

# Horizontal Electrophoresis

Protocol

Bulletin 6189

## Sub-Cell<sup>®</sup> Model 96 Comb Selection Guide

Catalog #	# of Wells	Height	Thickness, mm	Width of Well, mm	Length of Teeth, mm	Volume, $\mu$ l (in 5 mm deep gel)
<b>Adjustable-Height Combs for Sub-Cell Model 96 Systems</b>						
170-4528	51	—	0.75	3	15	11.25
170-4529	51	—	1.5	3	15	22.5
170-4526	26	—	0.75	6	15	22.5
170-4527	26	—	1.5	6	15	45
170-4530	2 or 4 preparative 2 reference	—	0.75	46 or 98 6	15	172.5 or 367.5 22.5
170-4531	2 or 4 preparative 2 reference	—	1.5	46 or 98 6	15	345 or 735 45

**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 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