

## ***READY GEL CELL***

### **INSTRUCTION MANUAL**

**Catalog Numbers**

**165-3125**

**165-3126**

**BIO-RAD**

## TABLE OF CONTENTS

---

### Safety

<b>Section 1.0 General Information</b> .....	1
1.1 Unpacking.....	1
1.2 Accessories .....	2
<b>Section 2.0 Set Up and Basic Operation</b> .....	3
2.1 Assembly of the Ready Gel Cell .....	3
2.2 Sample Loading .....	6
2.3 Running the Gel .....	7
2.4 Removing the Gel.....	7
<b>Section 3.0 Maintenance</b> .....	8
<b>Appendix A. Specifications</b> .....	9
<b>Appendix B. Warranty and Ordering Information</b> .....	10

### List of Figures

1. Ready Gel Cell Prepared for Use.....	1
2. Assembling the Ready Gel Cell.....	4
3. Sealing the Inner Chamber .....	5
4. Using the Sample Loading Guide.....	6

## SAFETY

---

Power to the Ready Gel Cell is supplied by an external DC voltage power supply. The output of this power supply must be isolated from external ground in such a way that the DC voltage output floats with respect to ground. All Bio-Rad power supplies meet this important safety requirement. Regardless of which power supply is used, the maximum specified operating parameters for the Ready Gel Cell are:

- 600 VDC            maximum voltage limit
- 15 Watts            maximum power limit
- 50° C            maximum ambient temperature limit

Current to the cell, provided from the external power supply, enters the unit through the lid assembly, which provides a safety interlock to the user. Current to the cell is broken when the lid is removed. Always turn off the power supply before removing the lid. **Do not attempt to use the cell without the safety lid.**

### Important:

This Bio-Rad product is designed and certified to meet EN61010-1 safety standards. Certified products are safe to use when operated in accordance with the instruction manual. This instrument should not be modified or altered in any way. Alteration of this instrument will:

- Void the manufacturer's warranty,
- Void the EN61010-1 certification, and
- Create a potential safety hazard.

Bio-Rad is not responsible for any injury or damage caused by use of this instrument for purposes other than for which it is intended or by modifications of the instrument not performed by Bio-Rad or an authorized agent.

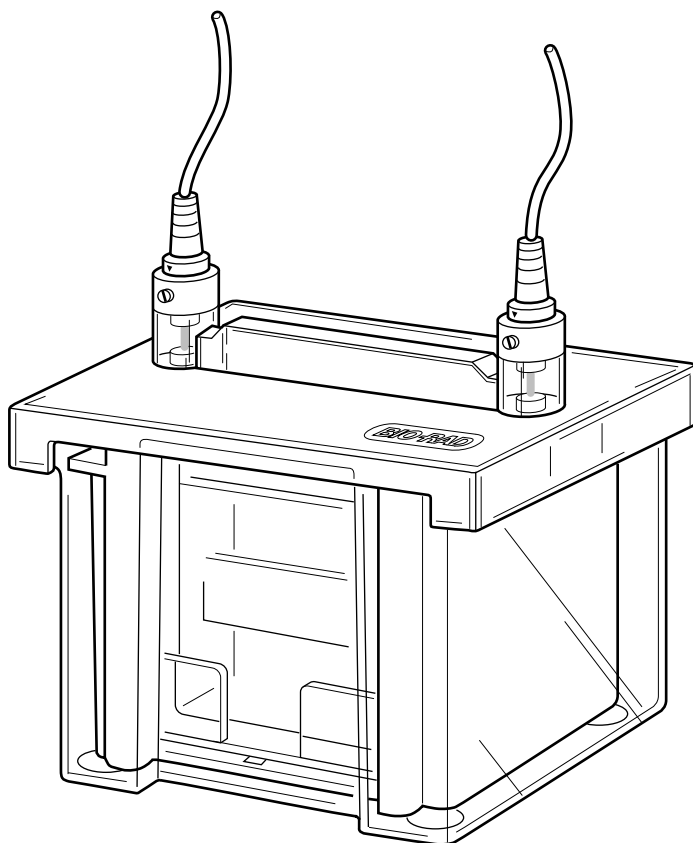
This product conforms to the class A standards for Electromagnetic Emissions, intended for laboratory equipment applications. It is possible that emissions from this product may interfere with some sensitive appliances when placed nearby or on the same circuit as those appliances. The user should be aware of this potential and take appropriate measures to avoid interference.

