Electrophoresis Power Supply

EC200XL Operating and Maintenance Manual 7217579 Rev. 0



Visit us online to register your warranty www.thermoscientific.com/warranty





MANUAL NUMBER 7217579

0		5/3/12	Transfer to Marietta (was LTM1595X2 8/28/07)	CCS
REV	ECR/ECN	DATE	DESCRIPTION	Ву



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. \blacktriangle

Caution All internal adjustments and maintenance must be performed by qualified service personnel. ▲

This equipment has been designed and tested to conform to CSA1010 safety standards, as applicable to laboratory instrumentation. This applies only to the EC200 when used as specified in the documentation, in its intended applications, with Thermo Fisher Scientific approved electrophoresis apparatus only. Usage in any other manor may not provide similar performance or safety protection.

This equipment is provided with a 3-conductor, grounded AC line cord. The protective earth ground is necessary for safe operation. Do not use any other AC line cord with this instrument.

The EC200 is a high voltage power supply capable of generating dangerous levels of voltage and current during operation. Exercise caution when working around and with the electrical connections of this equipment. Always check electrical connectors, wires, and associated apparatus for any signs of wear or damage before using with this equipment. Be sure to use only electrophoresis equipment that is suitably rated for the voltage and current capabilities of the EC200 power supply.

The output of the EC200 power supply is intended for connection to electrically isolated electrophoresis apparatus only. Use only with electrically isolated electrophoresis apparatus with minimum isolation of 600V. Do not connect any terminal of the EC200 output to earth ground. This may impair the safety protection provided by the equipment, or cause equipment damage.

The high voltage output of the EC200 power supply takes some amount of time to decay when unloaded or lightly loaded. Wait a minimum of 60 seconds after stopping a run before touching the power supply leads.

This equipment has a protective ground leakage current of 0.5mA at 120VAC, and 1mA at 230VAC. While significantly below the 3.5mA limitation for laboratory instrumentation, some applications and locations require a leakage current below 0.5mA. Check the specific requirements of your application before using this equipment.

This equipment is for indoor use only.

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

©2012 Thermo Fisher Scientific. All rights reserved.



Important operating and/or maintenance instructions. Read the accompanying text carefully.



Potential electrical hazards. Only qualified persons should perform procedures associated with this symbol.



Equipment being maintained or serviced must be turned off and locked off to prevent possible injury.



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



Marking of electrical and electronic equipment, which applies to electrical and electronic equipment falling under the Directive 2002/96/EC (WEEE) and the equipment that has been put on the market after 13 August 2005.

This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96/EC. It is marked with the WEEE symbol. Thermo Fisher Scientific has contracted with one or more recycling/disposal companies in each EU Member State European Country, and this product should be disposed of or recycled through them. Further information on Thermo's compliance with this directive, the recyclers in your country and information on Thermo products will be available at www.thermofisher.com.

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

Do You Need Information or Assistance on Thermo Scientific Products?

If you do, please contact us 8:00 a.m. to 6:00 p.m. (Eastern Time) at:

1-740-373-4763 1-800-438-4851 1-877-213-8051 http://www.thermoscientific.com service.led.marietta@thermofisher.com www.unitylabservices.com Direct Toll Free, U.S. and Canada FAX Internet Worldwide Web Home Page Tech Support Email Address Certified Service Web Page

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

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

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

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

Regardless of your needs, our professional telephone technicians are available to assist you Monday through Friday from 8:00 a.m. to 6:00 p.m. Eastern Time. Please contact us by telephone or fax. If you wish to write, our mailing address is:

> Thermo Fisher Scientific 401 Millcreek Road, Box 649 Marietta, OH 45750

International customers, please contact your local Thermo Scientific distributor.

Table of Contents

1-2 1-2
1-2 2-1
2_1
.2-1
2-2
2-3
2-3
2-4
2-4
.3-1
.4-1
4-1
4-1
4-1
4-1
-

Section 1 Introduction

Thank you for selecting a Thermo Scientific EC200 Electrophoresis power supply. This manual describes the operation of the EC200. The power supply that you have purchased is the most productive and easy-to-use unit available anywhere. This manual should answer any questions that might arise in operating your power supply; however, don't hesitate to call our Technical Services department if you need any assistance.

