



## Dodeca™ Stainers



Simplified high-throughput staining and improved handling for large format 2-D gels.

**BIO-RAD**

# The New Movement in Gel

The **shaker motor** provides variable settings for optimal staining efficiency. The patent-pending shaking motion is gentle to protect gels from breaking.

The **shaker control unit** includes the power switch and controls the oscillating speed of the shaking rack.

The **reagent access door** offers easy access for pouring in staining solutions.

The clear **staining trays** hold the gels during staining and facilitate large format gel handling. They can be used to transport gels throughout the laboratory. Gels easily slide off the tray onto an imager or spot cutter platform. The white staining tray holds the top gel to easily monitor the final development of the staining process.

The **tray attachments** connect to the staining trays to hold certain gel sizes in an appropriately sized area to prevent excessive movement and breakage.

The **solution tank** has two drain ports for rapid draining of staining solutions.



Dodeca stainers are high-throughput gel staining devices that accommodate up to 12 large format gels. Their design ensures high-quality, consistent results and eliminates gel breakage, allowing thorough analysis of results. Dodeca stainers match the capacity of the PROTEAN® IEF cell for first-dimension isoelectric focusing and the PROTEAN® Plus Dodeca cell for second-dimension electrophoresis to help streamline the 2-D gel electrophoresis workflow.

## Simplified Staining

With the convenient and compact design of the Dodeca stainers, staining multiple gels is no longer a chore. The shaking mechanism is integrated into the design, eliminating the need to purchase an additional shaker. This patent-pending shaking design allows gentle, consistent, and thorough staining. The Dodeca stainers include a shaking rack that holds up to 12 gels at once, allowing them to be easily managed as a single unit.

## Improved Gel Handling

The ease of handling large format gels is enhanced using Dodeca stainers. Features of the Dodeca stainers minimize direct hands-on gel manipulation and ensure that gels remain intact for downstream protein analysis.

# Staining

The **shaking rack** holds the staining trays at an angle to improve reagent flow and allow air bubbles to escape, ensuring uniform gel staining. The shaking rack also features built-in handles for easy placement in and out of the solution tank.



## Gel Clip

### Enhanced Large Format Gel Handling

The unique gel clip is a convenient tool that facilitates large format gel handling and eliminates gel breakage by minimizing direct hands-on gel manipulation. The gel clip gently but securely clamps along one edge of a gel, distributing the weight evenly so that the gel can easily be lifted without tearing. You can transfer a gel to a staining tray and then to an imaging or spot cutter platform without ever touching the gel.

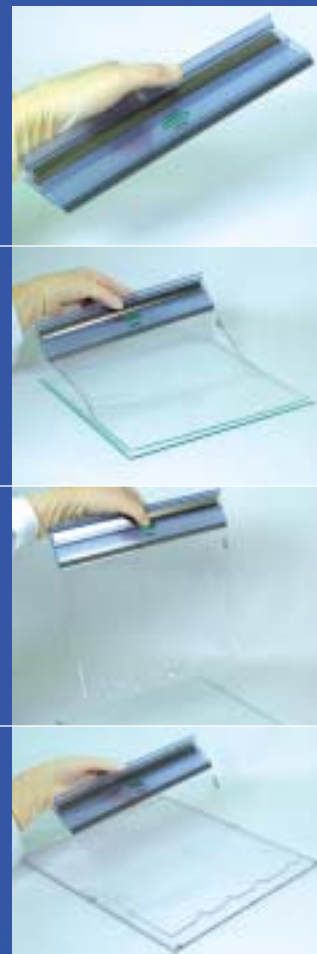
The gel clip can be used with any gel size.

The gel clip

Clamping onto a gel (PROTEAN Plus precast gel shown)

Lifting the gel off the glass plate

Transferring the gel to a large staining tray



### Available in Two Sizes

Dodeca stainers are available in two sizes to accommodate as many gel sizes as possible while maintaining high-quality staining. Both large and small Dodeca stainers provide optimized volumes for a particular gel size and eliminate gel breakage due to excessive or constricted movement.

### Stain Compatibility

The Dodeca stainers are compatible with Bio-Safe™ colloidal Coomassie Brilliant Blue G-250 stain, Coomassie Brilliant Blue R-250 stain, SYPRO Ruby protein gel stain, the new mass spectrometry compatible Dodeca silver stain kit, and the original silver stain kit (not mass spectrometry compatible). They are not compatible with the Silver Stain Plus™ kit.

