

28  
cm

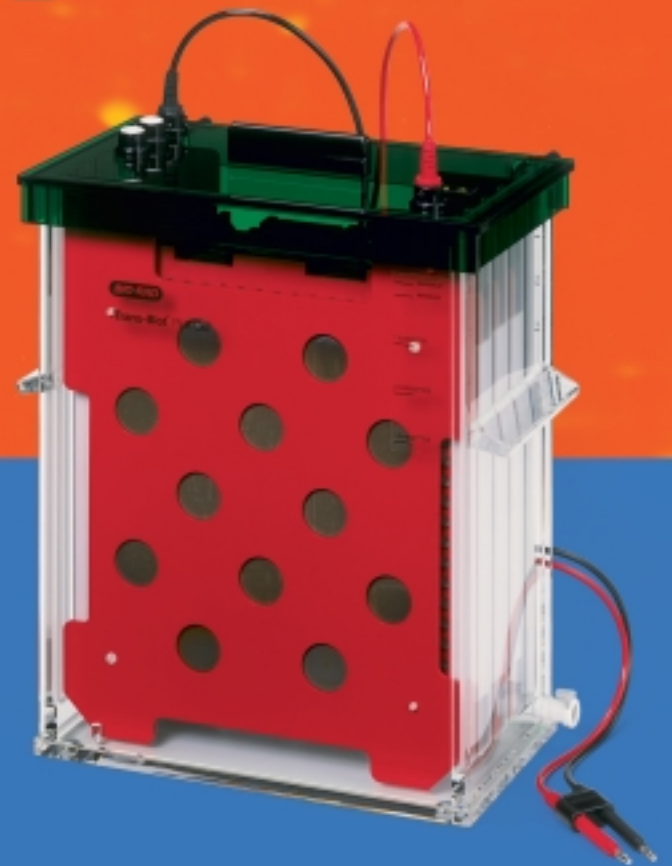
Now do blots in a

**BIG**

**way!**

26.5  
cm

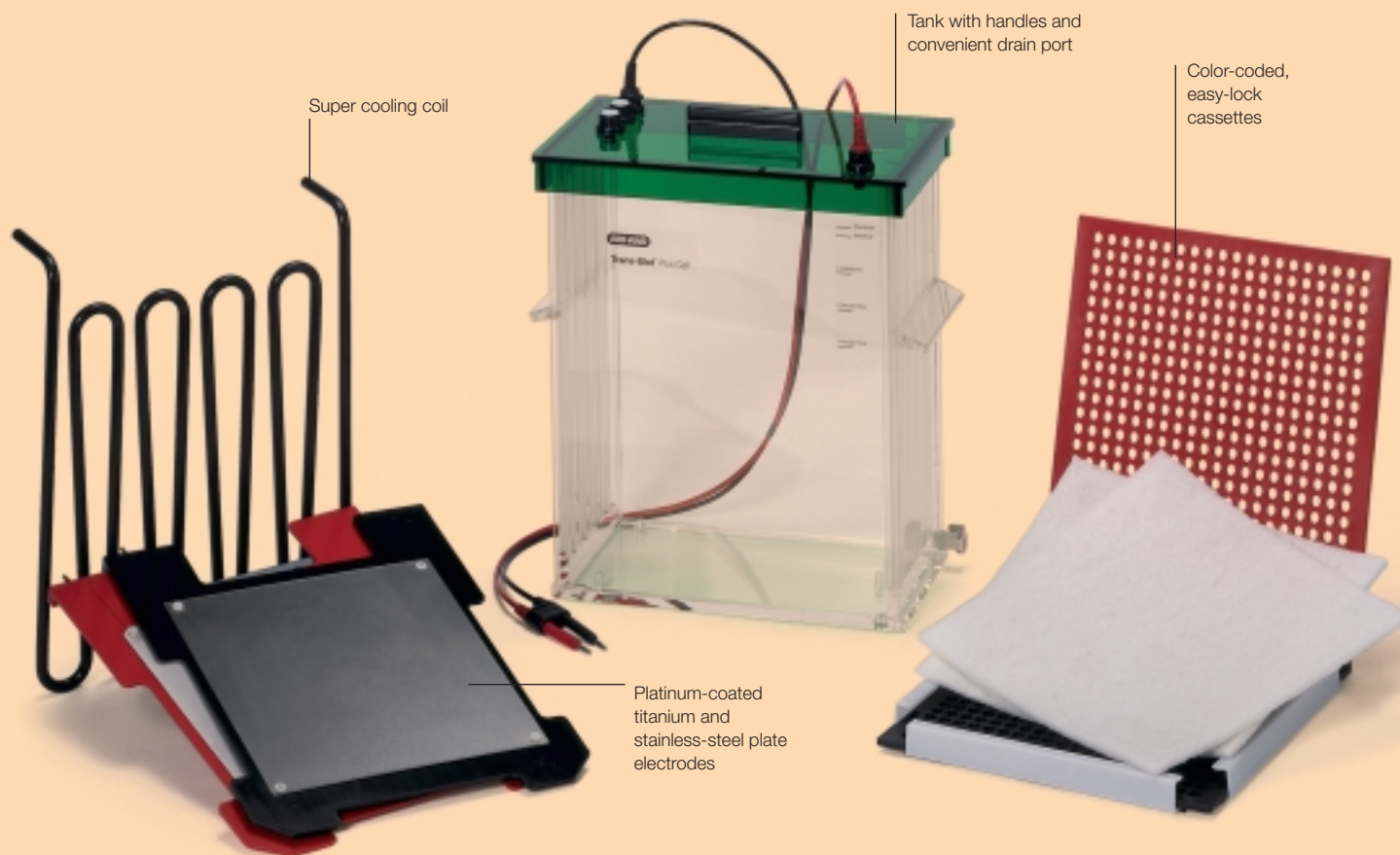
Now blot your large format gels with the efficiency of wet transfer and the speed of semi-dry transfer using the new Trans-Blot® Plus cell.



