

# It All Fits



Aurum™ Plasmid 96 Kit

## Vector Isolation

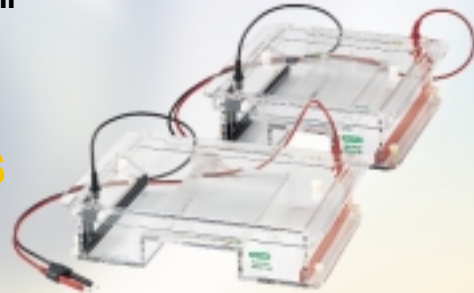


Sub-Cell® GT Cell



iCycler™ Thermal Cycler

## PCR\* Screening



Sub-Cell Model 96 and 192 Cells

## Gel-Based Analysis



Mini-Sub® and Wide Mini-Sub Cells

## Quantitation



SmartSpec™ 3000 Spectrophotometer

## Clone Identification



Gene Pulser® II Electroporator

Performing Nucleic Acid Electrophoresis?  
We Have the Cell for You.

Bio-Rad's extensive line of products and reagents meets all your needs. Whether you are screening amplified DNA, identifying clones, or using precast agarose gels for instant results, we offer the ideal Sub-Cell electrophoresis system to complete the job.

**BIO-RAD**

## Ordering Information

Catalog #	Description
<b>Sub-Cell GT Systems</b>	
170-4406	Mini-Sub Cell GT System, with 7 x 7 cm tray (tray included fits gates for casting gels in electrophoresis cell)
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 (tray included fits gates for casting gels in electrophoresis cell)
170-4467	Mini-Sub Cell GT System, with 7 x 10 cm tray, mini-gel caster
170-4405	Wide Mini-Sub Cell GT System, with 15 x 7 cm tray (tray included fits gates for casting gels in electrophoresis cell)
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 (tray included fits gates for casting gels in electrophoresis cell)
170-4469	Wide Mini-Sub Cell GT System, with 15 x 10 cm tray, mini-gel caster
170-4401	Sub-Cell GT System, with 15 x 10 cm tray
170-4402	Sub-Cell GT System, with 15 x 15 cm tray (tray included fits gates for casting gels in electrophoresis cell)
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 (tray included fits gates for casting gels in electrophoresis cell)
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

Each system includes Sub-Cell unit, UV-transparent gel tray, casting gates, two 1.5 mm fixed-height combs (15- and 20-well for the Sub-Cell and wide Mini-Sub cell, 8- and 15-well for the Mini-Sub cell).

### ReadySub-Cell™ GT Cells

170-4487	Mini ReadySub-Cell GT Cell, includes buffer tank, lid and electrodes, leveling bubble; accommodates 8-well and 12-well mini ReadyAgarose™ gels
170-4489	Wide Mini ReadySub-Cell GT Cell, includes buffer tank, lid and electrodes, leveling bubble; accommodates 20-well, 32-well, and 2 x 32-well wide mini ReadyAgarose gels

### Sub-Cell Model 96 and 192 Systems

170-4502	Sub-Cell Model 96 System, with 25 x 10 cm tray
170-4503	Sub-Cell Model 96 System, with 25 x 15 cm tray
170-4500	Sub-Cell Model 96 System, with 25 x 10 cm tray, gel caster
170-4501	Sub-Cell Model 96 System, with 25 x 15 cm tray, gel caster
170-4508	Sub-Cell Model 192 System, with 25 x 10 cm tray
170-4509	Sub-Cell Model 192 System, with 25 x 15 cm tray
170-4510	Sub-Cell Model 192 System, with 25 x 20 cm tray

### Sub-Cell Model 96 and 192 Systems (continued)

170-4511	Sub-Cell Model 192 System, with 25 x 25 cm tray
170-4504	Sub-Cell Model 192 System, with 25 x 10 cm tray, gel caster
170-4505	Sub-Cell Model 192 System, with 25 x 15 cm tray, gel caster
170-4506	Sub-Cell Model 192 System, with 25 x 20 cm tray, gel caster
170-4507	Sub-Cell Model 192 System, with 25 x 25 cm tray, gel caster

Each system includes Sub-Cell unit, UV-transparent gel tray, two 1.5 mm combs (26- and 51-well), comb holder, casting gates. For more information on power supplies compatible with Sub-Cells, visit us on the Web at [www.discover.bio-rad.com](http://www.discover.bio-rad.com)

### Plasmid Purification Systems

732-6400	Aurum Plasmid Mini Kit,** 100 preps
732-6440	Aurum Plasmid Midi Kit,** 20 preps
732-6450	Aurum Plasmid Maxi Kit,** 10 preps
732-6460	Aurum Plasmid 96 Kit, 2 x 96-well preps