### Compatibility of different gel sizes with the Dodeca stainers.

	Gel Size (W x L)	Gel Format
<b>Large Dodeca stainer</b>	25,6 x 23 cm	PROTEAN Plus precast gels
	25 x 20,5 cm	PROTEAN Plus handcast gels (require one attachment per tray)
<b>Small Dodeca stainer</b>	20 x 20,5 cm	PROTEAN Plus handcast gels
	18,5 x 20 cm	PROTEAN II XL handcast gels
	18,3 x 19,3 cm	PROTEAN II XL precast gels
	16 x 20 cm	PROTEAN II xi handcast gels
	16 x 16 cm	PROTEAN II xi handcast and precast gels
	13,3 x 8,7 cm	Criterion™ gels (up to 24 gels, require one attachment per tray)

# Dodeca Silver Stain Kit



The Dodeca silver stain kit was designed specifically for the Dodeca stainers; it has been formulated to minimize silver deposits in the stainers. The kit is conveniently packaged with stock solutions that allow easy preparation for high-throughput staining (up to 12 large format or 24 Criterion gels).

**Available in two sizes** — The small and large sizes of the kit are packaged to fill the appropriate Dodeca stainer to its maximum capacity. A minimum of 4 gels is recommended when silver staining in the Dodeca stainers.

**Reduced background** — The stain formulation ensures reduced background staining and artifacts to enhance the detection of low-abundance proteins.

**High sensitivity (0.25–0.5 ng BSA)** — High detection sensitivity allows improved protein detection. Sensitivity down to 0.25 ng BSA is possible in carefully controlled situations.

**Mass spectrometry compatible** — The stain does not modify protein residues.

For more information on the Dodeca silver stain kit, request bulletin 2971.

## Dodeca Stainer Specifications

Number of gels	1–12 large format gels 1–24 Criterion gels in the small Dodeca stainer (minimum of 4 gels recommended for silver staining)
Shaker device	Built-in shaker motor
Maximum staining solution volume	10 L for the large Dodeca stainer 7 L for the small Dodeca stainer
Compatible stains	Bio-Safe colloidal Coomassie Brilliant Blue G-250 stain, Coomassie Brilliant Blue R-250 stain, SYPRO Ruby protein gel stain, and Dodeca silver stain kit
Overall dimensions	41.3 x 46.2 x 38.9 cm (W x L x H)
Overall weight	9.1 kg for the large Dodeca stainer 7.5 kg for the small Dodeca stainer

## Ordering Information

Catalog #	Description
<b>Dodeca Stainers</b>	
165-3400	Dodeca Stainer, large, 100–240 V, includes 13 trays (12 clear and 1 white), 12 tray attachments, shaking rack, solution tank, lid with shaker, shaker control unit, gel clip, instructions
165-3401	Dodeca Stainer, small, 100–240 V, includes 13 trays (12 clear and 1 white), 12 Criterion tray attachments, shaking rack, solution tank, lid with shaker, shaker control unit, gel clip, instructions
165-3403	Dodeca Stainer and Dodeca Silver Stain Kit, large, 100–240 V, includes one large Dodeca stainer (165-3400) and one Dodeca silver stain kit for the large tank (161-0480)
165-3404	Dodeca Stainer and Dodeca Silver Stain Kit, small, 100–240 V, includes one small Dodeca stainer (165-3401) and one Dodeca silver stain kit for the small tank (161-0481)
<b>Accessories</b>	
165-3414	Gel Clip, holds any gel size, 1
165-3429	Storage Box, large, holds up to 4 gels on large staining trays, 1
165-3430	Storage Box, small, holds up to 4 gels on small staining trays, 1
<b>Dodeca Silver Stain Kits</b>	
161-0480	Dodeca Silver Stain Kit, for large Dodeca stainer tank, stains 12 PROTEAN Plus (25.6 x 23 cm or 25 x 20.5 cm) gels
161-0481	Dodeca Silver Stain Kit, for small Dodeca stainer tank, stains 12 PROTEAN (16 x 16 cm to 20 x 20.5 cm) or 24 Criterion (13.3 x 8.7 cm) gels

Coomassie is a trademark of Imperial Chemical Industries PLC. SYPRO is a trademark of Molecular Probes, Inc.



## Easy Gel Storage

Storage boxes are available in two sizes for convenient sealed storage of gels at 4°C. The storage boxes accommodate up to 4 gels (on the staining trays) plus one additional tray to hold the top gel in position.

**BIO-RAD**

**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 Czech Republic + 420 2 41 43 05 32 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84-177 Hong Kong 852-2789-3300 India (91-124)-6398112/113/114, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 55-52-00-05-20 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 7 095 721 1404 Singapore 65-6415 3188 South Africa 00 27 11 4428508 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717-9555 Taiwan (8862) 2578-7189/2578-7241 United Kingdom 020 8328 2000