

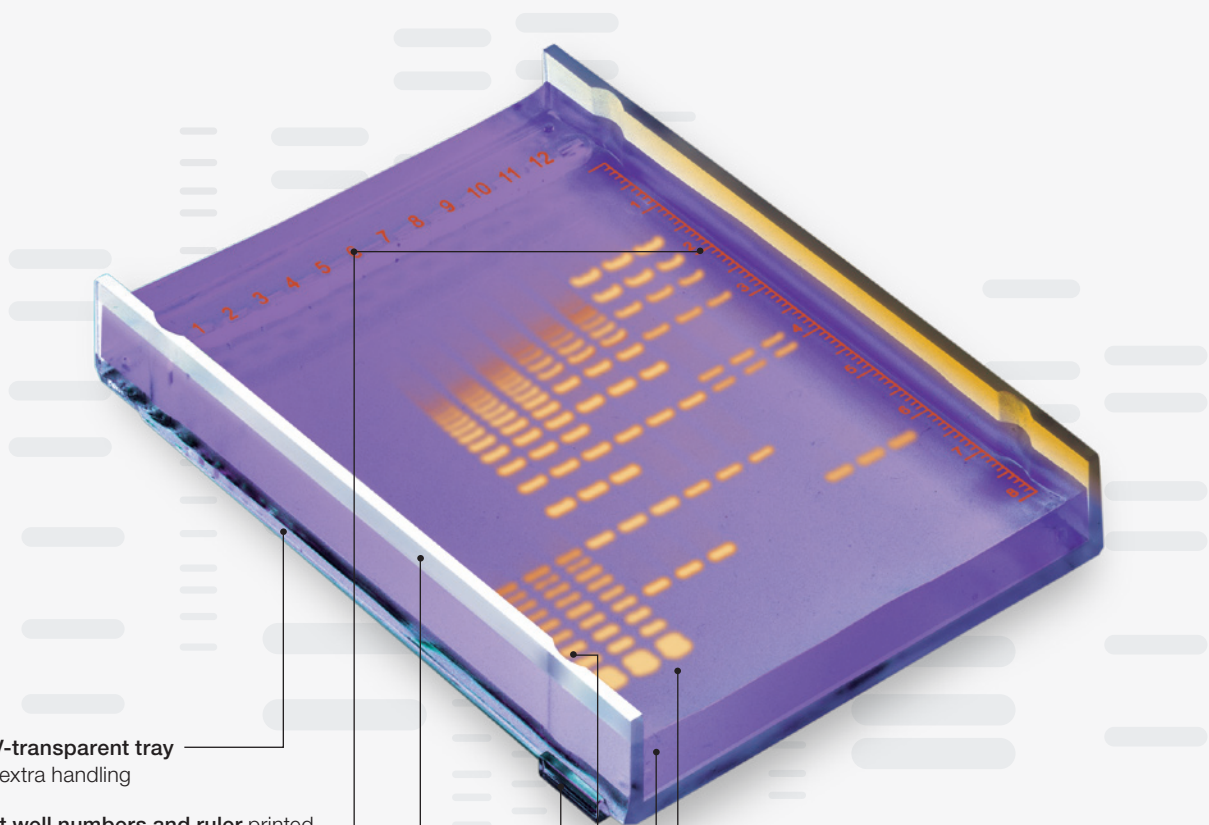
ReadyAgarose Precast Gels

Exemplary Performance. Diverse Formats.



Reload with ReadyAgarose Precast Gels

ReadyAgarose Precast Gels are made with Bio-Rad Certified Agarose and prepared in ultraviolet (UV)-transparent gel trays designed to fit securely in Mini-Sub Cell GT and Wide Mini-Sub Cell GT Systems as well as other manufacturers' electrophoresis cells. The UV-transparent trays are clearly marked with well numbers and a ruler for easy visual reference. The individually packaged gels come in a variety of formats, which resolve nucleic acids from 20 to 10,000 bp.



