

Instruction Manual

Savant[®] SPD121P SpeedVac[®] Concentrator



197-3002-00 Rev. C

Analyze • Detect • Measure • Control™

Thermo
ELECTRON CORPORATION

Contents

Section	Overview	Page
1.0	Overview of SPD121P Unit	1
1.1	Cover Lock Override	2
1.2	Installation of the SPD121P	3
1.3	How to Hook Up the SPD121P	3

Section	Control Panel	Page
2.0	The Control Panel	4
2.1	Description of the Control Panel	4

Section	Manual Run	Page
3.0	The Manual Run	5
3.1	The Auto Run	6-7

Section	Rotor Selection	Page
4.0	Rotor Selection Guide	8

Section	Guidelines	Page
5.0	Guideline for Solvent Choice	9

Section	Maintenance	Page
6.0	Maintenance & Service	10

Section	Specifications	Page
7.0	Specifications & Warranty	10
	Appendix 1 - Additional SPD121P System Set-ups	11

Disclaimer:

All statements, information and data given herein are believed to be accurate and reliable but are presented without guarantee, or responsibility of any kind, expressed or implied. Statements or suggestions concerning possible use of our products are made without representation that any such use is free of patent infringement and are not recommendations to infringe any patent. The user should not assume that all safety measures are indicated, or that other measures may not be required.

1.0 Overview of the SPD121P Unit

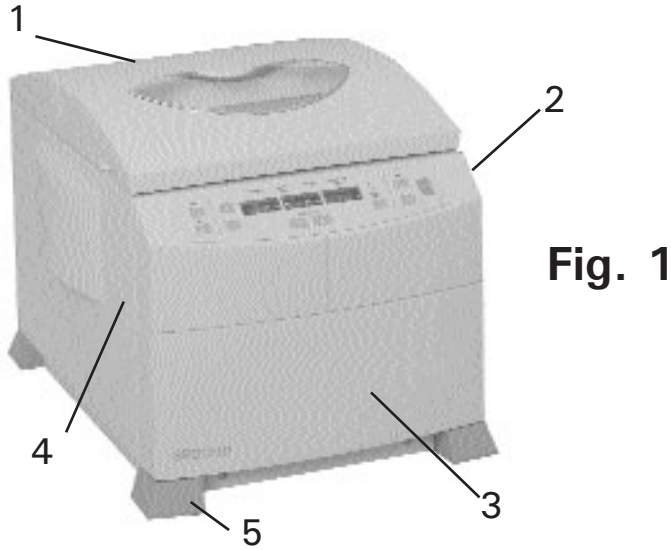
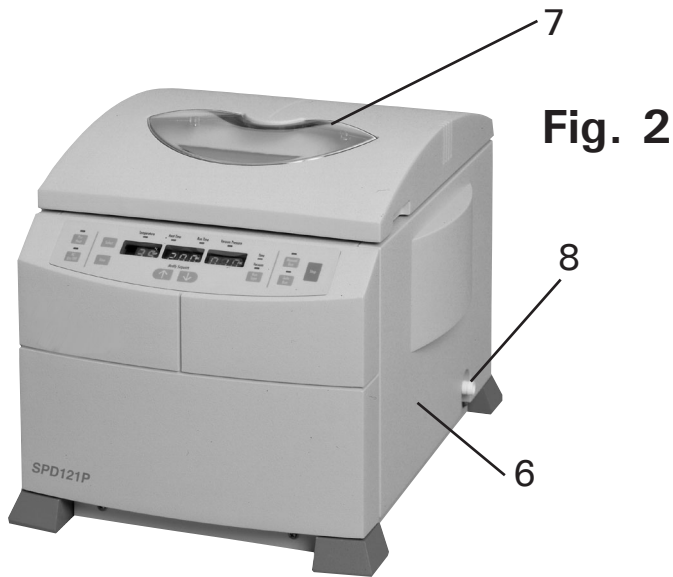


