



Molecular Imager[®] VersaDoc[™] MP 5000 System

Unsurpassed Sensitivity in Chemiluminescence Plus Multiplex Versatility

Product Feature

Benefit

Unique detection technology for highest sensitivity and widest dynamic range among molecular imaging products

- Reliable imaging with perfect data linearity and precision of the challenging chemiluminescence blots that emit the weakest signals

Red, green, blue, and ultraviolet (UV) sample excitation

- Widest range of fluorophores, including all variations of multiplex fluorescence, can be detected for unlimited flexibility in gel and blot applications

Unique dynamic flat fielding technology (patented*)

- Perfect uniformity through the entire image area for reliable data from every pixel in the image

Epi-white and transmitted light for sample illumination

- Unlimited range of fluorescent and colorimetric applications on gels and blots in addition to superior chemiluminescence detection

Integrated software controls

- 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)

- Application-driven workflow
- Mac and PC compatible
- Image reporting
- Comprehensive analysis feature set
- Options for CFR 21 Part 11

* US patent 5,951,838.



Molecular Imager **VersaDoc MP 5000** System

Ordering Information

Catalog # Description

Molecular Imager VersaDoc MP 5000 System

170-8650 **Molecular Imager VersaDoc MP 5000 System**, PC or Mac, 100/240 V, includes CCD camera, darkroom, power supply, cables, epi-illumination, transillumination, fluorescent reference plate, focusing target, Quantity One software, instructions

Accessories

170-8647 **Red Illumination Modules, 2**
170-8648 **Green Illumination Modules, 2**
170-7706 **Optional Fixed Lens, 105 mm**
170-8001 **White Light Conversion Screen**
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.

Mac is a trademark of Apple Computer, Inc. Mitsubishi is a trademark of Mitsubishi Corporation. Sony is a trademark of Sony Corporation.

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 905 364 3435 **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 6361 7000 **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