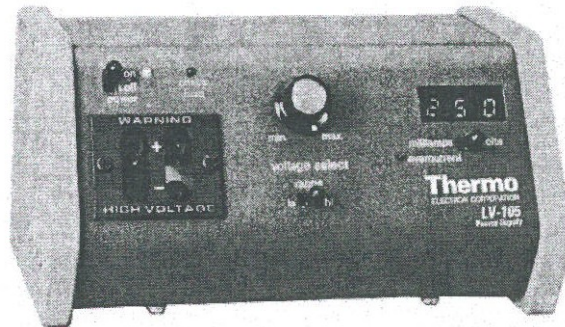
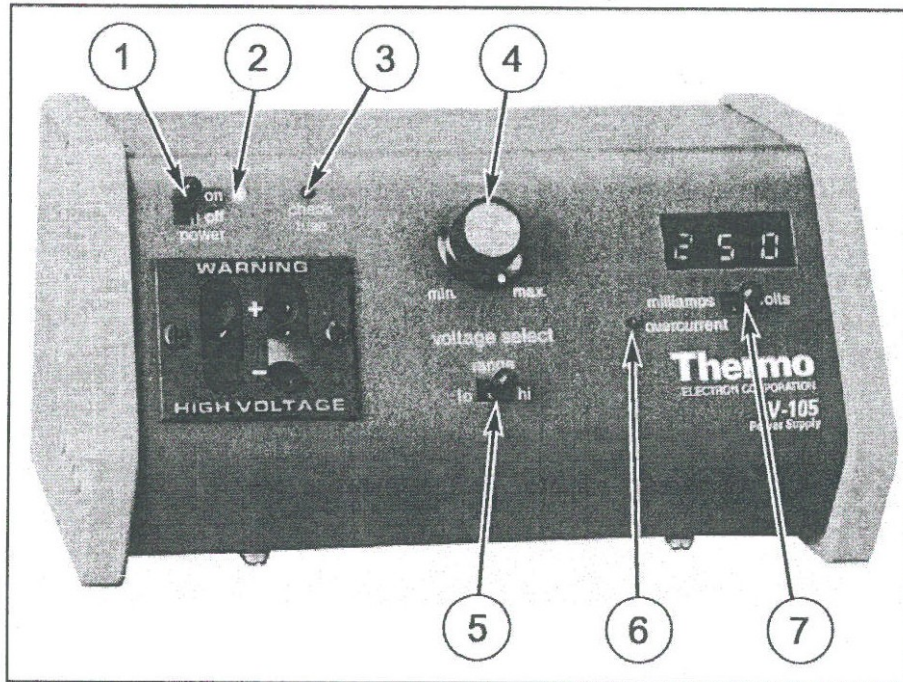


**Instructions**  
**EC-105**  
**Electrophoresis Power Supply**



**IMPORTANT**  
Please read these instructions carefully  
before using this apparatus

**Thermo**  
ELECTRON CORPORATION



- 1- Main Power Switch
- 2- Power On LED
- 3- Check Fuse LED
- 4- Voltage Select Control
- 5- Lo/Hi Range Switch
- 6- Overcurrent LED
- 7- Volts/Milliamp Display

## **INTENDED USES AND SPECIFICATIONS**

The EC-105 Power Supply is intended for use with any electrophoretic device designed to operate below 250 Volts and 500 milliAmps DC. Two sets of outputs operate in parallel to provide between 20 and 250 volts, up to 500 milliamps, and up to 75 watts in the constant voltage mode. A single LED display provides output voltage or current readings. Overcurrent protection is provided by audible and visual alarms and a single 1 amp slo blow fuse on the main circuit.

### **(1) SET UP**

#### **(a) Unpack and Check:**

Unpack and carefully examine the power supply. Report any damage to the distributor. Do not attempt to operate this device if physical damage is present. Save all packing material if damage is found.

#### **(b) Electricity Supply:**

Before connecting this device to the electricity supply, check the information given on the power supply rating label and confirm the following:

- Your electricity supply is single phase A.C. (alternating current) of the stated frequency with neutral nominally at earth potential.
- Your supply voltage is within the stated range.
- The current rating is within the capacity of your supply outlet.

### **WARNING: THIS EQUIPMENT MUST BE EARTHED.**

The wires in the electrical lead are colored in accordance with the following code:

Green	Earth (Ground)	White	Neutral (Common)	Black
			Live	(Phase)

Connect the wires to a non-reversible plug as follows:

Green wire to the terminal marked E (Earth), G (Ground), or colored Green.

White wire to the terminal marked N (Neutral) or Common.

Black wire to the terminal marked L (Live) or Phase.

### **IMPORTANT ELECTRICAL SAFETY NOTES:**

Consult a qualified electrician if any doubt or if your electricity supply system has any of the following:

- No Earth
- A Color Code Different from the Above
- Reversible Plugs
- Supply and Return Leads that are Both Above Earth Potential

#### **(c) Locate:**



Place the power supply in proximity to the electrophoretic cell(s) with which it is to be connected. Be sure to place the power supply in a safe, dry location with all the controls accessible.

**(d) Connect to Cell and Mains:**

Connect power supply to electrophoretic cell, then to the AC receptacle.

**(2) OPERATION**

**Power On:**

After the power supply is properly connected to the electrophoretic cell(s), select the lo range and set Voltage Select to the fully counter-clockwise (minimum) position. Set the display switch to Volts and turn on power using the Power switch. When "On," the red LED is illuminated.

