



# **Mini-PROTEAN® II Multiscreen Apparatus**

## **Instruction Manual**

**Catalog number  
170-4017**



# Table of Contents

|  | Page     |
|--|----------|
| <b>Section 1 Introduction .....</b>                                    | <b>1</b> |
| 1.1 Specifications .....   | 1        |
| <b>Section 2 Equipment and Reagents.....</b>                           | <b>1</b> |
| <b>Section 3 Special Handling Features.....</b>                        | <b>3</b> |
| 3.1 Chemical Stability .....   | 3        |
| <b>Section 4 Multiscreen Assembly Procedure .....</b>                  | <b>3</b> |
| 4.1 Preparation for the Immunoassay in the Multiscreen Apparatus ..... | 3        |
| 4.2 Assembly of the Multiscreen Apparatus.....                         | 3        |
| 4.3 Sample Loading and Washing .....                                   | 5        |
| 4.4 Color Development of Enzyme Conjugates .....                       | 7        |
| 4.5 Detection with Colloidal Gold Conjugates.....                      | 8        |
| <b>Section 5 Solution for Immunoassay Applications.....</b>            | <b>8</b> |
| <b>Section 6 References.....</b>                                       | <b>9</b> |

## Section 1 Introduction

With the Mini-PROTEAN II multiscreen apparatus you can quickly and easily screen up to 40 different antibody or serum samples on Western blots, without having to cut the membranes into strips. Only 600  $\mu$ l of sample is used per channel, eliminating waste of precious antibody. Two individual and independent sample templates allow screening of either one or two mini blots.

The multiscreen apparatus is simple to operate. Electrophorese an antigen sample on an SDS-PAGE mini gel using the Mini-PROTEAN II cell, and blot it onto nitrocellulose or Zeta-Probe<sup>®</sup> membrane with the Mini Trans-Blot<sup>®</sup>, Trans-Blot, or Trans-Blot SD cell. After blocking the unreacted sites, clamp the membrane between the gasket and sample template. The assembly is held together with four screws, and the rubber sealing gasket prevents any well-to-well leakage. Pipet serum or antibody samples into each of the channels for incubation with the antigen. Wash solutions can be easily introduced with the Eppendorf<sup>®</sup> Repeater<sup>™</sup> pipet, and are rapidly removed by vacuum aspiration. The Multiscreen apparatus is compatible with all common Western blotting procedures.

### 1.1 Specifications

#### Materials

|                              |                                    |
|------------------------------|------------------------------------|
| Multiscreen apparatus        | Acrylic plastic                    |
| Multiscreen gasket           | Silicone rubber                    |
| Shipping weight              | 1.4 kg                             |
| Overall size                 | 11 x 27 x 6 cm (W x L x H)         |
| Membrane size                | 7 x 8.4 cm                         |
| Channel dimensions           | 2.5 mm x 5.2 cm x 5 mm (W x L x H) |
| Channels per sample template | 20                                 |

## Section 2 Equipment and Reagents

| Catalog Number              | Product Description   |
|-----------------------------|---|
| 170-4017                    | Mini-PROTEAN II Multiscreen Apparatus, includes Multiscreen sample templates, 2 Gaskets and Base Plate  |
| 170-4018                    | Mini-PROTEAN II Multiscreen Gaskets, 2  |
| <b>Mini-Protean II Cell</b> |   |
| 165-2940                    | Mini-PROTEAN II Cell, includes 10 well combs, 0.75 mm spacers (4), Electrode Core with Gaskets, lower buffer chamber, lid with cables, 3 sets Glass Plates, Clamp Assemblies (2), Casting Stand with Gaskets, leveling bubble, and instructions |
| <b>Transfer Cells</b>       |   |
| 170-3930                    | Mini Trans-Blot Electrophoretic Transfer Cell   |
| 170-3935                    | Mini Trans-Blot Module  |
| 170-3946                    | Trans-Blot Electrophoretic Transfer Cell, with plate electrodes   |
| 170-3910                    | Trans-Blot Electrophoretic Transfer Cell, with standard electrodes  |
| 170-3940                    | Trans-Blot SD Semi-Dry Electrophoretic Transfer Cell  |