Cast in a **UV-transparent tray** to eliminate extra handling

Fluorescent well numbers and ruler printed on tray for easy reference

Two-sided tray for consistently straight lanes and familiar run conditions

Lock-tabs on patented tray prevent tray movement during loading and running

Supplied in **TBE** (Tris/Boric Acid/EDTA) or **TAE** (Tris/Acetic Acid/EDTA) Buffer, with or without ethidium bromide (EtBr)

Cast with **high-strength, high-resolution agaroses** that are qualified for molecular biology applications

Available in **1% and 3% agarose** to separate DNA from 20 to 10,000 bp

Notched tray holds gel in place

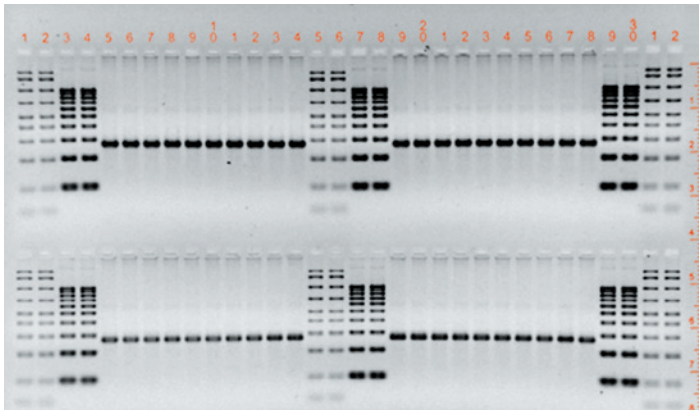
Offered in **mini and wide mini gel formats** with 8, 12, 2 x 8, 20, 32, 2 x 32, or 4 x 26 wells for all loading volume needs

Gel Selection Guide

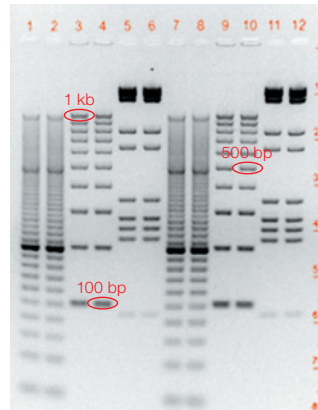
Gel, %	Resolution Range, bp
1	200–10,000
3*	20–1,000

*Our 3% gels are made with high-strength, high-resolution agarose, making them easier to handle while providing the same resolution as a standard 4% agarose gel.

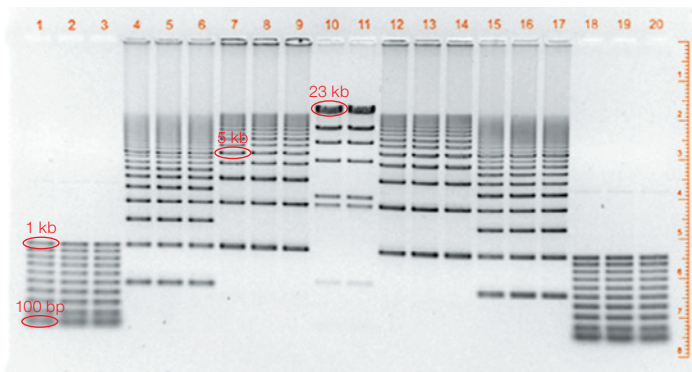
Obtain Publication-Quality Results



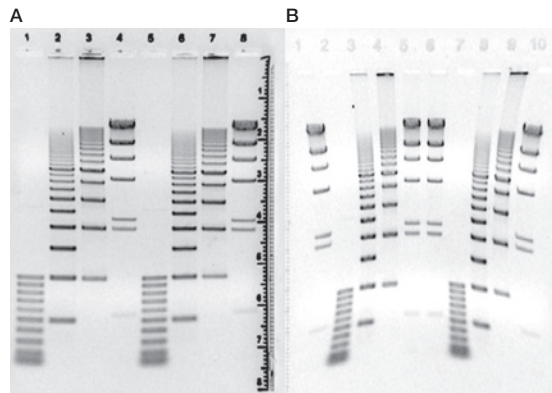
High-throughput capacity. A 3% Wide Mini Ready Agarose Precast Gel was repeatedly loaded with an Amplisize Molecular Ruler, an EZ Load 100 bp Molecular Ruler, and PCR samples containing a 280 bp product. The gel was run at 100 V for 40 min in TAE Buffer.