**(b) Select Voltage:**

- (a) Adjust the Voltage Select control to the desired output voltage. The Select Control will increase the output voltage approximately 10 volts for every incremental adjustment. The lo/hi range switch is set to the hi position whenever operating voltages above 140 volts are required.

PLEASE NOTE: The overcurrent alarm will sound whenever maximum output current (500 mA) is exceeded. Should this occur, reduce the operating voltage until the overcurrent alarm goes off.

The volts/milliamps display switch will display either volts to nearest volt or milliamps to the nearest milliamp, depending on the switch setting. This switch may be changed at any time without affecting the output.

**(c) Disconnection:**

When disconnecting the power supply from the electrophoretic cell, for any reason, first turn off the power switch and confirm that the red LED is no longer illuminated.

**(3) PRECAUTIONS AND SERVICE**

Use the same precautions as with any electrical device. Do not operate without the cover in place or with any possible short circuit. Do not operate in a damp, humid environment where condensing moisture may short out internal electrical components. When moving to or from a cold room, allow at least eight (8) hours for the unit to temperature-equilibrate.

This power supply is equipped with a slo blow 1 amp fuse in the mains circuit. Should this fuse open, the check fuse LED on the front panel will illuminate. USE ONLY 1 AMP (3/4 AMP FOR 240 VAC MAINS) SLO BLOW FUSES IN THIS EQUIPMENT.

Should this product become excessively warm, smoke, or display erratic information during operation, discontinue use immediately and contact a qualified service representative.

#### **(4) CLEANING**

Use only mild, non-abrasive detergents. DO NOT use solvents or immerse in water, as this may cause permanent damage.

#### **(5) TECHNICAL SPECIFICATIONS**

##### **EC-105 Minicell Power Supply (115V/60Hz or 230V/50Hz)**

**Output mode:** Constant Voltage  
**Voltage Steps:** Lo Range: 20,30,40,50,60,70,80,90,100,110,120,130  
Hi Range: 140,150,160,170,180,190,200,210,220,230,240,250  
**Maximum Voltage:** 250 Volts  
**Maximum Current:** 300 milliAmps @ 150–250 Volts  
500 milliAmps below 150 Volts  
**Maximum Power:** 75 Watts

**Number of Outputs:** 2

**Fuse Replacement:** 1 Amp slo blow

**Temperature Range:** 0°C – 40°C (non-condensing atmosphere)

**Dimensions:** 8" (W) x 6" (D) x 4.5" (H)  
20.3 cm x 13.3 cm x 11.4 cm

**Weight: (net)** 5 lbs 2.3 kg

**Weight: (shipping)** 7 lbs 3.2 kg

***Thank You***

***for purchasing our  
EC-105 Power Supply.***

**The world's most popular power supply.**

If you should ever have any concern about the operation of this,  
or any other Thermo Electron product, please feel free to contact us at

(508) 482-7000, (800) 327-2643(U.S. toll free) or online at [www.thermo.com](http://www.thermo.com).

### GUARANTEE

This laboratory equipment was produced by Thermo Electron with the highest practical standards of materials, workmanship, and design. The design and manufacture of parts have been conceived with one purpose — to produce a unit which will give superior service.

Thermo Electron guarantees this unit to be free from defects in materials or workmanship under normal use or service for four years from date of shipment. If, during this time, this unit proves defective in materials or workmanship, the Company will repair or replace it free of charge if returned to us prepaid. This guarantee does not cover damage in transit, damage caused by carelessness, misuse or neglect, or unsatisfactory performance as a result of conditions beyond our control; or consequential losses as a result of failure of our product.



## Thermo Electron Corporation

Bioscience Technologies

450 Fortune Boulevard Milford, MA  
01757 866.9.THERMO (866.984.3766)

Fax: 508.634.2199

**Laboratory Pipetting and Consumables**  
[info.pipettes@thermo.com](mailto:info.pipettes@thermo.com)

**Molecular Biology**  
[info.molbio@thermo.com](mailto:info.molbio@thermo.com)

**Microplate Instrumentation**  
[info.microplateinstruments@thermo.co](mailto:info.microplateinstruments@thermo.co)

**Sample Preparation**  
[info.sampleprep@thermo.com](mailto:info.sampleprep@thermo.com)

**Laboratory Automation & Integration**  
[info.labautomation@thermo.com](mailto:info.labautomation@thermo.com)

**New Labs**  
[newlabs@thermo.com](mailto:newlabs@thermo.com)

**Controlled Environment**  
[info.controlenv@thermo.com](mailto:info.controlenv@thermo.com)

**Services** [services.biosciencetech@thermo.com](mailto: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 69

### 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

### Madrid

+34 9165 74930  
Fax: +34 9165 74937

### 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  
+44 870 609 9203 Fax: +44  
870 609 9202

### 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

Navi Mumbai  
+91 22 2778 1101  
Fax: +91 22 2778 1103

### Japan

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

### Finland

Vantaa  
+358 9 329 100  
Fax: +358 9 3291 0414

**Thermo**  
ELECTRON CORPORATION

© 2003 Thermo Electron Corporation. All rights reserved. Thermo Electron Corporation, and Analyze, Detect, Measure, Control are trademarks of Thermo Electron Corporation. Windows is a registered trademark of Microsoft Corporation.