



PowerPac™ HV

high-voltage power supply



PROTEAN® II Ready Gel® Precast Gel System



Model 111 Mini IEF Cell



Rotofor® Cell and Mini Rotofor Cell

Plug in for great performance.

BIO-RAD

Powered Up!

The PowerPac HV high-voltage power supply supports an increased output of 5,000 V, 500 mA, or 400 W, which allows you to drive all high-voltage applications, including low-current applications in the microampere range. It is ideal for isoelectric focusing (IEF) and DNA sequencing. With 400 W output, the PowerPac HV offers enough power to run the most demanding IEF or DNA sequencing experiments.



Data Tracking via Easy-to-Use Software

With the optional PowerPac™ data transfer software, it's easy to organize, display, print, analyze, export, and annotate run data from the PowerPac HV power supply. Send run data first to a personal digital assistant (PDA) and then to a PC, or directly to a PC with a peripheral IR-receiving device. Once data are collected from the PowerPac HV, the data transfer software allows the operator to:

- View and track the programmed method
- See actual method and run conditions (voltage, current, power, method edits, error conditions, run date and time) recorded periodically during the run
- Annotate a run with notes
- Identify the serial number of the unit that generated the data
- Document all run-related information with reports
- Organize run data into projects
- Customize data presentation



IEF separation with voltage ramping.

Innovative Wireless Data Transfer

The PowerPac HV high-voltage power supply uses an integral IR port to transmit data to a PC either via a handheld PDA or directly via an IR-receiving device with the optional PowerPac data transfer software. With the IR receiver and PowerPac data transfer software, the PowerPac HV can be placed anywhere in the laboratory with respect to the PC. The recorded data include:

- The unit's unique serial number (to identify each data-generating PowerPac power supply)
- Programmed methods
- Actual run conditions recorded throughout the run
- User notes entered post-run
- A date/time stamp from the onboard real-time clock

Built for the Real World

The PowerPac HV power supply is designed for long-term reliability. Each power supply is manufactured in a state-of-the-art facility with automated testing equipment to ensure that it arrives at your laboratory ready to deliver top performance.

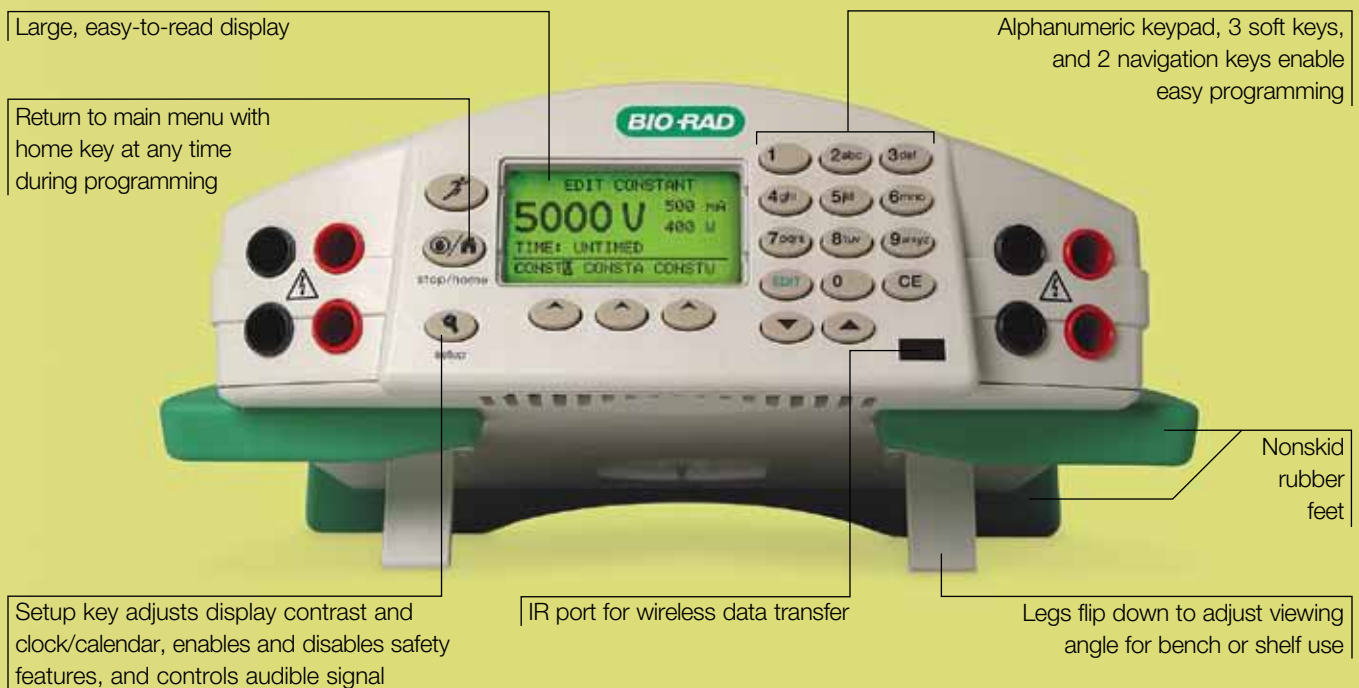
Extraordinary User Interface and Programmability