\* EN61010-1 is an internationally accepted electrical safety standard for laboratory instruments.

## 1.0 GENERAL INFORMATION

The Ready Gel Electrophoresis Cell will run all Bio-Rad precast Ready Gel products. The Ready Gel Cell is designed to make running precast gels very simple. Ready Gels are easily inserted into the cell and self align with the gasket to form a leak proof seal for the upper buffer. At the end of a run, the Ready Gels are easily removed.

The Ready Gel Cell can run two precast Ready Gels; one gel can be run when the Mini Cell Buffer Dam is used in place of the second gel. The Ready Gel mini tank and lid can be used with a variety of different electrode modules to run handcast gels, to blot mini gels, to run 2-D electrophoresis, or for electroelution.



*Figure 1. Ready Gel Cell Prepared for Use*

### 1.1 UNPACKING

When you receive the Ready Gel Cell, carefully inspect the shipping container for any damage which may have occurred in shipping. Severe damage to a container may indicate damage to its contents. If you suspect damage to the contents may have occurred in shipping, immediately file a claim with the carrier in accordance with their instructions before contacting Bio-Rad Laboratories.

Inspect the unit for external damage. If any part is missing or damaged, contact Bio-Rad Laboratories immediately.

## 1.2 ACCESSORIES

The following options are available with your Ready Gel Cell.

- **Sample Loading Guides, 10 and 15 Well:** The two sided sample loading guides match the well spacing of the 10 and 15 well Ready Gels. The sample guides serve to align a Hamilton Syringe or pipet tip with each of the sample wells in the gels.
- **Mini Cell Buffer Dam:** The Mini Cell Buffer Dam substitutes for a Ready Gel cassette when running only one gel. The buffer dam can also be used with the Mini-PROTEAN® II cell.

## 2.0 SET UP AND BASIC OPERATION

---

To get the best performance from your Ready Gel Cell, familiarize yourself with the unit's components by assembling and disassembling the cell before using it. Refer to Figure 2, which shows the assembling of the Ready Gel Cell, tank, and lid.

Feature	Description
<b>Clamping Frame</b>	This serves as the frame for the inner chamber. The Clamping Frame includes pressure plates and closure cams. When the cams are closed the Electrode Assembly is clamped snugly and evenly against the back of the Clamping Frame, sealing the inner chamber.
<b>Electrode Assembly</b>	The Electrode Assembly houses the molded U-shaped gasket, the upper and lower electrodes, and the connecting banana plugs. The anode (lower electrode) jack is identified with a red marker, and the cathode (upper electrode) jack with a black marker. The Electrode Assembly holds the Ready Gel cassettes.
<b>Inner Chamber</b>	The Electrode Assembly, two Ready Gel cassettes, and the Clamping Frame form the inner chamber.
<b>Mini Tank and Lid</b>	The mini tank and lid combine to fully enclose the inner chamber during electrophoresis. The lid cannot be removed without disconnecting the electrical circuit. The mini tank and lid can be used with a variety of different Mini-PROTEAN II electrode modules to run handcast gels, blot mini gels, to run 2-D electrophoresis or for electroelution.

### 2.1 ASSEMBLY OF THE READY GEL CELL

1. Remove the Electrode Assembly from the Clamping Frame. Rotate the cams outward to release the Electrode Assembly. Refer to Figure 2.
2. Prepare a Ready Gel by removing the pull tab from the bottom of the cassette and removing the comb. Rinse the sample wells with running buffer before use. For complete instructions on using the Ready Gel, see the Ready Gel box or request the Ready Gel Application Guide, catalog # 161-0993.
3. Repeat the procedure for a second Ready Gel cassette. If you will not be preparing a second Ready Gel cassette, you will need to use the mini cell buffer dam.

