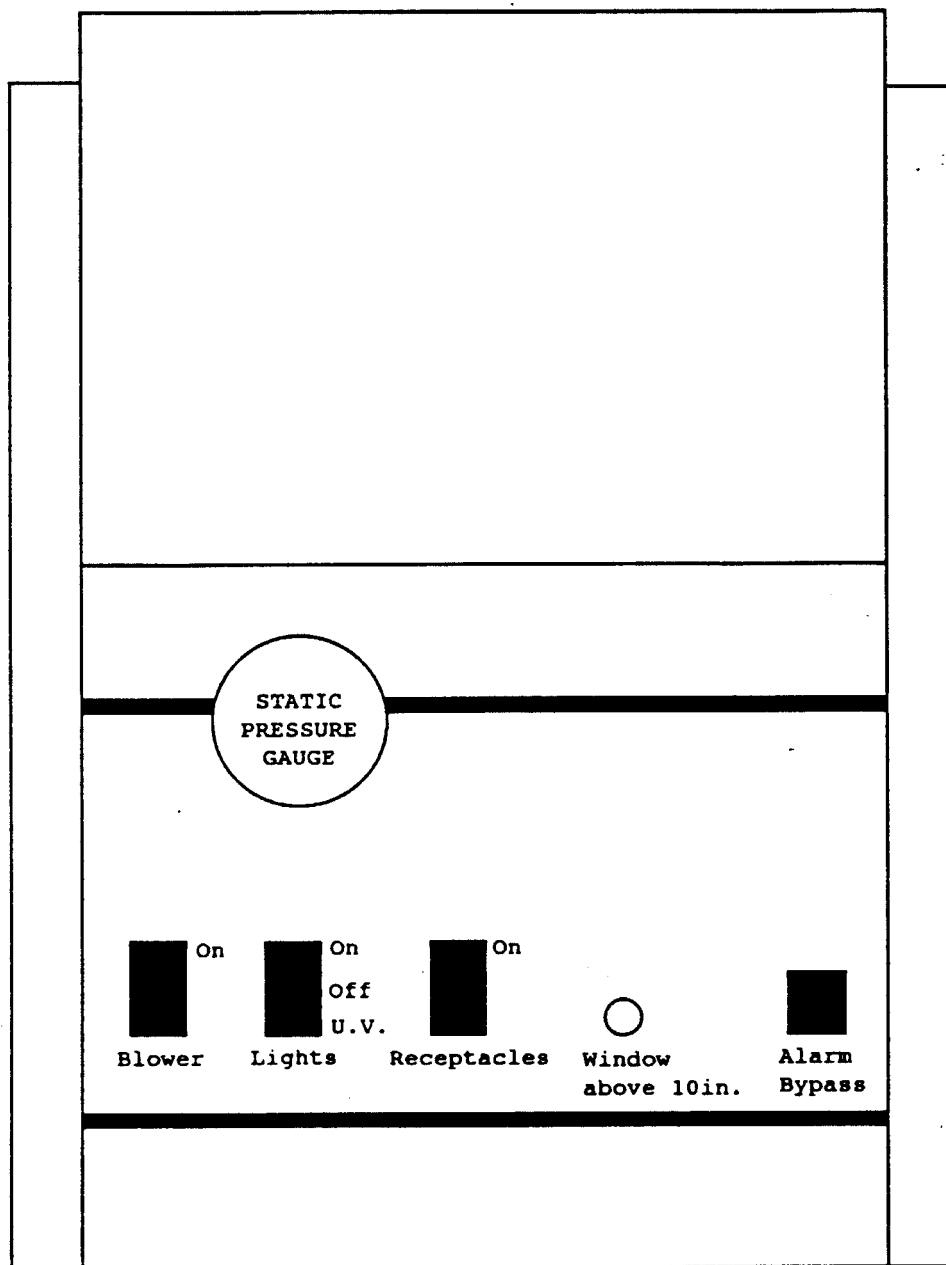
Note: If the sliding window is above the 10" level, a red indicating light and audible buzzer will warn that an unsafe condition exists.

