



# **GS Gene Linker<sup>®</sup> UV Chamber**

## **Instruction Manual**

For Technical Service  
Call Your Local Bio-Rad Office or  
in the U.S. Call **1-800-4BIORAD**  
(1-800-424-6723)

## Warranty Statement

Bio-Rad Laboratories GS Gene Linker chamber is warranted against defects in materials and workmanship for 1 year. If any defects occur in the instrument during this warranty period, Bio-Rad will repair or replace the defective parts free of charge. The following defects, however, are specifically excluded:

1. Defects caused by improper operation.
2. Damage caused by accident or misuse.
3. Damage caused by disaster.
4. Corrosion due to use of improper solvents or detergents.
5. Repair or modification performed by anyone other than Bio-Rad Laboratories or authorized agent.
6. Use of spare parts supplied by anyone other than Bio-Rad Laboratories or authorized agent.

The following components are not covered under warranty:

1. Fuses
2. Bulbs

**Note:** There are no user-serviceable parts or components inside the unit. Unauthorized service or repair may void the warranty.

For any inquiry regarding instrument operation, call Bio-Rad Technical Service at 1-(800) 4BIORAD, or contact your local representative.

For any request for repair service, determine the model number and serial number of your instrument, the purchase order number and invoice number, and in the U.S. call Bio-Rad Instrument Service at 1-(800) 876-7614.

## Table of Contents

<b>Section 1</b>	<b>Introduction</b> .....	1
<b>Section 2</b>	<b>Specifications</b> .....	1
2.1	Safety .....	1
2.2	Specifications .....	2
2.3	Components .....	3
<b>Section 3</b>	<b>Operating Instructions</b> .....	4
3.1	Set Up .....	4
3.2	Start/Stop .....	4
3.3	Modes .....	4
3.4	Monitoring .....	5
3.5	Quick Reference Chart .....	6
<b>Section 4</b>	<b>Methods</b> .....	7
4.1	DNA Crosslinking .....	7
4.2	Sterilization .....	8
4.3	DNA Nicking .....	8
<b>Section 5</b>	<b>Cleaning and Maintenance</b> .....	10
5.1	Cleaning .....	10
5.2	Bulb Replacement .....	10
<b>Section 6</b>	<b>Equipment and Accessories</b> .....	10
6.1	Gene Linker UV Chamber and Accessories .....	10
6.2	Electrophoresis Reagents and Equipment .....	10
6.3	Transfer Reagents and Equipment .....	11
<b>Section 7</b>	<b>References</b> .....	12
7.1	UV Crosslinking of DNA to Membrane .....	12
7.2	UV Crosslinking of DNA to Protein .....	12
7.3	UV Nicking of DNA .....	13
7.4	UV Footprinting of DNA .....	13
7.5	Repair of UV-Induced DNA Damage .....	13
7.6	UV Crosslinking of RNA to Membrane .....	14
7.7	UV Crosslinking of RNA to Protein .....	14
7.8	Cellular UV .....	15
7.9	UV-Induced Mutagenesis .....	15
7.10	UV Sterilization of PCR Products .....	15

## Section 1 Introduction

The GS Gene Linker UV chamber is useful for all molecular biology applications requiring UV irradiation. The GS Gene Linker chamber has three different operational modes: Program, Energy, and Time. The Program mode has preset energy or time parameters for specific applications. You can use one of these preset programs, or you can set your own desired energy level or time using the Energy mode or the Time mode. This manual provides information on how to operate the GS Gene Linker chamber along with valuable application guidelines.

## Section 2 Specifications

### 2.1 Safety



#### Definition of Symbols



Caution, risk of UV exposure



Caution (refer to accompanying document)

### Warning

This instrument is intended for laboratory use only.

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



#### Lamps

The lamps inside this instrument emit short wave UV radiation. Overexposure to direct or reflected UV light can cause severe damage to the eyes and skin. Never look into an illuminated UV lamp without proper eye protection.

Ozone may be formed near UV lamps. Excessive ozone exposure can cause eye irritation and discomfort in the respiratory tract. Operate the GS Gene Linker chamber in an adequately ventilated area if ozone is detected by measurement or odor.

#### Sensor

The UV emission sensor is located on the right-hand inside wall. The UV sensor detects only the energy output in the range of 200-400 nm, and will automatically turn off the bulbs when the desired amount of cumulative UV energy has been delivered. The UV sensor also measures maximum energy output from the bulbs, and the Replace Bulb light will illuminate when the energy output is too low (approximately 30%). For these reasons, the sensor should remain clear and clean for accurate energy readings (refer to Section 5 for maintenance).

## Safety Features

<i>Power Switch</i>	The power switch can be used at any time to stop any of the operations.
<i>Start/Stop</i>	The Start/Stop button will stop any of the on-going operations. If pressed again the operation will start at the beginning of the cycle.
<i>Open Door</i>	The GS Gene Linker chamber will not operate when the chamber door is open. If the door is open, the Door Open light will be on. When the door is properly closed the Door Open light will turn off. The chamber will stop operating when the door is opened, but will resume operation when the door is closed.
<i>Chamber</i>	The sealed chamber protects you from any UV radiation. The viewing window will protect your eyes and body from UV radiation.
<i>Maximum Operation</i>	The GS Gene Linker chamber is equipped with an automatic shut off after 999 seconds of energy emission if the detector is covered, or 24 minutes if the radiometer check is not turned off.

## 2.2 Specifications

<b>Dimensions</b>	<b>Width x Depth x Height</b>
<b>Outside</b>	42.7 x 30.5 x 26.4 cm (16.8 x 12 x 10.4 inches)
<b>Inside</b>	31.7 x 24.1 x 15.2 cm (12.5 x 9.5 x 6 inches)
<b>Weight</b>	10 kg (22 lb)

### Functional

<b>Input voltage range</b>	100 VAC/50 Hz/1 amps 120 VAC/60 Hz/1 amps 220 VAC/50 Hz/0.5 amps 240 VAC/50 Hz/0.5 amps
<b>Fuses</b>	2.0 amp Slow-Blow (100/120 V) or 1.0 amp Slow-Blow (220/240 V) Type T
<b>Environmental</b>	
Operating	50 ° F (10 °C) to 90 °F (32 °C) temperature 30–80% humidity
Storage	32 ° F (0 °C) to 140 °F (60 °C) temperature 10–90% humidity