Figure 1 = Front View
 Figure 2 = Side View
 Figure 3 = Back View

Fig. 1



- 1. Top Radiant Cover
- 2. Control Panel
- 3. Front Panel
- 4. Left Side Panel
- 5. Anti Skid/Vibration
- 6. Right Side Panel
- 7. Chamber View Window
- 8. Vacuum Port
- 9. Vacuum Hose
- 10. Main On/Off Switch
- 11. Manual Cover Lock Release
- 12. Rear Access Panel
- 13. AC Socket

Fig. 2

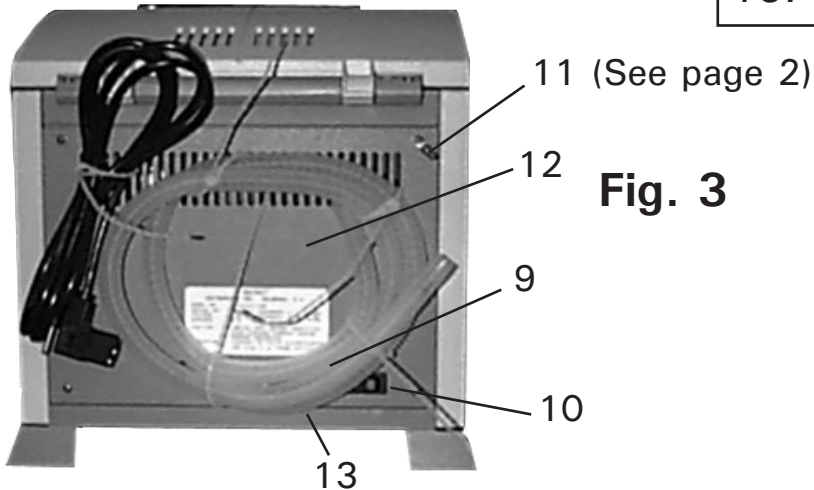
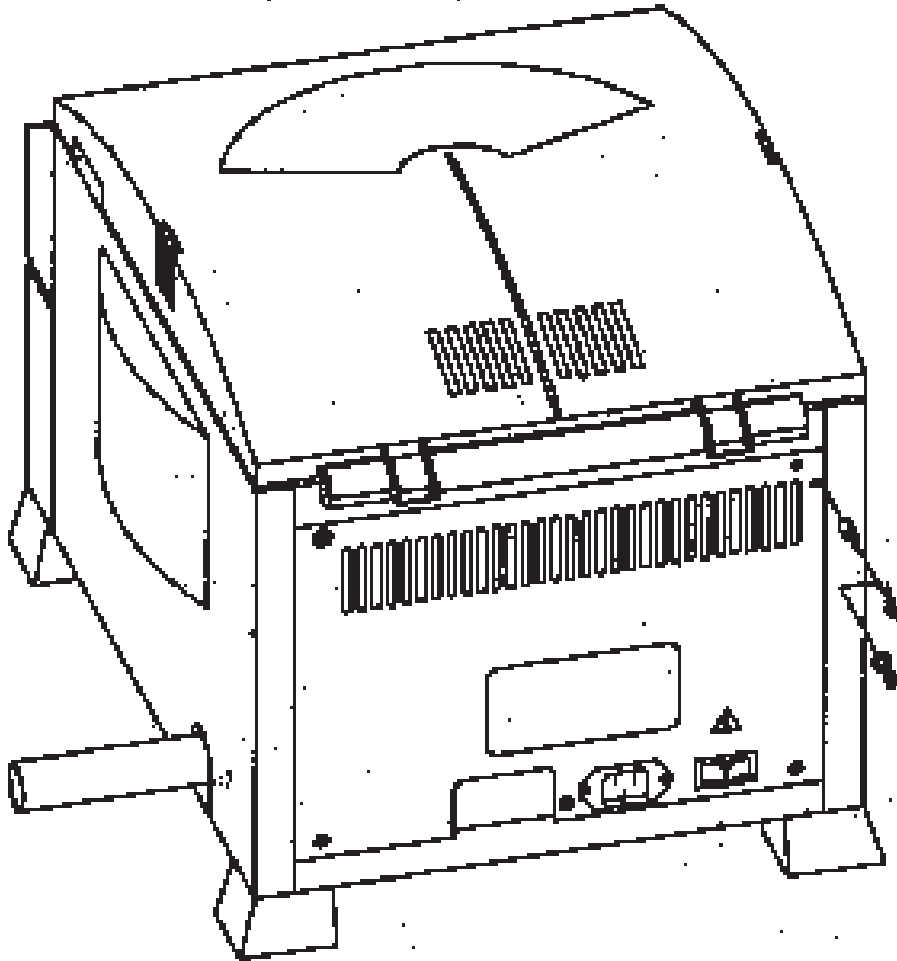