| <b>Catalog Number</b>  | <b>Product Description</b>   |
|--|--|
| <b>Power Supplies</b>  |  |
| 165-5052   | PowerPac 200 Power Supply, 100/120 V                               |
| 165-5053   | PowerPac 200 Power Supply, 220/240 V                               |
| <b>Related Instruments</b>   |  |
| 170-6545   | Bio-Dot® Microfiltration Apparatus                                 |
| 170-6542   | Bio-Dot SF Microfiltration Apparatus                               |
| <b>Blotting Media</b>  |  |
| <b>Nitrocellulose Membrane (0.45 micron)</b>   |  |
| 162-0115   | Nitrocellulose Membrane, roll, 33 cm x 3 m, 1                      |
| 162-0113   | Nitrocellulose Membrane, sheets, 20 x 20 cm, 5                     |
| 162-0116   | Nitrocellulose Membrane, sheets, 15 x 15 cm, 10                    |
| 162-0114   | Nitrocellulose Membrane, sheets, 15 x 9.2 cm, 10                   |
| 162-0117   | Nitrocellulose Membrane, sheets, 9 x 12 cm, 10                     |
| 162-0145   | Nitrocellulose Membrane, sheets, 7 x 8.4 cm, 10                    |
| <b>Nitrocellulose Membrane (0.2 micron)</b>  |  |
| 162-0112   | Nitrocellulose Membrane, roll, 33 cm x 3 cm, 1                     |
| 162-0146   | Nitrocellulose Membrane, sheets, 7 x 8.4 cm, 10                    |
| 162-0147   | Nitrocellulose Membrane, sheets, 13.5 x 16.5 cm, 10                |
| <b>Immun-Blot® Assay Kits</b>  |  |
| Bio-Rad's Immun-Blot assay kits contain the necessary components and instructions for performing immune detection assays on blotted membranes. |  |
| 170-6460   | Immun-Blot Assay Kit - Goat Anti-Rabbit IgG (H+L)<br>AP Conjugate  |
| 170-6461   | Immun-Blot Assay Kit - Goat Anti-Mouse IgG (H+L)<br>AP Conjugate   |
| 170-6462   | Immun-Blot Assay Kit - Goat Anti-Human IgG (H+L)<br>AP Conjugate   |
| 170-6463   | Immun-Blot Assay Kit - Goat Anti-Rabbit IgG (H+L)<br>HRP Conjugate |
| 170-6464   | Immun-Blot Assay Kit - Goat Anti-Mouse IgG (H+L)<br>HRP Conjugate  |
| 170-6465   | Immun-Blot Assay Kit - Goat Anti-Human IgG (H+L)<br>HRP Conjugate  |
| 170-6466   | Immun-Blot Assay Kit - Protein A HRP                               |
| 170-6467   | Immun-Blot Assay Kit - Protein G HRP                               |
| <b>Total Protein Detection Kits</b>  |  |
| 170-6512   | Biotin-Blot Protein Detection Kit                                  |
| 170-6517   | Enhanced Colloidal Gold Total Protein Detection Kit                |
| <b>Blotting Standards</b>  |  |
| 161-0305   | Prestained SDS-PAGE Standards, Low Range                           |
| 161-0309   | Prestained SDS-PAGE Standards, High Range                          |
| 161-0318   | Prestained SDS-PAGE Standards, Broad Range                         |
| 161-0307   | Biotinylated SDS-PAGE Standards Kit, Low Range, HRP                |
| 161-0308   | Biotinylated SDS-PAGE Standards Kit, Low Range, AP                 |
| 161-0312   | Biotinylated SDS-PAGE Standards Kit, High Range, HRP               |
| 161-0313   | Biotinylated SDS-PAGE Standards Kit, High Range, AP                |
| 161-0321   | Biotinylated SDS-PAGE Standards Kit, Broad Range, HRP              |
| 161-0322   | Biotinylated SDS-PAGE Standards Kit, Broad Range, AP               |

## Section 3 Special Handling Features

The multiscreen apparatus can be cleaned with a mild, non-abrasive detergent but not be autoclaved. Do not subject the unit to temperatures greater than 50 °C, as this will warp acrylic plates. If the unit becomes warped, it will no longer provide a proper seal. Heating the apparatus to >50 °C voids all warranties.