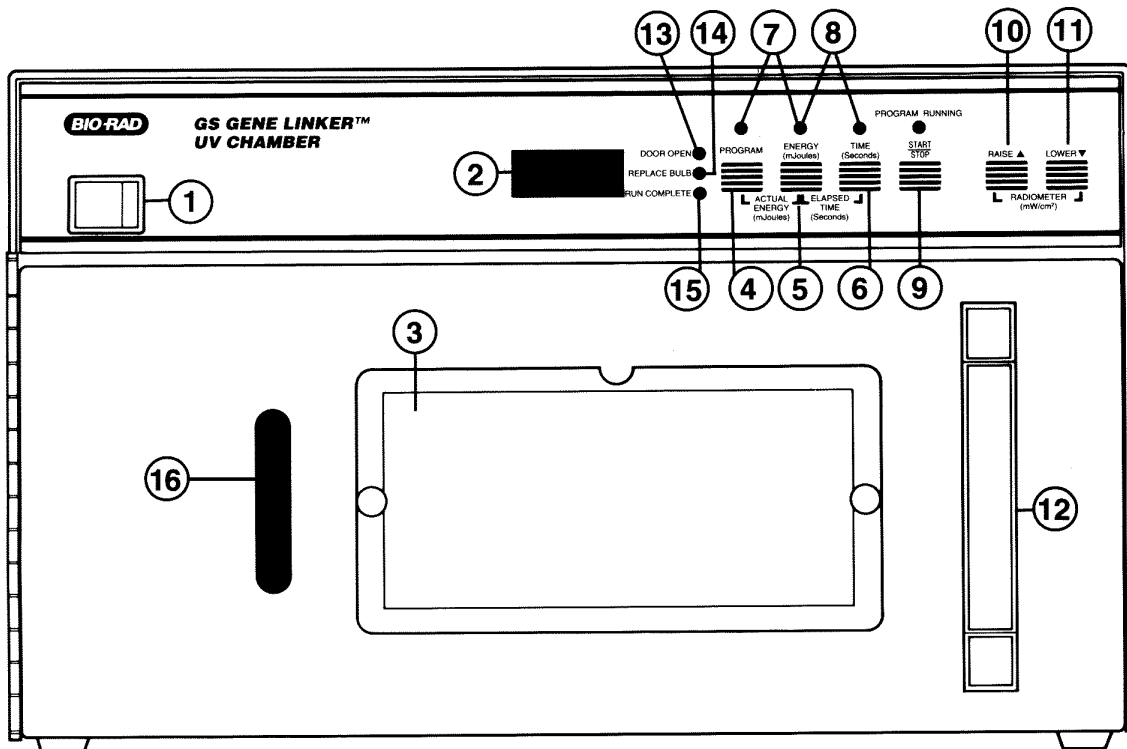
### UV Energy Source

<b>Germicidal bulbs</b>	(5) G8T5 format, minibipin
<b>Output</b>	253.7 nm energy maximum

**Note:** This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to Part 15 of the FCC rules. These limits provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference in which case the user will be required to correct the interference at his own expense.

## 2.3 Components

- |  |   |
|--|---|
| <ol style="list-style-type: none"> <li>1. Power switch</li> <li>2. LED display</li> <li>3. Reference chart</li> <li>4. Program</li> <li>5. Energy</li> <li>6. Time</li> <li>7. Actual energy</li> <li>8. Elapsed time</li> <li>9. Start/Stop</li> <li>10. Raise</li> <li>11. Lower</li> <li>12. Handle</li> <li>13. Door light</li> <li>14. Bulb light</li> <li>15. Run light</li> <li>16. Window</li> </ol> | <p>Power switch to turn the GS Gene Linker chamber On or Off.</p> <p>Selected Program, Energy, or Time is displayed here.</p> <p>Quick reference on preset programs.</p> <p>Program button is used to select the program mode.</p> <p>Energy button is used to select the energy mode. Energy is measured in milliJoules.</p> <p>Time button is used to select the time mode. Time is measured in seconds.</p> <p>Program light and Energy light are both on when the cumulative energy is displayed on the LED.</p> <p>Energy light and Time light are both on when the elapsed time is displayed on the LED.</p> <p>Start/Stop button is used to start the irradiation cycle or to stop a cycle.</p> <p>Increases selected mode.</p> <p>Decreases selected mode.</p> <p>To open and close chamber door.</p> <p>Door Open light indicates the chamber door is open.</p> <p>Replace Bulb light indicates the UV bulbs need replacement.</p> <p>Run Complete light is on when the irradiation cycle has ended.</p> <p>Viewing window</p> |
|--|---|



## **Section 3**

### **Operating Instructions**

#### **3.1 Set Up**

1. Plug the chamber power cord into the appropriate electrical outlet.
2. Press the power switch On. The digital display should be **C-L**. If the display shows **ERR**, the chamber is malfunctioning and needs repair. Contact your local Bio-Rad representative, or call Bio-Rad Instrument Service at 1-800 876-7614.

#### **3.2 Start/Stop**

The Start/Stop button is used to start a given irradiation cycle. When a cycle is started, the Start/Stop button can be used to stop the cycle. The cycle parameters will remain in the temporary memory and can be restarted at the beginning of the cycle by pressing the Start/Stop button again.

The light above the Start/Stop button will turn on when in use and turn off after the irradiation cycle has ended.

#### **3.3 Modes**

The GS Gene Linker chamber can be operated in 3 different modes: Program, Time, or Energy. The mode is indicated by the light above these buttons. The following instructions describe how to set and use each of these modes.

##### **Program**

Each Program is set in either energy or time. To see the preset program value, press the Energy button to display the energy for this program, or press the Time button to display the time for this program. The LED will display - - - when the Time or Energy parameter is not preset. Pressing the Raise or Lower buttons while viewing the set Time or Energy switches out of the Program mode and into the indicated mode.

1. Place the material to be irradiated inside the GS Gene Linker chamber. Use filter paper or plastic wrap to support membranes. Close the chamber door.
2. Press the Program button. Select the desired program by pressing the Raise or Lower button.
3. Start the selected program by pressing the Start/Stop button. The light above the Start/Stop button will be on during the irradiation. The LED will display Actual Energy or Elapsed Time with the two lights above the buttons indicating which units are being displayed.
4. The GS Gene Linker unit will automatically stop after reaching the set time or the set energy level. At the end of the program, the GS Gene Linker chamber will sound a tone for a few seconds. This tone can be stopped by pressing any button. The light above the Start/Stop button will be off and the Run Complete light will be on. The LED will display either energy or time indicated by the light above the button.

