



PROTEAN[®]
Plate
Washer/Holder

Instruction
Manual

Catalog Number

165-1991

165-1992

165-1993

BIO-RAD

Table of Contents

	Page
Section 1 General Information.....	1
1.1 Introduction	1
1.2 Specifications	1
1.3 Chemical Compatibility	1
1.4 Equipment and Accessories	2
Section 2 Description of Major Parts	3
2.1 Plate Washer/Holder Components.....	3
2.2 Tank and Lid	3
2.3 Plate Holder	3
Section 3 Operating Instructions	4
Section 4 Maintenance.....	5

Section 1

General Information

1.1 Introduction

The PROTEAN II xi plate washer/holder consists of two plate holders, a washing tank, and a lid. The plate holders minimize operator handling of glass plates, which reduces the likelihood of chipping and cracking the fragile plates. The washing tank accommodates all PROTEAN II xi, PROTEAN II, and Mini-PROTEAN® II glass plates for washing, and makes an ideal storage holder for clean glass plates.

1.2 Specifications

Tank and Lid

Construction	Polypropylene
Wall thickness	1.5 mm (minimum)
Size	38 cm (W) x 38 cm (L) x 37.0 cm (H)
Weight	1.9 kg

Plate Holder

Construction	
Plate holder	Polypropylene
Handle	Stainless steel
Nuts and screws	Nylon
Size	7.9 cm (W) x 22.9 cm (L) x 12.7 cm (H)
Weight	0.5 kg

1.3 Chemical Compatibility

The plate holder, tank, and lid are constructed of polypropylene. Polypropylene is not compatible with halogens, fuming nitric acid, or aromatic and chlorinated hydrocarbons at high temperatures. These chemicals will attack and destroy these parts.

Chromerge will weaken these parts after extended exposure. We do not recommend using the plate holder with Chromerge. If you must use these holders for Chromerge cleaning, do not store the plate holders in Chromerge solutions. NEVER USE THE TANK TO STORE CHROMERGE SOLUTION. USE A GLASS HOLDING TANK FOR ALL CLEANING WITH CHROMERGE.

1.4 Equipment and Accessories

Catalog

Number	Product Description
165-1991	PROTEAN II xi Plate Washer/Holder , includes tank, lid, plate holders (2), instructions, and a 400 g sample of Bio-Rad's Cleaning Concentrate
165-1992	PROTEAN II xi Plate Holder , and instructions
165-1993	PROTEAN II xi Plate Washer Tank and Lid , and instructions
165-0722	Bio-Rad Cleaning Concentrate , 1 kg bottle

PROTEAN II xi Cells

Catalog

Number	Product Description
165-1801	PROTEAN II xi Cell , 16 cm, without spacers and combs
165-1802	PROTEAN II xi Cell , 16 cm, 1.5 mm spacers (4) 15 well comb (2)
165-1803	PROTEAN II xi Cell , 16 cm, 1.0 mm spacers (4) 15 well comb (2)
165-1804	PROTEAN II xi Cell , 16 cm, 0.75 mm spacers (4) 15 well comb (2)
165-1811	PROTEAN II xi Cell , 20 cm, without spacers and combs
165-1812	PROTEAN II xi Cell , 20 cm, 1.5 mm spacers (4) 15 well comb (2)
165-1813	PROTEAN II xi Cell , 20 cm, 1.0 mm spacers (4) 15 well comb (2)
165-1814	PROTEAN II xi Cell , 20 cm, 0.75 mm spacers (4) 15 well comb (2)

PROTEAN II xi Glass Plates

Catalog

Number	Product Description
165-1821	Inner Plate , 16 cm cell, 16 x 20 cm (2)
165-1822	Outer Plate , 16 cm cell, 18.3 x 20 cm (2)
165-1823	Inner Plate , 20 cm cell, 20 x 20 cm (2)
165-1824	Outer Plate , 20 cm cell, 22.3 x 20 cm (2)
165-1825	Frosted Inner Plate (Agarose Gels) , 16 cm cell, 16 x 20 cm (2)
165-1826	Frosted Inner Plate (Agarose Gels) , 20 cm cell, 20 x 20 cm (2)
165-1827	Beveled Inner Plate (2-D procedures, bevel length = 16 cm) , 16 x 20 cm (2)
165-1828	Beveled Inner Plate (2-D procedures, bevel length = 16 cm) , 20 x 20 cm (2)
165-1832	Notched Inner Plate (double-up procedures, bevel length = 16 cm) , 16 cm cell, 16 x 20 cm (1)
165-1833	Notched Inner Plate (double-up procedures, bevel length = 16 cm) , 20 cm cell, 20 x 20 cm (1)