Fig. 3

1.1 COVER LOCK OVERRIDE



To Open The Top Cover During Power Failure:

Remove the screw holding the cover lock's manual release cord and lightly pull it as shown, while lifting the cover.

1.2 INSTALLATION OF THE SPD121P

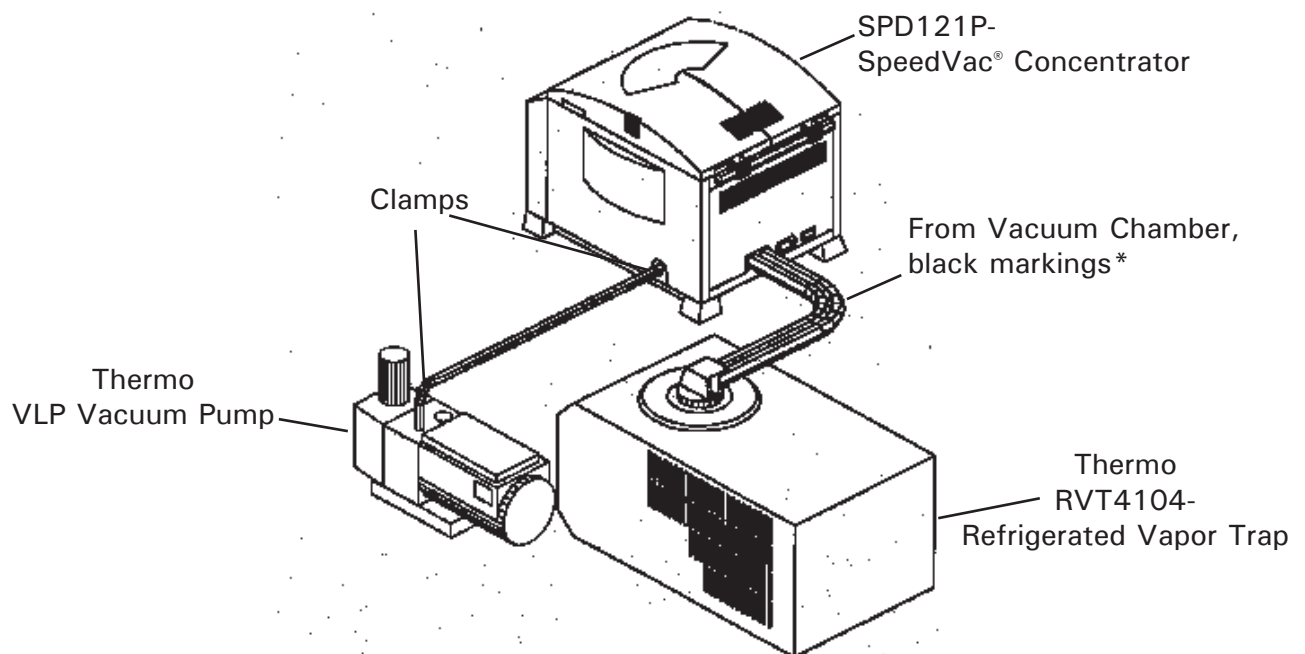


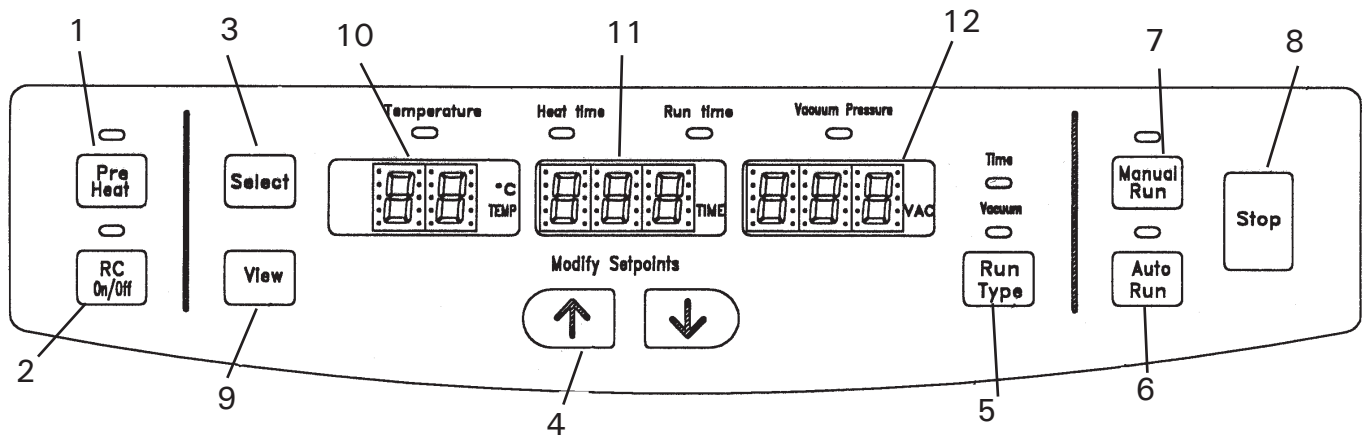
Figure 4-Rear View of System

1.3 How To Hook Up The SPD121P SpeedVac® System

1. Unpack unit from the box and verify that all the parts match packing list.
2. Read instruction manual carefully! If assistance is required, contact Thermo.
3. If the unit is part of a larger system, use the vacuum tubing supplied to hook up unit (See Figure 4).
4. If this SPD SpeedVac® is replacing an older single outlet vacuum tube SpeedVac, please see Appendix 1 on how to properly plumb the new SPD Concentrator.
5. If being hooked up to existing system, simply attach vacuum tubing to port on side of unit. (See item 8, figure 2-vacuum port).

CAUTION - Cold trap should be in line between the vacuum source and the SPD Unit.

2.0 SPD121P Control Panel



2.1 Description of SPD121P Control Panel

1. **PRE-HEAT** - use to pre-heat chamber to 45 °C prior to or between runs. Once run is initiated the pre-heat stops.