## **Energy**

1. Place the material to be irradiated inside the GS Gene Linker chamber. Close the chamber door.
2. Press the Energy button. Select the desired energy level by pressing the Raise or Lower button.
3. Start the selected energy level by pressing the Start/Stop button. The light above the Start/Stop button will be on during the irradiation. The LED will display the Actual Energy with the light above the Program and Energy buttons on.
4. The GS Gene Linker unit will automatically shut down after reaching the set energy level. At the end of the irradiation cycle, the GS Gene Linker chamber will sound a tone for a few seconds. This tone can be stopped by pressing any button. The light above the Start/Stop button will be off and the Run Complete light will be on. The LED will display cumulative energy.

## **Time**

1. Place the material to be irradiated inside the GS Gene Linker chamber. Close the chamber door.
2. Press the Time button. Select the desired irradiation time by pressing the Raise or Lower button.
3. Start the selected time by pressing the Start/Stop button. The light above the Start/Stop button will be on during the irradiation. The LED will display the Elapsed Time and the light above the Time and Energy buttons will be on.
4. The GS Gene Linker unit will automatically shut down after reaching the set time. At the end of the irradiation cycle the GS Gene Linker chamber will sound a tone for a few seconds. The tone can be stopped by pressing any button. The light above the Start/Stop button will be off and the Run Complete LED will be on. The LED will display elapsed time.

## **3.4 Monitoring**

Any time during or after an irradiation cycle, the elapsed time or cumulative energy can be determined.

1. To determine the elapsed time, press both the Energy button and the Time button simultaneously. The elapsed time in seconds will appear on the LED.
2. To determine the cumulative energy, press both the Energy button and the Program button simultaneously. The cumulative energy in milliJoules will appear on the LED.
3. Press the Energy or Time button alone and the LED will display the set values.



### 3.5 Quick Reference Chart

The GS Gene Linker chamber quick reference chart reproduced below, is located behind a plexiglass shield on the chamber door. This chart can be used as an applications guide. Bio-Rad will publish updated charts as new programs or protocols are developed. To receive these updates, it is important for you to be listed in the GS Gene Linker customer data base. Contact your Bio-Rad representative to be listed.

#### GS Gene Linker UV Chamber Quick Reference Chart

Make sure material is in the chamber and the door is properly closed. Turn ON power. The LED display should read C-L. Select operation mode by pressing the Program, Energy, or Time button. Press the Raise or Lower button to set the desired program, energy level or time. Press the Start/Stop button to start the irradiation cycle. At the completion of the cycle, a tone will sound and the UV light will automatically turn off.

Application	Conditions	Program* (LED Reading)	Setting
Crosslinking	Dot blot/NaOH NH <sub>4</sub> OAc dry Zeta-Probe	C-L	125 mJoule
Nicking	Pulsed field gels	nic	60 mJoule
Sterilization	UV resistant material	Str	90 sec
Crosslinking	Dot blot/damp Zeta-Probe	C1	30 mJoule
Crosslinking	Southern dry membrane	C2	50 mJoule
Crosslinking	Southern damp membrane	C3	150 mJoule
Crosslinking	Dot blot/NaOH NH <sub>4</sub> OAc dry membrane 312 nm	C4	250 mJoule

*Program (LED Reading)	Section
C-L	4.1
nic	4.2
Str	4.1
C1	4.1
C2	4.3
C3	4.3
C4	[312 nm bulbs (to be released)]

## Section 4 Methods

Bio-Rad Laboratories has developed protocols for UV crosslinking of nucleic acids to nylon membrane, UV-induced nicking of DNA prior to transfer to membrane, and UV sterilization. We have an ongoing research program to develop new protocols for molecular biology applications. Our studies indicate that the energy required for optimal crosslinking of nucleic acid to membrane is dependent upon several parameters. These include the type of membrane (charged vs. neutral nylon), the transfer buffer, whether the membrane is wet or dry, as well as the application (dot blot vs. genomic Southern). The following protocols are recommended, depending upon the application.

### 4.1 DNA Crosslinking

#### Slot or Dot Blot/damp Zeta-Probe membrane

1. Denature the DNA in 0.4 N NaOH final concentration.
2. Assemble the microfiltration apparatus according to the manufacturer's recommendations. Use a positively charged nylon membrane like Zeta-Probe® GT membrane.
3. Apply the DNA samples directly onto the membrane.
4. Remove the membrane from the apparatus and place the damp membrane inside the GS Gene Linker chamber. Use filter paper or plastic wrap to support the membrane.
5. Follow the instructions in Section 3.3, Program mode. Select the program with the LED reading **C1** (30 mJ). Start the irradiation cycle by pressing the Start/Stop button. Proceed with preferred hybridization protocol.

#### Slot or Dot Blot/NaOH NH<sub>4</sub>OAc/dry Zeta-Probe Membrane

1. Denature the DNA sample by adding NaOH and EDTA to a final concentration of 0.4 N NaOH and 10 mM EDTA. Heat the sample to 100 °C for 10 minutes.
2. Neutralize the DNA sample by adding an equal volume of cold 2 M ammonium acetate (NH<sub>4</sub>OAc), pH 7.0.
3. Assemble the microfiltration apparatus according to the manufacturer's recommendations. Use a positively charged nylon membrane like Zeta-Probe GT membrane.
4. Apply the DNA samples directly onto the membrane.
5. Remove the membrane from the apparatus and place the damp membrane between 2 filter papers. Allow the membrane to dry. Remove the top filter paper and place membrane inside the GS Gene Linker chamber.
6. Follow the instructions in Section 3.3, Program mode. Select the program with the LED reading **C-L** (125 mJ). Start the irradiation cycle by pressing the Start/Stop button. Proceed with preferred hybridization protocol.

**Note:** When using a neutral membrane select program **C3** (150 mJ).

#### Southern Transfer

1. Depurinate the DNA by soaking the gel in 0.25 N HCl for 10-15 minutes.
2. Denature the DNA by placing the gel in a bath of 0.5 N NaOH for 30 minutes.
3. Neutralize the gel by soaking it in 0.5 M Tris-HCl, pH 7.4, 1 M NaCl for 30 minutes.