High resolution. A 3% Mini Ready Agarose Precast Gel was loaded with an EZ Load 20 bp Molecular Ruler (lanes 1, 2, 7, and 8), an EZ Load 100 bp Molecular Ruler (lanes 3, 4, 9, and 10), and Avall/EcoRI-digested pBR322 (lanes 5, 6, 11, and 12). The gel was run at 100 V for 70 min in TAE Buffer.



Routine broad-range separations. A 1% TAE Wide Mini Ready Agarose Precast Gel was loaded with multiple replicates of 100 bp, 500 bp, and 1 kb molecular rulers and HindIII-digested lambda DNA. The gel was run at 75 V for 80 min in TAE Buffer.

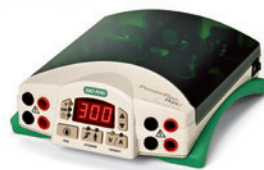


Distortion-free electrophoresis. **A**, a 1% TAE Ready Agarose Precast Gel in its 2-sided tray; **B**, a precast gel in a 4-sided tray. Only the 4-sided tray distorts bands during electrophoresis. Both gels were loaded with duplicate samples and run at 74 V for 80 min until the bromophenol blue ran 5.2 cm from the wells. Gels were stained with EtBr.



The **Sub-Cell GT System** makes running agarose gels simple and efficient, with a flexible design that offers optimal performance with handcast or precast gels.

Visit bio-rad.com/SubCells for more information.



Power your Sub-Cell System with the **PowerPac Basic Power Supply**. The unit is lightweight and stackable, and the vertical display is clearly visible on shelves and lab benches alike.

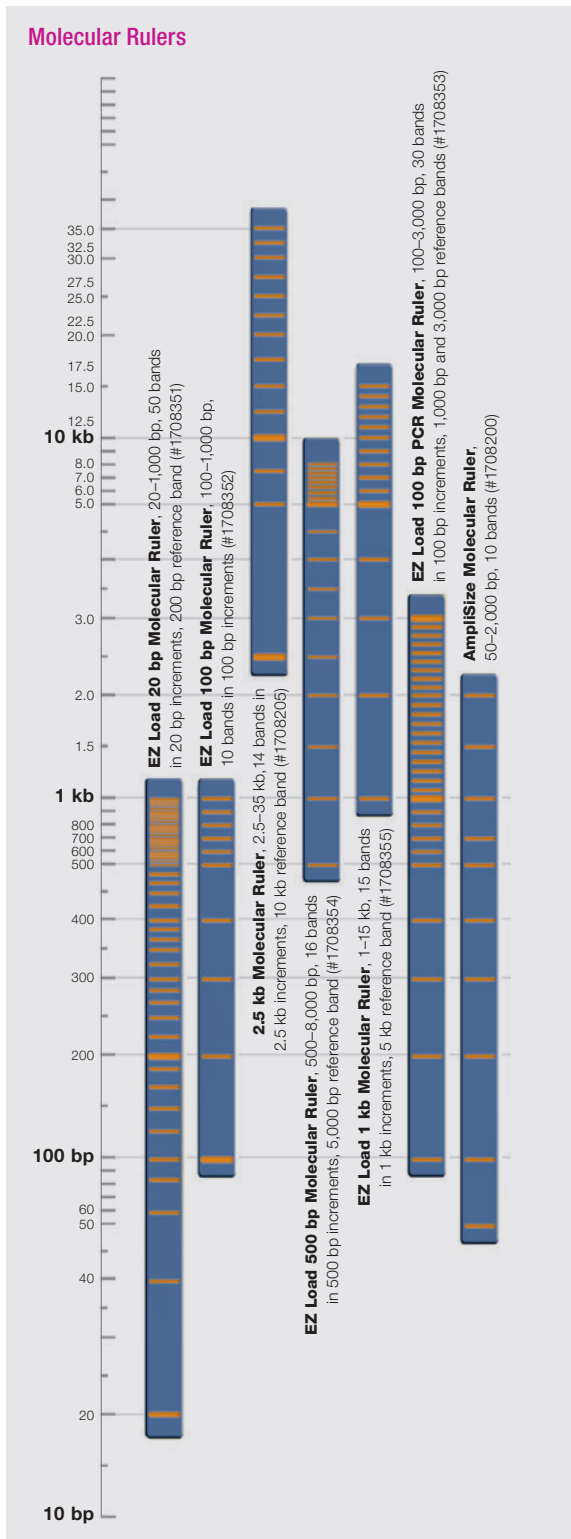
Visit bio-rad.com/PowerSupplies for more information.