SECTION 4.2



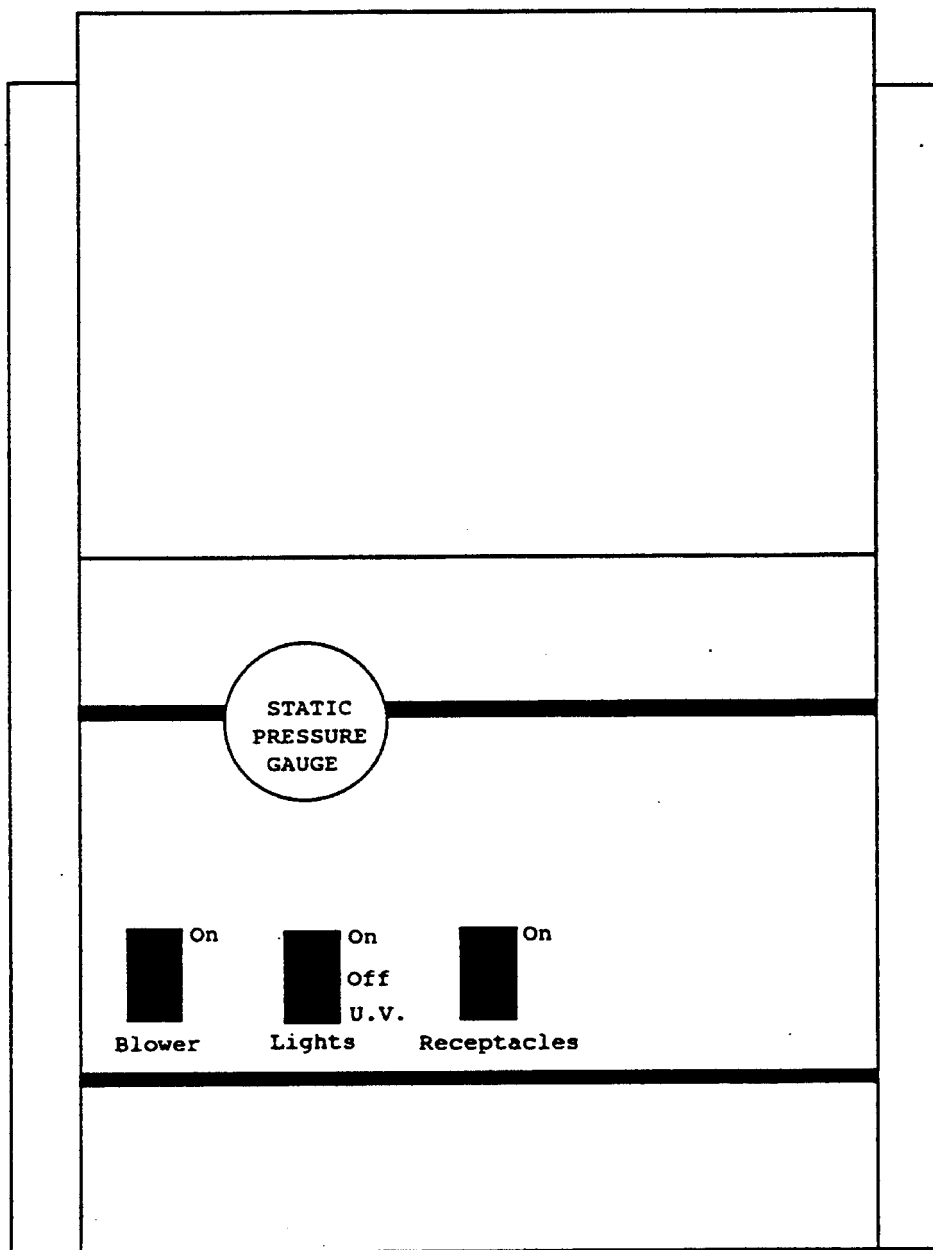
DRAWING #4.0
OVERVIEW OF CABINET

SECTION 4.3



DRAWING #4.1
OVERVIEW OF 1184/1186
CONTROL PANEL

SECTION 4.4



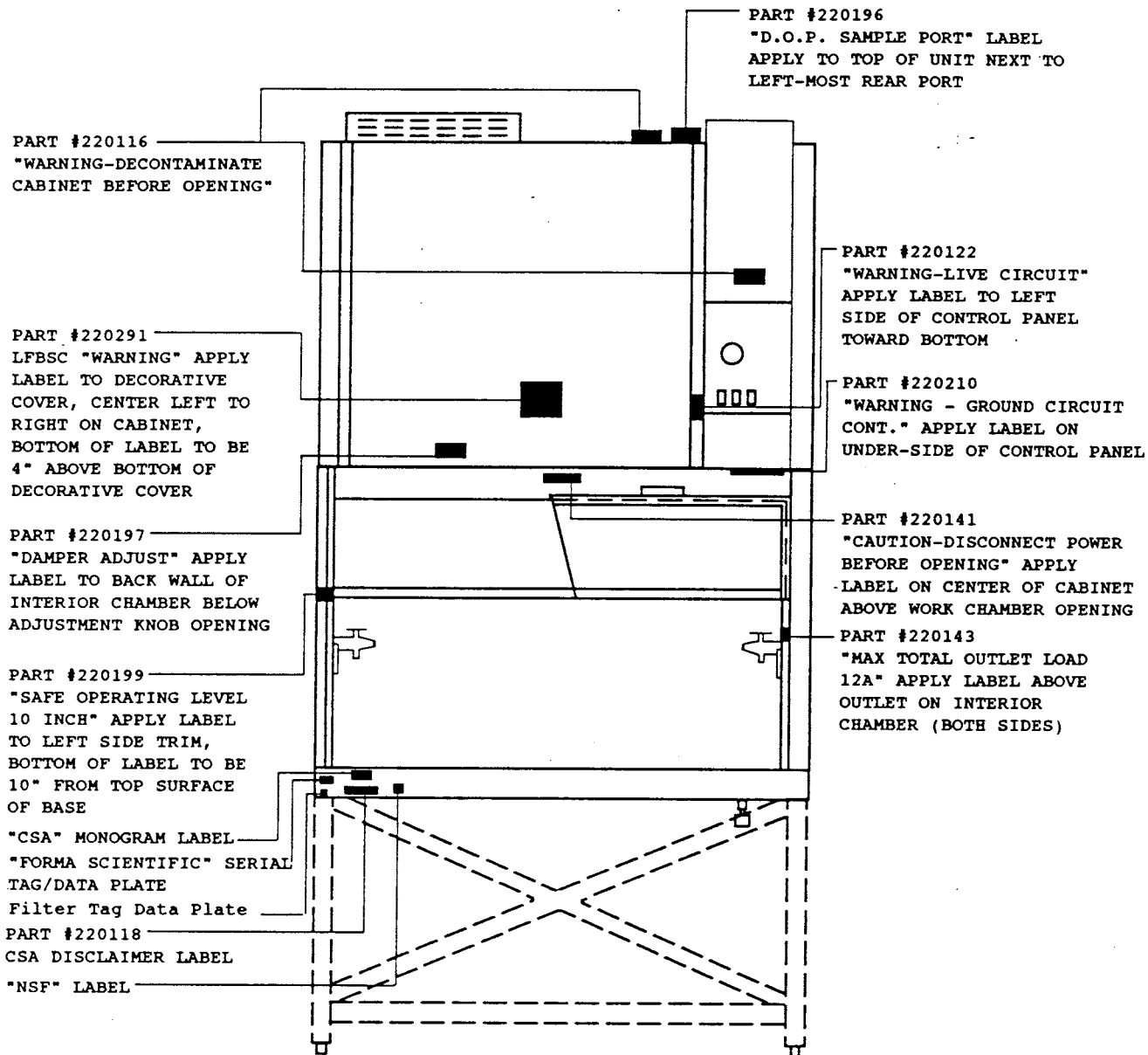
**DRAWING #4.2
OVERVIEW OF 1194/1196
CONTROL PANEL**

SECTION 5 - GENERAL CAUTIONS

5.1 Caution Notes

1. Following initial installation, the unit must be thoroughly tested and certified.
2. All activities to be performed within the cabinet should be approved by a biological safety officer or other qualified individual.
3. Since the HEPA filters remove particulates only (not gas), explosive/flammable substances should never be used in the cabinet, unless approved and monitored by a biological safety officer or other qualified individual.
4. Ultra-violet lighting should not be used while personnel are using the cabinet. If exposure cannot be avoided the proper safety gear/clothing must be worn. Consult a biological safety officer or other qualified individual for proper procedures.
5. If the cabinet is to be used for biological or toxicological applications, it must be monitored by a biological safety officer or other qualified individual.
6. If the unit needs to be serviced, it must be decontaminated to protect service personnel from contamination. The cabinet must then be recertified by a qualified certifying agency.
7. None of the perforations in the work area may be covered or blocked, as air flow will be disrupted and contamination may occur.

5.2 IMPORTANT LABEL LOCATION DRAWING



NOTE:

IF ANY OF THE ABOVE "LABELS" ARE LOST OR DAMAGED IN ANYWAY, PLEASE CONSULT THE
FORMA SCIENTIFIC SERVICE DEPARTMENT FOR A FREE REPLACEMENT!

PHONE 1-800-848-3080

FAX 614-373-4189

TELEX 29-8205

SECTION 6 - CABINET START-UP

6.1 General Recommendations

1. Keep the movement in the room to a minimum when the cabinet is in use.
2. Keep all laboratory doors closed to prevent drafts that may disturb critical air flow.
3. Pre-plan cabinet use and place everything needed in the cabinet so that nothing passes through the air barrier (in or out) during the procedure.
4. Practice good aseptic technique to insure safe use of the cabinet.
5. If a spill occurs, clean it up immediately. Decontaminate the work area and all affected equipment.
6. Do not cover or block the exhaust grille.
7. Do not cover or block any perforations (air holes) in the work area.

6.2 Use of Auxiliary Equipment in the Cabinet

Use auxiliary equipment in the cabinet only if proper precautions are taken. Any appliance used in the work area will cause turbulence and disturb the air flow. Use of such equipment should be carefully managed. The equipment should be placed at the rear of the work space where air turbulence will have a minimal effect.

Blender: A blender may be used in the cabinet. But because of the amount of aerosol it produces and the increased turbulence it causes, it is recommended that it be removed from the cabinet as soon as possible.

6.3 Cabinet Checklist

1. Check to see that the Drain Valve is closed (the handle horizontal), so that if a spill should occur, it will remain in the system and not drain onto the laboratory floor.
2. Check to see that All Service Valves are closed.

6.4 Start-Up Procedure

1. Turn the light switch on.
2. Check the intake and exhaust grills to ensure they are not blocked.
3. Turn the blower switch on.
4. Place everything needed into the cabinet.
5. Place viewing window at 10" (Sliding Type).
6. Close the viewing window (Hinged Type).

SECTION 7 - TROUBLESHOOTING

7.1 Troubleshooting Guide

The following is a guide to troubleshooting the system. If a contaminated area of the cabinet must be entered to determine and/or resolve the source of a particular problem, *the cabinet must first be decontaminated.*

Servicing of the unit must be performed by qualified service personnel.

Problem 1: Air flow in the cabinet work area and through the exhaust filter is inadequate.

Possible causes:

- a. Exhaust filter is blocked by laboratory materials or the protective shipping cover.
- b. If the biological safety cabinet is connected to an exhaust system, there may be inadequate exhaust suction or back pressure in the duct system. The system must be rebalanced to handle the correct air volume. A Biological Safety Officer should be consulted.
- c. Low voltage being applied to the blower motor.
- d. Blower motor or speed control is defective.
- e. Supply HEPA filter and Exhaust HEPA filter may be loaded. Decontaminate the unit and replace both HEPA filters.

Caution: Before any maintenance work is performed in the biological safety cabinet, the unit must first be decontaminated.

Problem 2: Ultra violet light malfunction

Possible causes:

- a. Check lamp pins and socket ends for contact.
- b. Defective starter for the U/V light.

Problem 3: Fluorescent light malfunction

Possible causes:

- a. Check lamp pins and socket ends for contact.
- b. Defective lamp.

Problem 4: Loud screeching noise

Possible causes:

- a. Bad bearings in motor blower unit.
- b. Blower scroll is rubbing against housing.

SECTION 8 - ROUTINE MAINTENANCE

8.1 Checking the Static Pressure Gauge "Zero"

Note: In order to provide an accurate reading, the indicating needle of the static pressure gauge should be precisely at zero when the cabinet is shut off. If the cabinet is connected to a central exhaust system, the exhaust system must also be shut off.

Following HEPA filter replacement, the static pressure gauge should be checked for zero when the cabinet is shut off. When the cabinet is started up and proper air flow balance has been reached, the reading on the gauge should be recorded. This initial reading will serve as a base line indication of subsequent filter loading. When the reading increases by approximately 50% , the air flow balance should again be checked. Replacement of the filters will probably be required. Refer to Section 8.2, Re-zeroing the Static Pressure Gauge.

8.2 Re-Zeroing the Static Pressure Gauge

1. The cabinet must be turned off.
2. Remove the front cover from the static pressure gauge by grasping the front cover and turning it counter-clockwise.
3. Locate the allen-type adjustment screw beside the gauge needle.
4. Turn the adjustment screw counter-clockwise to lower the reading or clockwise to raise it.

8.3 Adjusting the Damper

Since the HEPA filter resistance may vary considerably from filter to filter (even filters of the same size), a damper has been installed in the cabinet exhaust system for maintaining proper airflow balance. The purpose of the damper is to regulate the amount of exhaust air, intake velocity and supply velocity. The damper has been preset at the factory and should not be readjusted unless the proper velocities cannot be obtained.

