

Horizontal Electrophoresis



Bulletin 6186

Sub-Cell[®] GT Complete Systems

Catalog #	Casting Gates	Gel Caster	UVTP Tray (cm)				PowerPac Basic Power Supply
			15 x 10	15 x 15*	15 x 20	15 x 25	
170-4401			•				
170-4402	•			•			
170-4403					•		
170-4404						•	
170-4481		•	•				
170-4482	•	•		•			
170-4483		•			•		
170-4484		•				•	
164-0302	•	•			•		•

* 15 x 15 cm tray allows casting gels in the cell using casting gates. Other gel sizes can be cast with a gel caster.





**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 11 5044 5699
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** 03 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** 061 717 95 55
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000