



Molecular Imager[®] Gel Doc[™] XR System

The Sensible Solution for Uncompromised Gel Documentation

Product Feature

Benefit

Real-time imaging with short refresh interval	<ul style="list-style-type: none"> ▪ One-touch sample positioning and focusing for easy gel documentation
Megapixel resolution in a single snapshot with wide dynamic range and high sensitivity	<ul style="list-style-type: none"> ▪ Accurate data output suitable for publication and a variety of downstream analyses
Epi-white and transmitted light for sample illumination	<ul style="list-style-type: none"> ▪ Wide range of fluorescent and colorimetric applications on gels and blots
FireWire interface	<ul style="list-style-type: none"> ▪ Easy to install and use on Mac and PC computers
Integrated software controls	<ul style="list-style-type: none"> ▪ Simple to set up and begin image acquisition ▪ Easy to customize new applications
Image acquisition and analysis with Quantity One [®] 1-D analysis software (included)	<ul style="list-style-type: none"> ▪ Application-driven workflow ▪ Mac and PC compatible ▪ Image reporting ▪ Comprehensive analysis feature set ▪ Options for CFR 21 Part 11



BIO-RAD

Molecular Imager **Gel Doc XR** System

Ordering Information

Catalog # Description

Molecular Imager Gel Doc XR System

- 170-8170 **Molecular Imager Gel Doc XR System**, PC, includes darkroom, UV transilluminator, epi-white illumination, camera, FireWire card, cables, Quantity One software, instructions
- 170-8171 **Molecular Imager Gel Doc XR System**, Mac, includes darkroom, UV transilluminator, epi-white illumination, camera, FireWire card, cables, USB to serial converter, Quantity One software, instructions

Accessories

- 170-7950 **White Light Transilluminator**, universal hood
- 170-8001 **White Light Conversion Screen**
- 170-8074 **Filter**, 520DF30 62 mm for SYBR Green/GFP/SYBR Gold/fluorescein
- 170-8075 **Filter**, 560DF50 62 mm for Cy3/rhodamine
- 170-8076 **Filter**, 630BP30 62 mm for SYPRO Ruby/Texas Red
- 170-8077 **Filter**, 480BP70 62 mm for Hoechst/coumarin
- 170-8098 **254 nm UV Lamps**, 6
- 170-6887 **365 nm UV Lamps**, 6
- 170-8097 **Standard 302 nm UV Lamps**, 6
- 170-8089 **Mitsubishi P93DW Printer**
- 170-7581 **Mitsubishi Thermal Printer Paper**, 4 rolls
- 170-8067 **Sony UPD895 Printer Paper**, 5 rolls

Software

- 170-9600* **Quantity One 1-D Analysis Software**
- 170-9615 **Quantity One CFR Module**

* Included with imaging system.

Cy is a trademark of GE Healthcare. FireWire and Mac are trademarks of Apple Computer, Inc. Hoechst is a trademark of Hoechst Aktiengesellschaft. Mitsubishi is a trademark of Mitsubishi Corporation. Sony is a trademark of Sony Corporation. SYBR, SYPRO, and Texas Red are trademarks of Molecular Probes, Inc.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 905 712 2771 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000