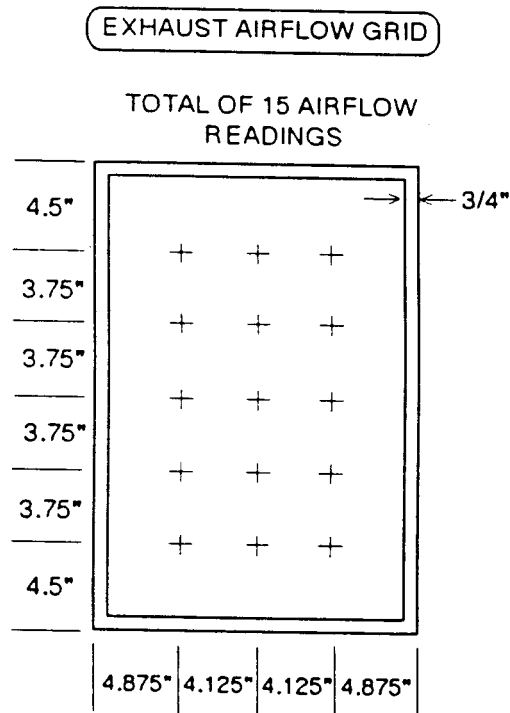
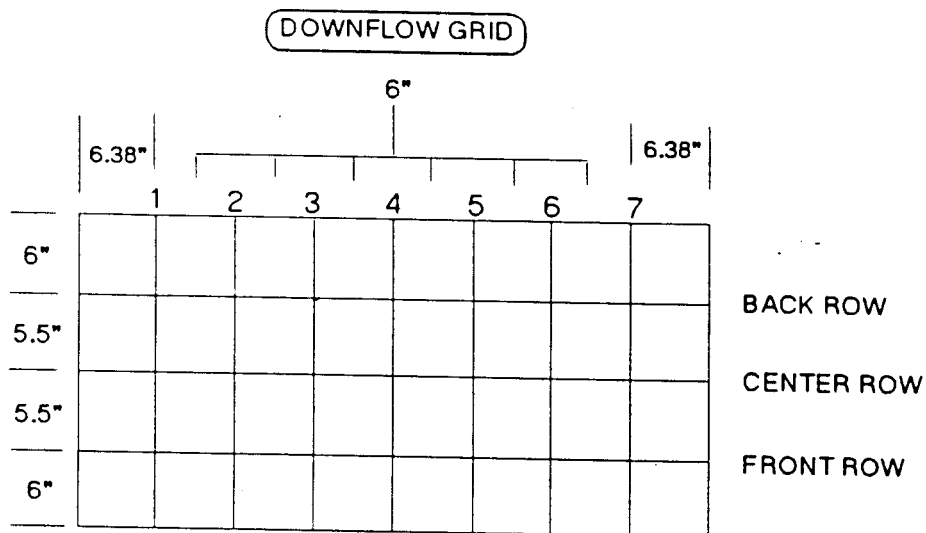
Caution! Adjustments must be made only by qualified personnel!

1. Layout test grids (refer to Sections 8.4 and 8.5).
2. Start-up the cabinet and allow it to run for at least twenty minutes.
3. Take airflow measurements. If airflow specifications are not sufficient, open the control panel and check the voltage on the power switch. **Note:** airflow measurements at the factory are recorded with the cabinet connected to a 115 volt power supply.
4. Locate the triac board in the upper left hand corner (behind control panel). Measure (using a true RMS Voltmeter) and record the voltage drop across the white and black wire leading to the terminal strip. The voltage pot is located on the lower right side and is labeled "Voltage Pot". Clockwise adjustment of this pot increases voltage supply to the blower motor. Adjust 2-3 volts accordingly (either up or down, depending upon the airflow required). Retake the airflow measurements.

If it is determined that the damper must be adjusted in order for the proper airflow balance to be maintained, adjust it as follows.

1. The damper adjustment is located above the supply diffuser screen (toward the left back wall) and is labeled "Damper Adjust Access".
2. The sprocket adjustment wheel is numbered from 0 to 9. 0 = is fully closed, cw, and 9 = is fully open, ccw.
3. Note the position (from 0-9) on the damper adjust sprocket. With a screwdriver or similiar tool, turn the adjustment wheel clockwise to fully close the damper. Damper adjustments must start from the fully closed position.
4. Example: If the present position is 4 and the supply velocity is High and the exhaust velocity is Low, adjust the damper sprocket fully closed (0) and then to approximately 5. Re-take airflow readings and repeat until the correct airflow balance is achieved.

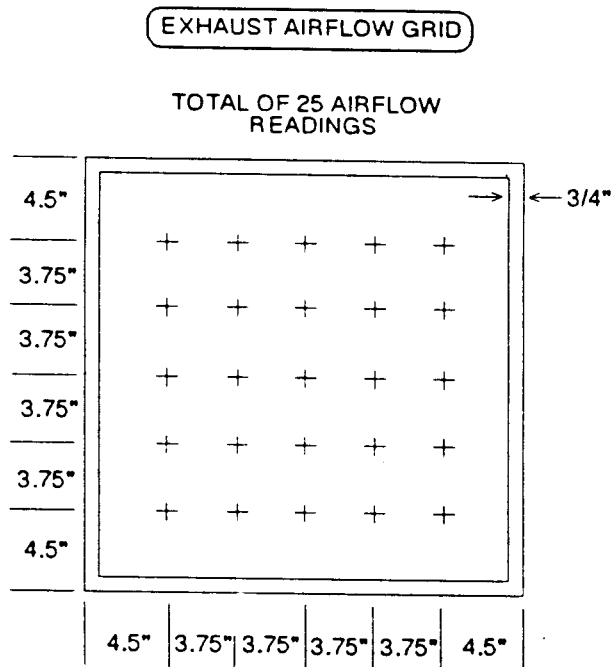
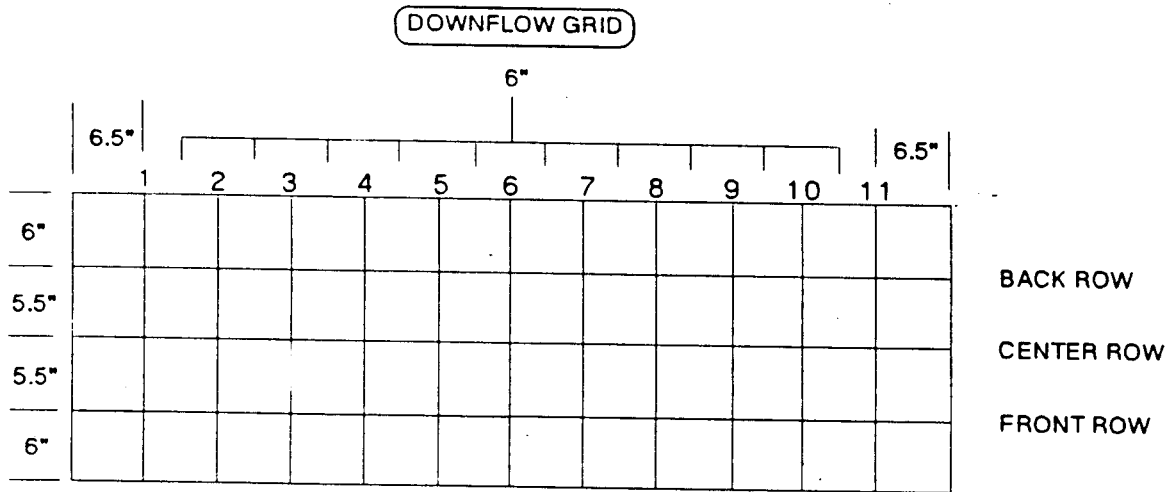
8.4 Biological Safety Cabinet Test Grids (Models 1184/1186)



WORK ACCESS OPENING FACE AREA= 3.385 (ft.sq.)
 OPEN EXHAUST AREA= 2.428 (ft.sq.)

• Total area minus obstructed area

8.5 Biological Safety Cabinet Test Grids (Models 1186/1196)



WORK ACCESS OPENING FACE AREA= 5.052 (ft.sq.)

OPEN EXHAUST AREA= 3.398 (ft.sq.)

* Total area minus obstructed area

SECTION 9 - SPECIFICATIONS

9.1 Model - 1184 (4' Cabinet With Sliding Window)

a. Construction

Work Surface: Type 304 Stainless Steel, #4 Finish
Cabinet: Cold Rolled Steel
Finish: Antique White with Windsor Blue Trim
Baked-on Powder Coat Epoxy Paint

b. Dimensions

Exterior: 54"W x 64"H x 32.5"F-B
Interior: 49"W x 28.33"H x 22.25"F-B

c. Electrical Requirements

Main: 100-125 VAC, 1 Phase, 2 Wire, 60 Hz, 10 FLA
Receptacle: 115 VAC, 1 Phase, 2 Wire, 60 Hz, 12 FLA
Circuit Breaker: 15 Amp

d. Filters

- (1) Supply HEPA Filter (18"W x 48"F-B x 5-7/8"H)
- (1) Exhaust HEPA Filter (18"W x 24"F-B x 11.5"H)

e. Lights

- (1) Fluorescent 60W, (F48T12/CW/HO)
- (1) Optional UV 30W, (G30T8) Germicidal Lamp

f. Blower Motor

3/4 HP, 1625 RPM

g. Drain Pan Capacity

21.14 Gallons

9.2 Model - 1185 (4' Cabinet With Sliding Window)

a. Construction

Work Surface: Type 304 Stainless Steel, #4 Finish
Cabinet: Cold Rolled Steel
Finish: Antique White with Windsor Blue Trim
Baked-on Powder Coat Epoxy Paint

b. Dimensions

Exterior: 54"W x 64"H x 32.5"F-B
Interior: 49"W x 28.33"H x 22.25"F-B

c. Electrical Requirements

Main: 220 VAC, 1 Phase, 2 Wire, 50 Hz, 6 FLA
Receptacle: 220 VAC, 1 Phase, 2 Wire, 50 Hz, 12 FLA
Circuit Breaker: 15 Amp

d. Filters

(2) Supply HEPA Filter (18"W x 48"F-B x 5-7/8"H)
(1) Exhaust HEPA Filter (18"W x 24"F-B x 11.5"H)

e. Lights

(2) Fluorescent 60W, (F48T12/VHO)
(1) Optional UV 30W, (G30T8) Germicidal Lamp

f. Blower Motor

3/4 HP, 1625 RPM

g. Drain Pan Capacity

21.14 Gallons

9.3 Model - 1186 (6' Cabinet With Sliding Window)

a. Construction

Work Surface: Type 304 Stainless Steel, #4 Finish

Cabinet: Cold Rolled Steel

Finish: Antique White with Windsor Blue Trim

Baked-on Powder Coat Epoxy Paint

b. Dimensions

Exterior: 78"W x 64"H x 32.5"F-B

Interior: 73"W x 28.33"H x 22.25"F-B

c. Electrical Requirements

Main: 100-125 VAC, 1 Phase, 2 Wire, 60 Hz, 12 FLA

Receptacle: 115 VAC, 1 Phase, 2 Wire, 60 Hz, 12 FLA

Circuit Breaker: 15 Amp

d. Filters

(1) Supply HEPA Filter (18"W x 72"F-B x 5-7/8"H)

