

Sub-Cell[®] GT 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)
Combs for Sub-Cell GT Systems						
170-4449	30	Fixed	1.5	2.7	14	20.25
170-4344	30	Adjustable	1.5	2.7	19.1	20.25
170-4447	20	Fixed	0.75	5	15	18.75
170-4448	20	Fixed	1.5	5	15	36.3
170-4321	20	Adjustable	0.75	4.8	19.1	17
170-4322	20	Adjustable	1.5	4.8	19.1	34
170-4445	15	Fixed	0.75	6	15	20.7
170-4446	15	Fixed	1.5	6	15	45
170-4323	15	Adjustable	0.75	5.5	19.1	21
170-4324	15	Adjustable	1.5	5.5	19.1	42
170-4443	10	Fixed	0.75	10	14	37.5
170-4444	10	Fixed	1.5	10	14	75
170-4325	10	Adjustable	0.75	9.9	19.1	37
170-4326	10	Adjustable	1.5	9.9	19.1	74
Preparative Combs for Sub-Cell GT Systems						
170-4442	4 preparative 2 reference	Fixed	1.5	27 3	14	202 22
170-4441	2 preparative 2 reference	Adjustable	1.5	50 4	14	375 30
170-4440	1 preparative 2 reference	Fixed	1.5	106 4	20	795 30
170-4328	1 preparative 2 reference	Adjustable	3	106 4	19.1	1,590 60
Multichannel Pipet-Compatible Combs for Sub-Cell GT Systems						
170-4456	26	Fixed	0.75	3	15	11.25
170-4457	26	Fixed	1.5	3	15	22.5
170-4454	18	Fixed	0.75	3	14	11.25
170-4455	18	Fixed	1.5	3	14	22.5
170-4452	14	Fixed	0.75	6	20	22.5
170-4453	14	Fixed	1.5	6	20	45
170-4450	10	Fixed	0.75	6	14	22.5
170-4451	10	Fixed	1.5	6	14	45



**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
