



Plug In for
Great Performance.

PowerPac™ Basic 300V Power Supply



Criterion™ Precast Gel System



Sub-Cell® GT Electrophoresis Cells



Mini-PROTEAN® 3 Electrophoresis Cell

BIO-RAD



Performance, Performance, Performance

For reliability and convenience, the PowerPac Basic power supply is your best choice. The unit is stackable and compact and has a small footprint. Its choice of constant voltage or current settings and its built-in timer with pause/resume run options let you stay focused on research and easily adjust programmed settings without having to reset the timer during an experiment.

PowerPac Basic power supply units are loaded with safety features and are available individually or as part of complete electrophoresis systems with mini vertical or submerged horizontal cells. When outstanding performance counts, the PowerPac Basic power supply drives research to a maximum. For more information, go to www.bio-rad.com/powersupplies/

Specifications

Output range (programmable)	10–300 V, fully adjustable in 1 V steps 4–400 mA, fully adjustable in 1 mA steps 75 W (maximum)
Type of output	Constant voltage or constant current with automatic crossover
Output terminals	4 pair recessed banana jacks in parallel
Timer	1–999 min, fully adjustable
Pause/resume function	Yes
Display	3-digit LED
Operating conditions	0–40°C; 0–95% humidity (noncondensing)
Regulatory	EN-61010, CE
Safety features	No-load detection; sudden load change detection, overload/short-circuit detection; overvoltage protection
Input protection	Fuse on both hot and neutral
Input power (actual)	90–120 or 198–264 VAC, 50/60 Hz, autoswitching
Dimensions (W x D x H)	21 x 24.5 x 6.5 cm
Weight	1.1 kg (2.4 lb)

Ordering Information

Catalog #	Description
164-5050	PowerPac Basic Power Supply , 100–120/220–240 V; includes power cord, instructions
165-5061	PowerPac™ Adaptor , 1
165-5066	PowerPac Adaptors , 2
Mini Vertical Electrophoresis and Blotting Systems, 100–120/220–240 V	
165-3301	Mini-PROTEAN 3 Electrophoresis Cell , 10-well, 0.75 mm thickness; complete system, includes 2 combs, 5 sets of glass plates, casting stand, casting clamp assembly, sample loading guide, electrophoresis module (electrode assembly, clamping frame, tank, lid with power cables, mini cell buffer dam)
165-3319	Mini-PROTEAN 3 Cell and PowerPac Basic Power Supply , includes 165-3301 and 164-5050
165-3323	Mini-PROTEAN 3 Electrophoresis Cell, Mini Trans-Blot® Module, and PowerPac Basic Power Supply
165-3324	Mini-PROTEAN 3 Electrophoresis Module, Mini Trans-Blot Module, and PowerPac Basic Power Supply
165-6019	Criterion Cell and PowerPac Basic Power Supply
170-3989	Mini Trans-Blot Cell and PowerPac Basic Power Supply
Submerged Horizontal Electrophoresis Systems, 100–120/220–240 V	
164-0300	Mini-Sub® Cell GT Cell and PowerPac Basic Power Supply
164-0301	Wide Mini-Sub Cell GT Cell and PowerPac Basic Power Supply
164-0302	Sub-Cell GT Cell and PowerPac Basic Power Supply
164-0303	Mini ReadySub-Cell™ GT Cell and PowerPac Basic Power Supply
164-0304	Wide Mini ReadySub-Cell GT Cell and PowerPac Basic Power Supply
164-0305	Sub-Cell Model 96 Cell and PowerPac Basic Power Supply
164-0306	Sub-Cell Model 192 Cell and PowerPac Basic Power Supply

Bright 3-digit LED display

Lights monitor output

Pause/resume capability



Easy-to-use rubber keys

Nonskid rubber feet

Legs flip down to adjust viewing angle for bench or shelf use

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
Canada 905 712 2771 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 4029300/5013478 Israel 03 963 6050
Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 55 5200 05 20 The Netherlands 0318 540666 New Zealand 64 9415 2280
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 095 721 14 04 Singapore 65 6415 3188 South Africa 27 0861 246 723
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189/2578 7241 United Kingdom 020 8328 2000