

AnyGel™ stands, for any electrophoresis laboratory.

Unique, Convenient Docking Stations for Any Size of Electrophoresis Gel.

AnyGel stands are stations for docking and prepping gels throughout an electrophoresis procedure. The stands safely hold any size of gel cassette. They're ideal for handcasting in empty Criterion™ or Ready Gel® cassettes, or for loading samples onto gels. The 6-row AnyGel stand offers an angled position for IPG strip loading onto Criterion or PROTEAN®-sized gel cassettes. The 6-row stand also provides an efficient transition from a PROTEAN IEF cell to a Dodeca™ cell for labs performing 2-D electrophoresis.

BIO-RAD

AnyGel Stands Are Great Tools when You're Preparing Gels.

Stable and Secure Gel Positioning

Both stands support the largest Bio-Rad gel size (handcast, 25 x 20.5 cm) with no danger of tipping on the lab bench. The stands are ideal for handcasting because they offer a large, level base to stabilize the empty gel cassettes while supporting them in a secure vertical position, without squeezing the cassettes, for level gel casting.

Complete Gel Visibility and Accessibility

The rows in the 6-row stand are spaced 3.8 cm apart with each row 1.9 cm higher than the row in front of it, allowing easy access for handling of individual gels. In addition to the stair-step design, the clear clamps make all gels clearly visible for casting and loading.

Ideal Design for IPG Strip Loading

The 6-row stand includes a mechanism to angle a gel cassette in an ideal position to facilitate IPG strip loading. The 6-row AnyGel stand is the only stand that provides both vertical and angled gel positioning in one stand!

Quality Built In

All bases and clamps are molded out of durable polycarbonate.

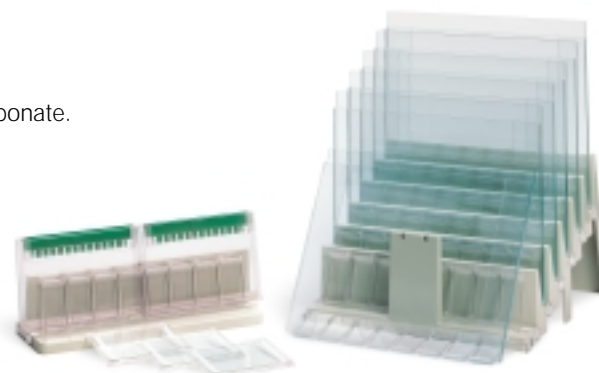
Specifications

AnyGel Stand, Single-Row

Overall dimensions	27.9 x 10.2 x 7.9 cm (W x D x H)
Capacity	1 vertical PROTEAN-sized gel 2 vertical Criterion gels 3 vertical mini gels

AnyGel Stand, 6-Row

Overall dimensions	27.9 x 25.4 x 17.8 cm (W x D x H)
Capacity	6 vertical PROTEAN-sized gels, or 5 vertical plus 1 angled in front 12 vertical Criterion gels plus 1 angled in front 18 vertical mini gels*



* Small mini gels and 7 cm strips can be easily handled without a specialized stand for loading. Mini-PROTEAN 3 gels are not positioned appropriately for IPG strip loading in the 6-row AnyGel stand.

Ordering Information

Catalog #	Description
165-4131	AnyGel Stand, single-row, includes single-row base, 1 clamp, instructions
165-5131	AnyGel Stand, 6-row, includes 6-row base, 7 clamps, instructions
165-6020	Criterion Cell and Single-Row AnyGel Stand (includes catalog #165-6001 and 165-4131)
165-3321	Mini-PROTEAN 3 Cell and Single-Row AnyGel Stand (includes catalog #165-3301 and 165-4131)
165-4132	Replacement Clamps (2)

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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 86-10-8201-1366/68 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 4124 Italy 34 91 590 5200 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 7 095 979 98 00 Singapore 65-2729877 Spain 34-91-590-5200 Sweden 46 (0)8-55 51 27 00 Switzerland 061-717-9555 United Kingdom 0800-181134