

Mini-Sub[®] Cell Comb Selection Guide

Catalog #	# of Wells	Height*	Thickness, mm	Width of Well, mm	Length of Teeth, mm	Volume, μ l (in 5 mm deep gel)
170-4464	15	Fixed	0.75	3	10	11.25
170-4465**	15	Fixed	1.5	3	10	22.5
170-4332	15	Adjustable	1.0	2.6	10.2	13
170-4462	8	Fixed	0.75	5.5	11	20
170-4463**	8	Fixed	1.5	5.5	11	40
170-4333	8	Adjustable	1.0	5.5	10.2	27.5
170-4461	2 preparative 2 reference	Fixed	1.5	20 4	10	150 30
170-4460	1 preparative 2 reference	Fixed	1.5	44 3	10	330 22.5
170-4342	1 preparative 2 reference	Adjustable	3	43.3 3	10.2	645 45

* Fixed-height combs must be used with Mini-Sub cell GT system gel trays. Adjustable-height combs require comb holder, catalog #170-4331.

** Combs included in systems.



**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