

NUCLEIC ACID ELECTROPHORESIS

ReadyAgarose™ 96 Plus Precast Gel System

ReadyAgarose 96 Plus Gels

- High-throughput capacity — 96 samples and 8 reference lanes per gel
- Highly reproducible results
- Wide selection of gels to separate a variety of DNA fragments

Convenience and Capacity for High-Throughput Runs

Integrated, Timesaving Products

The ReadyAgarose 96 Plus system delivers a complete solution for high-throughput nucleic acid electrophoresis, using convenient, quality products designed to integrate with each other:

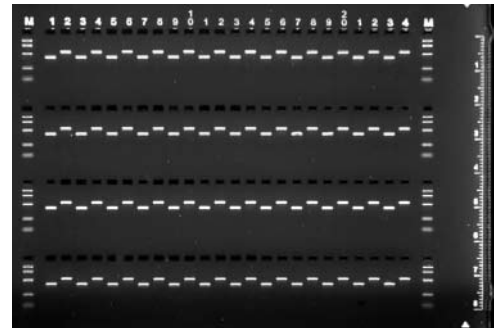
- ReadyAgarose 96 Plus precast gels
- Wide mini ReadySub-Cell™ GT cell
- ReadyAgarose 96 Plus software wizard

The ReadyAgarose 96 Plus system is suited for analysis of PCR and RT-PCR products, reaction digests, plasmid preparations, DNA libraries and arrays, and SNPs.

ReadyAgarose 96 Plus Gels

ReadyAgarose 96 Plus gels are wide mini precast agarose gels designed to accommodate 96 samples and 8 reference lanes in just one run. The wells are defined for easy loading of samples directly from microplates and are compatible with 4- and 12-channel pipets.

Each lane offers a 2 cm resolution distance, delivering up to 20% more resolving power than other commercially available 96-well precast gels. ReadyAgarose 96 Plus gels are available in 1% and 3% agarose concentrations in TAE and TBE buffers, and contain ethidium bromide. The optimal resolution ranges of the gels are 500 bp–10 kb for 1% gels and 20 bp–2 kb for 3% gels. ReadyAgarose 96 Plus gels can be stored at room temperature for up to 12 months.

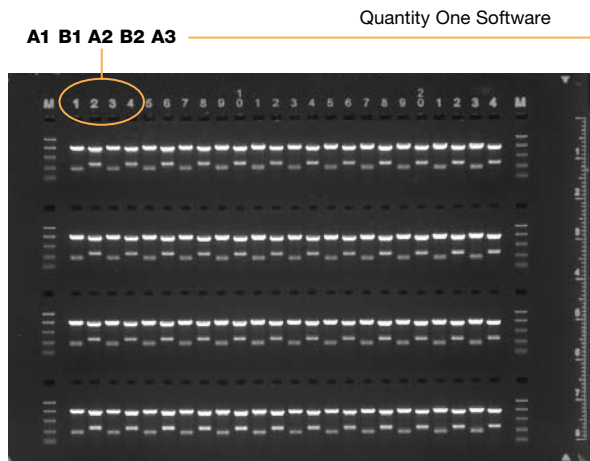


Separation of DNA fragments on ReadyAgarose 96 Plus gels. A 3% gel with 375 bp and 530 bp fragments in alternating numbered lanes and molecular marker, 100 bp–2 kb, in the "M" marker lanes.

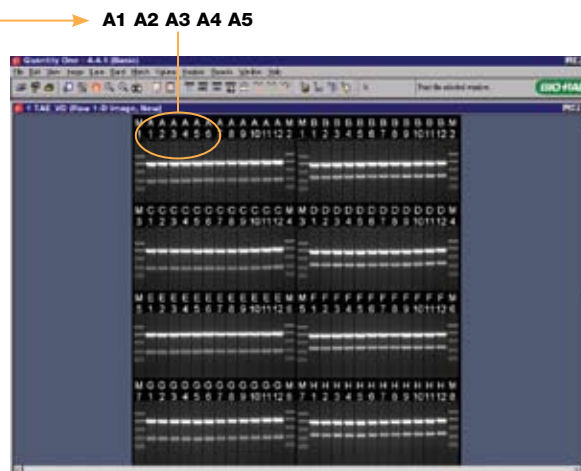
To ensure the highest quality and reproducibility, ReadyAgarose 96 Plus gels are manufactured with genetic quality tested (GQT) Certified™ agarose powders. In addition, they have all the features that have made the ReadyAgarose line of precast gels so popular, including the UV-transparent tray with fluorescent well numbers and a ruler.



BIO-RAD



Lanes on a ReadyAgarose 96 Plus gel loaded with a multichannel pipet.



Lanes on a ReadyAgarose 96 Plus gel image after rearrangement with the ReadyAgarose 96 Plus wizard in Quantity One Basic software.

Wide Mini ReadySub-Cell GT Cell

ReadyAgarose 96 Plus gels fit Bio-Rad's wide Mini-Sub[®] Cell GT cell and wide mini ReadySub-Cell GT cells. The patent-pending lock tabs on the ReadyAgarose 96 Plus gel tray lock into the dedicated Bio-Rad subcells to prevent errors due to gel movement during sample loading.

ReadyAgarose 96 Plus Wizard

The ReadyAgarose 96 Plus wizard is a feature in the free Quantity One Basic software available for download. ReadyAgarose 96 Plus gels are multichannel pipet compatible. However, due to the tip-to-tip spacing of multichannel pipets, samples from adjacent wells in a 96-well microplate are loaded in alternating wells on the gel. The ReadyAgarose 96 Plus wizard rearranges the gel image and displays it in the same format as the original 96-well plate.

You can take advantage of this feature with **free** Quantity One[®] Basic software. To download this software, go to www.bio-rad.com/software/

For more information on the ReadyAgarose 96 Plus wizard in Quantity One, request bulletin 2980.

Ordering Information

Catalog #	Description
ReadyAgarose 96 Plus TBE Gels*	
161-3060	1% ReadyAgarose 96 Plus Gel , with ethidium bromide
161-3062	3% ReadyAgarose 96 Plus Gel , with ethidium bromide
ReadyAgarose 96 Plus TAE Gels*	
161-3063	1% ReadyAgarose 96 Plus Gel , with ethidium bromide
161-3065	3% ReadyAgarose 96 Plus Gel , with ethidium bromide
ReadySub-Cell GT Cell	
170-4489	Wide Mini ReadySub-Cell GT Cell , includes buffer tank, lid and electrodes, leveling bubble

* Dimensions are 15 x 10 cm (W x L) and fit Bio-Rad's wide mini subcells as well as other manufacturers' subcells of similar size.

The polymerase chain reaction (PCR) process is covered by patents owned by Hoffman-LaRoche. Use of the PCR process requires a license.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 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 0508 805 500
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 United Kingdom 020 8328 2000