2. **RC ON/OFF** - Use to add radiant heat to chamber. Manual activation ON/OFF at any point during run. Also will turn off at end of run.
3. **SELECT** - Press this button to select the parameters to be modified. Selection is from left to right. GREEN light indicates the parameter to be modified.
NOTE: To select "VACUUM PRESURE", "RUN TYPE" light should light under "VACUUM".
4. **MODIFY SETPOINTS UP/DOWN** - Modifies selected parameter.
5. **RUN TYPE** - Selects a "Time" or "Vacuum" run.
6. **AUTO RUN** - Starts an automated "TIME" or "VACUUM" run (See page 5).
7. **MANUAL RUN** - Starts a "Manual" run (See page 4).
8. **STOP** - Terminates "Manual" or "Auto" run.
9. **VIEW** - Press to view. Displays preset parameters all at once.
10. **TEMPERATURE DISPLAY (RED DISPLAY)** - Indicates set temperature or actual temperature during a run in °C.
11. **TIME DISPLAYS (GREEN DISPLAY)** - Indicates heat timer or run timer set-points. In the process of a run it indicates elapsed time or time left.
12. **VACUUM PRESSURE DISPLAY (AMBER DISPLAY)** - Displays chamber vacuum in torrs or microns by shifting a decimal point. "Hpr." represents atmospheric pressure. No decimal point indicates microns.

