

The ProteomeWorks™ System



1st dimension just went  
**extreme.**



The **Cup Loading Tray** for the PROTEAN® IEF Cell

**BIO-RAD**

# Cup loading is ...

an effective method for improved resolution and separation by IEF at **extreme** pH ranges.

## Ultimate Resolving Power

- Combine the new cup loading tray for the PROTEAN IEF cell with the new micro-range ReadyStrip™ IPG strips for ultimate resolving power and unbeatable performance for first-dimension runs!

## Versatility and Flexibility

- Cup loading expands the versatility and range of applications for first-dimension isoelectric focusing using the PROTEAN IEF cell
- The cup loading tray accommodates all ReadyStrip sizes, from 7 to 24 cm
- The movable electrodes provide the flexibility to run any ReadyStrip length, from 7 to 24 cm, in one universal focusing tray

## High Capacity and Throughput

- The cup loading tray accommodates up to 12 ReadyStrip IPG strips in a single tray per run
- To further increase throughput, rehydrate a second set of strips in the appropriate rehydration/equilibration tray while running another set in the cup loading tray

## Easy to Use

- The disposable sample cups easily snap into the cup loading tray on top of the strips. The cups are available in two sizes, 150 µl and 100 µl, and come in a convenient chain of 12, so the number of cups needed for each experiment can be easily cut off
- The durable polycarbonate cup loading tray ensures sufficient heat transfer for accurate and reproducible pI determination, and is easy to clean

## Ordering Information

Catalog # Description

### PROTEAN IEF Cell

- 165-4000 PROTEAN IEF System, complete, includes basic unit, 17, 11, and 7 cm focusing trays with lids, 1 pack each of 17, 11, and 7 cm rehydration/equilibration trays with lids, 2 pairs of forceps, pack of electrode wicks, mineral oil, cleaning brushes, instructions
- 165-4001 PROTEAN IEF Cell, 90–240 VAC, basic unit, includes cell, instructions

### Cup Loading Tray

- 165-4050 Cup Loading Tray, includes 1 pair movable electrodes, 1 pack each of large (150 µl) and small (100 µl) replacement cups
- 165-4055 Cup Loading Tray With Forceps
- 165-4051 Large Replacement Cups (150 µl), 1 pack (120 count)
- 165-4052 Small Replacement Cups (100 µl), 1 pack (120 count)
- 165-4053 Replacement Movable Electrodes, 1 pair
- 165-4054 Replacement Tray Base, 1

SYPRO is a trademark of Molecular Probes, Inc. Bio-Rad is licensed by Molecular Probes to sell SYPRO products for research use only, under US patent 5,616,502.

**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) 2-4141 0532 **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, 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** 64-9-4152280 **Norway** 47-23-38-41-30 **Poland** (48) 22-8126 672 **Portugal** 351-21-472-7700 **Russia** 7 095 721 1404 **Singapore** 65-2729877 **South Africa** 00 27 11 4428508 **Spain** 590 5200 **Sweden** 46 (0)8-55 51 27 00 **Switzerland** 061 717-9555 **United Kingdom** 0800-181134

A. Conventional in-gel sample loading



B. Sample loading with cup loading tray



*E. coli* sample (300 µg) from Bio-Rad's ReadyPrep™ 2-D starter kit was loaded onto 11 cm ReadyStrip IPG strips, pH 7–10. A, results using the conventional in-gel sample loading protocol. B, results using the cup loading method. The sample cup was placed close to the anode. The strips were focused for a total of 30 kV-hr on the PROTEAN IEF cell. The second-dimension separation was carried out using 8–16% Criterion™ Tris-HCl precast gels and the Criterion cell. Gels were stained with SYPRO Ruby protein gel stain and imaged using Bio-Rad's Personal Molecular Imager FX™ system.



### ReadyStrip IPG Strips, 12 per package\*

pH Range	7 cm	11 cm	17 cm	18 cm
3–6	163-2003	163-2017	163-2010	163-2035
7–10	163-2005	163-2019	163-2012	163-2037
3.9–5.1	163-2028	163-2024	163-2020	163-2038
4.7–5.9	163-2029	163-2025	163-2021	163-2039
5.5–6.7	163-2030	163-2026	163-2022	163-2040
6.3–8.3	163-2031	163-2027	163-2023	163-2041

\* Refer to the Bio-Rad Life Science catalog or visit us on the Web at [discover.bio-rad.com](http://discover.bio-rad.com) for complete information on ReadyStrip IPG strips.