(1) Exhaust HEPA Filter (24"W x 24"F-B x 11.5"H)

e. Lights

(1) Fluorescent 85W, (F72T12/CW/HO)

(1) Optional UV 30W, (G30T8) Germicidal Lamp

f. Blower Motor

3/4 HP, 1050 RPM

g. Drain Pan Capacity

30.38 Gallons

9.4 Model - 1194 (4' Cabinet With Hinged Window)

a. Construction

Work Surface: Type 304 Stainless Steel, #4 Finish

Cabinet: Cold Rolled Steel

Finish: Antique White with Windsor Blue Trim

Baked-on Powder Coat Epoxy Paint

b. Dimensions

Exterior: 54"W x 64"H x 32.5"F-B

Interior: 49"W x 28.33"H x 22.25"F-B

c. Electrical Requirements

Main: 100-125 VAC, 1 Phase, 2 Wire, 60 Hz, 10 FLA

Receptacle: 115 VAC, 1 Phase, 2 Wire, 60 Hz, 12 FLA

Circuit Breaker: 15 Amp

d. Filters

(1) Supply HEPA Filter (18"W x 48"F-B x 5-7/8"H)

(1) Exhaust HEPA Filter (18"W x 24"F-B x 11.5"H)

e. Lights

(1) Fluorescent 60W, (F48T12/CW/HO)

(1) Optional UV 30W, (G30T8) Germicidal Lamp

f. Blower Motor

3/4 HP, 1625 RPM

g. Drain Pan Capacity

21.14 Gallons

9.4 Model - 1196 (6' Cabinet With Hinged Window)

a. Construction

Work Surface: Type 304 Stainless Steel, #4 Finish

Cabinet: Cold Rolled Steel

Finish: Antique White with Windsor Blue Trim

Baked-on Powder Coat Epoxy Paint

b. Dimensions

Exterior: 78"W x 64"H x 32.5"F-B

Interior: 73"W x 28.33"H x 22.25"F-B

c. Electrical Requirements

Main: 100-125 VAC, 1 Phase, 2 Wire, 60 Hz, 12 FLA

Receptacle: 115 VAC, 1 Phase, 2 Wire, 60 Hz, 12 FLA

Circuit Breaker: 15 Amp

d. Filters

(1) Supply HEPA Filter (18"W x 72"F-B x 5-7/8"H)

(1) Exhaust HEPA Filter (24"W x 24"F-B x 11.5"H)

e. Lights

(1) Fluorescent 85W, (F72T12/CW/HO)

(1) Optional UV 30W, (G30T8) Germicidal Lamp

f. Blower Motor

3/4 HP, 1050 RPM

g. Drain Pan Capacity

30.38 Gallons

SECTION 10 - ACCESSORIES

10.1 ACCESSORIES - TYPE A/B₃ CABINETS

DESCRIPTION:	ORDER NO.	1184	1186	1194	1196
Service Valve, Vacuum	191141	-	-	-	-
Service Valve, Air	191142	-	-	-	-
Service Valve, Nitrogen	191143	-	-	-	-
Service Valve, Gas	191144	-	-	-	-
UV Light, 30W	191419	-	-	-	-
UV Light, 30W Portable	191070	-	-	-	-
Ground Fault, Duplex Outlet (R.S.)	191540	-	-	-	-
IV Rod, Stainless Steel 4' Cabinet	191522	-	-	-	-
IV Rod, Stainless Steel 6' Cabinet	191523	-	-	-	-
(* See note below)					
*30" Modular Stand, 4' Cabinet	191552	-	-	-	-
*30" Modular Stand, 6' Cabinet	191553	-	-	-	-
*36" Modular Stand, 4' Cabinet	191556	-	-	-	-
*36" Modular Stand, 6' Cabinet	191557	-	-	-	-
Locking Caster for modular stand	191416	-	-	-	-
Adjustable Foot Rest 4' Cabinet	191127	-	-	-	-
Adjustable Foot Rest 6' Cabinet	191129	-	-	-	-
Low Air Flow Alarm	191168	-	-	-	-
Exhaust Transition	191570	-	-	-	-

NOTE:

* APPROXIMATE HEIGHT FROM FLOOR TO WORK STATION.

ACCESSORIES ARE CUSTOMER INSTALLED!

SECTION 11 - PARTS LIST

11.1 Model 1184/1185

Stock #	Description
191510	1184, 4' Sliding Window
156079	3/4 HP Blower Motor (1625 RPM)
170045	Capacitor, Motor 25MFD, 370V
190396	Motor Speed Control
225250	Ballast (Fluorescent Lighting)
225414	48" Fluorescent Lamp (60W, 120V)
230054	Circuit Breaker, 15A SP
430200	Line cord Assembly, 15A, 120V, Hospital Grade
760063	Filter, Exhaust HEPA 18x24x11-1/2
760154	Filter, Supply HEPA 18x48x6
500009	Ballast, (U.V. Lighting)
141014	30W Germicidal Lamp
250013	Relay, DPDT. 10A. 120V
280005	Pilot Light, #312, Red
300055	Delay Relay, 3 to 300 Seconds
360095	Rocker Switch, SPST, Flat Black
360096	Pushbutton Switch, SPDT
360105	Rocker Switch, SPDT
249025	Valve Body w/Tip
104008	Gauge, Static Pressure
7011186	1186/1196 Technical Manual

Model 1185

Stock #	Description
275012	Transformer, 1.5 KVA, 240/120V
460052	Plug, European, 220V, 16A
500023	Ballast, F48T12, 220V, 50 Hz
500022	Ballast, 1 F30T8, 220V, 50 Hz

11.2 Model 1186 - Replacement Parts

Stock #	Description
191511	1186, 6' Sliding Window
156039	3/4 HP Blower Motor (1050 RPM)
170039	Capacitor, Motor 15MFD, 370V
190396	Motor Speed Control
225250	Ballast (Fluorescent Lighting)
225418	72" Fluorescent Lamp (85W, 120V)
230054	Circuit Breaker, 15A SP
430200	Line cord Assembly, 15A, 120V, Hospital Grade
285812	Plug 15A, 120V Hospital Grade
360157	Rocker Switch (5A)
249025	Valve Body w/Tip
104008	Gauge, Static Pressure
760064	Filter, Exhaust HEPA 24x24x11-1/2
760153	Filter, Supply HEPA 18x72x6
500009	Ballast, (U.V. Lighting)
141014	30W Germicidal Lamp
420057	25VA Transformer
7011186	1186/1196 Technical Manual