3.0 THE MANUAL RUN SPD121P

1. Connect the unit to its required voltage.
2. Turn the power switch to ON, located on the back of the unit. The cover lock, disengages, allowing the top cover to be opened.

The display lights up, showing the following default values:

Temperature:	35 °C in RED
Run Time:	2.00 HRS. in GREEN
Vacuum Pressure:	01.0 in AMBER

Press "SELECT" to "HEAT TIME": 1.00 HR default in GREEN.

3. Place sample tubes in rotor so load is balanced. Secure rotor with the supplied knob (hand tight). Close cover.
4. Using the "SELECT" button and the up/down keys, set temperature between 35 °C-80 °C, or "no", for no heat.
5. Using the "SELECT" button and the up/down keys select and modify "TEMPERATURE", and "HEAT TIME" to between 0.01 and 9.59 hours or CCC (for continual heating).

When the heat timer expires, the heater will shut off no matter what the temperature setpoint reads (except if CCC).

6. Close cover.
7. Press the "MANUAL RUN" button. The cover locks and rotor starts turning. The decimal point blinks and the time display is counting up. The temperature rises to the set temperature. The "HEAT TIME" will count down and the SAV valves will both actuate, applying vacuum to the chamber.

NOTE: If the cover is not closed, the display will show "Lid" and the run will not start.

8. Press R/C for radiant cover heat. Press at any time to turn OFF and ON.
9. To end the manual run, press "STOP" button. The display will show "End", the SAV valves will click, isolating the chamber from the vacuum pump and also allowing air to bleed into the chamber.
10. After the rotor stops spinning, the cover lock disengages and the display reverts to last set parameters.
11. Open the cover and remove samples.

3.1 THE AUTO RUN TIME OR VACUUM SPD121P TIME



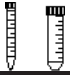



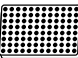
1. Refer to the "MANUAL RUN" section for start up.
2. Place the sample tubes in rotor so load is balanced. Secure rotor with supplied knob (hand tight). Close cover.
3. To execute an AUTO "TIMED" RUN, execute the following steps:
 - a) Press "RUN TYPE" to select "TIME" run
 - b) Use "SELECT" button and the up/down keys to select and modify "TEMPERATURE, "HEAT TIME, "RUN TIME" parameters ONLY, (The unit will not allow setting of vacuum pressure). RUN and HEAT TIME can be set from 0.01 to 9.59 hours (HEAT TIME also has "CCC").
 - c) Press "AUTO RUN" button to start the run. The cover locks and the rotor starts spinning. The time display is counting down in 1 minute intervals and the decimal point blinks. The heat time is counting down (use select button to view "HEAT TIME"). The temperature rises in 1 °C increments to set temperature. The vacuum pressure begins decrementing down from "HPr" (atmospheric pressure), after both SAV valves actuate, applying vacuum to the chamber.
 - d) The vacuum display will indicate vacuum pressure in the chamber.
NOTE: If the cover is not closed, the display will show "Lid" and the run will not start.
 - e) Press R/C for radiant cover heat. Press at any time to turn OFF and ON.
 - f) Once the time decrements to 0.00 HRS. the run will automatically stop, the display will show "End", the SAV valves will click, also allowing air to bleed into the chamber.
 - g) After the rotor stops spinning, the cover unlocks and the display reverts to last set parameters.
 - h) Open the cover and remove samples.

