



Plug In for
Great Performance.

PowerPac™ HC High-Current Power Supply



Trans-Blot® Plus Cell



Trans-Blot® SD Cell



PROTEAN® Plus Dodeca™ Cell





Amped Up

If you want to bring new power to your separations, plug into the PowerPac HC power supply. With 3.0 A of current and 300 W of power, no other power supply available today matches the output of the PowerPac HC power supply. It easily handles all your blotting applications and drives high-throughput SDS-PAGE to maximum performance with the PROTEAN plus Dodeca cells. To help you optimize every run, the PowerPac HC power supply provides fully adjustable voltage, current, and power control.

A two-line, 16-character LCD shows all parameters during running and gives you sophisticated programming options, including three-step methods.

The PowerPac HC power supply will energize your high-current applications, and with its streamlined design, stackable case, and vibrant, translucent green lid, it'll turn a lot of heads in your laboratory.

The PowerPac HC power supply is available as part of complete systems for electrophoresis. See below for available combinations.

Specifications

Output specifications	250 V, 3.0 A, 300 W
Output range (programmable)	5–250 V, fully adjustable in 1 V increments 0.01–3.0 A, fully adjustable in 0.01 A increments
Type of output	1–300 W, fully adjustable in 1 W increments Constant voltage, current, or power with automatic crossover
Output terminals	4 pair recessed banana jacks in parallel
Timer	up to 99 hr, 59 min
Pause/resume function	Yes
Display	2-line, 16-character backlit LCD
Operating conditions	0–40°C; 0–95% humidity in absence of condensation
Safety compliance	EN61010
EMI	Conforms to CE standards for emissions and immunity class A, tested only at 230 V. TÜV EMC certification
Safety features	No-load detection; rapid resistance change detection; ground leak detection; overload/short circuit detection; overvoltage protection; over-temperature protection
Input protection	Fuse on hot and neutral
Input power (nominal)	100–120/220–240 VAC, autoswitching
Stackable	Yes
Dimensions	25 x 28.5 x 8 cm (W x D x H)
Weight	2.0 kg

Ordering Information

Catalog #	Description
PowerPac HC Power Supply, 100–120/220–240 V	
164-5052	PowerPac HC Power Supply, includes power cord, instructions
165-5051	PowerPac™ Adaptor, 1 each
165-5066	PowerPac Adaptor, 2 each
Blotting Systems, 100–120/220–240 V	
170-3872	Criterion™ Blotter With Plate Electrodes and PowerPac HC Power Supply
170-3874	Criterion Blotter With Wire Electrodes and PowerPac HC Power Supply
170-3836	Mini Trans-Blot® Cell and PowerPac HC Power Supply
170-3853	Trans-Blot System With Plate Electrodes, Super Cooling Coil, and PowerPac HC Power Supply
170-3850	Trans-Blot System With Plate Electrodes and PowerPac HC Power Supply
170-3825	Trans-Blot System With Wire Electrodes and PowerPac HC Power Supply
170-3848	Trans-Blot SD System and PowerPac HC Power Supply
170-3991	Trans-Blot Plus Cell and PowerPac HC Power Supply
High-Throughput Electrophoresis Systems, 100–120/220–240 V	
165-4140	PROTEAN Plus Dodeca Cell (100/120 V) and PowerPac HC Power Supply
165-4141	PROTEAN Plus Dodeca Cell (220/240 V) and PowerPac HC Power Supply
165-4138	Criterion Dodeca Cell and PowerPac HC Power Supply



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050
Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500
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 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000