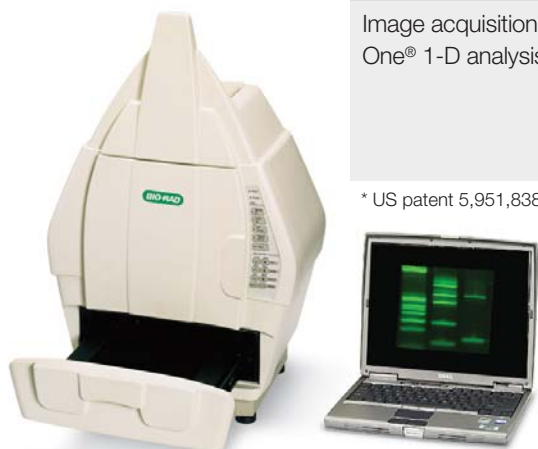


# Molecular Imager® ChemiDoc™ XRS System

## Value-Added Sensitivity for Uncompromised Gel and Blot Imaging

Product Feature	Benefit
Extended integration time for detection of most faint and bright light signals	<ul style="list-style-type: none"> <li>Chemiluminescence imaging with perfect linearity and sensitivity to detect low picogram substrate quantities</li> </ul>
Unique dynamic flat fielding technology (patented*)	<ul style="list-style-type: none"> <li>Perfect uniformity through the entire image area for reliable data from every pixel in the image</li> </ul>
Real-time imaging with short refresh interval	<ul style="list-style-type: none"> <li>One-touch sample positioning and focusing for easy gel documentation</li> </ul>
Megapixel resolution in a single snapshot with wide dynamic range and superior sensitivity	<ul style="list-style-type: none"> <li>Accurate data output suitable for publication and a variety of downstream analyses</li> </ul>
Epi-white and transmitted light for sample illumination	<ul style="list-style-type: none"> <li>Wide range of fluorescent and colorimetric applications on gels and blots in addition to superior chemiluminescence detection</li> </ul>
FireWire interface	<ul style="list-style-type: none"> <li>Easy to install and use on Mac and PC computers</li> </ul>
Integrated software controls	<ul style="list-style-type: none"> <li>Simple to set up and begin image acquisition</li> <li>Easy to customize new applications</li> </ul>
Image acquisition and analysis with Quantity One® 1-D analysis software (included)	<ul style="list-style-type: none"> <li>Application-driven workflow</li> <li>Mac and PC compatible</li> <li>Image reporting</li> <li>Comprehensive analysis feature set</li> <li>Options for CFR 21 Part 11</li> </ul>

\* US patent 5,951,838.



**BIO-RAD**

# Molecular Imager **ChemiDoc XRS** System

## Ordering Information

Catalog # Description

### Molecular Imager ChemiDoc XRS System

- 170-8070 **Molecular Imager ChemiDoc XRS System**, PC, includes darkroom, UV transilluminator, epi-white illumination, fluorescent reference plate, camera, power supply, PCI digitizing card, cables, Quantity One software, instructions
- 170-8071 **Molecular Imager ChemiDoc XRS System**, Mac, includes darkroom, UV transilluminator, epi-white illumination, fluorescent reference plate, camera, power supply, PCI digitizing card, cables, USB to serial converter, Quantity One software, instructions

### Accessories

- 170-7950 **White Light Transilluminator**, universal hood
- 170-8001 **White Light Conversion Screen**
- 170-8072 **f/0.95 25 mm Wide-Angle Lens**
- 170-8073 **f/0.95 17 mm Wide-Angle Lens**
- 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](http://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