### 3.1 Chemical Stability

#### Chemicals compatible with acrylic plastic:

|                   |               |
|-------------------|---------------|
| hydrochloric acid | < 50% ethanol |
| sodium hydroxide  | < 50% ethanol |

#### Chemicals that will attack acrylic plastic:

all polar aromatic solvents or chlorinated hydrocarbons, esters, and ketones  
glacial acetic acid  
chromic acid  
trichloroacetic acid  
> 50% ethanol  
> 50% methanol

## Section 4 Multiscreen Operating Procedure

### 4.1 Preparation for the Immunoassay in the Multiscreen Apparatus

1. Electrophorese the antigen sample into a mini gel following the instructions provided with the Mini-PROTEAN II cell. The maximum length of the separating gel should not exceed 5.2 cm, the length of the channels on the multiscreen sample template. The stacking gel should be cast with the Mini-PROTEAN II preparative comb. This comb contains one large sample well and one reference well.
2. Blot the gel to nitrocellulose or Zeta-Probe membrane, using a 7 x 8.4 cm membrane size. Refer to the Mini Trans-Blot, Trans-Blot or Trans-Blot SD cell instruction manual for electrophoretic transfer procedures.
3. Following the transfer, mark the outline of the gel on the membrane using a pen or pencil. This will aid in aligning the blot with the sample template. Block the unreacted sites on the blot with a blocking solution, 1 hour at room temperature for nitrocellulose, 2 hours at room temperature for Zeta-Probe membrane. Rinse the blocked blot in TBS before applying to the multiscreen apparatus. See Section 5 for buffer formulations.

**Note:** Always use forceps or wear gloves when handling membranes.

### 4.2 Assembly of the Multiscreen Apparatus

1. Clean and dry the multiscreen apparatus and gaskets prior to assembly.

**Note:** Do not heat the apparatus to temperatures greater than 50 °C. This will cause the unit to warp.

2. Place the sealing gasket onto the base plate with the raised surface down, using the guide pins to help align the gasket.

