


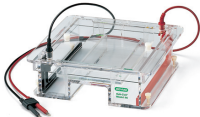
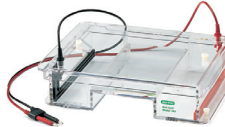


Horizontal Electrophoresis

Protocol

Bulletin 6181

Sub-Cell Selection Guide

	Mini-Sub Cell GT*	Wide Mini-Sub Cell GT**	Sub-Cell GT	Sub-Cell Model 96	Sub-Cell Model 192
					
Cell size (W × L × H)	9.2 × 25.5 × 5.6 cm	17.8 × 25.5 × 6.8 cm	18 × 40.5 × 9.4 cm	29 × 30 × 9 cm	29 × 40 × 9 cm
Gel tray sizes (outside dimensions, W × L)	7 × 7 cm 7 × 10 cm	15 × 7 cm 15 × 10 cm	15 × 10 cm 15 × 15 cm 15 × 20 cm 15 × 25 cm	25 × 10 cm 25 × 15 cm	25 × 10 cm 25 × 15 cm 25 × 20 cm 25 × 25 cm
ReadyAgarose™ gels accommodated	Yes (mini format 8-, 12-, 2 × 8-well)	Yes (wide mini and 96 Plus formats)	No	No	No
Sample throughput	8–30***	10–60***	1–120†	24–96***	24–192†
Base buffer volume	~270 ml	~650 ml	~1 L	~2 L	~3 L
Buffer recirculation	No	No	No	Yes	Yes
Bromophenol blue migration	~4.5 cm/hr (at 75 V)	~4.5 cm/hr (at 75 V)	~3.0 cm/hr (at 75 V)	~6.2 cm/hr (at 200 V)	~5.2 cm/hr (at 200 V)

* The mini ReadySub-Cell™ GT cell is a Mini-Sub cell GT cell dedicated to running ReadyAgarose precast gels, gel size 7 × 10 cm; sample throughput is 8, 12, or 2 × 8. This cell does not include casting gates, tray, or combs.

** The wide mini ReadySub-Cell GT cell is a wide Mini-Sub cell GT cell dedicated to running ReadyAgarose precast gels, gel size 15 × 10 cm; sample throughput is 20, 32, 2 × 32, or 4 × 26. This cell does not include casting gates, tray, or combs.

*** Sample throughput value assumes 1–2 combs per gel.

† Sample throughput value assumes 1–4 combs per gel.

BIO-RAD



**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