VACUUM

1. Executing an AUTO "VACUUM" RUN type:
(GREEN light ON above vacuum).
 - a) Press "RUN TYPE" to select "VACUUM" RUN.
 - b) Using "SELECT" button and the up/down keys set "VACUUM" pressure at which the run should end. (Ex. 500 Microns), would be entered as 00.5 torr.
 - c) Close cover.
 - d) Press "AUTO RUN" button to start the run. The cover locks and the rotor starts turning. The display will count up in 1 minute intervals and the decimal point blinks. The heat time will count down in 1 minute intervals. The temperature rises in 1 °C increments from 35 °C to set temperature. The vacuum pressure is declining from Hpr. (Atmospheric Pressure), after both SAV valves actuate, allowing vacuum into the chamber.
 - e) Press R/C for radiant cover heat. Press at any time to turn OFF and ON.
 - f) Once the unit senses the set vacuum pressure in the chamber has been achieved (< 500 Microns) it ends the run and the display will show "End", the SAV valves will click, also allowing air to bleed into the chamber.
 - g) After the rotor stops spinning, the cover unlocks and the display reverts to last parameters.
 - h) Open the cover and remove samples.

GENERAL: During the run, display shows actual parameters.
To check set parameters press "VIEW" button and "SELECT".
The display will revert temporarily to set points.

4.0 Rotor Selection Guide

	Working Volume (ml)	Number of Tubes	Description	Rotor Model
MICROCENTRIFUGE TUBES 	1.2 - 1.6	40	1.5 - 2.0 ml	RH40-11
	1.2 - 1.6	64	1.5 - 2.0 ml	RH64-11
	1.2 - 1.6	120	1.5 - 2.0 ml	RH120-11
GLASS AND PLASTIC TUBES 	0.3	100	0.5ml (8 x 29mm)	RH100-8
	0.3	40	0.4ml (6x 50mm)	RH40-6
	0.3	100	0.4ml (6 x 50mm)	RH100-6
	4	20	12 x 75mm	RH20-12
	4	40	1.5 - 2.0 (12 x 75mm)	RH40-12
	4	72	12 x 75mm	RH72-12
	8	10	13 x 100mm	RH20-12
	8	32	13 x 100mm	RH32-13
	10	8	18 x 100mm, 17 x 95, 16 x 100	RH8-18
CENTRIFUGE TUBES 	12	10	15ml conical (16 x 120mm)	RH10-15
	40	6	50ml conical (28 x 115mm)	RH6-50
FLASKS 	35	8	50ml pear shaped flask	RH8-50
	80	4	100ml pear shaped flask	RH4-100
VIALS 	2	60	12 x 32mm	RH60-12-40
	2.0	60	12 x 40mm vials	RH60-12-40
	2.4	12	20 x 47mm v-vials	RH12-20
	3	24	1 dram vials (15 x 45mm), 4ml	RH24-15
	4	12	20 x 60mm v-vials	RH12-20
	5.6	24	18 x 52mm mini-scintillation vials	RH24-18
	16	12	28 x 60mm scintillation vials	RH12-28
MICROTITER PLATES (Shallow) 	0.3	2 plates	Microplates	RH2MP
	0.3	6 plates	Microplates	RHSW6M6
DEEPWELL PLATES 	2.0	2 plates	Deepwell Microplates	RHDW2MP

5.0 GUIDELINES FOR SOLVENT CHOICE

Part I Solvents suited for the SPD121P unit

- Ethanol
- Methanol
- Formic Acid
- Water
- Acetonitrile
- Methylene Chloride
- Chloroform
- Ethyl Acetate
- Hydrochloric Acid
- Trifluoroacetic Acid
- Dimethyl Sulfoxide (DMSO)

Part II All other solvents not listed above

Contact a Thermo Application Scientist
to see if your solvent can be used with the SPD121P concentrator.

6.0 MAINTENANCE/SERVICE

1. Maintenance: The SpeedVac® SPD121P requires no scheduled maintenance.
2. Cleaning: The SpeedVac® SPD121P should be cleaned if solvents spill on or inside the unit. Always clean up any spills immediately using absorbent towels.

Always wear gloves when cleaning and dispose of paper towel in appropriate designated refuse containers.

3. Replace chamber seal if cracked (Part number is 197-6020-00).
4. Outside of unit can be cleaned with dilute solution of soap and water.
5. For any other maintenance or service issues or service problems, please contact Thermo.

