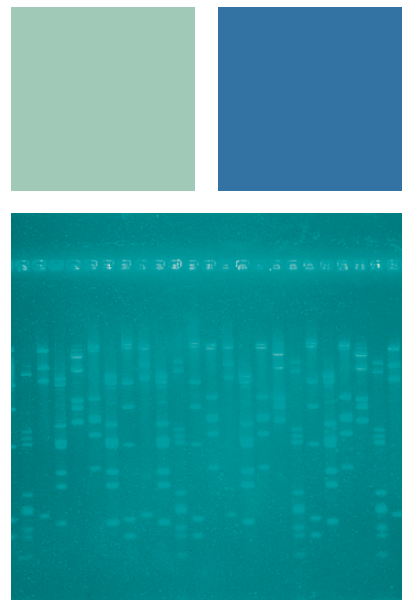
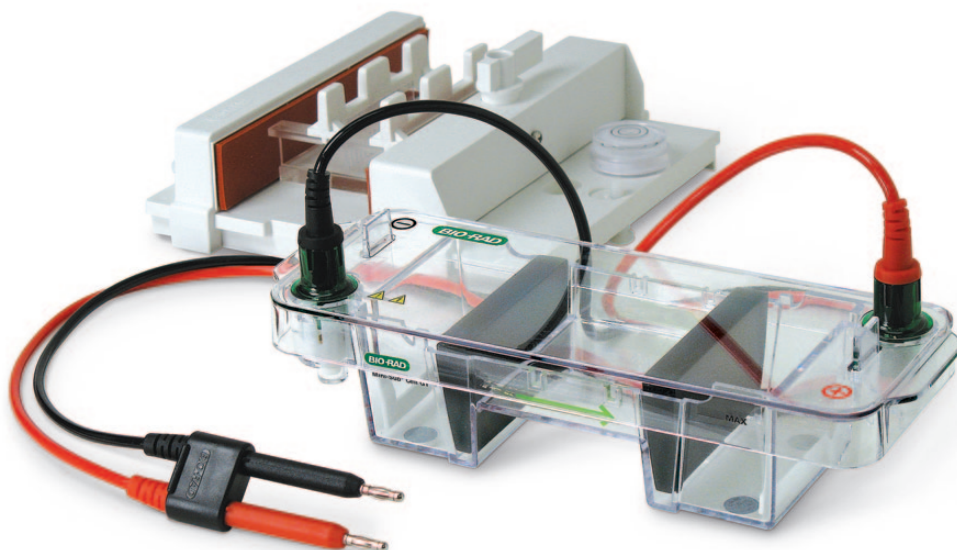


Submerged Horizontal Electrophoresis: Sub-Cell Systems



New and Improved Sub-Cell® Systems

Whether you are running precious PCR samples, or performing RFLP analysis or any other application requiring nucleic acid separation, Bio-Rad has a solution. The Mini-Sub® cell GT, wide Mini-Sub cell GT, and Sub-Cell GT cell models have been redesigned for even easier electrophoresis and more intuitive experiment setup.

- **Easy to clean** — patent-pending QuickSnap electrodes are easy to remove, simplifying cleaning
- **Leakproof** — new electrode design prevents buffer leaks from the base
- **Flexible design** — UV-transparent gel trays, combs, and other accessories are compatible with both new and old models
- **Easy lid removal** — longer tabs on the base prevent incorrect lid positioning and also enable easy removal of the lid, reducing buffer spillage
- **Intuitive setup** — a printed arrow on the side of the base indicates the direction of the run and ensures proper gel orientation
- **Gel casting** — multiple options are available for hand casting of different gel sizes
- **Time savings and gel reproducibility** — the new Mini-Sub cell GT and wide Mini-Sub cell GT cells accommodate ReadyAgarose™ precast gels

Bio-Rad is your one-stop shop for everything electrophoresis.
For more information, visit us at www.bio-rad.com/horizontalelectro.