The alphanumeric keypad and graphics display combine for straightforward programming of single- or multi-step methods. The PowerPac HV user interface provides:

- Simple methods programming for starting a run
- Repeat of the previous run with a single keystroke
- Continuous, timed, or volt-hour operation
- A running screen displaying large numbers that can be seen from a distance, as well as helpful run details such as the constant variable, elapsed time, and volt-hours

A Helping Hand With IQ/OQ

To support installation qualification and operational qualification (IQ/OQ) within GLP- and FDA-regulated environments, a PowerPac IQ/OQ protocol binder and test box are available for use exclusively with the PowerPac HV and PowerPac™ Universal power supplies. The first part of the user-customizable PowerPac IQ/OQ protocol provides an installation checklist and procedure to check performance of a newly installed PowerPac HV. The second part of the protocol checks actual performance against specified performance over time. The test box houses a factory-calibrated resistive load. When used with the appropriate meters to monitor output voltage and current, the test box determines the power supply's performance against its specifications.



PowerPac HV high-voltage power supply



Specifications

Output (programmable)	
Voltage	20–5,000 V
Current	0.01–500 mA
Power	1–400 W
Type of output (with automatic crossover)	Constant voltage, constant current, constant power, or constant temperature
Timer control	1 min to 99 hr, 59 min
Volt-hour control	Yes, 99,999 V-hr
Pause/resume function	Yes
Display functions	128 x 64 pixel, yellow-green backlit graphics LCD
Programmable methods	Stores up to 9 basic and 9 IEF methods, each with up to 9 steps
Real-time editing	Yes
Real-time clock	Yes
Automatic recovery after power failure	Yes, user-selectable; setup values maintained
Data transfer/archiving	Yes
Temperature control	Yes, via temperature probe; 30–90°C ± 2°C
Microampere readout and control	Yes
Safety features	No-load detection, sudden load change detection, ground leak detection, overload/short circuit protection, overvoltage protection
Operating conditions	0–40°C, 0–95% humidity
Stackable	Yes
Number of output jacks	4 sets in parallel
Regulatory	EN-61010, CE
IQ/OQ protocols	Yes
Input power (actual)	90–120 or 198–264 VAC, 50 or 60 Hz, autoswitching
Dimensions (W x D x H)	27.5 x 34 x 10 cm
Weight	2.85 kg (6.3 lb)

Minimum Hardware Requirements for PowerPac Data Transfer Software

PC requirements	Windows XP or 2000 operating system 400 MHz processor 256 MB RAM minimum, 512 MB recommended 1,024 x 768 pixel screen resolution with true-color mode (24 or 32 bits) 6 GB hard drive CD-ROM drive IR port
PDA requirements	Palm OS software version 4.0 or 5.0 8 MB memory

Ordering Information

Catalog # Description

PowerPac HV Power Supply and Accessories, 100–120/220–240 V	
164-5056	PowerPac HV Power Supply, includes power cord, instructions
164-5059	PowerPac HV Power Supply With Temperature Probe
164-5097	PowerPac HV Data Transfer Software
164-5098	PowerPac HV IQ/OQ Protocol Binder and Test Box
164-5099	PowerPac HV IQ/OQ Protocol Binder

Palm OS is a trademark of Palm Computing, Inc. Windows and Windows 2000 and XP are trademarks of Microsoft Corp.



BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Malaysia** 60 3 2117 5260 **Mexico** 52 555 488 7670
The Netherlands 0318 540666 **New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700
Russia 7 495 721 14 04 **Singapore** 65 6415 3170 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700
Switzerland 061 717 95 55 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 6518311 **United Kingdom** 020 8328 2000