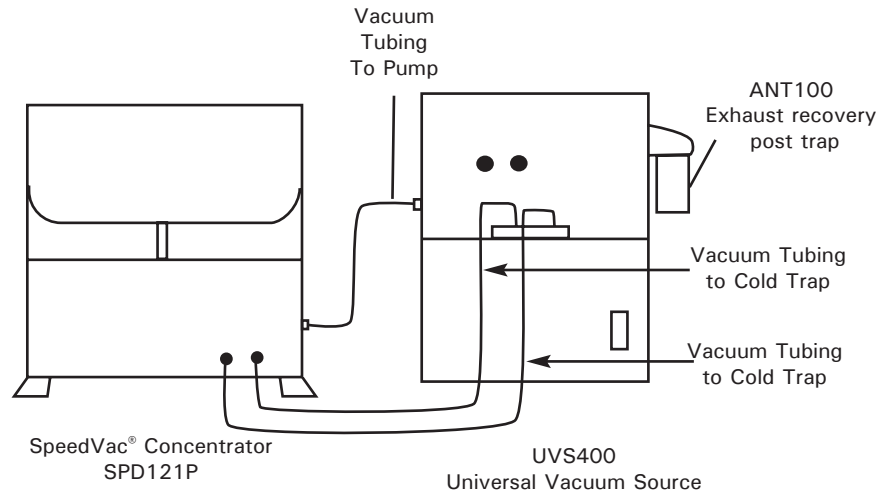
7.0 SPECIFICATIONS/WARRANTY

Model:	SPD121P
Bleeder Valve:	Integrated Automatic Bleeder Valve
Temperature Range:	35 °C-80 °C
Maximum Tube Size:	18 x 150 mm
Maximum Tube Capacity:	6
Maximum Carrier Capacity:	2 (96-deepwell plates) see chart on page 6
Dimensions: (W x D x H) in.:	14 x 18 x 13
cm:	36 x 45 x 33
Weight: lbs. (kg)	31 (14)
Power requirements:	115 VAC/60Hz, 5.0 A (Part number SPD121P-115) or 230 VAC/50Hz, 3.0 A (Part number SPD121P-230)

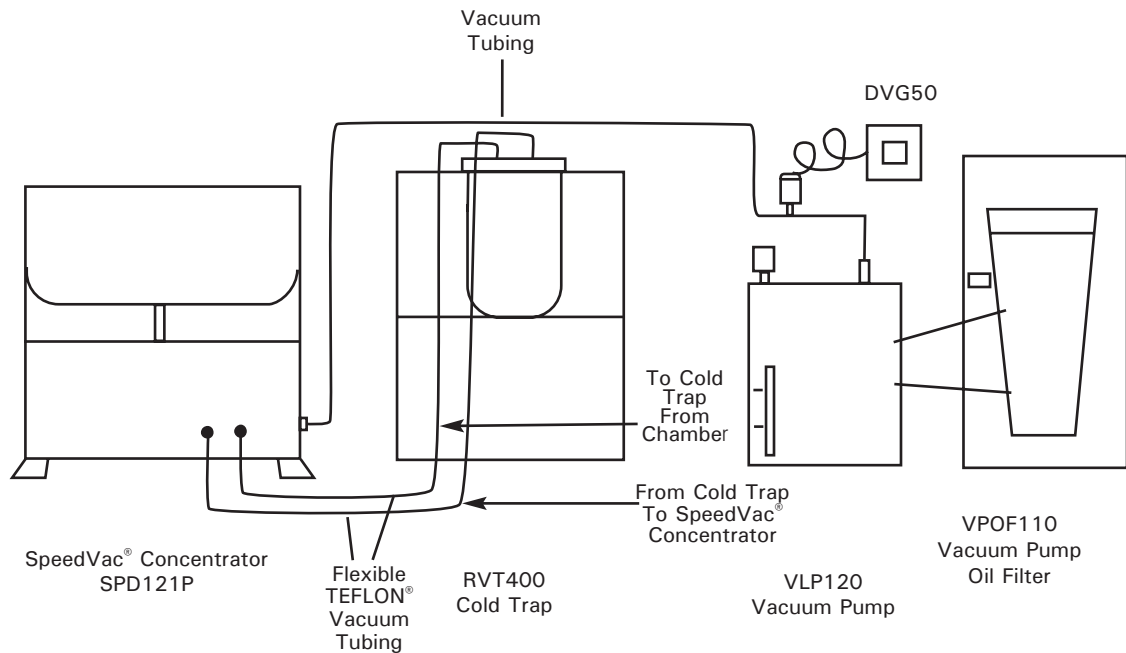
All Thermo products mentioned in this manual (except glassware) are warranted against defects in material and workmanship for one year after the date of delivery to the original purchaser. Thermo's warranty is limited to defective materials and workmanship, and does not cover incidental or consequential damages. Warranty work is subject to our inspection of the unit. No instruments, equipment, or accessories will be accepted without a Return Material Authorization (RMA) number issued by Thermo. Costs of shipping the unit are not covered under warranty. The warranty obliges you to follow the precautions in this manual. It is the responsibility of the user to dispose of ALL materials in a manner in accordance with all federal, state and local regulations. ALL RETURNED UNITS MUST BE DECONTAMINATED AND FREE OF RADIOACTIVITY AND SHOULD BE ACCOMPANIED WITH A DECONTAMINATION FORM. PLEASE CALL THERMO TO HAVE THIS FORM FAXED TO YOU! Under no circumstances shall Thermo be liable for damages due to the improper handling or use of its products. Thermo assumes no liability, express or implied, for your use of this equipment.