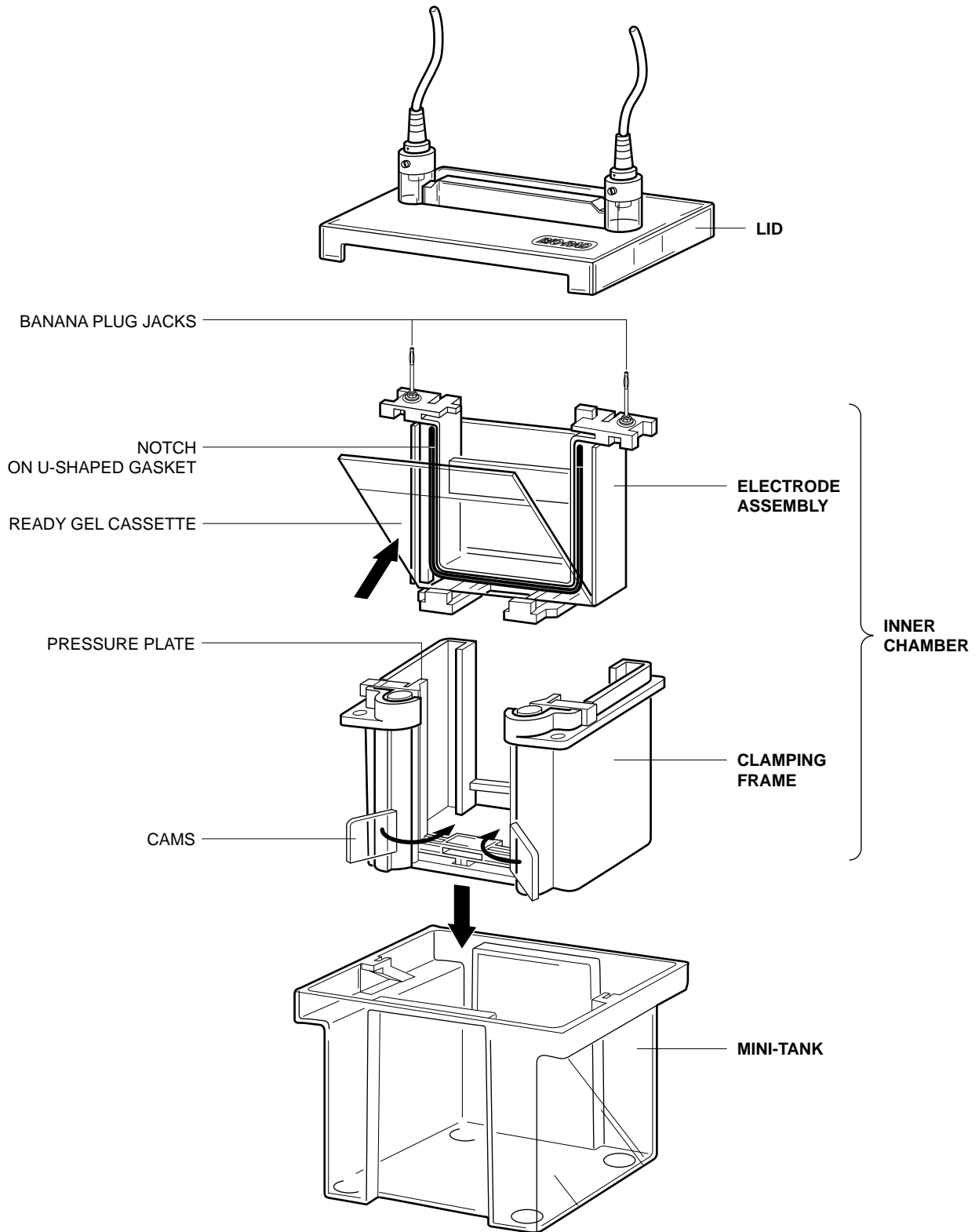
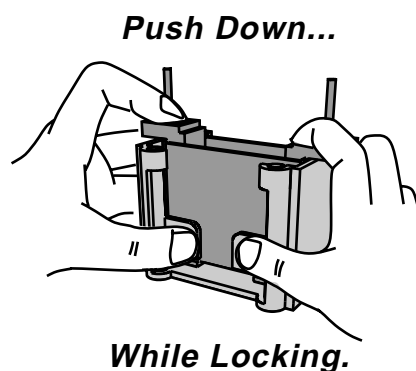


Figure 2. Assembling the Ready Gel Cell

4. Place the Ready Gel cassettes into the slots at the bottom of each side of the Electrode Assembly. Be sure that the short glass plate of the Ready Gel cassette faces inward toward the notches on the green U-shaped gaskets.
5. Press the Ready Gel cassettes up against the gaskets. The Ready Gel cassettes and the Electrode Assembly fit together to form the inner chamber.
6. Transfer the Electrode Assembly and gels into the Clamping Frame.
7. Press down on the Electrode Assembly while closing the two cam levers of the Clamping Frame. See Figure 3.

