



Plug In for  
Great Performance.

# PowerPac™ Basic 300V Power Supply



Criterion™ Precast Gel System



Sub-Cell® GT Electrophoresis Cells



Mini-PROTEAN® 3 Electrophoresis Cell





## 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/](http://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	<b>PowerPac Basic Power Supply</b> , 100–120/220–240 V; includes power cord, instructions
165-5061	<b>PowerPac™ Adaptor</b> , 1
165-5066	<b>PowerPac Adaptors</b> , 2
<b>Mini Vertical Electrophoresis and Blotting Systems, 100–120/220–240 V</b>	
165-3301	<b>Mini-PROTEAN 3 Electrophoresis Cell</b> , 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	<b>Mini-PROTEAN 3 Cell and PowerPac Basic Power Supply</b> , includes 165-3301 and 164-5050
165-3323	<b>Mini-PROTEAN 3 Electrophoresis Cell, Mini Trans-Blot® Module, and PowerPac Basic Power Supply</b>
165-3324	<b>Mini-PROTEAN 3 Electrophoresis Module, Mini Trans-Blot Module, and PowerPac Basic Power Supply</b>
165-6019	<b>Criterion Cell and PowerPac Basic Power Supply</b>
170-3989	<b>Mini Trans-Blot Cell and PowerPac Basic Power Supply</b>
<b>Submerged Horizontal Electrophoresis Systems, 100–120/220–240 V</b>	
164-0300	<b>Mini-Sub® Cell GT Cell and PowerPac Basic Power Supply</b>
164-0301	<b>Wide Mini-Sub Cell GT Cell and PowerPac Basic Power Supply</b>
164-0302	<b>Sub-Cell GT Cell and PowerPac Basic Power Supply</b>
164-0303	<b>Mini ReadySub-Cell™ GT Cell and PowerPac Basic Power Supply</b>
164-0304	<b>Wide Mini ReadySub-Cell GT Cell and PowerPac Basic Power Supply</b>
164-0305	<b>Sub-Cell Model 96 Cell and PowerPac Basic Power Supply</b>
164-0306	<b>Sub-Cell Model 192 Cell and PowerPac Basic Power Supply</b>

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](http://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