



Molecular Imager® VersaDoc™ MP 4000 System

Multiplex Imaging for Ultimate Versatility

Product Feature	Benefit
Red, green, blue, and ultraviolet (UV) sample excitation	<ul style="list-style-type: none"> Widest range of fluorophores, including all variations of multiplex fluorescence, can be detected for unlimited flexibility in gel and blot applications
Extended integration time for detection of most faint and bright light signals	<ul style="list-style-type: none"> Chemiluminescence imaging with perfect linearity and sensitivity to detect low picogram substrate quantities
Unique dynamic flat fielding technology (patented*)	<ul style="list-style-type: none"> Perfect uniformity through the entire image area for reliable data from every pixel in the image
Multimillion pixel resolution in a single snapshot with widest dynamic range	<ul style="list-style-type: none"> Superior resolution and high precision data for the most advanced gel and blot applications including 2-D proteomics and arrays
Epi-white and transmitted light for sample illumination	<ul style="list-style-type: none"> Unlimited range of fluorescent and colorimetric applications on gels and blots in addition to superior chemiluminescence detection
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



* U.S. patent 5,951,838.



Molecular Imager **VersaDoc MP 4000** System

Ordering Information

Catalog # Description

Molecular Imager VersaDoc MP 4000 System

170-8640 **Molecular Imager VersaDoc MP 4000 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

Software

170-9600* **Quantity One 1-D Analysis Software**
170-9615 **Quantity One CFR Module**
170-9620 **PDQuest™ Basic 2-D Analysis Software**
170-9630 **PDQuest Advanced 2-D Analysis Software**
170-9645 **PDQuest Advanced CFR Module**

* Included with imaging system.

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