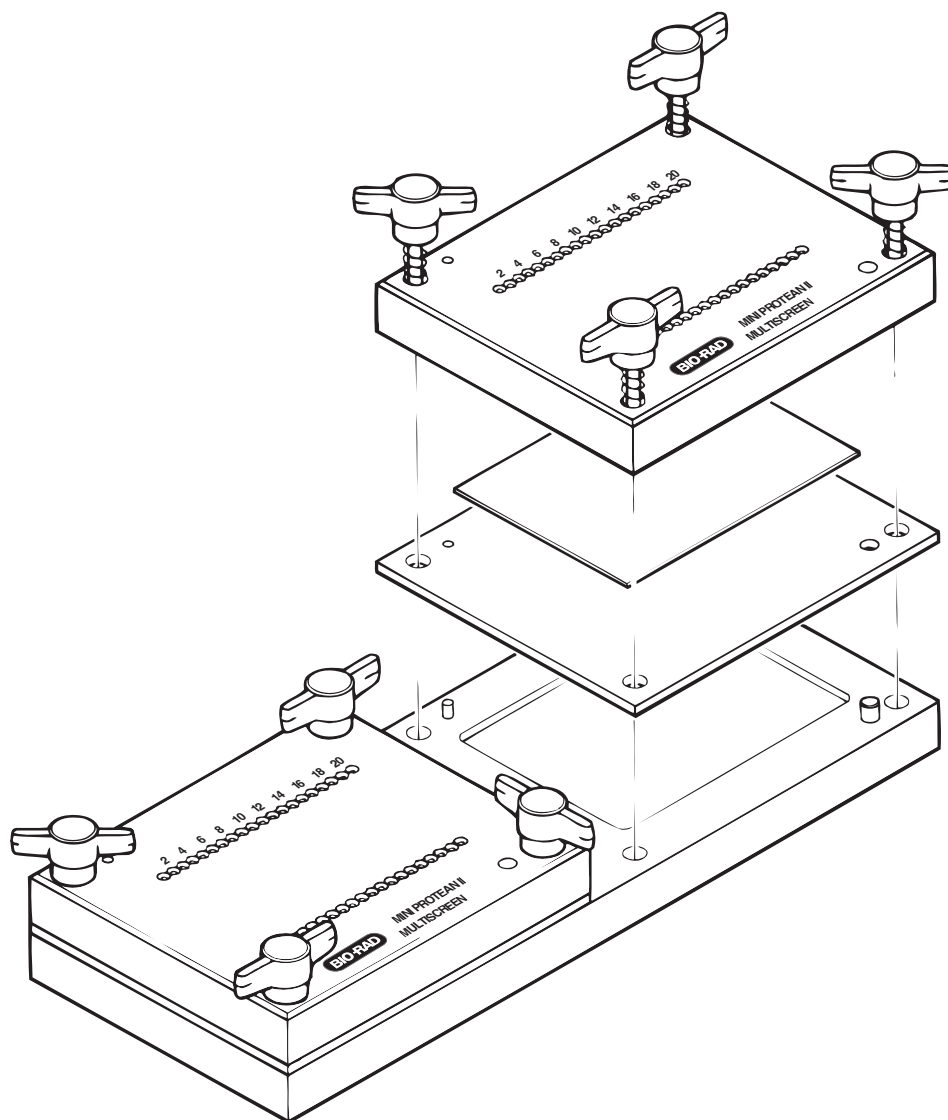


Fig. 1. Assembly of the multiscreen apparatus.

3. Lay the blocked blot on the gasket with the antigen side facing up. Center the membrane so that the channels of the sample template cover the length of the blotted sample.
4. Place the sample template on top of the membrane. The guide pins insure that the template will be properly aligned. Finger tighten the four screws. When tightening the screws, use a diagonal crossing pattern to insure even pressure on the membrane surface (see Figure 2). The multiscreen apparatus is ready for sample application.

**Note:** Use of excessive force when tightening the screws is not necessary to prevent well-to-well leakage. Finger tightening is sufficient to obtain a good seal. Overtightening can cause the channels to cut into the membrane.

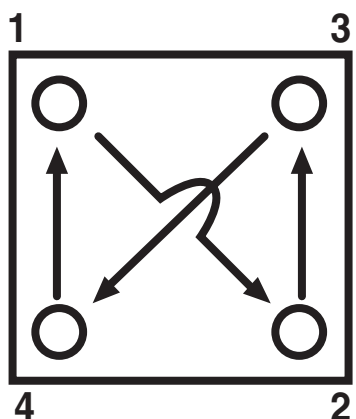


Fig. 2. Diagonal crossing pattern for tightening screws in the multiscreen apparatus.

### 4.3 Sample Loading and Washing

Detailed instructions for performing immunoassays, including a comprehensive troubleshooting guide, can be found in any of the Immun-Blot assay instruction manuals. See page 2 for a complete listing of Immun-Blot assay kits available from Bio-Rad.

1. To load an antibody or serum sample, tilt the multiscreen apparatus toward you so that the back of the unit is tilted up  $\sim 30^\circ$  (see Figure 3). Using a syringe or Eppendorf pipet, load the solution into the bottom unmarked holes of the channels. Slow, careful delivery of sample is necessary to avoid trapping bubbles inside the channels. Titrating the apparatus helps the bubbles rise to the top, towards the numbered holes of the channels. Fill the channel with 600  $\mu\text{l}$  antibody solution.

**Note:** Antibody buffers containing BSA or BLOTTO\* are recommended for use with the multiscreen apparatus. Do not use antibody buffers with gelatin, as this may cause coagulation of gelatin within the channels of the unit.