11.3 Model 1194 - Replacement Parts

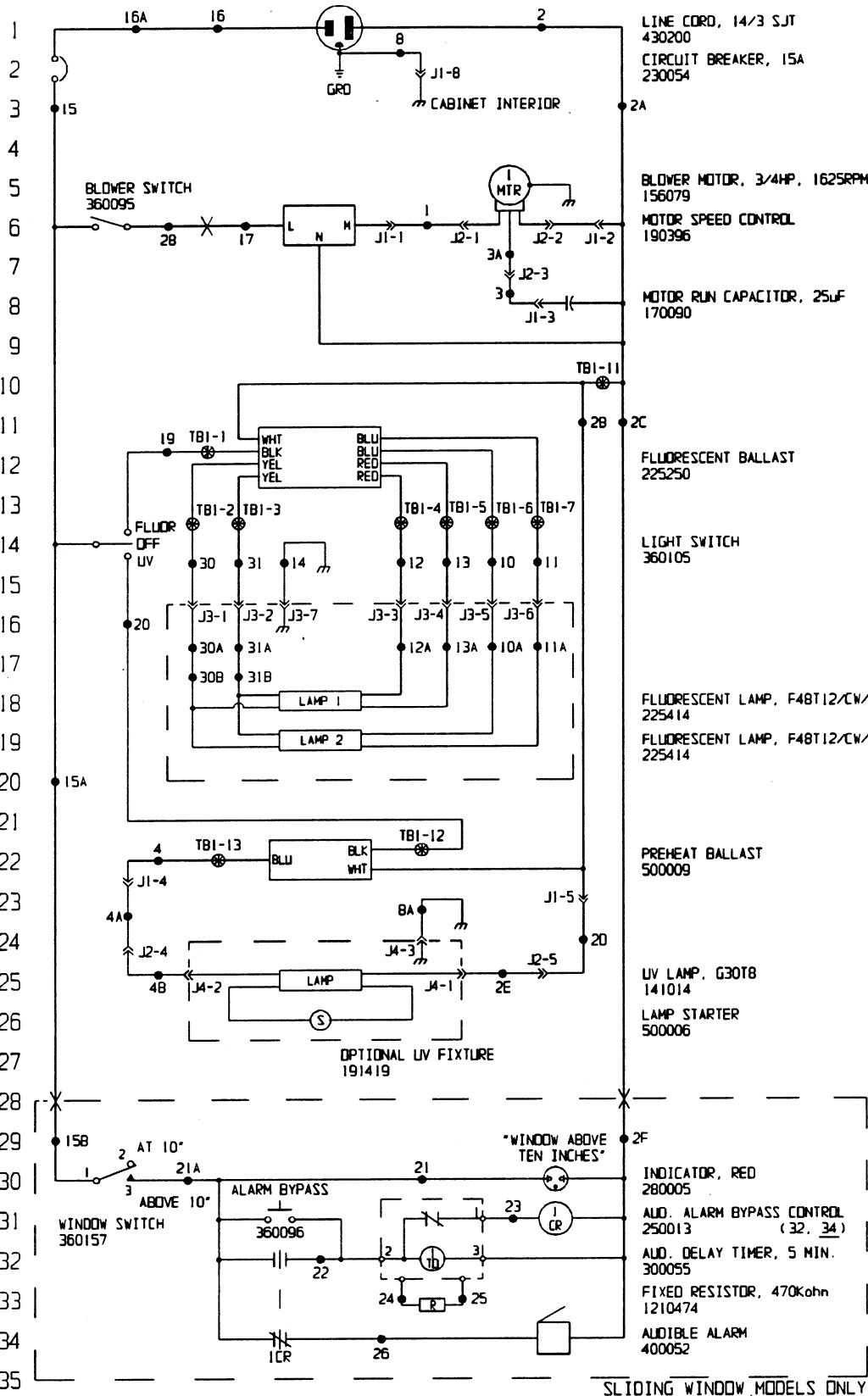
Stock #	Description
191530	1194, 4' Hinged Window
156079	3/4 HP Blower Motor (1625 RPM)
170045	Capacitor, Motor 25MFD, 370V
190396	Motor Speed Control
225250	Ballast (Fluorescent Lighting)
225414	48" Fluorescent Lamp (60W, 120V)
230054	Circuit Breaker, 15A SP
430200	Line cord Assembly, 15A, 120V, Hospital Grade
760063	Filter, Exhaust HEPA 12x24x11-1/2
760154	Filter, Supply HEPA 18x48x6
500009	Ballast, (U.V. Lighting)
141014	30W Germicidal Lamp
250013	Relay, DPDT. 10A. 120V
280005	Pilot Light, #312, Red
300055	Delay Relay, 3 to 300 Seconds
360095	Rocker Switch, SPST, Flat Black
360096	Pushbutton Switch, SPDT
360105	Rocker Switch, SPDT
249025	Valve Body w/Tip
104008	Gauge, Static Pressure
7011186	1186/1196 Technical Manual

11.4 Model 1196 - Replacement Parts

Stock #	Description
191531	1196, 6' Sliding Window
156039	3/4 HP Blower Motor (1050 RPM)
170039	Capacitor, Motor 15MFD, 370V
190396	Motor Speed Control
230054	Circuit Breaker, 15A SP
430200	Line cord Assembly, 15A, 120V, Hospital Grade
285812	Plug, 15A, 120V, Hospital Grade
225250	Ballast (Fluorescent Lighting)
225418	72" Fluorescent Lamp (85W, 120V)
360157	Rocker Switch (5A)
249025	Valve Body w/Tip
104008	Gauge, Static Pressure
760064	Filter, Exhaust HEPA 24x24x11-1/2
760153	Filter, Supply HEPA 18x72x6
420057	25VA Transformer
500009	Ballast, (U.V. Lighting)
141014	30W Germicidal Lamp
7011186	1186/1196 Tech Manual

SECTION 12 – ELECTRICAL SCHEMATICS

POWER CONNECTION
100-130 VAC, 1PH, 2W, 60HZ, 10FLA

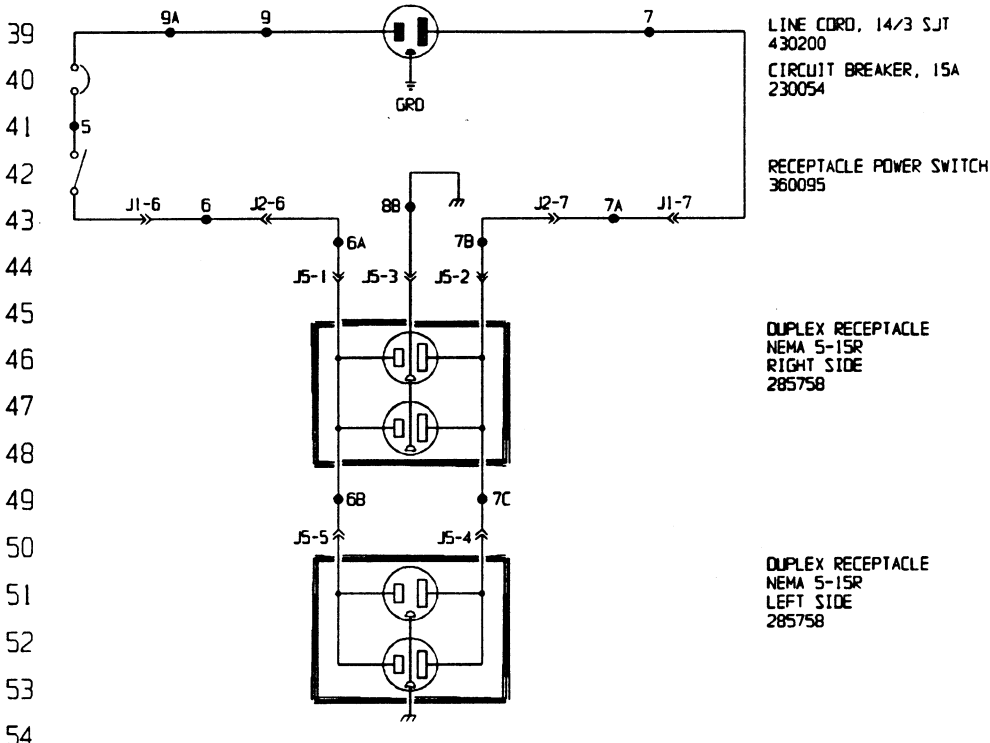


- LINE CORD, 14/3 SJT
430200
- CIRCUIT BREAKER, 15A
230054
- BLOWER MOTOR, 3/4HP, 1625RPM
156079
- MOTOR SPEED CONTROL
190396
- MOTOR RUN CAPACITOR, 25 μ F
170090
- FLUORESCENT BALLAST
225250
- LIGHT SWITCH
360105
- FLUORESCENT LAMP, F48T12/CW/HO
225414
- FLUORESCENT LAMP, F48T12/CW/HO
225414
- PREHEAT BALLAST
500009
- UV LAMP, G30T8
141014
- LAMP STARTER
500006
- INDICATOR, RED
280005
- AUD. ALARM BYPASS CONTROL
250013 (32, 34)
- AUD. DELAY TIMER, 5 MIN.
300055
- FIXED RESISTOR, 470Kohm
1210474
- AUDIBLE ALARM
400052

SLIDING WINDOW MODELS ONLY

Electrical Schematic
Forma Models:
1184 and 1194

POWER CONNECTION - STANDARD DUPLEX RECEPTACLES
 115 VAC NOMINAL, 60HZ, 12FLA MAX.



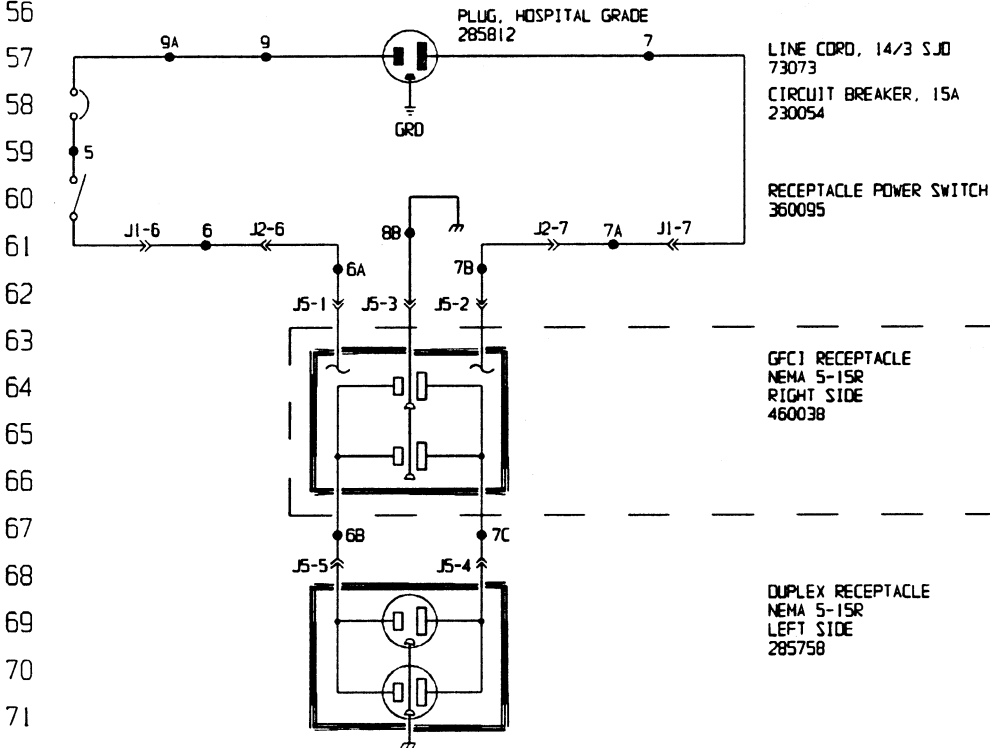
LINE CORD, 14/3 SJT
 430200
 CIRCUIT BREAKER, 15A
 230054

RECEPTACLE POWER SWITCH
 360095

DUPLEX RECEPTACLE
 NEMA 5-15R
 RIGHT SIDE
 285758

DUPLEX RECEPTACLE
 NEMA 5-15R
 LEFT SIDE
 285758

POWER CONNECTION - OPTIONAL GFCI RECEPTACLES
 115 VAC NOMINAL, 60HZ, 12FLA MAX.