4. Transfer the DNA onto a nylon membrane using 10x SSC or 10x SSPE as the transfer buffer. The membrane can be a neutral nylon or a positively charged nylon membrane like Zeta-Probe GT membrane.
5. After transfer, rinse the membrane in 2x SSC for 5 minutes. The membrane can be damp or dry.
6. Follow the instructions in Section 3.3 Program mode. Select the program with the LED reading **C2** (50 mJ) if the membrane is dry, or **C3** (150 mJ) if the membrane is damp. Start the irradiation cycle by pressing the Start/Stop button. Proceed with preferred hybridization protocol.

## 4.2 Sterilization

The GS Gene Linker chamber has a sterilization cycle preset. Follow the instructions in Section 3.3 Program mode. Select the third program with the LED reading **Str** (90 sec). Start the irradiation cycle by pressing the Start/Stop button. The GS Gene Linker chamber will sterilize only the exposed areas on any object in the chamber.

## 4.3 DNA Nicking<sup>25</sup>

The transfer efficiency of large DNA (>20 kb) from agarose gels is poor unless the DNA is nicked prior to transfer to nylon membrane. The DNA can be cleaved either by HCl depurination or by UV-irradiation. The depurination reaction is harder to control and is extremely sensitive to temperature. Exposure to short wavelength UV light (254 nm) is a more reliable method for nicking DNA in pulsed field gels before transfer. For optimal results, the following protocol must be followed rigorously.

1. Stain the gel with 1.0 µg/ml ethidium bromide (EtBr) for exactly 30 minutes with constant agitation. Do not destain the gel prior to nicking. Use a fresh dilution of EtBr for each gel.
2. Place the gel on a glass tray for transporting the gel to the GS Gene Linker chamber.
3. Place the gel inside the chamber and close the door.
4. Follow the instructions in Section 3.3 Program mode. Select the second program with the LED reading **nic** (60 mJ).
5. Start the irradiation cycle by pressing the Start/Stop button. The gel can be photographed, but exposure to UV radiation must be minimized (< 10 seconds). The gel can be destained prior to photography if desired.
6. Soak the gel in 0.4 N NaOH, 1.5 M NaCl for 15 minutes.
7. Transfer the DNA onto Zeta-Probe GT membrane using two liters of 0.4 N NaOH, 1.5 M NaCl as the transfer solvent.
8. Set up the capillary transfer as follows, from bottom to top:
  - A. Corning Pyrex glass dish (28 x 18 x 4 cm).
  - B. A plexiglass or plastic box for support, about 3 cm high and small enough to fit in the glass dish (*e.g.*, Eppendorf yellow pipette tip rack).
  - C. Glass Plate (16 x 20 cm).
  - D. Three sheets of blotting paper as wick (18 x 30 cm) (S&S, GB002).

- E. Agarose gel (well side down).
  - F. Zeta-Probe membrane cut to the same size as the gel and pre-wetted with distilled water.
  - G. Three sheets of blotting paper (18 x 15 cm) (S&S, GB002).
  - H. A stack of paper towels 10 cm thick.
9. Transfer the DNA for 24–48 hours.
  10. Carefully remove the paper towel and blotting papers. Remove the membrane together with the gel, turn over the membrane and gel, lay them gel side up, and mark the location of the wells and the orientation marker on the top of the gel. The position of the wells can be accurately marked on the membrane by using a fine point permanent alcohol marker pen, cutting through the bottoms of the wells.
  11. Neutralize the membrane in 0.5 M Tris, pH 7.0 (neutralization buffer) for 5 minutes followed by rinsing briefly in 2x SSC. Transferred DNA can be visualized on the membrane by placing the damp blot on a transilluminator.
  12. Dry the membrane by blotting on 3MM or other adsorbent paper and proceed to hybridization. UV crosslinking of the DNA to the membrane is not recommended with this alkaline transfer method.

### **Discussion**

1. The procedure is based on gels approximately 6 mm thick. If thicker gels are used, the staining period may be prolonged to allow diffusion of EtBr into the middle of the gels. DNA that is not stained with EtBr will not be nicked by the UV light and thus will not be transferred from the gel.
2. Presoaking the gel in NaOH prior to transfer decreases background and increases transfer efficiency.
3. Pulsed field gels can also be blotted onto membranes using 10x SSC as the transfer buffer with standard alkaline denaturation followed by neutralization. Alkaline transfer onto nylon membranes gives as good or better sensitivity as standard transfers onto nitrocellulose filters. The alkaline procedure is much simpler and faster. In addition, nylon membranes can be reused many more times than nitrocellulose filters. Some blots may be reused as many as twenty times.
4. DNA separated on the CHEF-DR<sup>®</sup> II or CHEF Mapper<sup>®</sup> system can also be vacuum transferred onto nylon membrane in 4 hours using the Model 785 Vacuum Blotter (165-5001, 120 V) and NaOH as the transfer buffer.
5. The DNA is transferred from the back of the gel (the side opposite the wells) onto the membrane because irregularities in the surface of the gel frequently occur during solidification of these high percentage gels (1%). These surface artifacts will interfere with the transfer of the DNAs from the gel. Transfer from the other side of the gel insures smooth surface contact between the gel and the membrane.
6. It is essential to neutralize the membrane after transfer to prevent changing the pH of the hybridization buffer during the hybridization.
7. It is not necessary to bake nylon membranes after alkaline transfer since the DNA should be fixed onto the membrane by NaOH.

- To monitor the efficiency of the transfer, stain the gel in neutralization buffer for 30 minutes with 1.0  $\mu\text{g/ml}$  EtBr. Photograph the post-transferred gel and compare with the original picture.

## **Section 5**

### **Cleaning and Maintenance**

#### **5.1 Cleaning**

*Outside* - Clean the outside of the GS Gene Linker UV chamber with a damp towel. Do not use solvents or strong detergents.

*Inside* - Clean the inside aluminum walls of the chamber with ethanol (reagent grade) and a soft towel. The sensor is located on the right-hand inside wall. Clean the sensor with a soft towel and ethanol. Be careful not to scratch the sensor with an abrasive towel.

#### **5.2 Bulb Replacement**

The GS Gene Linker chamber is equipped with a radiometer mode. This mode is used to determine the bulb brightness. Enter into the radiometer mode by pressing both the Raise and Lower buttons at the same time. The UV lights will turn on and the LED will display an 'L' along with the UV illuminance measured in  $\text{mW/cm}^2$ . After 1 minute, the reading should stabilize. If the UV illuminance is  $\leq 2 \text{ mW/cm}^2$  ALL 5 bulbs should be replaced. Stop the radiometer mode by pressing any button.