APPENDIX 1

Additional SPD121P System Set-ups



SEMI-INTEGRATED SYSTEM



COMPONENT HIGH VACUUM SYSTEM

SPD-SERIES SPEEDVAC® FAMILY



SPD101B



SPD111V



SPD121P



Thermo Electron Corporation

Bioscience Technologies

450 Fortune Boulevard
Milford, MA 01757
866.9.THERMO (866.984.3766) • Fax: 508.634.2199
www.thermo.com

Laboratory Pipetting and Consumables

info.pipettes@thermo.com

Molecular Biology

info.molbio@thermo.com

Microplate Instrumentation

info.microplateinstruments@thermo.com

Sample Preparation

info.sampleprep@thermo.com

Laboratory Automation & Integration

info.labautomation@thermo.com

New Labs

newlabs@thermo.com

Controlled Environment

info.controlenv@thermo.com

Services

services.biosciencetech@thermo.com

Bioscience
Technologies

International Sales Office Locations

Belgium

Brussels
+32 2 482 30 30
Fax: +32 2 482 30 31

France

Cergy Pontoise Cedex
+33 1 34 32 51 51
Fax: +33 1 34 32 51 59

Germany

Dreieich
+49 6103 408 0
Fax: +49 6103 408 1222

Netherlands

Breda
+31 76 571 4440
Fax: +31 76 587 9757

Russia

Saint-Petersburg
+7 812 325 8045
Fax: +7 812 186 1194

Moscow

+7 095 755 9045
Fax: +7 095 755 9046

Spain

Barcelona
+34 93 2233154
Fax: +34 93 2230857

Sweden

Stockholm
+46 8 742 03 90
Fax: +46 8 742 09 47

Lund

+46 46 90 96 60
Fax: +46 46 32 87 70

United Kingdom

Basingstoke, Hampshire
+44 01256 81782
Fax: +44 01256 81792

China

Beijing
+8610 5850 3588
Fax: +8610 6621 0847

Shanghai

+8621 5465 7588
Fax: +8621 6445 7830

Hong Kong

Wanchai
+852 2885 4613
Fax: +852 2567 4447

India

Bangalore
+91 22 2778 1101
Fax: +91 22 2778 1103

Japan

Yokohama-City
+81 45 453 9122
Fax: +81 45 453 9222

Thermo
ELECTRON CORPORATION

© 2001-2004 Thermo Electron Corporation, All rights reserved.
Thermo Electron Corporation, SpeedVac, and Analyze, Detect, Measure,
Control, are trademarks of Thermo Electron Corporation.