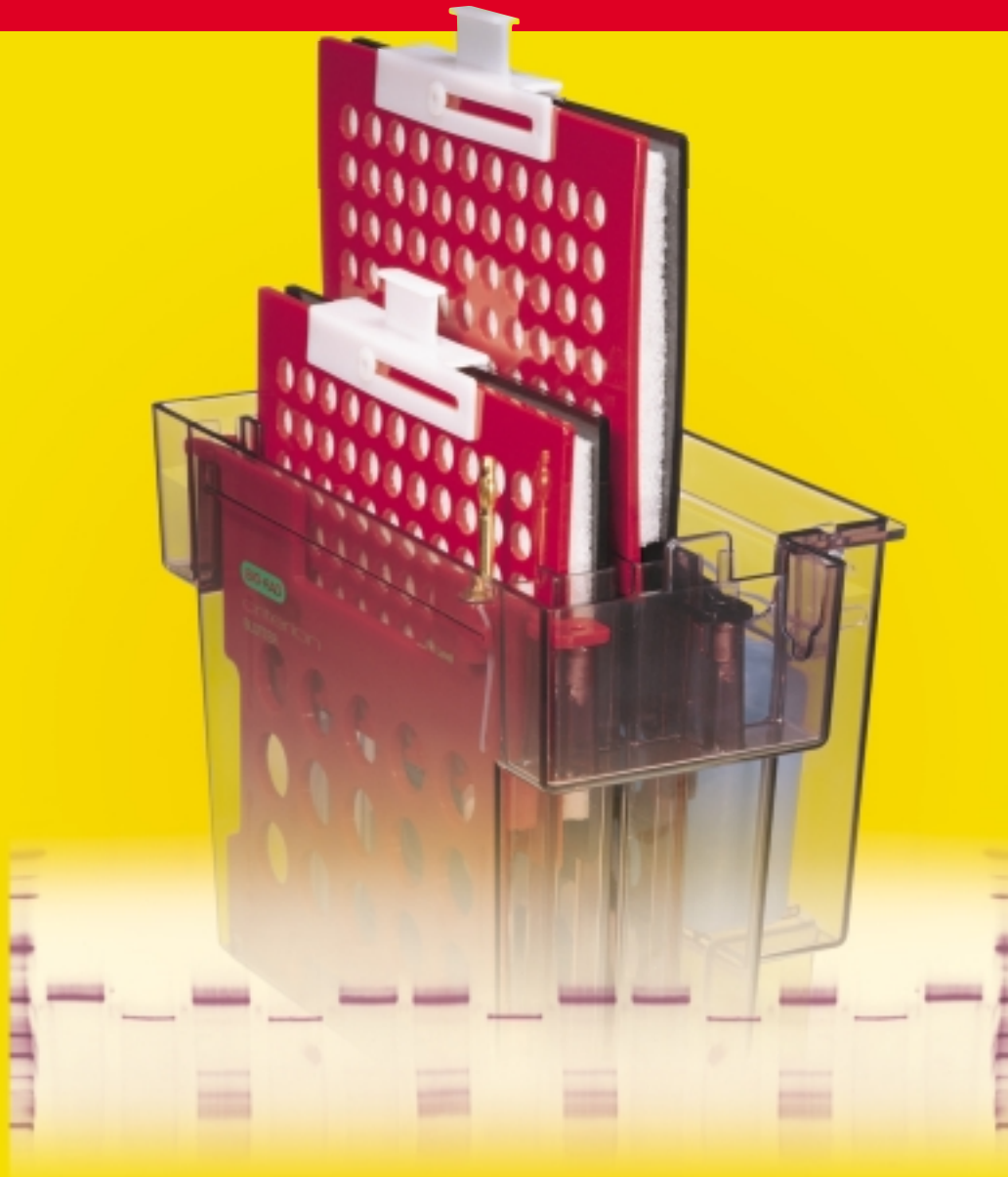


THE BLOTTER FOR THE NEXT PRECAST GEL SYSTEM



the  
**crit**erion™  
**blotter**

The Perfect Blotting Complement  
to the Criterion Precast Gel System

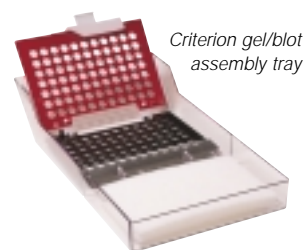
A blotter fit for use with the incredible new Criterion precast gel system, the Criterion blotter combines excellent blotting efficiency with the flexibility of the Trans-Blot™ system and the ease-of-use of the Mini Trans-Blot system.

**BIO-RAD**

# The Criterion Blotter



The Criterion blotter as assembled with wire electrodes (#170-4071)



Criterion gel/blot assembly tray

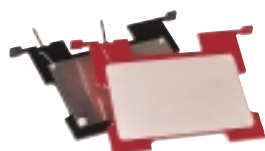


Plate electrodes (installed in #170-4070)



Fiber pads and paper



Sealed ice cooling unit



Optional Criterion blotter cooling coil

## Blots 2 Criterion Gels

Matches the sample throughput efficiency of the wide Criterion precast gels.

## Blots 4 Mini-PROTEAN 3 Gels

Allows mini-gel transfer with Criterion performance features.

## Close Electrode Proximity

Efficient transfers due to high field strength, resulting in transfer times of 30 min to 1 hr for most proteins.