BIO-RAD

Sub-Cell System Selection Guide

	New, Convenient Sub-Cell Systems			High-Throughput Sub-Cell Models	
	Mini-Sub Cell GT System	Wide Mini-Sub Cell GT System	Sub-Cell GT System	Sub-Cell Model 96 Cell	Sub-Cell Model 192 Cell
Gel tray sizes, cm (W x L)	7 x 7 7 x 10	15 x 7 15 x 10	15 x 10 15 x 15 15 x 20 15 x 25	25 x 10 25 x 15	25 x 10 25 x 15 25 x 20 25 x 25
ReadyAgarose gels accommodated	Yes 8-, 12-, 2 x 8-well	Yes 20-, 32-, 2 x 32-, 4 x 26-well	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
Cell size, cm (W x L x H)	9.2 x 25.5 x 5.6	17.8 x 25.5 x 6.8	18 x 40.5 x 9.4	29 x 30 x 9	29 x 40 x 9
Bromophenol blue dye migration rate	~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)

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

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

Ordering Information

Catalog # Description

Sub-Cell GT Systems*

170-4406	Mini-Sub Cell GT System, with 7 x 7 cm tray
170-4466	Mini-Sub Cell GT System, with 7 x 10 cm tray
170-4486	Mini-Sub Cell GT System, with 7 x 7 cm tray, mini-gel caster
170-4467	Mini-Sub Cell GT System, with 7 x 10 cm tray, mini-gel caster
170-4487	Mini ReadySub-Cell™ GT System, includes Sub-Cell GT tank, lid, leveling bubble
170-4405	Wide Mini-Sub Cell GT System, with 15 x 7 cm tray
170-4468	Wide Mini-Sub Cell GT System, with 15 x 10 cm tray
170-4485	Wide Mini-Sub Cell GT System, with 15 x 7 cm tray, mini-gel caster
170-4469	Wide Mini-Sub Cell GT System, with 15 x 10 cm tray, mini-gel caster
170-4489	Wide Mini ReadySub-Cell GT System, includes Sub-Cell GT tank, lid, leveling bubble
170-4401	Sub-Cell GT System, with 15 x 10 cm tray
170-4402	Sub-Cell GT System, with 15 x 15 cm tray
170-4403	Sub-Cell GT System, with 15 x 20 cm tray
170-4404	Sub-Cell GT System, with 15 x 25 cm tray
170-4481	Sub-Cell GT System, with 15 x 10 cm tray, gel caster

170-4482	Sub-Cell GT System, with 15 x 15 cm tray, gel caster
170-4483	Sub-Cell GT System, with 15 x 20 cm tray, gel caster
170-4484	Sub-Cell GT System, with 15 x 25 cm tray, gel caster

Sub-Cell GT/PowerPac™ Basic Power Supply Systems**

164-0300	Mini-Sub Cell GT/PowerPac Basic System, 100/120 V
164-0301	Wide Mini-Sub Cell GT/PowerPac Basic System, 100/120 V
164-0302	Sub-Cell GT/PowerPac Basic System, 100/120 V

Upgrade Kits***

170-4491	Mini Handcasting Kit
170-4497	Wide Mini Handcasting Kit

* All Sub-Cell GT systems include a buffer tank, a safety lid with cables, a leveling bubble, casting gates, and two 1.5 mm thick fixed-height combs: 15- and 20-well combs for Sub-Cell and wide Mini-Sub cell, 8- and 15-well combs for Mini-Sub cell.

** All Sub-Cell GT/PowerPac Basic systems include the same accessories listed above and a gel caster.

*** Handcasting kits can be purchased to upgrade ReadySub-Cell systems to Mini-Sub or wide Mini-Sub cell systems. The kits include casting gates and 1.5 mm thick 8- and 15-well combs for mini ReadySub-Cell models or 15- and 20-well combs for wide mini ReadySub-Cell models.

Practice of the polymerase chain reaction (PCR) may require a license.



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 31 3689 6600 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 30 210 777 4396 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 Malaysia 60 3 2117 5260 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 3170 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 66 2 6518311 United Kingdom 020 8328 2000 Vietnam 84 8 38131140