The EC200 power supply is designed to provide constant voltage or constant current output to apparatus used in electrophoresis applications. One to four sets of electrophoresis cells can be connected in parallel and run simultaneously. The EC200 can deliver up to 200W of total output power. When operating in constant voltage or constant current mode, the power supply automatically limits the other parameter to either the power supply maximum, or a lower limit if set by the user. If this non-constant limit is reached, the power supply will automatically switch control modes, from constant voltage to constant current, or vice versa. In this way, the EC200 protects your electrophoresis cells from damaging over power conditions. The EC200 power supply also provides for timed operation in either voltage or current modes, and allows an automatic completion in the event of a power loss if enabled by the user.

- 5-200V, adjustable in 1 volt steps
- 0.01-2.00A, adjustable in 0.01A steps
- 200W maximum output
- Automatic control mode crossover
- 0-99 hour 59 minute timed run
- Automatic restart if loss of AC power (if enabled)
- Large LCD display

Unpacking the Power Supply

When unpacking your EC200 power supply, be sure you have received the EC200 unit plus the line cord.

Inspect your equipment and packaging material for signs of damage. Damage to the shipping container may indicate rough handling which could cause internal damage to the power supply. If you suspect shipping damage to the power supply, contact your carrier for instructions on filing a claim. If you are missing any of the above items, contact your supplier for instructions.

Specifications

AC Input Power 95-265VAC, 50-60Hz, 250VA Max

Environmental

Operating temperature: 0-40°C, 0-95% R.H. non-condensing Altitude: 2000m Overvoltage category II, IEC664 Pollution degree 2, IEC664

DC Output Power

5 - 200VDC, 200W Max 0.01 - 2.00A, 200W Max Ripple: ±1% Drift: ±1%, after 30-minute warm-up

Compliance

CSA 1010

This equipment has been designed and tested to conform to CSA1010 safety standards, as applicable to laboratory instrumentation. This applies only to the EC200 when used as specified in the documentation, in its intended applications, with Thermo Scientific approved electrophoresis apparatus only. Usage in any other manner may not provide suitable protection.

Section 2 Set-Up and Operation

Select a location that allows for 3" clearance behind the power supply, and comfortable reach of the front panel controls and cell connections. Do not block the vented area of the case - on the front bottom of the unit, or the fan area at the rear. Connect the power supply to a 3-prong grounded AC outlet, using the AC cord provided with the unit only. Connect the electrophoresis apparatus to the power supply, making sure to match the red positive lead to the red positive jack, and the black negative lead to the black negative jack. Similarly, connect one or more electrophoresis apparatus to the power supply if you plan to run more than one in parallel. Power the unit on using the AC power switch located at the rear of the unit, next to the AC line cord entry.



Figure 2-1. Control Panel

Blue On/Off key

Using the Power Supply

Press the blue key on the front of the unit to enable/disable the control logic. The LCD display will illuminate and show the values of the last saved run setup. The EC200 preserves the run settings each time you start a run. One set of conditions is saved for each mode, constant voltage or constant current. The setpoint value, limit parameter, time duration, and power-fail restart settings are all saved. This allows convenient setup for repetitive runs.

To change the control mode, press the Exit key. From the resulting screen, select either constant volts or constant amps using the function keys to the right of the display.

After selecting a control mode, either constant voltage or current, and setting the limit and timed run parameters if desired (see Constant Voltage/Current Operations), press the Run key to energize the power supply output. The power supply output will ramp up to the appropriate setpoint, while not allowing the limit parameter to be exceeded. If the limit parameter should be exceeded, then the power supply will crossover control modes, making the limit parameter the new control setpoint. The run mode display shows the controlled parameter highlighted, and will automatically change if a crossover has occurred.

To stop a run in progress, press the Run key, the Stop key, or the blue ON/OFF key. The display will change to "STOP" signifying that the output is no longer energized. To pause a run in progress, press the Pause key. The display will change to "PAUSE" signifying that the output is no longer energized, and that the run can be resumed. Pressing Run or Pause resumes the run from the previous point (i.e. for a timed run). Only a run that has been paused can be restarted. After Stopping a run, or to change parameters of a (paused) run, you must cycle the control logic off and then back on by pressing the blue ON/OFF key.

During a run, the display will show the actual values of voltage and current. The time value displays elapsed time for an untimed run, or time left for a timed run.

