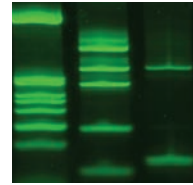
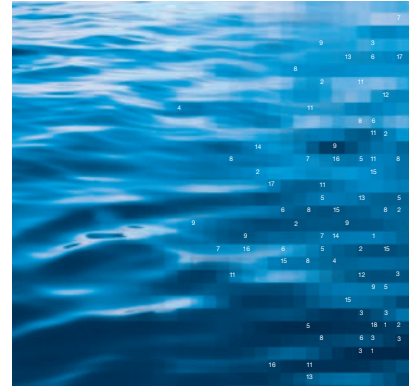
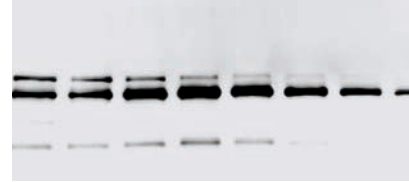


## Imaging



Molecular Imager® **ChemiDoc™ XRS System**



**Extra Resolution and Sensitivity**



**BIO-RAD**

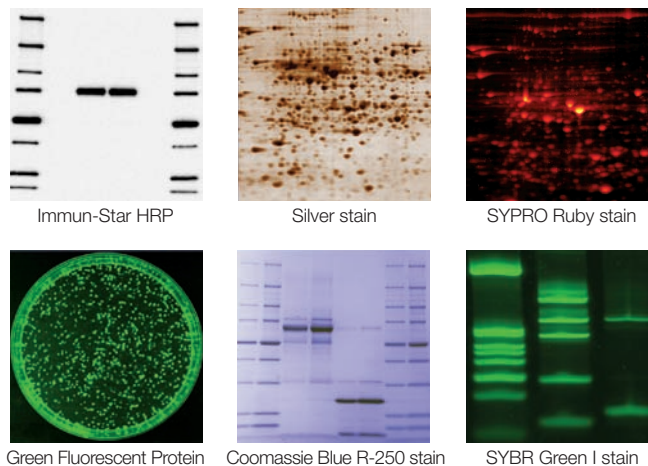
# Molecular Imager ChemiDoc XRS System

The ChemiDoc XRS system is the most advanced chemiluminescence detection system in Bio-Rad's gel documentation line. The system features flat fielding technology for superior uniformity and quantitation, and the flexibility to image chemiluminescent, fluorescent, and colorimetric samples. The system includes a supersensitive CCD, a light-tight compact darkroom, a slide-out transilluminator with preparative mode for added safety when manually visualizing samples, a software-driven motorized lens, and powerful, user-friendly Quantity One® 1-D analysis software. A touch pad gives you total access and control right at the instrument.

The ChemiDoc XRS system can be used for imaging in a wide variety of applications (western blotting, nucleic acid detection, 2-D gel electrophoresis, dot blotting, densitometry, and colony counting) using a large portfolio of detection methods (Immun-Star™ HRP and AP chemiluminescent detection, and ethidium bromide, SYBR Green I, SYPRO Ruby, Cy3, rhodamine, Green Fluorescent Protein, Hoechst, silver stain, copper stain, and Coomassie Brilliant Blue staining).

## Optimized for High-Resolution Chemiluminescence and Fluorescence Imaging

- Supercooled high-resolution CCD camera with 1.44 megapixel resolution
- Real-time imaging
- Software-driven motorized control of zoom, focus, and iris (standard on every system)
- True dynamic flat fielding
- User control of camera bin and gain settings
- Excellent collection efficiency; f/1.2 motorized zoom lens or f/0.95 wide-angle lens
- Exceptional performance with -50°C Peltier-cooled supersensitive camera
- Choice of 3 UV transillumination wavelengths
- Integrated software controls



## Specifications

CCD resolution (H x V)	1,392 x 1,040 µm
Pixel density	65,536 gray levels
Imaging area	25 x 26 cm
Camera cooling temperature	-50°C
Camera cooling system	Peltier
Dynamic range	4 orders of magnitude
Motorized zoom lens	C-mount, f/1.2, 12-75 mm
Operating system compatibility	Windows 2000 and XP, Mac OS X
Illumination modes	Epi-white, trans-UV (302 nm), and white
Dimensions (W x D x H)	60 x 36 x 96 cm
Weight	32 kg (70.4 lb)

## Ordering Information

Catalog #	Description
<b>Molecular Imager ChemiDoc XRS System</b>	
170-8070	Molecular Imager ChemiDoc XRS System, PC, includes darkroom, UV transilluminator, epi-white illumination, fluorescent reference plate, camera, power supply, FireWire cable, 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, FireWire cable, 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
170-8098	254 nm UV Lamps, 6
170-6887	365 nm UV Lamps, 6
170-8097	Standard 302 nm UV Lamps, 6

For optional emission filters or computer systems, please inquire.

Coomassie is a trademark of BASF Aktiengesellschaft. Cy is a trademark of GE Healthcare. Hoechst is a trademark of Hoechst Aktiengesellschaft. FireWire, Mac, and Mac OS are trademarks of Apple Computer, Inc. SYBR and SYPRO are trademarks of Molecular Probes, Inc. Windows and Windows 2000 and XP are trademarks of Microsoft Corp.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 4BIORAD Australia 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 4029300/5013478 Israel 03 963 6050  
Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 55 5200 05 20 The Netherlands 0318 540666 New Zealand 64 9415 2280  
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 095 721 14 04 Singapore 65 6415 3188 South Africa 27 0861 246 723  
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189/2578 7241 United Kingdom 020 8328 2000