## Choice of Electrodes

Durable plate electrodes consisting of a platinum-coated titanium anode and stainless steel cathode pair, or an economical platinum wire electrode pair.

## Tidy-Lab Features

Cassettes are set low, with handles for easy removal, and overflow chambers make spillover impossible.

## Minimal Buffer Requirement

Only 1.3 L of buffer is required for 2 Criterion blots.

## Quality Built In

Tank, cassettes and electrode cards are molded from durable polysulfone.

## Sealed Ice Block

Ice block will provide sufficient cooling for most normal blotting applications. Sealed block is held submerged in transfer buffer. Ice block will not spill during refreezing.

## Optional Cooling Coil

For temperature-sensitive procedures, cooling coil is available for precise temperature control.

## Assembly Tray Bonus Feature

Assembly tray is the perfect size for gel soaking and sandwich assembly, and provides for correct orientation of gel and blot.

## Ordering Information

| Catalog # | Description  |
|-----------|--|
| 170-4070  | Criterion Blotter with Plate Electrodes, includes cell assembled with plate electrodes, lid with cables, 2 Criterion gel holder cassettes, filter paper pack, fiber pad pack, gel blot assembly tray, roller, sealed ice cooling unit and holder, instructions |
| 170-4071  | Criterion Blotter with Wire Electrodes, includes cell assembled with wire electrodes, lid with cables, 2 Criterion gel holder cassettes, filter paper pack, fiber pad pack, gel blot assembly tray, sealed ice cooling unit, instructions                      |
| 165-6024  | Criterion Gel/Plate Blot System, includes Criterion cell for running Criterion precast gels and Criterion blotter with plate electrodes  |
| 165-6025  | Criterion Gel/Wire Blot System, includes Criterion cell for running Criterion precast gels and Criterion blotter with wire electrodes  |
| 170-4072  | Criterion Blotter with Plate Electrodes (170-4070) and PowerPac 200 Power Supply, 110/120 V  |
| 170-4073  | Criterion Blotter Plate Electrodes (170-4070) and PowerPac 200 Power Supply, 220/240 V   |
| 170-4074  | Criterion Blotter with Wire Electrodes (170-4071) and PowerPac 200 Power Supply, 110/120 V   |
| 170-4075  | Criterion Blotter with Wire Electrodes (170-4071) and PowerPac 200 Power Supply, 220/240 V   |
| 170-4076  | Optional Criterion Blotter Cooling Coil  |

## CRITERION BLOTTER ACCESSORIES

|          |   |
|----------|---|
| 170-4080 | Criterion Gel Holder Cassettes, 2                         |
| 170-4081 | Criterion Blotter Platinum Anode Plate Electrode          |
| 170-4082 | Criterion Blotter Stainless Steel Cathode Plate Electrode |
| 170-4083 | Criterion Blotter Wire Electrode Card, anode              |
| 170-4084 | Criterion Blotter Wire Electrode Card, cathode            |
| 170-4085 | Filter Paper, 9.5 x 15.2 cm, 50                           |
| 170-4086 | Fiber Pads, 9.5 x 15.2 cm, 4                              |
| 170-4087 | Sealed Ice Cooling Unit, 2                                |
| 170-4089 | Criterion Gel/Blot Assembly Tray                          |
| 165-1279 | Roller  |

## CRITERION MEMBRANE FILTER PAPER SANDWICHES

Each sandwich consists of one membrane (nitrocellulose or PVDF) and two sheets of thick filter paper cut to fit Criterion gels, 8.5 x 13.5 cm. Supplied in packs of 20 or 50 sandwiches.

|          |   |
|----------|---|
| 162-0232 | 0.2 µm Nitrocellulose/Filter Paper Sandwich, 8.5 x 13.5cm, 20 pack  |
| 162-0233 | 0.2 µm Nitrocellulose/Filter Paper Sandwich, 8.5 x 13.5cm, 50 pack  |
| 162-0234 | 0.45 µm Nitrocellulose/Filter Paper Sandwich, 8.5 x 13.5cm, 20 pack |
| 162-0235 | 0.45 µm Nitrocellulose/Filter Paper Sandwich, 8.5 x 13.5cm, 50 pack |
| 162-0236 | 0.2 µm PVDF/Filter Paper Sandwich, 8.5 x 13.5cm, 20 pack            |
| 162-0237 | 0.2 µm PVDF/Filter Paper Sandwich, 8.5 x 13.5cm, 50 pack            |

**BIO-RAD**

**Bio-Rad  
Laboratories**

Life Science  
Group

Web Site [www.bio-rad.com](http://www.bio-rad.com) U.S. (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771  
China 86-10-62051850/51 Denmark 45 39 17 99 47 Finland 358 (0)9 804 2200 France 01 43 90 46 90 Germany 089 318 84-0  
Hong Kong 852-2789-3300 India (91-11) 461-0103 Israel 03 951 4127 Italy 39-02-216091 Japan 03-5811-6270 Korea 82-2-3473-4460  
Latin America 305-894-5950 Mexico 514-2210 The Netherlands 0318-540666 New Zealand 64-9-4152280 Norway 22-74-18-70 Russia 7 095 979 98 00  
Singapore 65-2729877 Spain 34-91-661-7085 Sweden 46 (0)8-55 51 27 00 Switzerland 01-809 55 55 United Kingdom 0800-181134