LINE CORD, 14/3 SJT
 73073
 CIRCUIT BREAKER, 15A
 230054

RECEPTACLE POWER SWITCH
 360095

GFCI RECEPTACLE
 NEMA 5-15R
 RIGHT SIDE
 460038

DUPLEX RECEPTACLE
 NEMA 5-15R
 LEFT SIDE
 285758

Electrical Schematic
 Forma Models:
 1184 and 1194

WIRE REFERENCE CHART

	WIRE NO.	GAUGE	COLOR
77			
78	1	16	BLACK
	2	14	WHITE
79	2A	16	WHITE
	2B	20	WHITE
80	2C	20	WHITE
	2D	16	WHITE/BLACK
81	2E	16	WHITE
	2F	18	WHITE
82	3	16	YELLOW
	3A	16	RED
83	4	16	ORANGE
	4A	16	ORANGE
84	4B	16	BLACK
	5	16	BLUE
85	6	16	BROWN
	6A	14	ORANGE
86	6B	16	BLACK
	7	14	WHITE
87	7A	16	GRAY
	7B	14	RED
	7C	16	WHITE
88	8	16	GREEN
	8A	16	GREEN
89	8B	14	GREEN
	9	14	BLACK
90	9A	16	BLACK
	10	20	BLUE
91	10A	22	BLUE
	11	20	BLUE
92	11A	22	BLACK
	12	20	RED
93	12A	22	RED
	13	20	RED
94	13A	22	BROWN
	14	16	GREEN
95	15	16	BROWN
	15A	20	BROWN
96	15B	18	BROWN
	16	14	BLACK
97	16A	16	BLACK
	17	16	RED
	18	NOT USED	
98	19	16	PURPLE
	19A	18	BLACK
99	20	20	YELLOW
	21	20	RED
100	21A	18	RED
	22	20	BLUE
101	23	20	PURPLE
	24	RESISTOR	LEAD
102	25	RESISTOR	LEAD
	26	18	BLACK
103	27	NOT USED	
	28	16	ORANGE
104	29	NOT USED	
	30	20	YELLOW
	30A	22	YELLOW
105	30B	18	YELLOW
	31	20	YELLOW
106	31A	22	ORANGE
107	31B	18	YELLOW

NOTES:

<input checked="" type="checkbox"/> Denotes Terminal Strip Connection	Parts List Reference Number
1CR Lost Relay Number	<input type="checkbox"/> Assembly
13 Lost Terminal Number	<input type="checkbox"/> Panel
31 Lost Wire Number	<input type="checkbox"/> Refrigeration
	<input type="checkbox"/> 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



Forma Scientific

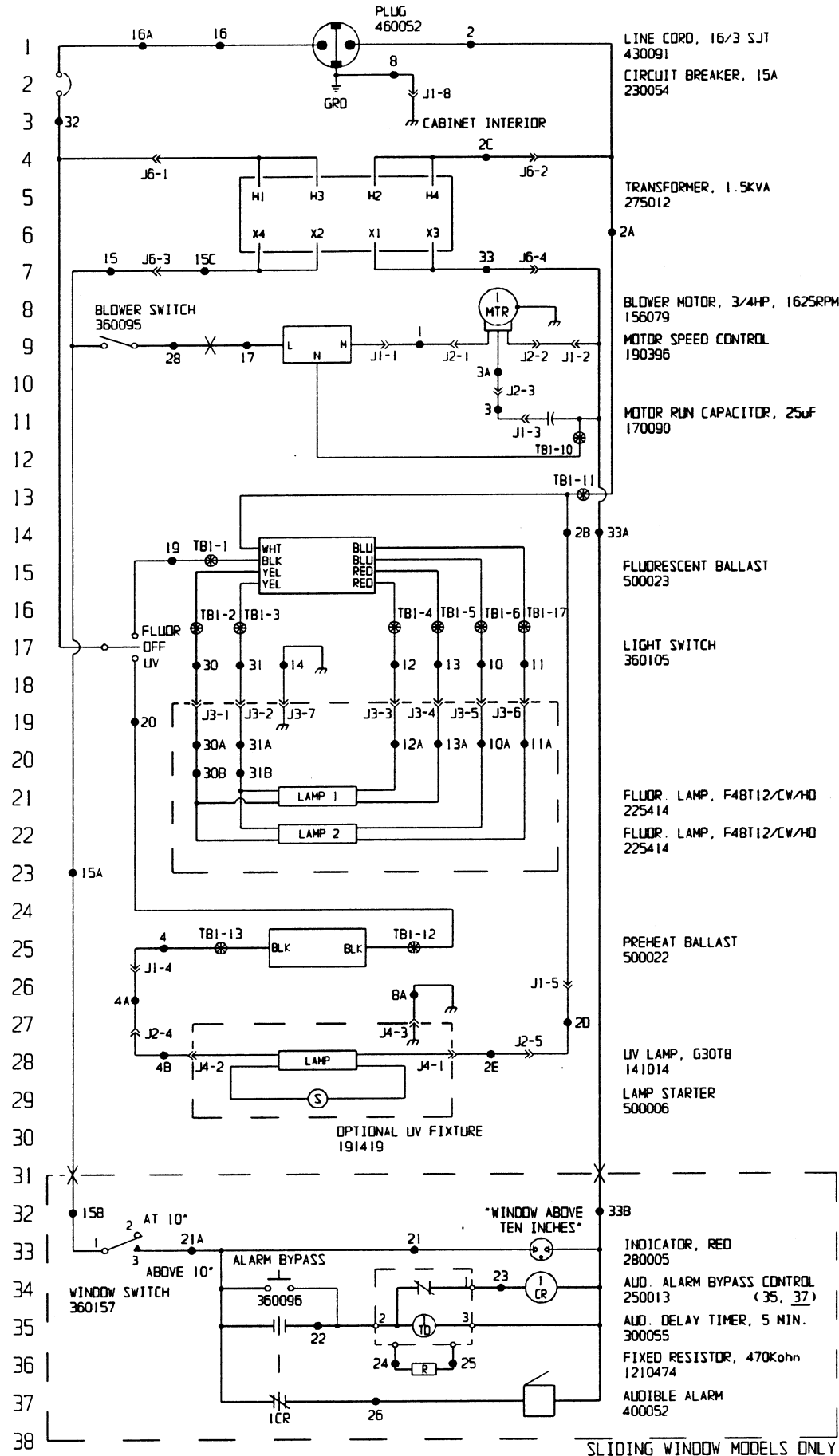
REV. 6/98 BAP/ETL, 0140 4570 1845 24-530
TELL FREE USA 800-848-3300, 0408 614-373-4183

6	HD-738	08-01-95	JAS	KDG		CORRECT RELAY LINE NUMBERS
5	HD-679	5-11-93	JAS	JAS	LON	REVISED SPEED CONTROL
4	HD-638	9-09-92	JAS	KDG	LON	REV. PLUG & CORD
3	SI-2991	4-28-92	JAS	JAS	LON	REV. ALARM WIRE GAUGES
2	HD-515	1-30-91	JAS	JAS	LON	ADD 2ND FLUOR. LAMP
REV ECR NO.	DATE	BY	CAO	APPD	DESCRIPTION OF REVISION	
DATE	3-3-90	DWN	GLM	CAO	JAS	APPD LON
CUSTOMER						
JOB TITLE	1184, 1194 LFBS					
DWG TITLE	ELECTRICAL SCHEMATIC					
LOCATION	HOODSO1		JOB NUMBER	DRAWING NUMBER		
				1184-70-0-D		

**Electrical Schematic
Forma Models:
1184 and 1194**

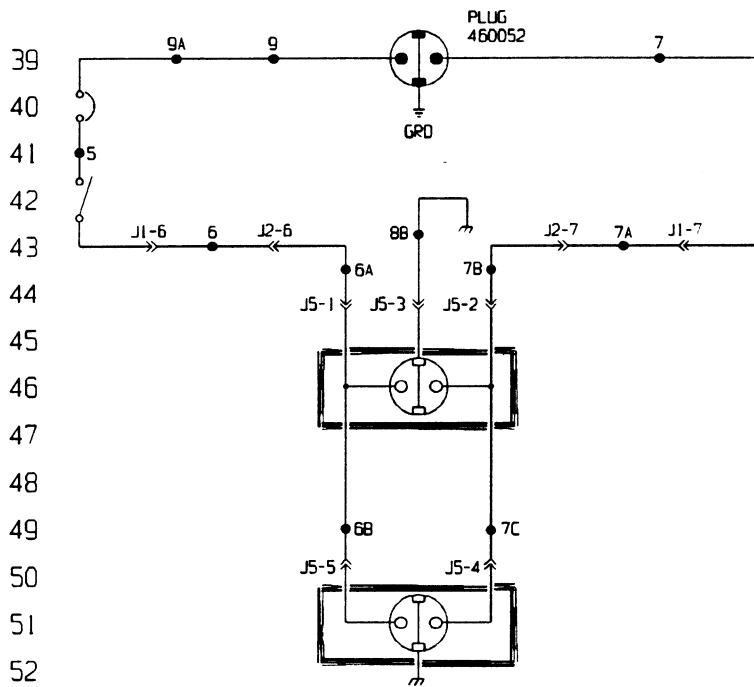
**1184-70-0-D Rev. 6
Page 3 of 3**

POWER CONNECTION
220VAC, 1PH, 2W, 50HZ, 6FLA



Electrical Schematic
Forma Models:
1185 and 1195

POWER CONNECTION - RECEPTACLES
 220VAC NOMINAL, 1PH, 2W, 50HZ, 12FLA MAX.