Turn off the GS Gene Linker chamber power and unplug the power cord. Remove the bulb by holding either metal end and turn the bulb 1/4 turn. The bulb will slip out of the socket through the slots on the bottom. Insert the new bulb through the slots and turn the bulb 1/4 turn until the bulb snaps into place.

## **Section 6**

### **Equipment and Accessories**

#### **6.1 GS Gene Linker UV Chamber and Accessories**

165-5031	<b>GS Gene Linker UV Chamber</b> , 120 VAC/60 Hz; includes five 254 nm bulbs and instruction manual
165-5032	<b>GS Gene Linker UV Chamber</b> , 220 VAC/50 Hz; includes five 254 nm bulbs and instruction manual
165-5033	<b>GS Gene Linker UV Chamber</b> , 240 VAC/50 Hz; includes five 254 nm bulbs and instruction manual
165-5034	<b>GS Gene Linker UV Chamber</b> , 100 VAC/50 Hz; includes five 254 nm bulbs and instruction manual
165-5035	<b>GS Gene Linker UV Chamber Replacement Bulbs</b> , 254 nm, 5 bulbs

#### **6.2 Electrophoresis Reagents and Equipment**

170-4300	<b>Sub-Cell DNA Electrophoresis Cell without Casting Tray</b>
170-4304	<b>Sub-Cell DNA Electrophoresis Cell with 15x20 cm Casting Tray</b>
170-4343	<b>Wide Mini-Sub Cell DNA Electrophoresis Cell Basic Unit</b>
170-4307	<b>Mini-Sub Cell DNA Electrophoresis Cell Basic Unit</b>

## **Electrophoresis Reagents and Equipment (Continued)**

165-5052      **PowerPac 200 Power Supply, 100/120V**

165-5053      **PowerPac 200 Power Supply, 220/240V**

### **Ultra Pure DNA Agaroses**

162-0017      **Low Melt Preparative Grade Agarose, 25 gm**

162-0020      **Low Melt Preparative Grade Agarose, 250 gm**

162-0125      **High Strength Analytical Grade Agarose, 100 gm**

162-0126      **High Strength Analytical Grade Agarose, 500 gm**

162-0133      **Molecular Biology Certified Agarose, 100 gm**

162-0134      **Molecular Biology Certified Agarose, 500 gm**

162-0135      **Chromosomal Grade Agarose, 25 gm**

162-0136      **Chromosomal Grade Agarose, 100 gm**

161-0733      **Premix 10x TBE Buffer, 1 liter**

## **6.3 Transfer Reagents and Equipment**

165-5000      **Model 785 Vacuum Blotter with Regulator**, includes: Vacuum Regulator; Base unit with Vacuum Stage; Porous Vacuum Plate; Reservoir Seal O-ring; Sealing Frame; Assortment Window Gaskets; Vacuum Blotter lid; Instruction manual

165-5001      **Model 785 Vacuum Blotter System**: same as 165-5000 except with vacuum pump (120 V)

165-5002      **Model 785 Vacuum Blotter System**: same as 165-5000 except with vacuum pump (220/240 V)

165-5003      **Model 785 Vacuum Blotter Basic Unit**

170-6545      **Bio-Dot Apparatus**

170-6542      **Bio-Dot Slot Format**

170-3910      **Trans-Blot Electrophoretic Transfer Cell**

162-0190      **Zeta-Probe GT Membrane**, sheets, 9 x 12 cm, 15

162-0191      **Zeta-Probe GT Membrane**, sheets, 10 x 15 cm, 15

162-0192      **Zeta-Probe GT Membrane**, sheets, 15 x 15 cm, 15

162-0193      **Zeta-Probe GT Membrane**, sheets, 15 x 20 cm, 15

162-0194      **Zeta-Probe GT Membrane**, sheets, 20 x 20 cm, 15

162-0195      **Zeta-Probe GT Membrane**, sheets, 20 x 25 cm, 3

162-0196      **Zeta-Probe GT Membrane**, roll 30 cm x 3.3 m, 1

162-0197      **Zeta-Probe GT Membrane**, roll 20 cm x 3.3 m, 1

162-0198      **Zeta-Probe GT Membrane**, roll 30 cm x 30 m, 1

170-3557      **Random Primer DNA Labeling Kit**, 25 reactions

165-0962      **Filter Paper Backing**, 35 x 45 cm, 25 sheets

165-0921      **Filter Paper Backing**, 18 x 34 cm, 25 sheets

## Section 7

### References

#### 7.1 UV Crosslinking of DNA to Membrane

1. Allefs, J.J.H.M., Salentijn, E.M.J., Krens, F.A. and Rouwendal, G.J.A., Optimization of Non-radioactive Southern Blot Hybridization: Single Copy Detection and Reuse of Blots. *Nucleic Acids Research*, **18**, (10) 3099-3100. (1990).
2. Cannon, Gordon, Heinhorst, Sabine, and Weissbach, Arthur, Quantitative Molecular Hybridization on Nylon Membranes. *Anal. Biochem.* **149**, 229-237. (1985).
3. Church, George M. and Gilbert, Walter, Genomic Sequencing. *Proc. Natl. Acad. Sci.* **81**, 1991-1995. (1984).
4. Khandjian, E.W., Optimized Hybridization of DNA Blotted and Fixed to Nitrocellulose and Nylon Membranes. *BioTech.*, **5**, 165-167 (1987).
5. Knight, Pamela., Nucleic Acid and Protein Blotting. *BioTech.* **8**, 166-167 (1990).
6. Nierzwicki-Bauer, Sandra A., Gebhardt, Joan S., Linkkila, Leslie and Walsh, Kieron, A Comparison of UV Cross-linking and Vacuum Baking for Nucleic Acid Immobilization and Retention. *BioTech.* **9**, 472-478 (1990).
7. Razin, S.V., Yarovaya, O.V. and Georgiev, G.P. Low Ionic Strength Extraction of Nuclease-Treated Nuclei Destroys the Attachment of Transcriptionally Active DNA to the Nuclear Skeleton. *Nucleic Acids Research*, **13**, 7427-7444 (1985).
8. Twomey, Tara A. and Krawetz, S.A. Parameters Affecting Hybridization of Nucleic Acids Blotted onto Nylon or Nitrocellulose Membranes. *BioTech* **8**, 478-481 (1990).
9. Saluz, Hanspeter and Jost, Jean-Pierre, Optimized Genomic Sequencing as a Tool for the Study of Cytosine Methylation in the Regulatory Region of the Chicken Vitellogenin II Gene. *Gene* **42**, 151-157 (1986).