\* BLOTTO is an acronym for Bovine Lacto Transfer Technique Optimizer, and refers to non-fat dry milk.<sup>1,2</sup>



Fig. 3. Tilt the multiscreen apparatus toward you during sample application.

2. Wash solutions can be applied in the same manner as the antibody samples, or with the Eppendorf repeater pipet. Use 600  $\mu$ l per channel. The number and stringency of washes may vary and should be determined separately for each experiment. However, a minimum of three washes with a buffer containing a detergent such as Tween-20 is recommended after each antibody incubation. See Section 5 for buffer formulations.



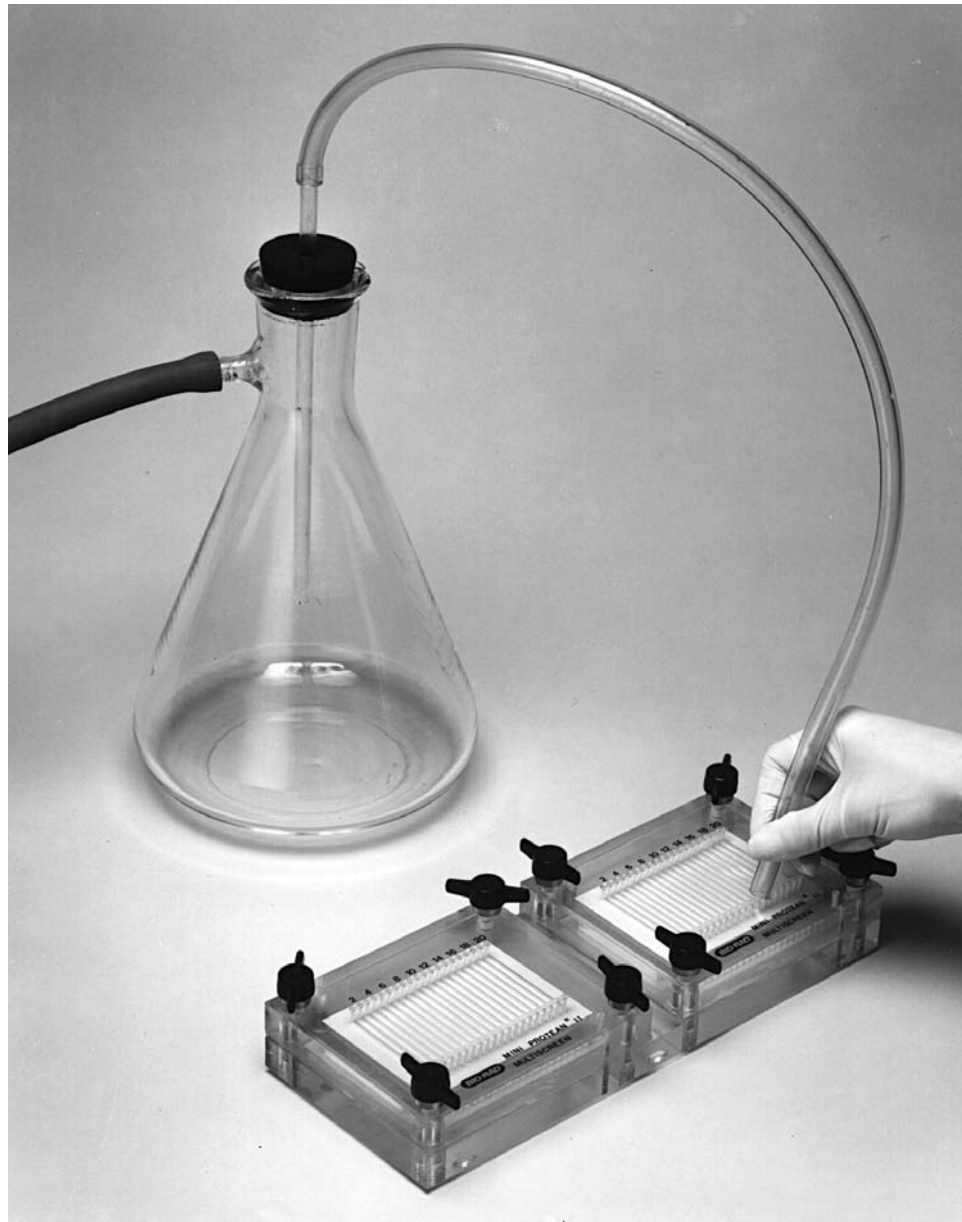


Fig. 4. Vacuum aspiration of sample and wash solutions.