**Note:** Gently pressing the top of the Electrode Assembly forces the top of the short glass plate on each Ready Gel cassette to seat against the rubber gasket. If you do not push down, the upper buffer may leak.



**Figure 3. Sealing the Inner Chamber**

8. Lower the Electrode Assembly and Clamping Frame into the mini tank. Fill the inner chamber with approximately 125 ml of running buffer, so that the buffer reaches a level between the tops of the short and long plates of the Ready Gel.

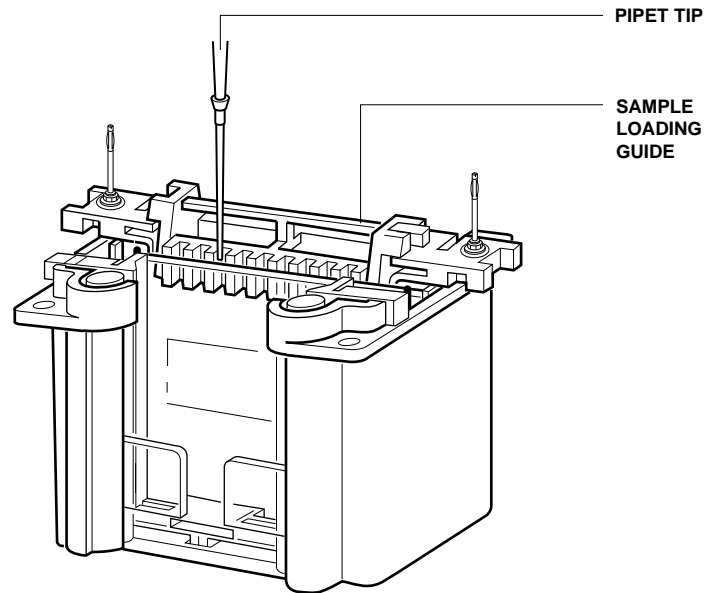
**Note:** Do **not** overfill the inner chamber so that buffer spills over the top of the long plate. Overfilling may cause siphoning of the buffer, resulting in buffer loss and interruption of the electrophoresis.

9. Add approximately 200 ml of running buffer to the mini tank.



## 2.2 SAMPLE LOADING

1. Load the samples into the wells with a Hamilton syringe or a pipet using gel loading tips.
2. If you are using a Sample Loading Guide, place the guide between the two gels in the Electrode Assembly. The guide rests across the top of the Electrode Assembly.



**Figure 4. Using the Sample Loading Guide**

3. Use the Sample Loading Guide to locate the sample wells. Insert the Hamilton syringe or pipet tip into the slots of the guide and fill the corresponding wells.

**Note:** Load samples slowly and allow them to settle evenly on the bottom of the well. Be careful not to puncture the bottom of the well with the syringe needle or pipet tip.

## **2.3 RUNNING THE GEL**

1. Align the electrode plugs and jacks and place the lid on top of the mini tank. The correct orientation is made by matching the colors of the plugs on the lid with the jacks on the electrode assembly. A stop on the lid prevents it from being orientated incorrectly.
2. Attach the electrical leads to a suitable power supply (150 V minimum ) with the proper polarity.
3. Apply power to the Ready Gel Cell and begin electrophoresis; 200 volts, constant is recommended. No adjustment of the setting is necessary for the number of gels. The usual run time is approximately 35 minutes.

## **2.4 REMOVING THE GEL**

1. After electrophoresis is complete, turn off the power supply and disconnect the electrical leads.
2. Remove the cell lid and carefully lift out the Electrode Assembly and Clamping Frame. Pour off and discard the buffer in the inner chamber.