LINE CORD, 16/3 SJT
 430091
 CIRCUIT BREAKER, 15A
 230054

RECEPTACLE POWER SWITCH
 360095

RECEPTACLE WITH COVER
 RIGHT SIDE
 460128


RECEPTACLE WITH COVER
 LEFT SIDE
 460128

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

Electrical Schematic
 Forma Models:
 1185 and 1195

WIRE REFERENCE CHART

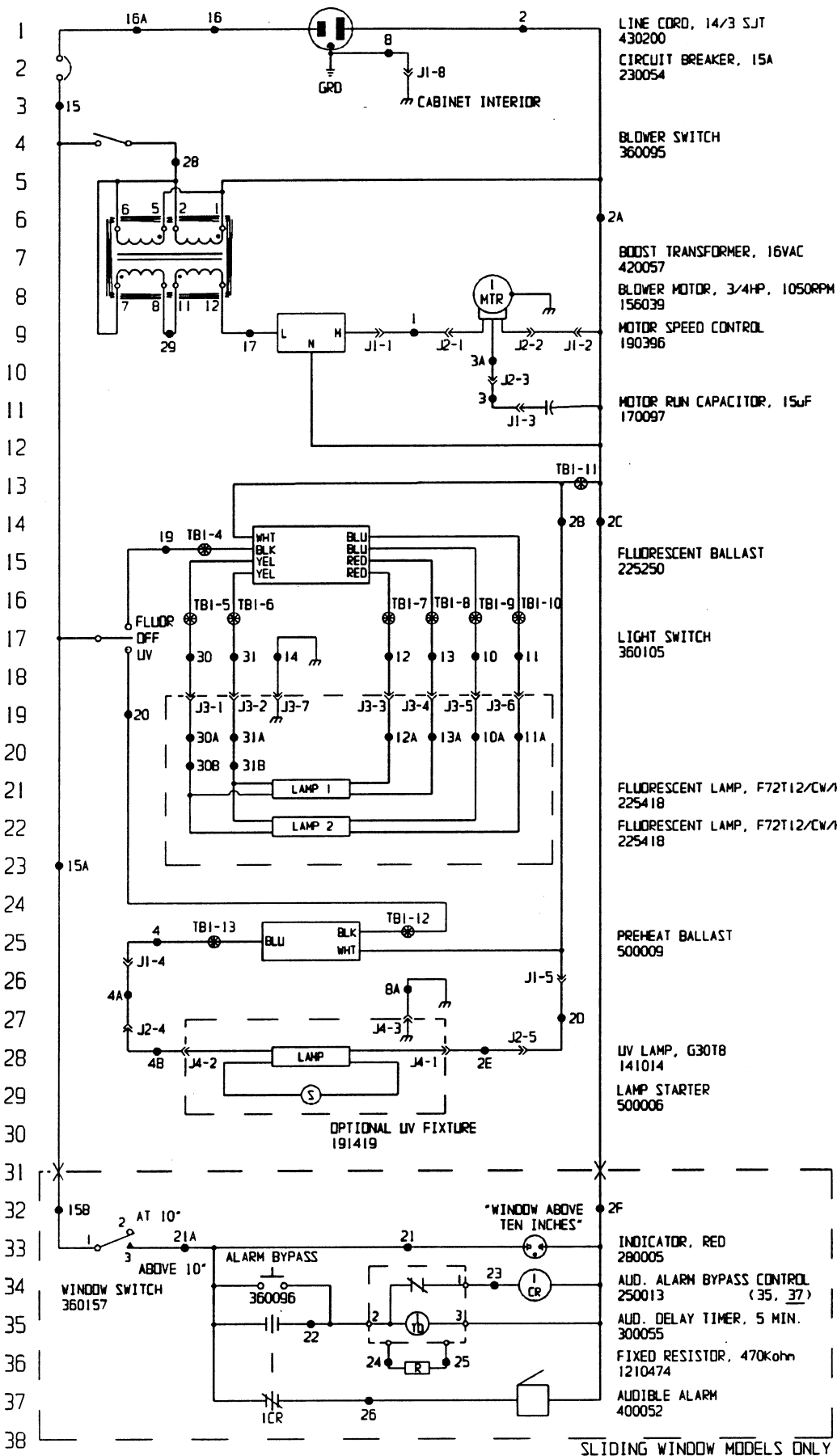
	<u>WIRE NO.</u>	<u>GAUGE</u>	<u>COLOR</u>
77			
78	1	16	BLACK
	2	16	BLUE
79	2A	16	WHITE
	2B	20	WHITE
80	2C	16	BLUE
	2D	16	WHITE/BLACK
81	2E	16	WHITE
	3	16	YELLOW
82	3A	16	RED
	4	16	ORANGE
83	4A	16	ORANGE
	4B	16	BLACK
84	5	16	BLUE
	6	16	BROWN
85	6A	14	ORANGE
	6B	16	BLACK
	7	16	BLUE
86	7A	16	GRAY
	7B	14	RED
87	7C	16	WHITE
	8	16	GREEN
88	8A	16	GREEN
	8B	14	GREEN
89	9	16	BROWN
	9A	16	BLACK
90	10	20	BLUE
	10A	22	BLUE
91	11	20	BLUE
	11A	22	BLACK
92	12	20	RED
	12A	22	RED
	13	20	RED
93	13A	22	BROWN
	14	16	GREEN
94	15	16	BROWN
	15A	20	BROWN
95	15B	18	BROWN
	15C	16	BLACK
96	16	16	BROWN
	16A	16	BLACK
	17	16	RED
97	18	NOT USED	
	19	16	PURPLE
98	19A	18	BLACK
	20	20	YELLOW
99	21	20	RED
	21A	18	RED
100	22	20	BLUE
	23	20	PURPLE
101	24	RESISTOR	LEAD
	25	RESISTOR	LEAD
	26	18	BLACK
102	27	NOT USED	
	28	16	ORANGE
103	29	NOT USED	
	30	20	YELLOW
104	30A	22	YELLOW
	30B	18	YELLOW
105	31	20	YELLOW
	31A	22	ORANGE
106	31B	18	YELLOW
	32	16	BROWN
	33	16	WHITE
107	33A	20	WHITE
	33B	18	WHITE

NOTES:		CUSTOMER APPROVAL/REFERENCE		5	HO-753	8-15-94	JAS	KDG		ADD TBI-10
<input checked="" type="checkbox"/> Denotes Terminal Strip Connection	Parts List Reference Number	APPROVED BY	DATE OF APPROVAL	4	HO-715	1-27-94	JAS	PKK	LON	ADDED MODEL 1195
ICR Last Relay Number	<input type="checkbox"/> Assembly	THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION AND SUCH INFORMATION IS NOT TO BE DISCLOSED TO OTHERS FOR ANY PURPOSE NOR USED FOR MANUFACTURING PURPOSES WITHOUT WRITTEN PERMISSION FROM FORMA SCIENTIFIC								
13 Last Terminal Number	<input type="checkbox"/> Panel	REV	ECR NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION		
33 Last Wire Number	<input type="checkbox"/> Refrigeration	DATE	12-5-91	DWN	CBG	CAD	CBG	APPD	LON	SCALE NTS
	<input type="checkbox"/> Wiring	 <p>Forma Scientific</p> <p>824 642 HANOVER, NH 03042 TEL: 603 241-5200 TOLL FREE USA 800-849-3880, 8410 844-373-4763</p>								
		CUSTOMER								
		JOB TITLE		1185/1195 LFBSC, 220V/50HZ						
		DWG TITLE		ELECTRICAL SCHEMATIC						
		LOCATION	JOB NUMBER	DRAWING NUMBER						
		HOODS01		1185-70-0-D						