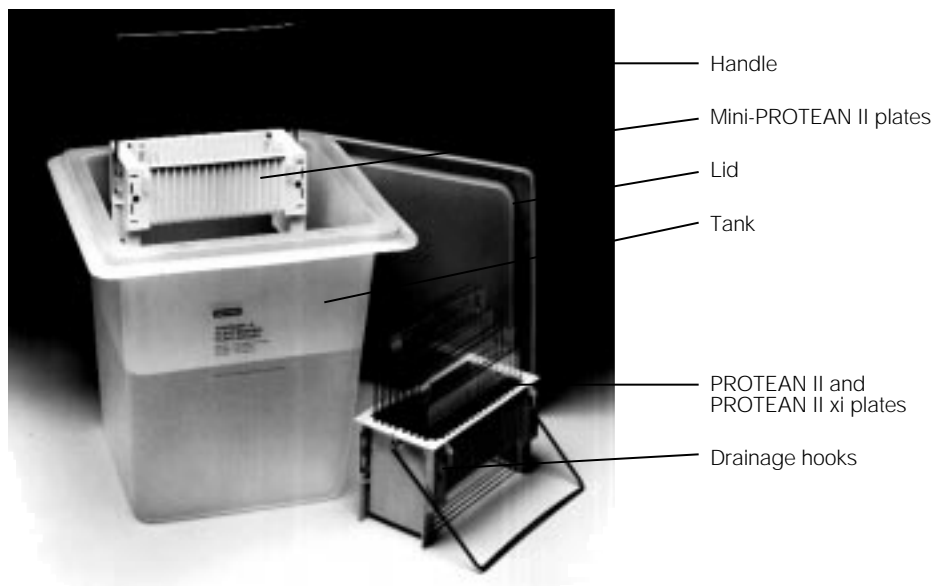
Mini-PROTEAN II Cell and Accessories

Catalog

Number	Product Description
165-2940	Mini-PROTEAN II Cell , 10 well combs, 0.75 mm spacers
165-2941	Mini-PROTEAN II Cell , same as 165-2940 without combs and spacers
165-2912	Glass Plates , 10 pairs of inner and outer plates
165-2907	Inner Glass Plates , 7.3 x 10.2 cm, 10
165-2908	Outer Glass Plates , 8.3 x 10.2 cm, 10

Section 2 Description of Major Parts

2.1 Plate Washer/Holder Components



2.2 Tank and Lid

The plate washer tank and lid can be used for both washing and plate storage. The polypropylene lid prevents accidental spillage of cleaning solution from the tank, and also provides a dust-free environment for plate storage.

Polypropylene plastic construction makes the tank resistant to the strong acid and base solutions commonly used in cleaning solutions. Any standard laboratory cleaning detergent will work well as a cleaning solution. See Section 1.3 for chemical compatibility before adding solutions other than detergents to the tank. Do not use the tank as a storage container for Chromerge.

Note: We recommend Bio-Rad's Cleaning Concentrate (catalog number 161-0722) for washing glass plates with the PROTEAN II xi plate washer/holder. The moderately alkaline cleaning concentrate is harmless to skin and clothing, and cleans by solubilizing and emulsifying contaminants in the glass surface. A 400 g sample bottle is provided with each plate washer/holder.

2.3 Plate Holder

The plate holder accommodates all types and sizes of PROTEAN II xi, PROTEAN II, and Mini-PROTEAN II glass plates. Lengthwise, each plate holder will accommodate 8 PROTEAN II xi and PROTEAN II plates; widthwise they can hold 18 Mini-PROTEAN II plates.

A unique plate holder feature is its drainage hooks. These hooks are used to hang the holder on the edge of the tank to drain the wet plates following cleaning and soaking. This eliminates messy transfer of plates.

The plate holder also has a convenient fold-down stainless steel handle for raising and lowering the holder without immersing your hands in the cleaning solution. The handle allows safe transportation of the plates to and from the tank or a drying oven.

Note: The handle is shipped separately from the plate holder and must be installed prior to use.

Section 3 Operating Instructions

1. Fill the tank with enough diluted Cleaning Concentrate (1:50 dilution with water) to completely submerge the glass plates. The approximate volume of cleaning solution and amount of cleaning concentrate required for each type of glass plate is given below. The maximum amount of solution that should be added is 15 liters. This will cover the 22.3 cm long PROTEAN II xi and PROTEAN II plates and still leave room to reach the handle without having to touch the cleaning solution.

<u>Type of glass plate</u>	<u>Quantity of cleaning solution required</u>
Mini-PROTEAN II plates	8 liters (use 160 ml of cleaning concentrate)
PROTEAN II xi and PROTEAN II 16 cm cell	12 liters (use 240 ml of cleaning concentrate)
PROTEAN II xi and PROTEAN II 20 cm cell	15 liters (use 300 ml of cleaning concentrate)

Note: We do not recommend the use of the strongly caustic Chromerge cleaning solution with the plate washer/holder due to the hazards associated with this highly corrosive material.

2. Insert up to 8 PROTEAN II xi or PROTEAN II (20 cm width) glass plates, or up to 18 Mini-PROTEAN II (10 cm width) plates into each holder.