#### 7.2 UV Crosslinking of DNA to Protein

10. Molitor, J. A., Walker, W., Doerre, S., Ballard, D.W., and Greene, W.C., NF-kB: A Family of Inducible and Differentially Expressed Enhancer-Binding Proteins in Human T Cells. *Proc. Natl. Acad. Sci.*, **87**, 10028-10032 (1990).
11. Gilmour, David S., Deitz, Thomas J. and Elgin, Sara C.R, UV Cross-linking Identifies Four Polypeptides That Require the TATA Box to Bind to the Drosophila hsp70 Promoter. *Mol. and Cell. Bio.*, **10**, 4233-4238 (1990).
12. Dooley, Steven, Welter, Cornelius, Theisinger, Brigit and Blin, Nikolaus, Generating Highly Labeled Oligonucleotides for DNA-Protein Interaction. *GATA*, **7**(5), 133-137 (1990).
13. Meffert, Rainer, Rathgeber, Gabriele, Schader, Hans-Jochen and Dose, Klaus, UV-Induced Cross-linking of Tet Repressor to DNA Containing Tet Operator Sequences and 8-Azidoadenines. *Nucleic Acids Research*, **18** (22), 6633-6636 (1990).
14. Becker, Michael M., Lesser, David, Kurpiewski, Michael, Baranger, Anne and Jen-Jacobson, Linda, Ultraviolet Footprinting Accurately Maps Sequence-Specific Contacts and DNA Kinking in the EcoRI Endonuclease - DNA Complex. *Proc. Natl. Acad. Sci.*, **85**, 6247-6251 (1988).
15. Schnapp, Andreas, Clos, Joachim, Hadelt, Wolfgang, Schreck, Ralf., Cvekl, Ales and Grummt, Ingrid, Isolation and Functional Characterization of TIF-IB, a Factor that Confers Promoter Specificity to Mouse RNA Polymerase I. *Nucleic Acids Research*, **18**(6), 1385-1393 (1990).
16. Schneider, Harald R., Waldschmidt, Rainer and Seifart, Klaus H., Human Transcription Factor III C Contains a Polypeptide of 55 kDA Specifically Binding to Pol III Genes. *Nucleic Acids Research*, **18**(16), 4743-4750 (1990).
17. Pfeleiderer, Christa, Smid, Amke, Bartsch, Ingrid and Grummt, Ingrid, An Undecamer DNA Sequence Directs Termination of Human Ribosomal Gene Transcription. *Nucleic Acids Research*, **18**(16), 4727-4736 (1990).

18. Kaufman, J.D., Valanda, G., Rodriguez, G., Bushar, G., Giri, C., Norcross, MA., *Mol. Cell Biol.*, **7**, 3759-3766, (1987).
19. Nabel, G. and Baltimore, D., *Nature*, **326**, 711-713, (1987).
20. Tong-Starksen, S.E., Luciw, P.C., Peterlin, B.M., *Proc. Natl. Acad. Sci. USA*, **84**, 6845-6849, (1987).
21. Rosen, C.A., Sodroski, J.G., Haseltine, W.A., *Cell*, **41**, 813-823, (1985).
22. Spandidos, D.A., Yiagnisis, M., Pintzas, A., *Anticancer Res.*, **9**, 383-386, (1989).
23. Meffert, R. and Dose, K., *FEBS Letters*, **239**, 190-194 (1988).

### 7.3 UV Nicking of DNA

24. Harding, Fiona A., Cohen, Nicholas and Litman, Gary W., Immunoglobulin Heavy Chain Gene Organization and Complexity in the Skate, *Raja erinaces*. *Nucleic Acids Research* **18**(4), 1015-1020 (1990).
25. Personal communication with Eric Lai University of North Carolina, Chapel Hill.

### 7.4 UV Footprinting of DNA

26. Uchida, Kiyoshi, Pyle, Anna Marie, Morii, Takashi and Barton, Jacqueline K., High Resolution Footprinting of EcoRI and Distamycin with Rh(phi)<sub>2</sub> (bpy)<sup>3+</sup>, a New Photofootprinting Reagent. *Nucleic Acids Research*, **17**(24), 10259-10279 (1989).

### 7.5 Repair of UV-Induced DNA Damage

27. Freeman, Ph.D., Steven, E., Gange, M.D., Richard W., Sutherland, Ph.D., John C., Matzinger, B.A., Ezra A. and Sutherland, Ph.D., Betsy M., Production of Pyrimidine Dimers in DNA of Human Skin Exposed *in Situ* to UVA Radiation. *The Journal for Investigative Dermatology*, **88**, 430-433 (1987).
28. Gekik, Catherine M. and Collins, Andrew R., Comparison of Effects of Fostriecin, Novobiocin, and Camptothecin, Inhibitors of DNA Topoisomerases, on DNA Replication and Repair in Human Cells. *Nucleic Acids Research*, **18**(4), 1007-1013 (1990).
29. Lyamichev, Victor I., Frank-Kamenetskii, Maxim D. and Soyfer, Valery N., Protection Against UV-Induced Pyrimidine Dimerization in DNA by Triplex Formation. *Nature*, **344**, 568-570 (1990).
30. Smerdon, Michael J., Bedoyan, Jirair and Thoma, Fritz., DNA Repair in a Small Yeast Plasmid Folded into Chromatin. *Nucleic Acids Research*, **18**(8), 2045-2051 (1990).
31. Chu, Gilbert and Chang, Elaine., Cisplatin-Resistant Cells Express Increased Levels of a Factor That Recognizes Damaged DNA. *Natl. Acad. Sci.*, **87**, 3324-3327 (1990).
32. Freeman, Steven E. and Thompson, Bryan D., Quantitation of Ultraviolet Radiation-Induced Cyclobutyl Pyrimidine Dimers in DNA by Video and Photographic Densitometry. *Analytical Biochemistry*, **186**, 222-228 (1990).
33. Tomkinson, Alan E., Bonk, R. Thomas., Kim, Joon., Bartfeld, Neil and Linn, Stuart., Mammalian Mitochondrial Endonuclease Activities Specific for Ultraviolet-Irradiated DNA., *Nucleic Acids Research*, **18**(4), 929-935 (1989).
34. Madura, Kiran, Prakash, Satya and Prakash, Louise, Expression of the *Saccharomyces cerevisiae* DNA Repair Gene RAD6 that Encodes a Ubiquitin Conjugating Enzyme, Increases in Response to DNA Damage and in Meiosis but Remains Constant During the Mitotic Cell Cycle, *Nucleic Acids Research*, **18**(4), 771-778 (1990).
35. Tsujimura, Tohru, Maher, Veronica M., Godwin, Alan R., Liskay, Michael and McCormick, J. Justin, Frequency of Intrachromosomal Homologous Recombination Induced by UV Radiation in Normally Repairing and Excision Repair-Deficient Human Cells. *Proc. Natl. Acad. Sci.*, **87**, 1566-1570 (1990).
36. Venema, J., Mullenders, L.H.F., Natarajan, A.T., van Zeeland, A.A. and Mayne, L.V., The Genetic Defect in Cockayne Syndrome is Associated With a Defect in Repair of UV-Induced DNA Damage in Transcriptionally Active DNA. *Proc. Natl. Acad. Sci.*, **87**, 4707-4711 (1990).