Constant Voltage Operations

From the Mode Selection screen, select Constant Voltage operation by pressing the appropriate function key located to the right of the display. To get to the selection screen from an operating screen, press Exit, or cycle the control logic off and then back on by pressing the blue ON/OFF key. The EC200 will display the last saved values for constant voltage mode. To change a parameter, press the function key located to the right of the display, next to the parameter you wish to change. The EC200 will allow a new value to be entered using the keypad numeric keys. Press the function key labeled "Enter" to complete the new setting. Press Clear to erase previous keystrokes. Press Exit to cancel entering a new value. The EC200 will not accept a voltage setpoint outside of the allowed range. Press Clear to re-enter an appropriate value.

The limit parameter, in this case current, is normally set to the power supply maximum value of 2.00 A. The EC200 will not accept a current limit outside of the allowed range.

In addition, the EC200 will never exceed the maximum power output specification of 200W. When starting a run, the power supply calculates a maximum limit parameter, above which the power supply maximum power output rating would be exceeded. If this value is lower than the limit setpoint entered, the EC200 will automatically lower the limit parameter.

Constant Current Operation

From the mode selection screen, select constant current operation by pressing the appropriate function key located to the right of the display. To get to the selection screen from an operating screen, press Exit, or cycle the control logic off and then back on by pressing the blue ON/OFF key. The EC200 will display the last saved values for constant current mode. To change a parameter, press the function key located to the right of the display, next to the parameter you wish to change. The EC200 will allow a new value to be entered using the keypad numeric keys. Press the function key labeled "Enter" to complete the new setting. Press Clear to erase previous keystrokes. Press Exit to cancel entering a new value. The EC200 will not accept a current setpoint outside of the allowed range. Press Clear to re-enter an appropriate value.

The limit parameter, in this case voltage, is normally set to the power supply maximum value of 200 volts. The EC200 will not accept a voltage limit outside of the allowed range.

In addition, the EC200 will never exceed maximum power output specification of 200W. When starting a run, the power supply calculates a maximum limit parameter, above which power supply maximum power output rating would be exceeded. If this value is lower than the limit setpoint entered, the EC200 automatically lowers the limit parameter.

Timed Operations It is possible to enter an amount of time for the power supply to provide power, after which it will automatically shut off. Use this feature for timed runs. Enter a time parameter as described above for setting a voltage or current setpoint. Enter a time duration of 0-99 hours, 59 minutes. A time duration of zero effectively disables timed run mode. When running a timed run, the time parameter displays the time remaining in the run. When running a non-timed run, the time parameter displays the elapsed time during the run **Automatic Power-fail** Automatic power-fail restart capability allows a timed run which is interrupted by loss of AC power to be restarted automatically, so that the Restart total time programmed for the time parameter will be met. **Note** Loss of AC power includes switching off the unit using the power switch located at the rear of the unit. When performing timed runs with power-fail restart enabled, always use the front panel controls to stop a run in progress. To enable power-fail restart, press the top function key when in the operating mode screen (constant voltage or current). The state of power-fail restart will toggle from off to on or on to off with each keypress. When AC power is restored during a timed run in which power-fail restart is enabled, the display will show "POWER-FAIL RESTART" for approximately 10 seconds indicating power-fail restart pending. During this time the output is not energized, to allow safely stopping the run (by pressing Run). After the power-fail restart pending delay is completed, the power supply output will ramp up to the setpoint value and the time will continue from that point. Any number of power interruptions can occur during the completion of a single run.

Section 3 Troubleshooting

The EC200 detects and reports several events and conditions that are considered errors. The EC200 will stop any run in process and display "ERROR XX" where XX is one of the below listed errors. When the EC200 is displaying an error indication, press the blue ON/OFF key to clear the error and return to setup mode, or press the Run key to resume the run (after correcting the cause of the error). Note that loss of AC power during a run is considered an error condition.

Always stop a run before turning off AC power to the EC200.

<u>'XX'</u>

Condition and Possible Remedy

00

Minimum load current detected. The output is not connected, or the electrophoresis apparatus is not set up properly. Check your setup and connections.

01

Maximum load current exceeded. The output is short circuit, or the electrophoresis apparatus is not set up properly. Check your setup and connections

02

Step load current change. The EC200 has detected a large change in load current. Check your setup and connections

04

Loss of AC power; PF disabled and not completed. The EC200 detected loss of AC power, the run was not completed since PF was not enabled.

05

Loss of AC power; run not timed. The EC200 detected loss of AC power.

06-99

The EC200 detected an internal error condition. Contact Technical Services for instructions.

