

WHITE LIGHT TRANSILLUMINATOR 2000

Instruction Manual

Catalog numbers

170-8124

and

170-7943

WARRANTY

The White Light transilluminator 2000 is warranted against defects in materials and workmanship for 1 year. If any defect occurs in the instrument or accessories during this warranty period, Bio-Rad Laboratories will repair or replace the defective parts at its own discretion without charge. The following defects, however, are specifically excluded:

1. Defects caused by improper operation.
2. Repair or modification done by anyone other than Bio-Rad Laboratories or an authorised agent.
3. Damage caused by substituting alternative parts.
4. Use of fittings or spare parts supplied by anyone other than Bio-Rad Laboratories.
5. Damage caused by accident or misuse.
6. Damage caused by disaster.
7. Corrosion caused by improper solvent or sample.

This warranty does not apply to parts listed below:

- Fuses
- Lamps
- Starters

The warranty general conditions are indicated in the attached warranty card that has to be filled in and sent back.

For any inquiry or request for repair service, contact your local Bio-Rad Laboratories office. Inform Bio-Rad of the model and serial number of your instrument.

REGULATORY NOTICE

IMPORTANT: This Bio-Rad instrument is designed and certified to meet EN 61010* safety standards and EMC regulations. 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 safety and EMC certification
- Create a potential safety hazard

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

*EN 61010 is an internationally accepted electrical safety standard for laboratory instruments.

IMPORTANT NOTICE

Before installing this unit, please read carefully the installation instructions.

This instrument is intended for clinical and research laboratory use and must be operated only by specialised personnel aware of the potential risks connected to the chemical - biological agents normally used with this unit.

Contact your local Bio-Rad Laboratories representative if service is required.

INSTALLATION

Unpacking

Carefully unpack the transilluminator as follows:

- First remove the cardboard inside the box.
- Remove the transilluminator from the two cardboard shells and place it on a stable, horizontal surface.
- Remove the blue protection film from the WL Filter.

Stand alone installation

- The instrument must be placed on a bench leaving at least 10 cm of space all around in order to avoid any obstacle that may reduce the ventilation.
- Connect the instrument to the power using the annexed cable. The power font must be able to deliver at least 250 VA with a voltage from 100 to 240 Vac. The plug must have a ground connection.
- Position the acrylic face shield sliding it below the transilluminator.

**The instrument is supplied in two versions: 230 Vac or 115 Vac.
Before connecting the cable to the power, verify that the voltage selectr is set to the right power.**

Installation with the Bio-Rad Gel Doc 2000 :

This instrument has been designed in order to be lodged inside the Gel Doc 2000 – Gel Documentation System and with the Instadoc system.

In this case follow the instructions included in the Gel Doc 2000 or Instadoc manuals or any other instrument that will be released in the future.

SWITCHING ON

To switch ON the instrument press the switch «Power ON/OFF».

The instrument is also provided with a thermal protection in order to avoid to exceed 50°C inside the instrument.

If the green key of the «Power ON/OFF» is lighted without that the lamps light it could be activated the thermal protection, this happens after an uninterrupted use of at least 1 hour . Switch Off the key and wait for cooling of the instrument.

LAMPS AND/OR STARTER REPLACEMENT

Attention: if substances dangerous for the health are used the instrument must be cleaned and treated for a perfect decontamination. When the instrument is open it would be better to use protection gloves.

All the internal circuits are fed at 230Vac independently by the line voltage through an autotransformer.

To ensure a good instrument running replace the lamps every 500 hours of working.

- Decontaminate the instrument as per the instruction given in the chapter “General precaution”;
- Remove the power cable;
- Unscrew out the screws on both sides of the instrument;
- Remove the lid putting it down on the back side making attention to the grounding cable now connected (do not put the filter directly on the bench);
- Remove the lamp rotating it until the lateral pins be in a vertical position;
- In case of replacement of a starter, remove it by rotating anticlockwise;
- Insert the new lamp into the lamp holders and rotate it until the pins be in a horizontal position;
- To replace the starter insert it in the holder and rotate clockwise;
- Reassemble the lid so that the grounding cable slides under the reflecting panel and retighten the screws on both sides.

It is essential to assemble the lid before making any test.
Follow in any case the instructions in the chapter SWITCHING ON.

FUSES REPLACEMENT

The fuses are on the side of the power module entry on the right lateral panel. It's very easy to replace them following this procedure:

- Remove the power cable;
- Unscrew the plastic screws of the fuse holder with a screw driver;
- Replace the fuses with others of suitable power (0,8A for 230-240Vac e 1,6A for 100 and 115 Vac).
- Retighten the fuse holder screws.

GENERAL PRECAUTIONS

- Connect the instrument **ONLY** with electric line connected to the ground or to the Gel Doc 2000.
- The instrument is equipped with a thermal protection for its overheating. During the normal use (5-15 min.) the instrument doesn't reach the temperature to switch off the thermal protection; if the instrument is left switched on for at least 1 hour the thermal protection automatically switch off the lamps (internal temperature around 48 °C) and the green light of the main switch remains lighted. When the temperature come back to 38°C the lamps turn on again.
- Do not pour out liquids directly on the instrument.
- Do not blank the aeration slit.
- Switch off the instrument immediately after its use.
- **Clean well the filter after its use. Avoid alcohol or organic solvents that can damage the plastic filter. Use only aqueous solutions of soaps and mild detergents Always wear disposable gloves.**

SPECIFICATIONS

Electrical data

Voltage 100/230 Vac 50/60 Hz +/- 10%

Fuses 0.8/1.6 Amp

Consumption 120 VA max

Physical data

Dimensions

Width 28 cm

Length 36 cm

Height 10 cm

Weight 9,5 kg

Operating conditions

Temperature 5-40°C

Relative Humidity max 80

Manufactured by :

Ordering Information

170-7943 WL Transilluminator 2000 (100-115 Vac/60 Hz)

Including : Instruction Manual

Power Cable

Spare Fuses 1.6 Amps

170-8124 WL Transilluminator 2000 (220-240 Vac/50 Hz)

Including : Instruction Manual

Power Cable

Spare Fuses 0.8 Amps

Bio-Rad Laboratories Srl

Via Cellini 18/A

20090 SEGRATE (Mi) - Italy

Tel. +39-2-21609.1

Fax. +39-2-21609.399

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
Australia, Bio-Rad Laboratories Pty Limited, Unit 11, 112-118 Talavera Rd P.O. Box 371, North Ryde, NSW 2113 • Phone 02-805-5000 • Fax 02-805-1920
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 Laboratones 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 Laboratones, Symbion Science Park, Fruebjergvej 3, DK-2100 Copenhagen • Phone 39 17 9947 • Fax 39 27 1698
Finland, Bio-Rad Laboratones, 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, 7-16, 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., 464 Siglap Road, #01-02 Flamingo Valley, Singapore 1545 • Phone (65) 4432529 • Fax (65) 4421667
Spain, Bio-Rad Laboratones, S. A. Avda Valdeiaparra 3, Pol. Ind. Alcobendas, E-28100 Alcobendas, Madrid • Phone (91) 661 70 65 • Fax (91) 661 96 98
Sweden, Bio-Rad Laboratones 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 Laboratones AG, Kanalstrasse 17, Postfach, CH-8152 Glattpfugg • Phone 01-809 55 55 • Fax 01-809 55 00
United Kingdom, Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemei Hempstead, Herts HP2 7TD • Free Phone 0800 181134 • Fax 01442 259118