37. Madura, Kiran and Prakash, Satya, Transcript Levels of *Saccharomyces cerevisiae* DNA Repair Gene RAD23 Increase in Response to UV Light and in Meiosis but Remain Constant in the Mitotic Cell Cycle. *Nucleic Acids Research*, **18**(16), 4737-4742 (1990).
38. Mitchell, David, L., Brash, Douglas E. and Nairn, Rodney S., Rapid Repair Kinetics of Pyrimidine(6-4)Pyrimidine Photoproducts in Human Cells are Due to Excision Rather Than Conformational Change. *Nucleic Acids Research*, **18**(4), 963-971 (1990).
39. Ohnishi, Takeo, Yuba, Shunsuke, Date, Takayasu, Utsumi, Hiroshi and Matsukage, Akio, Rat DNA Polymerase B Gene Can Join in Excision Repair of *Escherichia coli*. *Nucleic Acids Research*, **18**(19), 5673-5676 (1990).

## 7.6 UV Crosslinking of RNA to Membrane

40. Fingerle, J., Johnson, R., Clowes, A.W., Majesky, M.W. and Reidy, M.A. Role of uin Smooth Muscle Cell Proliferation and Migration After Vascular Injury in Rat Carotid artery. *Proc. Natl. Acad. Sci.*, **86**, 8412-8416 (1989).
41. Herrin, D.L. and Schmidt, G.W., Rapid, Reversible Staining of Northern Blots Prior to Hybridization. *Bio/Tech*, **6**, 196-198 (1988).
42. Khandjian, Edouard W. and Meric, Claude A., Procedure for Northern Blot Analysis of Native RNA. *Anal. Biochem.*, **159**, 227-232 (1986).
43. Kroczek, R.A., Mediate Visualization of Blotted RNA in Northern Analysis. *Nucleic Acids Research.*, **17**, 9497 (1989).
44. Lin, Leu-Fen, H., Mismar, Drzislav, Lile, Jack D., Armes, Lyman G., Butler, Eugene T., Vannice, James L., Collins, Frank, Purification, Cloning and Expression of Ciliary Neurotrophic Factor (CNTF). *Science*, 1023-1025 (1989).
45. Williams, Simon C., Bruckheimer, Sally M., Lusic, Aldons, J., LeBoeuf, R.C. and Kinniburgh, Alan J., Mouse Apolipoprotein A-IV Gene: Nucleotide Sequence and Induction by a High-Lipid Diet. *Mol. Cell. Bio.* **6**, 3807-3814 (1986).
46. Kroczek, Richard A. and Siebert, Erwin, Optimization of Northern Analysis by Vacuum-Blotting, RNA-Transfer Visualization and Ultraviolet Fixation. *Analytical Biochemistry*, **184**, 90-95 (1990).
47. Farrell, Jr., Robert E., Methodologies for RNA Characterization II: Quantitation by Northern Blot Analysis and the S1 Nuclease Assay. *Clinical Biotechnology*, **2**(2), 107-119 (1990).

## 7.7 UV Crosslinking of RNA to Protein

48. Garcia-Blanco, Mariano A., Anderson, Gordon J., Beggs, Jean and Sharp, Phillip A., A Mammalian Protein of 220 kDa Binds Pre-mRNAs in the Spliceosome: A Potential Homologue of the Yeast PRP8 Protein. *Proc. Natl. Acad. Sci.*, **87**, 3082-3086 (1990).
49. Marciniak, Robert A., Garcia-Blanco, Mariano A. and Sharp, Phillip A., Identification and Characterization of a HeLa Nuclear Protein that Specifically Binds to the Trans-Activation-Response (TAR) Element of Human Immunodeficiency Virus. *Proc. Natl. Acad. Sci.*, **87**, 3624-3628 (1990).
50. Leibold, Elizabeth, A., Laudano, Andrew and Yu, Yang., Structural Requirements of Iron-Responsive Elements for Binding of the Protein Involved in Both Transferring Receptor and Ferritin mRNA Post-Transcriptional Regulation. *Nucleic Acids Research*, **18**(7), 1819-1824 (1990).
51. Siebel, Christian W. and Rio, Donald C., Regulated Splicing of the *Drosophila P* Transposable Element Third Intron *in Vitro*: Somatic Repression. *Science*, **248**, 1200-1208 (1990).

## 7.8 Cellular UV

52. Satokata, Ichiro, Tanaka, Kiyoji, Miura, Naoyuki, *et al.*, Characterization of a Splicing Mutation in Group A Xeroderma Pigmentosum. *Proc. Natl. Acad. Sci.*, **87**, 9908-9912 (1990).
53. Lawrence, C.W., Borden, A., Banerjee, S.K. and LeCler, J.C., Mutation Frequency and Spectrum Resulting From a Single Abasic Site in a Single-Stranded Vector. *Nucleic Acids Research*, **18**(8), 2153-2157 (1990).