**Electrical Schematic
Forma Models:
1185 and 1195**

**1185-70-0-D Rev. 5
Page 3 of 3**

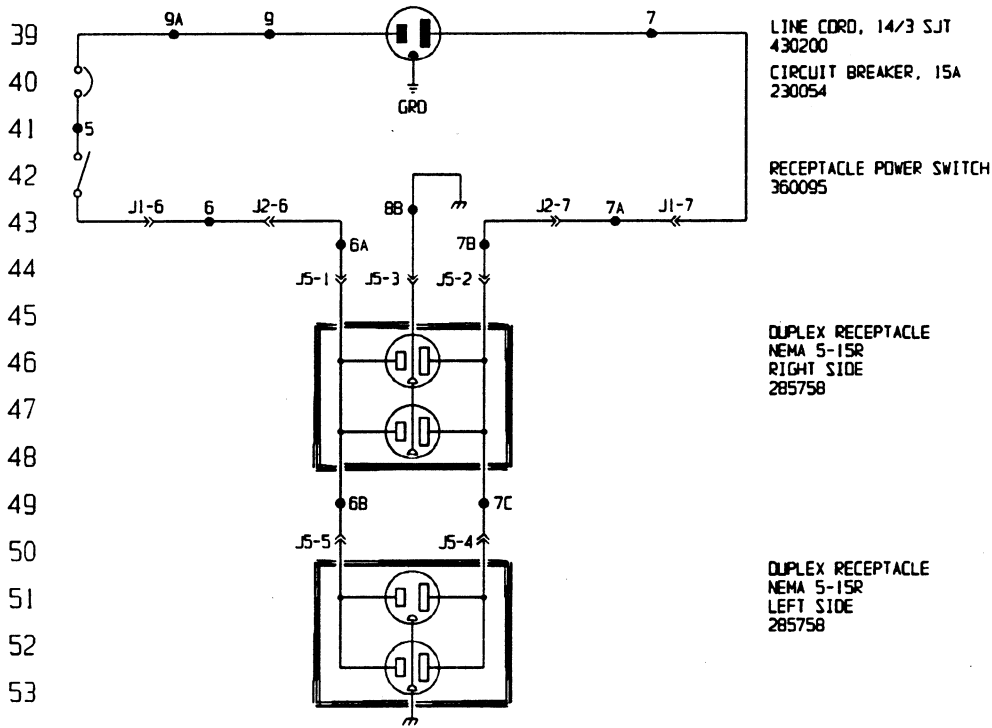
POWER CONNECTION
100-130 VAC, 1PH, 2W, 60HZ, 11FLA



- LINE CORD, 14/3 SJT
430200
- CIRCUIT BREAKER, 15A
230054
- BLOWER SWITCH
360095
- BOOST TRANSFORMER, 16VAC
420057
- BLOWER MOTOR, 3/4HP, 1050RPM
156039
- MOTOR SPEED CONTROL
190396
- MOTOR RUN CAPACITOR, 15uF
170097
- FLUORESCENT BALLAST
225250
- LIGHT SWITCH
360105
- FLUORESCENT LAMP, F72T12/CW/HO
225418
- FLUORESCENT LAMP, F72T12/CW/HO
225418
- PREHEAT BALLAST
500009
- UV LAMP, G30T8
141014
- LAMP STARTER
500006
- INDICATOR, RED
280005
- AUD. ALARM BYPASS CONTROL
(35, 37)
250013
- AUD. DELAY TIMER, 5 MIN.
300055
- FIXED RESISTOR, 470kOhm
1210474
- AUDIBLE ALARM
400052

Electrical Schematic
Forma Models:
1186 and 1196

POWER CONNECTION - STANDARD DUPLEX RECEPTACLES
 115 VAC NOMINAL, 60HZ, 12FLA MAX.



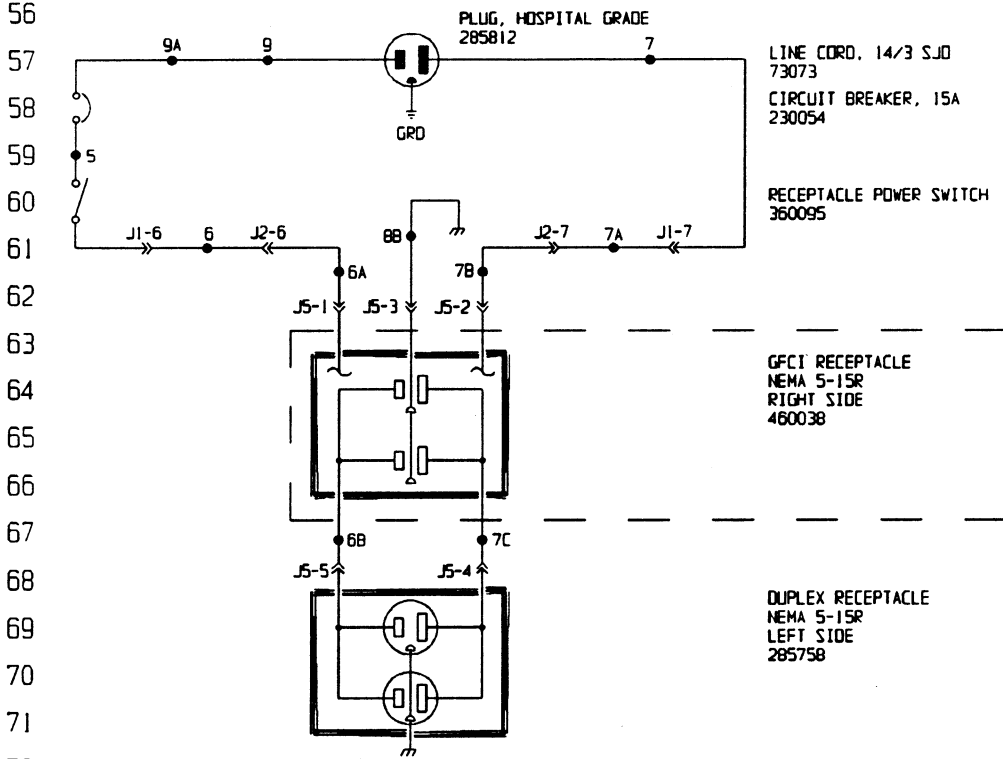
LINE CORD, 14/3 SJT
 430200
 CIRCUIT BREAKER, 15A
 230054

RECEPTACLE POWER SWITCH
 360095

DUPLEX RECEPTACLE
 NEMA 5-15R
 RIGHT SIDE
 285758

DUPLEX RECEPTACLE
 NEMA 5-15R
 LEFT SIDE
 285758

POWER CONNECTION - OPTIONAL GFCI RECEPTACLES
 115 VAC NOMINAL, 60HZ, 12FLA MAX.



PLUG, HOSPITAL GRADE
 285812
 LINE CORD, 14/3 SJO
 73073
 CIRCUIT BREAKER, 15A
 230054

RECEPTACLE POWER SWITCH
 360095

GFCI RECEPTACLE
 NEMA 5-15R
 RIGHT SIDE
 460038

DUPLEX RECEPTACLE
 NEMA 5-15R
 LEFT SIDE
 285758

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

WIRE REFERENCE CHART

	<u>WIRE NO.</u>	<u>GAUGE</u>	<u>COLOR</u>
77			
78	1	16	BLACK
	2	14	WHITE
79	2A	16	WHITE
	2B	20	WHITE
80	2C	20	WHITE
	2D	16	WHITE/BLACK
81	2E	16	WHITE
	2F	18	WHITE
82	3	16	YELLOW
	3A	16	RED
83	4	16	ORANGE
	4A	16	ORANGE
84	4B	16	BLACK
	5	16	BLUE
85	6	16	BROWN
	6A	14	ORANGE
86	6B	16	BLACK
	7	14	WHITE
87	7A	16	GRAY
	7B	14	RED
88	7C	16	WHITE
	8	16	GREEN
89	8A	16	GREEN
	8B	14	GREEN
90	9	14	BLACK
	9A	16	BLACK
91	10	20	BLUE
	10A	22	BLUE
92	11	20	BLUE
	11A	22	BLACK
93	12	20	RED
	12A	22	RED
94	13	20	RED
	13A	22	BROWN
95	14	16	GREEN
	15	16	BROWN
96	15A	20	BROWN
	15B	18	BROWN
97	16	14	BLACK
	16A	16	BLACK
98	17	16	RED
	18	NOT USED	
99	19	16	PURPLE
	19A	18	BLACK
	20	20	YELLOW
100	21	20	RED
	21A	18	RED
101	22	20	BLUE
	23	20	PURPLE
102	24	RESISTOR	LEAD
	25	RESISTOR	LEAD
103	26	18	BLACK
	27	NOT USED	
	28	16	ORANGE
104	29	16	YELLOW
	30	20	YELLOW
105	30A	22	YELLOW
	30B	18	YELLOW
106	31	20	YELLOW
	31A	22	ORANGE
107	31B	18	YELLOW

NOTES:	
⊗ Denotes Terminal Strip Connection	Parts List Reference Number
ICR Lost Relay Number	○ Assembly
13 Lost Terminal Number	○ Panel
31 Lost Wire Number	○ Refrigeration
	□ Wiring

CUSTOMER APPROVAL/REFERENCE	
APPROVED BY	_____
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	



Forma Scientific

800 640 4467 FAX 800 427 7811 TEL 640 427 7811 FAX 640 427 7811

7	HO-738	08-01-95	JAS	KDG		CORRECT RELAY LINE NUMBERS
6	HO-679	5-11-93	JAS	JAS	LON	REVISED SPEED CONTROL
5	HO-638	9-9-92	JAS	KDG	LON	REV. PLUG & CORD
4	SI-2991	5-1-92	JAS	JAS	LON	REV. ALARM WIRE GAUGES
3	HO-597	11-12-91	JAS	JAS	LON	CHANGE CAP. FROM 170039
REV	ECR NO.	DATE	BY	CAD	APPD	DESCRIPTION OF REVISION
	DATE	3-3-90	OWN	GLM	CAD	JAS APPD LON SCALE NTS
CUSTOMER						
JOB TITLE 1186, 1196 LF8SC						
DWG TITLE ELECTRICAL SCHEMATIC						
LOCATION			JOB NUMBER		DRAWING NUMBER	
HOODSOJ					1186-70-0-D	

**Electrical Schematic
Forma Models:
1186 and 1196**