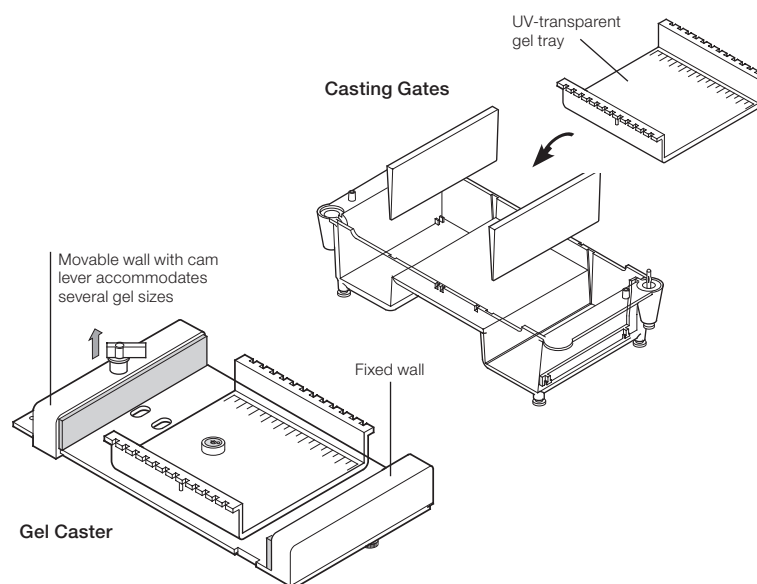


Hand Casting Options for a Variety of Gel Sizes

Two accessories are available for hand casting gels with the Sub-Cell family of electrophoresis cells. Gel casters offer the flexibility of accommodating several different-size UV-transparent (UVTP) gel trays and are available for the entire Sub-Cell product line. Casting gates offer the convenience of casting directly in the electrophoresis cells and are available in one specific size for each cell (see Casting Guide below).



Gel Caster Guide

| Tray Size | Mini Gel Caster | Full-Size Gel Caster | Model 96 Gel Caster | Model 192 Gel Caster |
|------------|-----------------|----------------------|---------------------|----------------------|
| 7 × 7 cm | • | • | • | • |
| 7 × 10 cm | • | • | • | • |
| 15 × 7 cm | • | • | • | • |
| 15 × 10 cm | • | • | • | • |
| 15 × 15 cm | | • | • | • |
| 15 × 20 cm | | • | • | • |
| 15 × 25 cm | | • | • | • |
| 25 × 10 cm | | | • | • |
| 25 × 15 cm | | | • | • |
| 25 × 20 cm | | | | • |
| 25 × 25 cm | | | | • |

Casting Guide

| Electrophoresis Cell | Tray Size | Casting Gates | Gel Caster |
|-----------------------|------------|---------------|------------|
| Mini-Sub cell GT | 7 × 7 cm | • | • |
| | 7 × 10 cm | | • |
| Wide Mini-Sub cell GT | 15 × 7 cm | • | • |
| | 15 × 10 cm | | • |
| Sub-Cell GT | 15 × 10 cm | | • |
| | 15 × 15 cm | • | • |
| | 15 × 20 cm | | • |
| | 15 × 25 cm | | • |
| Sub-Cell Model 96 | 25 × 10 cm | • | • |
| | 25 × 15 cm | | • |
| Sub-Cell Model 192 | 25 × 10 cm | | • |
| | 25 × 15 cm | • | • |
| | 25 × 20 cm | | • |
| | 25 × 25 cm | | • |



**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