Market-leading and easy-to-use, Bio-Rad gel documentation systems such as the **GelDoc Go System** are reliable for all your DNA and protein gel documentation needs.

Visit bio-rad.com/Imaging for more information.

Molecular Rulers



	Mini ReadyAgarose Precast Gels			Wide Mini ReadyAgarose Precast Gels			ReadyAgarose 96 Plus Precast 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)	4 x 26-Well (15 µl)
TBE Gels							
1% plus EtBr	1613004	1613010	—	1613028	1613034	—	—
3% plus EtBr	—	—	—	1613030	1613036	—	1613062
TAE Gels							
1%	1613015	—	1613057	—	—	—	—
1% plus EtBr	1613016	1613022	—	1613044	1613050	1613054	1613063
3%	1613017	—	—	—	—	—	—
3% plus EtBr	—	—	—	—	—	—	1613065

Mini ReadyAgarose Precast Gels fit in the Mini-Sub Cell GT Cell. Wide Mini and 96 Plus ReadyAgarose Precast Gels fit in the Wide Mini-Sub Cell GT Cell. Gel shelf life at room temperature: 12 months without EtBr, 6 months with EtBr.

Ordering Information

Catalog # Description

Horizontal Electrophoresis Systems and Power Supplies

- 1704487 **Mini ReadySub-Cell GT Horizontal Electrophoresis System**, includes lid and electrodes and accommodates Mini ReadyAgarose Gels
- 1704489 **Wide Mini ReadySub-Cell GT Horizontal Electrophoresis System**, includes lid and electrodes and accommodates Wide Mini ReadyAgarose Gels
- 1640300 **Mini-Sub Cell GT Horizontal Electrophoresis System and PowerPac Basic Power Supply**, includes 8- and 15-well combs, gel caster, 7 x 10 cm (W x L) UV-transparent tray, basic power supply
- 1640301 **Wide Mini-Sub Cell GT Horizontal Electrophoresis System and PowerPac Basic Power Supply**, includes 15- and 20-well combs, gel caster, 15 x 10 cm (W x L) UV-transparent tray, basic power supply
- 1645070 **PowerPac Universal Power Supply**, 100–120/220–240 V, power supply for a broad range of electrophoresis and blotting applications, includes power cord
- 1645050 **PowerPac Basic Power Supply**, 100–120/220–240 V, power supply for basic applications such as submerged horizontal gel electrophoresis, includes power cord

Buffers and Solutions

- 1610770 **10x TBE (Tris/Boric Acid/EDTA) Buffer**, 5 L
- 1610773 **50x TAE (Tris/Acetic Acid/EDTA) Buffer**, 5 L
- 1610433 **Ethidium Bromide Solution**, 10 mg/ml

Imaging

- 12009077 **GelDoc Go Gel Imaging System with Image Lab Touch Software**, includes UV/Stain-Free Imaging Tray

Visit bio-rad.com/ReadyAgarose for more information.

BIO-RAD, AMPLISIZE, MINI-SUB, and SUB-CELL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.



Bio-Rad PCR reagents and analytical instruments are manufactured under an ISO 13485:2016 certified Quality Management System and are quality control tested to ensure consistent product performance you can trust.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