### Quantum Prep® Plasmid Miniprep Kits

732-6100	Quantum Prep Plasmid Miniprep Kit, 100 preps
732-6120	Quantum Prep Midiprep Kit, 20 preps

### Gene Transfer Products

165-2105	Gene Pulser II Main Unit, 100/120 V, 50/60 Hz, includes shocking chamber, 15 sterile sample cuvettes (5 each of 0.1, 0.2, and 0.4 cm gap), cuvette rack
165-2106	Gene Pulser II Main Unit, 220/240 V, 50/60 Hz
165-2111	Gene Pulser II Shocking Chamber
165-2100	MicroPulser™ Electroporator, universal voltage, includes chamber with leads, 10 sterile cuvettes (5 each of 0.1 and 0.2 cm gap)
165-2088	Gene Pulser/MicroPulser Cuvette, 0.4 cm gap, 50, sterile
165-2086	Gene Pulser/MicroPulser Cuvette, 0.2 cm gap, 50, sterile
165-2089	Gene Pulser/MicroPulser Cuvette, 0.1 cm gap, 50, sterile

### Quantitation Products

170-2501	SmartSpec 3000 Spectrophotometer
170-2402	VersaFluor™ Fluorometer, 100/120/220 V, includes standard cuvette holder, 100 standard cuvettes, excitation filter and emission filter

### Amplification Products

170-8720	iCycler Thermal Cycler with 96 x 0.2 ml Reaction Module, includes iCycler base with 96 x 0.2 ml iCycler reaction module, in-sample temperature probe (0.2 ml tube size), PCR tubes, power cord, quick-reference card, instructions
170-8703	iCycler Thermal Cycler 96 x 0.2 ml Reaction Module, includes in-sample temperature probe (0.2 ml tube size)
170-8707	iCycler Thermal Cycler 60 x 0.5 ml Reaction Module, includes in-sample temperature probe (0.5 ml tube size)

### ReadyAgarose Mini Gels

### ReadyAgarose Wide Mini Gels

TBE Gels	8-well (20 µl)	12-well (10 µl)	2 x 8-well (20 µl)	20-well (20 µl)	32-well** (15 µl)	2 x 32-well** (15 µl)
0.8%	161-3001	161-3007	—	161-3025	161-3031	—
0.8% plus ethidium bromide	161-3002	161-3008	—	161-3026	161-3032	—
1%	161-3003	161-3009	—	161-3027	161-3033	161-3037
1% plus ethidium bromide	161-3004	161-3010	—	161-3028	161-3034	161-3038
3%	161-3005	161-3011	—	161-3029	161-3035	161-3039
3% plus ethidium bromide	161-3006	161-3012	—	161-3030	161-3036	161-3040
<b>TAE Gels</b>						
0.8%	161-3013	161-3019	—	161-3041	161-3047	—
0.8% plus ethidium bromide	161-3014	161-3020	—	161-3042	161-3048	—
1%	161-3015	161-3021	161-3057	161-3043	161-3049	161-3053
1% plus ethidium bromide	161-3016	161-3022	—	161-3044	161-3050	161-3054
3%	161-3017	161-3023	—	161-3045	161-3051	161-3055
3% plus ethidium bromide	161-3018	161-3024	—	161-3046	161-3052	161-3056

Gel shelf life at room temperature: 12 months without ethidium bromide, 6 months with ethidium bromide.

\* The polymerase chain reaction (PCR) process is covered by patents owned by Hoffman-LaRoche. Use of the PCR process requires a license. The iCycler system includes a licensed thermal cycler and may be used with PCR licenses available from Applied. Its use with authorized reagents also provides a limited PCR license in accordance with the label rights accompanying such reagents. Some applications may require licenses from other parties.

\*\* Check for availability.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191  
Canada (905) 712-2771 China (86-21) 63052255 Denmark 45 44 52-1000 Finland 358 (0)9 804 2200 France 01 47 95 69 65 Germany 089 318 84-177  
Hong Kong 852-2789-3300 India (91-124) 6398112/113/114 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460  
Latin America 305-894-5950 Mexico 52 5 534 2552 to 54 The Netherlands 0318-540666 New Zealand 64-9-4152280 Norway 47-23-38-41-30  
Russia 7 095 721 1404 Singapore 65-2729877 Spain 34 91 661 7085 Sweden 46 (0)8-55 51 27 00 Switzerland 061-717-9555 United Kingdom 0800-181134