



## Electrophoresis Power Supplies

Open a World of Powerful Possibilities

**BIO-RAD**

## PowerPac™ Basic Power Supply



- Recommended for basic horizontal and vertical electrophoresis applications
- Compact, stackable (as shown)
- Constant voltage or constant current output

## PowerPac™ HC High-Current Power Supply



- Recommended for high-current applications, such as western blotting
- Output of 250 V, 3.0 A, 300 W
- 2-line, 16-character LCD for programming
- Constant voltage, constant power, or constant current output

## PowerPac™ HV High-Voltage Power Supply



- Ideal for IEF and DNA sequencing
- Output of 5,000 V, 500 mA, and 400 W
- Protocol binder and text box to support IQ/OQ within GLP- and FDA-regulated environments

## PowerPac™ Universal Power Supply



- For all applications from mini vertical and high-throughput electrophoresis to blotting
- Protocol binder and text box to support IQ/OQ within GLP- and FDA-regulated environments

## PowerPac Power Supply Specifications



	PowerPac Basic	PowerPac HC	PowerPac HV	PowerPac Universal
Output range (programmable)				
Volts	10–300 V	5–250 V	20–5,000 V	10–500 V
Current	4–400 mA	0.01–3.0 A	0.01–500 mA	0.01–2.5 A
Power	75 W (maximum)	1–300 W	1–400 W	1–500 W
Type of output (with automatic crossover)	Constant voltage or constant current	Constant voltage, constant current, or constant power	Constant voltage, constant current, constant power, or constant temperature	Constant voltage, constant current, or constant power
Timer	1–999 min	1 min–99 hr, 59 min	1 min–99 hr, 59 min	1 min–99 hr, 59 min
Volt-hour control	—	—	• (99,000 V-hr)	• (99,000 V-hr)
Pause/resume function	•	•	•	•
Display	3-digit LED	16-character x 2-line LCD	128 x 64 pixel, backlit graphics LCD	128 x 64 pixel, backlit graphics LCD
Programmable methods	—	1 method up to 3 steps, no storage capability	Stores 9 methods, each with up to 9 steps	Stores 9 methods, each with up to 9 steps
Real-time clock	—	—	•	•
Automatic recovery after power failure	•	•	•	•
Data transfer/archiving	—	—	•	• (optional)
Temperature control	—	—	Via temperature probe; 30–90°C ± 2°C	—
Microampere readout	—	—	•	—
Safety features	Overload/short-circuit detection; overvoltage protection; no-load detection; sudden load change detection	Overload/short-circuit detection; overvoltage protection; no-load detection; sudden load change detection; ground leak detection	Overload/short-circuit detection; overvoltage protection; no-load detection; sudden load change detection; ground leak detection; arc detection	Overload/short-circuit detection; overvoltage protection; no-load detection; sudden load change detection; ground leak detection
Operating conditions	0–40°C; 0–95% humidity	0–40°C; 0–95% humidity	0–40°C; 0–95% humidity	0–40°C; 0–95% humidity
Stackable	•	•	•	•
Number of output jacks	4 sets in parallel	4 sets in parallel	4 sets in parallel	4 sets in parallel
Regulatory	EN-61010, CE	EN-61010, CE	EN-61010, CE	EN-61010, CE
IQ/OQ protocols	—	—	• (optional)	• (optional)
Input power (actual)	90–120 or 198–264 VAC, 50/60 Hz, autoswitching	90–120 or 198–264 VAC, 50/60 Hz, autoswitching	90–120 or 198–264 VAC, 50/60 Hz, autoswitching	90–120 or 198–264 VAC, 50/60 Hz, autoswitching
Dimensions (W x D x H)	21.0 x 24.5 x 6.5 cm	25.0 x 28.5 x 8.0 cm	27.5 x 34.0 x 10.0 cm	27.5 x 34.0 x 10.0 cm
Weight	1.1 kg (2.4 lb)	2.0 kg (4.4 lb)	2.85 kg (6.3 lb)	2.5 kg (5.5 lb)

## Ordering Information

Catalog # Description

### Power Supplies and Adaptors

164-5050	PowerPac Basic Power Supply, 100–120/220–240 V
164-5052	PowerPac HC Power Supply, 100–120/220–240 V
164-5056	PowerPac HV Power Supply, 100–120/220–240 V
164-5059	PowerPac HV Power Supply with Temperature Probe, 100–120/220–240 V
164-5097	PowerPac Data Transfer Software, version 2.0
164-5099	PowerPac HV IQ/OQ Protocol Binder
164-5098	PowerPac HV IQ/OQ Protocol Binder and Test Box
165-5058	PowerPac Temperature Probe
164-5070	PowerPac Universal Power Supply, 100–120/220–240 V
164-5069	PowerPac Universal IQ/OQ Protocol Binder and Test Box
164-5062	PowerPac Adaptor, 2 mm
164-5064	PowerPac Adaptor, 4 mm

### Mini-PROTEAN Tetra Systems

165-8025	Mini-PROTEAN® Tetra Cell and PowerPac Basic Power Supply, includes #165-8001 and #164-5050
165-8026	Mini-PROTEAN Tetra Cell and PowerPac Universal Power Supply, includes #165-8001 and #164-5070
165-8027	Mini-PROTEAN Tetra Cell and PowerPac HC Power Supply, includes #165-8001 and #164-5052
165-8028	Mini-PROTEAN Tetra Cell and PowerPac HV Power Supply, includes #165-8001 and #164-5056
165-8033	Mini-PROTEAN Tetra Cell, Mini Trans-Blot® Module, and PowerPac Basic Power Supply includes #165-8001, #170-3935, and #164-5050
165-8034	Mini-PROTEAN Tetra Cell for Mini Precast Gels, Mini Trans-Blot Module, and PowerPac Basic Power Supply, includes #165-8004, #170-3935, and #164-5050
165-8035	Mini-PROTEAN Tetra Cell, Mini Trans-Blot Module, and PowerPac HC Power Supply, includes #165-8001, #170-3935, and #164-5052
165-8036	Mini-PROTEAN Tetra Cell for Mini Precast Gels, Mini Trans-Blot Module, and PowerPac HC Power Supply, includes #165-8004, #170-3935, and #164-5052

### Criterion Systems

165-6019	Criterion™ Cell and PowerPac Basic Power Supply, 100–120/220–240 V, includes #165-6001 and #164-5050
----------	--

### Criterion Dodeca Systems

165-4138	Criterion™ Dodeca™ Cell and PowerPac HC Power Supply, includes #165-4130 and #164-5052
165-4139	Criterion Dodeca Cell and PowerPac Universal Power Supply, includes #165-4130 and #164-5070

For additional information, please visit [www.bio-rad.com/powersupplies](http://www.bio-rad.com/powersupplies).



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://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 11 3065 7550  
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 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300  
Israel 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 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 3188 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 026 674 55 05  
Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000