Note: Each plate holder can be filled with either PROTEAN II and PROTEAN II xi, or Mini-PROTEAN II plates; both types of plates cannot be held simultaneously in a single plate holder.

3. Lower the holders into the cleaning solution and soak the plates overnight. Two holders can be washed at a time. For more rapid washing, or to help remove traces of polymerized acrylamide from glass surfaces, repeatedly raise and lower a single plate holder into and out of the cleaning solution.
4. After soaking has been completed, withdraw the plate holder(s) and hook them onto the upper edge of the tank using drainage hooks. This allows the plates to drain completely before they are transferred to a sink for rinsing.
5. Rinse the plates thoroughly, first with tap water, then with distilled water. A final ethanol rinse is optional.
6. Air dry plates or place the plate holder(s) into a drying oven for rapid drying.
7. The clean glass plates are conveniently held upright in the plate holder. To protect the plates from dust or accidental breakage, we recommend that you store them in a clean, dry tank with lid (catalog number 165-1993).

Section 4 Maintenance

Bio-Rad's Cleaning Concentrate is a moderately alkaline solution. When diluted it has a pH of approximately 10.5. Replace the cleaning solution when the time required for cleaning begins to increase or when the pH of the solution drops. No general maintenance of the plate washer/holder is required with normal use.

BIO-RAD**Bio-Rad
Laboratories**

**Life Science
Group**

2000 Alfred Nobel Drive
Hercules, California 94547
Telephone (510) 741-1000
Fax: (510) 741-1060

Eastern Regional Office, 85A Marcus Dr., Melville, New York 11747 • Phone (516) 756-2575 • Fax (516) 756-2594
European Headquarters, Bio-Rad Laboratories, Inc., Drève du Sénéchal, 19, B-1180 Brussels • Phone 02 375 59 70 • Fax 02 374 61 62
Australia, Bio-Rad Laboratories Pty Limited, Unit 11, 112-118 Talavera Rd P.O. Box 371, North Ryde, N.S.W. 2113 • Phone 02-805-5000 • Fax 02-805-1920
Austria, Bio-Rad Laboratories Ges.m.b.H., Auhofstrasse 78D, A-1130 Wien • Phone 0222-877 89 01 • Fax 0222-876 56 29
Belgium, Bio-Rad Laboratories S.A./N.V., Begoniastraat 5, B-9810 Nazareth Eke • Phone 091-85 55 11 • Fax 091-85 65 54
Canada, Bio-Rad Laboratories (Canada) Ltd., 5149 Bradco Boulevard, Mississauga, Ontario L4W 2A6 • Phone (416) 624-0713 • Fax (416) 624-3019
China, Bio-Rad Laboratories, Yanshan Hotel Office Tower, #1307, 138A Haidian Road, Beijing • Phone 2563146 • Fax 2564308
France, Bio-Rad S.A., 94/96 rue Victor Hugo, B.P. 220, 94203 Ivry Sur Seine Cedex • Phone 01-49 60 68 34 • Fax 01-46 71 24 67
Germany, Bio-Rad Laboratories GmbH, Heidemannstraße 164, 80939 München/Postfach 450133, 80901 München • Phone 089-318 84-0 • Fax 089-318 84 100
Italy, Bio-Rad Laboratories S.r.l., Via Cellini, 18/A, 20090 Segrate Milano • Phone 02-21609.1 • Fax 02-21609-399
Japan, Nippon Bio-Rad Laboratories, K. K., Sumitomo Seimei Kachidoki Bldg 5-3-6 Kachidoki, Chuo-Ku, Tokyo 104 • Phone 03-3534-7515 • Fax 03-3534-8027
The Netherlands, Bio-Rad Laboratories B. V., Fokkerstraat 10, 3905 KV Veenendaal • Phone 08385-40666 • Fax 08385-42216
New Zealand, Bio-Rad Laboratories Pty Ltd., P. O. Box 100-051, North Shore Mail Centre, Auckland 10 • Phone 09-443 3099 • Fax 09-443 3097
Pacific, Bio-Rad Laboratories, Unit 1111, 11/F., New Kowloon Plaza, 38, Tai Kok Tsui Road, Tai Kok Tsui, Kowloon, Hong Kong • Phone 7893300 • Fax 7891257
Spain, Bio-Rad Laboratories, S. A. Avda Valdelaparra 3, Pol. Ind. Alcobendas, E-28100 Alcobendas, Madrid • Phone (91) 661 70 85 • Fax (91) 661 96 98
Sweden, Bio-Rad Laboratories AB, Gärdsvägen 7D, Box 1276, S-171 24 Solna • Phone 46-(0)8-735 83 00 • Fax 46-(0)8-735 54 60
Switzerland, Bio-Rad Laboratories AG, Kanalstrasse 17, 8152 Glattbrugg • Phone 01-810 16 77 • Fax 01-810 19 33
United Kingdom, Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead, Herts HP2 7TD • Free Phone 0800 181134 • Fax 0442 259118