**BIO-RAD**

# Get Ready for Big Blotting!

When it's time to blot your large format gels, the Trans-Blot Plus cell offers a huge 26.5 x 28 cm blotting area. This versatile tank transfer system provides more effective and uniform protein transfer of a broad molecular weight range than any semi-dry blotting system — without sacrificing efficiency. Blot up to three large format gels or up to 12 smaller format Criterion gels in as little as 15–30 minutes!



## Tremendous Performance

### Large Capacity

- Transfer up to three 26.5 x 28 cm gels
- Compatible with multiple gel sizes from a variety of vertical electrophoresis systems: simultaneous transfer of up to 3 PROTEAN® Plus or PROTEAN® II XL gels, 12 Criterion™ precast gels, or 27 Mini-PROTEAN® 3 or Ready Gel® precast gels

### Outstanding Flexibility

- Flexible electrode placement and variable power settings allow overnight standard-field transfers or rapid high-intensity field transfers

### Fast and Efficient

- Plate electrodes create a high-strength electrical field with higher current densities than wire electrodes to produce efficient transfers in 15–30 minutes

### Durable, Easy-to-Assemble Cassettes

- Sturdy construction ensures even contact across the entire gel and membrane surface for uniform transfers
- Detachable hinge mechanism eliminates slippage of gel sandwiches during assembly
- An optional tray for gel soaking and cassette assembly makes working with even the largest gels easy

### Cooling

- The cooling coil provides temperature regulation during native enzyme or high-intensity transfers

## Specifications

### Trans-Blot Plus Cell Tank

Material	Hand-fabricated acrylic
Buffer requirement	10–12 L (depends on number of cassettes)
Dimensions (W x D x H)	30 x 17.3 x 39.4 cm

### Electrodes

Material:	Support card	Red and black molded polycarbonate
	Anode plate	Platinum-coated titanium
	Cathode plate	Stainless steel
Distance:	3 cassettes	10 cm
	2 cassettes	7 cm
	1 cassette	4 cm

Dimensions (W x L)	28.3 x 36.2 cm
--------------------	----------------

### Cassettes (3 Cassettes per Tank)

Material:	Cassette	Molded polyphthalamide PPA
	Clamps	Delrin (acetal plastic)

Maximum gel size	26.5 x 28 cm
------------------	--------------

### Gel capacity (per cassette)

PROTEAN II XL	1 gel
PROTEAN Plus	1 gel
Criterion	4 gels
Mini-PROTEAN 3	9 gels
Ready Gel	9 gels

Dimensions (W x L)	28 x 30.7 cm
--------------------	--------------

### Cooling Coil

Material	Aluminum coil coated with KYNAR
----------	---------------------------------

### Recommended Power Supply

PowerPac™ 200 power supply	
PowerPac™ HC power supply (available soon — ask for availability)	

## Ordering Configurations

### Trans-Blot Plus Cell

170-3990	Trans-Blot Plus Cell With Plate Electrodes and Super Cooling Coil, includes 3 gel holder cassettes, cell with lid and power cables, 6 fiber pads, 1 pack blot absorbent paper (26.5 x 28 cm; pack of 30), roller, stirbar
----------	---

### Accessories

170-3994	Gel/Cassette Assembly Tray
170-3995	Fiber Pads, 27 x 28.5 cm, 2
170-3996	Blot Absorbent Paper, 26.5 x 28 cm, 60 sheets
170-3997	Stirbar
170-3998	Trans-Blot Plus Roller, 6" wide
170-3999	Trans-Blot Plus Gel Holder Cassette With Clamps, 1
170-4990	Trans-Blot Plus Super Cooling Coil
170-4991	Trans-Blot Plus Platinum Anode Plate Electrode
170-4992	Trans-Blot Plus Stainless-Steel Cathode Plate Electrode
170-4995	Trans-Blot Plus Cell Buffer Tank
170-4996	Trans-Blot Plus Cell Lid With Cables
170-4997	Trans-Blot Plus Gel Holder Cassette Clamps, 3

### Membranes

162-0251	Nitrocellulose Membrane, 0.45 μm, 26.5 x 28 cm, 10 sheets
162-0252	Nitrocellulose Membrane, 0.2 μm, 26.5 x 28 cm, 10 sheets
162-0253	Supported Nitrocellulose Membrane, 0.2 μm, 26.5 x 28 cm, 10 sheets
162-0254	Supported Nitrocellulose Membrane, 0.45 μm, 26.5 x 28 cm, 10 sheets
162-0255	Immun-Blot® PVDF Membrane, 26.5 x 28 cm, 10 sheets
162-0256	Sequi-Blot™ PVDF Membrane, 26.5 x 28 cm, 10 sheets

Delrin is a trademark of E.I. du Pont de Nemours and Co. KYNAR is a trademark of Atofina Chemicals, Inc.

**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 Czech Republic + 420 2 41 43 05 32 China (86-21) 63052255 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 52 5 534 2552 to 54 The Netherlands 0318-540666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland + 48 22 8126 672 Portugal 351-21-472-7700 Russia 7 095 721 1404 Singapore 65-62729877 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