3. Antibody samples and wash solutions can be rapidly removed by vacuum aspiration. The tubing from the vacuum source should be attached to the bottom, unmarked row of holes. Move the tubing back and forth along the row of holes until all the channels are dry (see Figure 4). To save an antibody or serum sample after incubation, remove the solution individually from the channel with a pipet or syringe.

#### 4.4 Color Development of Enzyme Conjugates

Incubation of the blot with antibodies conjugated to enzymes, such as horseradish peroxidase or alkaline phosphatase, can be conducted either in the multiscreen apparatus or in a separate vessel. If this step is carried out in the unit, wash the blot in the apparatus after the second antibody incubation as outlined in Section 4.3.

1. Color development of enzyme conjugated antibodies should be performed in a separate container to prevent permanent discoloration of the multiscreen apparatus. Remove the membrane by loosening the four screws. Lift out the sample template and move the membrane to a color development vessel.
2. Wash the membrane once with TBS for 5 minutes. After the color development solution has been prepared, incubate the membrane in the solution. Gently agitate the solution until development is complete. Remove the solution and rinse the membrane several times in distilled water to stop the reaction. Air dry the blot on filter paper.

#### 4.5 Detection with Colloidal Gold Conjugates

Incubation of the blot with colloidal gold conjugated antibodies, protein A, or protein G, should be conducted in a separate vessel to prevent discoloration of the multiscreen apparatus.

1. Remove the blot from the multiscreen apparatus after washing to remove excess first antibody. Place the membrane in a color development vessel.
2. Wash the membrane once with TBS for 5 minutes. Add the gold solution to the vessel until the membrane is completely covered. Gently agitate the solution. Red bands identifying antigen will appear on the membrane surface within 10–15 minutes at the sites of highest antigen concentrations. Allow the incubation to continue until the desired sensitivity is achieved.

## Section 5 Solutions for Immunoassay Applications

Tris Buffered Saline, 1 x TBS, 2 L  
20 mM Tris-HCl, 500 mM NaCl, pH 7.5  
Dissolve 4.84 g Tris, 58.48 g NaCl in ~1.5 L distilled, deionized H<sub>2</sub>O. Adjust to pH 7.5 with HCl. Adjust the volume to 2 L with dd H<sub>2</sub>O.

**Note:** Bio-Rad's Premixed Tris-Buffered Saline (catalog number 170-6430) eliminates weighing of buffer components. One bottle produces 1 L of 10 x TBS.

Tween-20 Wash Solution, 1X TTBS, 1 L  
20 mM Tris-HCl, 500 mM NaCl, 0.05% Tween-20, pH 7.5  
Add 0.5 ml Tween-20 to 1 L of TBS.  
Blocking Solution, 100 ml

Both of the following blocking solutions can be used with nitrocellulose. The solutions containing BLOTTO should be used with Zeta-Probe membrane. Incubate nitrocellulose blots for 1 hour at room temperature. Zeta-Probe membrane should be blocked for 2 hours at room temperature.

3% Gelatin - TBS  
Add 3.0 g gelatin to 100 ml TBS. Heat to 50 °C, stirring until dissolved. A microwave oven will quickly solubilize the gelatin, but do not heat above 65 °C.

OR:

5% BLOTTO in 100 ml of TBS

Add 5% g of BLOTTO to 100 ml of TBS.

Antibody Buffer, 200 ml

1% BSA - TTBS

Add 2 g BSA to 200 ml TTBS. Stir to dissolve.

OR

1% BLOTTO in TTBS.