## 7.9 UV-Induced Mutagenesis

53. Armstrong, John D. and Kunz, Bernard A., Site and Strand Specificity of UVB Mutagenesis in the SUP4-0 Gene of Yeast. *Proc. Natl. Acad. Sci.*, **87**, 9005-9009 (1990).

## 7.10 UV Sterilization of PCR Products

54. Rich, Jessica J., and Willis, David, A Single Oligonucleotide Can be Used to Rapidly Isolate DNA Sequences Flanking a Transposon Tn5 Insertion by the Polymerase Chain Reaction, *Nucleic Acids Research*, **18**(22), 6673-6676 (1990).
55. Sarkar, Gobinda, and Sommer, Steve S., Parameters Affecting Susceptibility of PCR Contamination to UV Inactivation, *Biotechniques* **10**(5), 591-593 (1991).

The PCR process is covered by U.S. patent numbers 4,683,195, 4,683,202, and 4,899,818 which are owned by Hoffmann-La Roche, Inc. and F. Hoffmann-La Roche, Ltd. The purchase of this product does not convey a license to use the process covered by these patents. The user of this product to perform PCR must obtain a license from Hoffmann-La Roche, Inc.



## Bio-Rad Laboratories

### Life Science Group

2000 Alfred Nobel Drive  
Hercules, California 94547  
Telephone (510) 741-1000  
Fax: (510) 741-5800  
[www.bio-rad.com](http://www.bio-rad.com)

**Australia**, Bio-Rad Laboratories Pty. Ltd., Block Y, Unit 1, Regents Park Industrial Estate, 391 Park Road, Regents Park, NSW 2143  
Phone 02 9914 2800 • Fax 02 9914 2889

**Austria**, Bio-Rad Laboratories Ges.m.b.H., Auhofstraße 78D, A-1130 Wien • Phone (01)-877 89 01 • Fax (01)-876 56 29

**Belgium**, Bio-Rad Laboratories S.A.-N.V., Begoniastraat 5, B-9810 Nazareth • Phone 09-385 55 11 • Free Phone 0800/97032 • Fax 09-385 65 54

**Brazil**, Bio-Rad Laboratories (Brazil), Rua dos Invalidos 212 - 5 andar, Lapa - Rio de Janeiro - RJ, CEP 20331-020 • Phone 55 21 507 6191

**Canada**, Bio-Rad Laboratories (Canada) Ltd., 5671 McAdam Road, Mississauga, Ontario L4Z 1N9 • Phone (905) 712-2771 • Fax (905) 712-2990

**China**, Bio-Rad China (Beijing), Rm 615, Shang Fang Plaza, No. 27, North Third Round Center Road, West District, Beijing 100029  
Phone 86-10-8201-1366/68 • Fax 86-10-8201-1367

**Denmark**, Bio-Rad Laboratories, Generatorvej 8 C, 2730 Herlev • Phone 45 44 52-1000 • Fax 45 44 52-1001

**Finland**, Bio-Rad Laboratories, Pihatörmä 1A, FIN-02240 Espoo • Phone 358 (0)9 804 2200 • Fax 358 (0)9 804 1110

**France**, Bio-Rad Laboratories, 3, Boulevard Raymond Poincaré, 92430 Marnes-la-Coquette • Phone 01 47 95 69 65 • Fax 01 47 41 9133

**Germany**, Bio-Rad Laboratories GmbH, Heidemannstraße 164, D-80939 München, Postfach 45 01 33, D-80901 München  
Phone 089 318 84-177 • Fax 089 318 84-123

**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

**India**, Bio-Rad Laboratories (India) Pvt. Ltd., B&B1, Enkay Towers Vanijyanikunj, Udhvy Vihar Phase V, Gurgaon, Haryana 122016  
Phone (91-124)-6398112/113/114 • Fax (91-124)-6398115

**Israel**, Bio-Rad Laboratories, Ltd., 14 Homa Street, P.O. Box 5044, Rishon Le Zion 75150 • Phone 03 951 4124 • Fax 03 951 4129

**Italy**, Bio-Rad Laboratories S.r.l., Via M. Peroglio 23, 00144 Rome • Phone 34 91 590 5200 • Fax 34 91 590 5211

**Japan**, Nippon Bio-Rad Laboratories KK, 7-18 Higashi-Nippori 5-chome, Arakawa-ku Tokyo 116-0014 • Phone 03-5811-6270 • Fax 03-5811-6272

**Korea**, Bio-Rad Korea Ltd., Cambridge Building, 1461-15 Seocho-Dong Seocho-Ku, Seoul 137-070 • Phone 82-2-3473-4460 • Fax 82-2-3472-7003

**Latin America**, Bio-Rad Latin America, 14100 Palmetto Frontage Road, Suite 101, Miami Lakes, Florida USA 33016 • Phone 305-894-5950 • Fax 305-894-5960

**Mexico**, Bio-Rad Laboratorios Mexico, Adolfo Prieto No. 1653, Col. De Valle, CP. 03100, Mexico D.F. • Phone 52 5 534 2552 to 54 • Fax 52 5 524 5971

**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., PO Box 300-571, Albany, Auckland • Phone 64-9-4152280 • Fax 64-9-443 3097

**Norway**, Bio-Rad Laboratories, Johan Scharffenbergs vei 91, N-0694 Oslo • Phone 47-23-38-41-30 • Fax 47-23-38-41-39

**Russia**, Bio-Rad Laboratorii, ul. Butirskaya 79 "B", office 156 RF-125015 Moscow • Phone 7 095 979 98 00 • Fax 7 095 979 98 56

**Singapore**, Bio-Rad Laboratories, Singapore, 211 Henderson Rd. #03-02, Henderson Industrial Park, 159552 • Phone 65-2729877 • Fax 65-2734835

**Spain**, Bio-Rad Laboratories, S.A., Lopez de Hoyos, 245-247, 28043 Madrid • Phone 34-91-590-5200 • Fax 34-91-590-5211

**Sweden**, Bio-Rad Laboratories AB, Vintergatan 1, Box 1097, S-172 22 Sundbyberg • Phone 46 (0)8-55 51 27 00 • Fax 46 (0)8-55 51 27 80

**Switzerland**, Bio-Rad Laboratories AG, Nenzlingerweg 2, CH-4153 Reinach • Phone 061-717-9555 • Fax 061-717-9550

**United Kingdom**, Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead, Hertfordshire HP2 7TD  
Phone 0181 328 2000 • Free Phone 0800-181134 • Fax 01442-259118