Section 4 Maintenance

Cleaning

Before cleaning the unit, be sure to always turn off AC power using the switch at the rear of the unit, and then disconnecting the AC line cord. The front of the unit is sealed and can be wiped clean with any mild detergent solution. Avoid harsh cleaners or agents as they may deteriorate the surface of the tactile membrane keys.

Replacing a Fuse

Before servicing the unit, be sure to always turn off AC power using the switch at the rear of the unit, and then disconnecting the AC line cord. The EC200 power supply requires no periodic servicing and should provide years of trouble free operation. Should you need to replace the fuses proceed as follows:

Caution The EC200 uses double pole neutral fusing. ▲

Turn off AC power by using the switch at the rear of the unit, and then disconnecting the AC line cord. Remove the fuse holder assembly with a small flatblade screwdriver.

Always replace both fuses with appropriate replacement fuses: 3A, 250V, 5X20mm, type T fuse (T3A,250V - Thermo catalog number FB-FUSE-2).

Replacement Parts

AC line cord	FB-CORD-1
Fuse, T3A,250V	FB-FUSE-2
For replacement parts, contact Technic	cal Services.

Accessories

Adapter for Cells		FBAD-1
Under Counter Bracket	FB	UB3000

The Warranty Period starts two weeks from the date your equipment is shipped from our facility. This allows for shipping time so the warranty will go
Warranty period. Warranty period. During the first year, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo's expense, labor included. Installation and calibration are not covered by this warranty agreement. The Technical Services Department must be contacted for warranty determination and direction prior to performance of any repairs. Expendable items, class, filters and paskets are excluded from this warranty
Replacement or repair of components parts or equipment under this warranty shall not extend the warranty to either the equipment or to the component or to the component part beyond the original warranty period. The Technical Services Department must give prior approval for return of any components or equipment. At Thermo's option, all non-conforming parts must be returned to Thermo Fisher Scientific postage paid and replacement parts are shipped FOB destination.
THIS WARRANTY IS EXCLUSIVE AND IN LIEU OF ALL OTHER WARRANTIES, WHETHER WRITTEN, ORAL OR IMPLIED. NO WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE SHALL APPLY. Thermo shall not be liable for any indirect or consequential damages including, without limitation, damages relating to lost profits or loss of products.
Your local Thermo Sales Office is ready to help with comprehensive site preparation information before your equipment arrives. Printed instruction man- uals carefully detail equipment installation, operation and preventive maintenance.
If equipment service is required, please call your Technical Services Department at 1-800-438-4851 (USA and Canada) or 1-740-373-4763. We're ready to answer your questions on equipment warranty, operation, maintenance, service and special application. Outside the USA, contact your local distributor for warranty information.
Rev. 4.4/09

Rep ner At ⁻ nat

EC200XL

5-2

The Warranty Period starts two months from the date your equipment is shipped from our facility. This allows for shipping Replacement or repair of components parts or equipment under this warranty shall not extend the warranty to either the approval for return of any components or equipment. At Thermo's option, all non-conforming parts must be returned to time so the warranty will go into effect at approximately the same time your equipment is delivered. The warranty protection extends to any subsequent owner during the first year warranty period. Dealers who stock our equipment are allowed During the first year, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo's expense, labor excluded. Installation and calibration are not covered by this warranty agreement. The Technical equipment or to the component part beyond the original warranty period. The Technical Services Department must give prior This warranty is exclusive and in Lieu of All other warranties, whether written, oral or Your local Thermo Sales Office is ready to help with comprehensive site preparation information before your equipment Contact your localdistributor for warranty information. We're ready to answer your questions on equipment warranty, operan additional six months for delivery and installation, provided the warranty card is completed and returned to the Technical Services Department must be contacted for warranty determination and direction prior to performance of any repairs. IMPLIED. NO WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE SHALL APPLY Thermo shall not be liable for any indirect or consequential damages including, without limitation, damages relating to lost THERMO FISHER SCIENTIFIC INTERNATIONAL DEALER WARRANTY arrives. Printed instruction manuals carefully detail equipment installation, operation and preventive maintenance. 2/09 Rev. 4 Expendable items, glass, filters, reagents, tubing, and gaskets are excluded from this warranty. Thermo postage paid and replacement parts are shipped FOB destination. ation, maintenance, service and special application profits or loss of products. Services Department.

Thermo Fisher Scientific 401 Millcreek Road Marietta, Ohio 45750 United States

www.thermofisher.com