Add 2 g of BLOTTO to 200 ml of TTBS.

### **First Antibody Solution**

Dilute antigen specific primary antibody to the appropriate titer in antibody buffer.

### **Second Antibody Solution, 100 ml**

Dilute species specific Bio-Rad second antibody enzyme conjugate, 1:3,000 by adding 33  $\mu$ l of conjugate to 100 ml of antibody buffer.

Consult the Immun-Blot assay kit instruction manual for dilution protocols of the colloidal gold conjugates.

### **Color Development Solution**

The specific chemicals and buffers are dependent on the enzyme conjugate being used. See the Immun-Blot assay kit instruction manual for details on how to make the appropriate solution.

## **Section 6 References**

1. Jerome, J. F. and Jaehning, J. A., *Mol. and Cell Bio.*, **6**, 1633 (1986).
2. Johnson, D. A., et. al., *Gene Anal. Tech.*, **1**, 3 (1984).

Eppendorf is a trademark of Eppendorf Geratebau Netheler & Hinz, GmbH



**Bio-Rad  
Laboratories**

---

**Life Science  
Group**

2000 Alfred Nobel Drive  
Hercules, California 94547  
Telephone (510) 741-1000  
Fax: (510) 741-5800

**Australia**, Bio-Rad Laboratories Pty Limited, Block Y Unit 1, Regents Park Industrial Estate, 391 Park Road, Regents Park, NSW 2143 • Phone 02-9414-2800 • Fax 02-9914-2888  
**Austria**, Bio-Rad Laboratories Ges.m.b.H., Auhofstrasse 78D, 1130 Wien • Phone (1) 877 89 01 • Fax (1) 876 56 29  
**Belgium**, Bio-Rad Laboratories S.A./N.V., Begoniastraat 5, 9810 Nazareth Eke • Phone 09-385 55 11 • Fax 09-385 65 54  
**Canada**, Bio-Rad Laboratories (Canada) Ltd., 5671 McAdam Road, Mississauga, Ontario L4Z 1N9 • Phone (905) 712-2771 • Fax (905) 712-2990  
**China**, Bio-Rad Laboratories, 14, Zhi Chun Road, Hai Dian District, Beijing 100088 • Phone (01) 2046622 • Fax (01) 2051876  
**Denmark**, Bio-Rad Laboratories, Symbion Science Park, Fruebjergvej 3, DK-2100 Copenhagen • Phone 39 17 9947 • Fax 39 27 1698  
**Finland**, Bio-Rad Laboratories, Business Center Länsikeskus, Pihatörmä 1A SF-02240, Espoo, • Phone 90 804 2200 • Fax 90 804 1100  
**France**, Bio-Rad S.A., 94/96 rue Victor Hugo, B.P. 220, 94 203 Ivry Sur Seine Cedex • Phone (1) 49 60 68 34 • Fax (1) 46 71 24 67  
**Germany**, Bio-Rad Laboratories GmbH, Heidemannstraße 164, D-80939 München/Postfach 450133, D-80901 München • Phone 089 31884-0 • Fax 089 31884-100  
**India**, Bio-Rad Laboratories, C-248 Defence Colony, New Delhi 110 024 • Phone 91-11-461-0103 • Fax 91-11-461-0765  
**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, 7-18, Higashi-Nippori 5-Chome, Arakawa-ku, Tokyo 116 • Phone 03-5811-6270 • Fax 03-5811-6272  
**The Netherlands**, Bio-Rad Laboratories B. V., Fokkerstraat 10, 3905 KV Veenendaal • Phone 0318-540666 • Fax 0318-542216  
**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  
**Singapore**, Bio-Rad Laboratories (Singapore) Ltd., 221 Henderson Rd #05-19, Henderson Building, Singapore 0315 • Phone (65) 272-9877 • Fax (65) 273-4835  
**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, Postfach, CH-8152 Glattbrugg • Phone 01-809 55 55 • Fax 01-809 55 00  
**United Kingdom**, Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead, Herts HP2 7TD • Free Phone 0800 181134 • Fax 01442 259118