**Note:** Always pour off the upper buffer before opening the cams. Otherwise the buffer will flow out of the chamber.

3. Open the cams of the clamping box. Pull the Electrode Assembly out of the Clamping Frame and remove the gel cassettes.
4. To remove the gel from the Ready Gel cassette, slice the tape along the sides of the Ready Gel cassette where the inner glass plate meets the outer plastic plate.

Separate the plates of the cassette and float the gel off the glass plate by inverting the gel and plate under fixative solution, agitating gently until the gel separates from the plate.

5. The Ready Gel Cell's electrode assembly, Clamping Frame, and mini tank should be rinsed thoroughly with distilled, deionized water after every use.

## 3.0 MAINTENANCE

---

Bio-Rad recommends that all Ready Gel Cell components and accessories be cleaned with suitable laboratory cleaner such as Bio-Rad Cleaning Concentrate, and rinsed thoroughly with distilled, deionized water before use.

**Note:** Ready Gel Cell components are not compatible with acetone or ethanol. Use of organic solvents voids all warranties. Call 1-800-4BIORAD or your local Bio-Rad representative, for technical information regarding chemical compatibility of the Ready Gel Cell components with various laboratory reagents.

## APPENDIX A. SPECIFICATIONS

---

<b>Clamping Frame</b>	Glass-filled liquid crystal polymer
<b>Pressure Plate and Cams</b>	Polycarbonate
<b>Electrode Assembly</b>	Glass-filled liquid crystal polymer
<b>Mini Tank and Lid</b>	Molded polycarbonate
<b>Electrodes</b>	Platinum wire, 0.010 inches diameter
<b>Gasket, electrode inner core</b>	Silicone Rubber
<b>Shipping Weight</b>	2.0 kg
<b>Overall Size</b>	16 cm (L) x 12 cm (W) x 18 cm (H)
<b>Ready Gel Size</b>	7 cm x 8 cm
<b>Voltage Limit</b>	600 VDC and 15 watts

## APPENDIX B. WARRANTY AND ORDERING INFORMATION

---

The Ready Gel Cell is warranted for 1 year against defects in materials and workmanship. If any defects should occur during this warranty period, Bio-Rad Laboratories will replace the defective parts without charge. However the following defects are specifically excluded:

1. Defects caused by improper operation.
2. Repairs or modifications done by anyone other than a Bio-Rad Laboratories or their authorized agent.
3. Damaged caused by accidental misuse.
4. Damage caused by disaster.
5. Common replacement parts including platinum wire and the rubber gasket.
6. Damage caused by the use of organic solvents.

For inquiry or request for repair service, contact your local Bio-Rad office.

### **Bio-Rad Laboratories**

2000 Alfred Nobel Drive  
Hercules, California 94547  
Phone: (510) 741-1000  
1-(800) 4-BIORAD  
1-(800) 424-6723  
Fax: (510) 741-5800 or  
1-(800) 879-2289  
Telex: 335-358

### **Canada**

Bio-Rad Laboratories Ltd.  
5671 McAdam Road  
Mississauga, Ontario L4Z 1N9  
Phone: (905) 712-2771  
1 (800) 268-0213  
Fax: (905) 712-2990

### **India**

Bio-Rad Laboratories Ltd.  
C-248 Defence Colony  
New Delhi 110 024  
Phone: (91-11) 461 0103  
Fax: (91-11) 461 0765  
(91-11) 4662 1863

### **Singapore**

Bio-Rad Laboratories Pte Ltd. (Singapore)  
221 Henderson Road  
05-19 Henderson Building  
Singapore 0315  
Phone: 65-2729877  
Fax: 65-2734835

### **Australia**

Bio-Rad Laboratories Pty., Ltd.  
Block Y, Unit 1  
Regents Park Industrial Estate  
391 Park Road  
Regents Park  
New South Wales 2143  
Phone: 61-2-9914-2800  
Fax: 61-2-9914-2889  
Telex: 79070166

### **Peoples Republic of China**

Bio-Rad China (Beijing Office)  
14 Zhi Chun Road  
HaiDian District  
Beijing 100 088  
Phone: 86-10-2051875  
Fax: 86-10-2051876

### **Italy**

Bio-Rad Laboratories S.r.l  
Via Cellini, 18A  
20090 Segrate - Milano  
Phone: 02/21609.1  
Fax: 02/21609.399

### **Spain**

Bio-Rad Laboratories, S.A.  
Avda, Valdelaparra,  
3 Poligono Industrial de Alcobendas  
E-28100 Alcobendas (Madrid)  
Phone: 34-1-6617085  
(900) 100 204  
Fax: 34-1-6619698

### **Austria**

Bio-Rad Laboratories Ges.m.b.H.  
Auhofstrasse 78D  
A-1130 Wien  
Phone: (1)-877 89 01  
Fax: (1)-876-56-29  
Telex: 13-6565

### **France**

Bio-Rad S.A.  
94/96 rue Victor Hugo  
B.P. 220  
94203 Ivry Sur Seine  
Cedex  
Phone: 01-4960-6834  
Fax: 01-4671-2467

### **Japan**

Nippon Bio-Rad Laboratories KK  
7-18 Higashi-Nippori 5 chome  
Arakawa-ku  
Tokyo 116  
Phone: 03-5811-6271  
Fax: 03-5811-6272

### **Sweden**

Bio-Rad Laboratories AB  
Box 1267  
S-171 24 Solna  
Gardsvagen 7D  
Phone: 46 8 735 83 00  
020-660 660  
Fax: 46 8 735 54 60

### **Belgium**

Bio-Rad Laboratories S.A.-N.V.  
Begoniastraat 5  
B-9810 Nazareth Eke  
Phone: 09-385-55-11  
Fax: 09-385-65-54

### **Germany**

Bio-Rad Laboratories GmbH  
Heidemannstrasse 164  
D-80939 Munchen  
Postfach 45 01 33  
D-80901 Munchen  
Phone: 089 318 84-0  
Fax: 089 318 84-100

### **Netherlands**

Bio-Rad Laboratories B.V.  
Fokkerstraat 10  
3905 KV Veenendaal  
Phone: 0318-540666  
Fax: 0318-542216

### **Switzerland**

Bio-Rad Laboratories, A.G.  
Kanalstrasse 17 Postfach  
CH-8152 Glattbrugg  
Phone: 41-1-8095555  
Fax: 41-1-8095500

### **Hong Kong**

Bio-Rad Pacific Ltd.  
Unit 1111, 11/F., New Kowloon Plaza  
38 Tai Kok Tsui Road  
Tai Kok Tsui, Kowloon  
Phone: 852-2789-3300  
Fax: 852-2789-1257

### **New Zealand**

Bio-Rad Laboratories Pty., Ltd.  
Unit 15 Poland Court  
21 Poland Road  
P.O. Box 100-051  
Glenfield, Auckland  
Phone: 09-443 3099  
0508-805 500  
Fax: 09-443 3097

### **United Kingdom**

Bio-Rad Laboratories, Ltd.  
Bio-Rad House  
Maylands Avenue  
Hemel Hempstead  
Hertfordshire HP2 7TD  
Phone: 01442-232552  
0800-181134  
Fax: 01442-259118

## WARRANTY INFORMATION

Model \_\_\_\_\_

Catalog Number \_\_\_\_\_

Date of Delivery \_\_\_\_\_

Serial Number \_\_\_\_\_

Invoice Number \_\_\_\_\_

Purchase Order No \_\_\_\_\_

## ORDERING INFORMATION

---

Catalog Number	Description
165-3125	<b>Ready Gel Cell</b> , includes Ready Gel Cell Electrode Assembly, Clamping Frame, mini tank, lid with power cables, mini cell buffer dam and instructions.
165-3126	<b>Ready Gel Cell Module</b> , includes Electrode Assembly, Clamping Frame, mini cell buffer dam and instructions.
165-3130	<b>Mini Cell Buffer Dam</b>
165-3146	<b>Sample Loading Guide, 10 well</b>
165-3132	<b>Sample Loading Guide, 15 well</b>
165-3149	<b>Ready Gel Cell Replacement Gaskets</b>
165-2975	<b>Mini Tank and Lid with Power Cables</b>
900-7680	<b>Platinum Wire</b> (Cathode = 8", Anode =11")

---



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