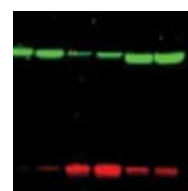
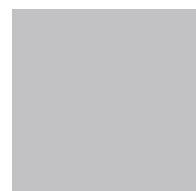


Imaging



Clear Advantages
in High-Resolution
Multiplex Images

Molecular Imager® **VersaDoc™ MP 4000** System



BIO-RAD

The Molecular Imager VersaDoc MP 4000 system — with red, green, and blue excitation, and 3.2 million pixels — offers clear advantages in high-resolution multiplex imaging.

The differences are striking. Multiple illumination and detection modes provide exceptional application versatility. Detect Flamingo™ fluorescent and SYPRO Ruby protein gel stains, Immun-Star™ enzyme-conjugated antibodies, quantum dots, ECL reagents, Cy, Pro-Q, and SYBR dyes, and many more. The multi-megapixel resolution of the VersaDoc MP 4000 system enables you to visualize samples that most imagers cannot resolve.

You can also combine pixels for greater sensitivity, and optimize lens and illumination characteristics for image uniformity and accurate quantitation. All this, along with outstanding ease of use and a clear upgrade path, make the VersaDoc MP 4000 system a definite standout.

Illumination Modes

- Red, green, blue, and white epi-illumination
- Transillumination (UV and white)

Multi-Megapixel Resolution

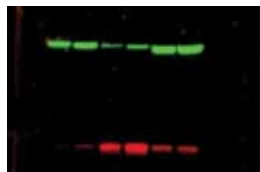
- 3.2 million pixels, with microlens technology and 53 µm resolution

Versatile Detection

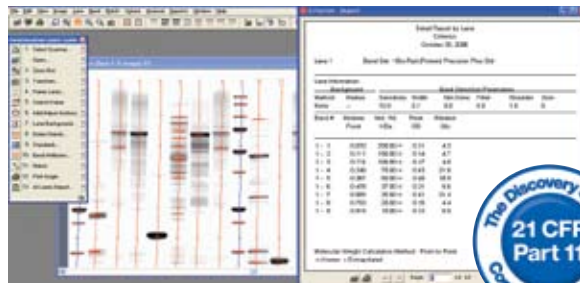
- Fluorescent
- Colorimetric
- Chemiluminescent
- Chemifluorescent

A Wide Range of Applications

- Proteomics and genomics
- 1-D nucleic acid and protein gels
- 2-D protein gels
- Northern, Southern, and western blots
- Dot and slot blots
- Colony counting
- Arrays



Cy3-labeled human β -actin and Cy5-labeled phosphorylated cofilin, imaged with green LED/605BP filter and red LED/695BP filter on VersaDoc MP 4000 system.



Imaging Software

Quantity One® 1-D analysis software and PDQuest™ 2-D analysis software offer sophisticated ways to visualize and analyze your electrophoretic data. Whether quantitating a DNA gel, looking for the absence or presence of a band, determining purity of a protein sample, calculating molecular weight of bands in an unknown sample, or analyzing protein expression, you can rely on comprehensive and sophisticated analyses of 1-D and 2-D samples.

For more information on Bio-Rad's line of gel analysis software, visit us on the Web at

www.bio-rad.com/software/

Ordering Information

Catalog # Description

Molecular Imager VersaDoc MP 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

Imaging Software

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

Bio-Rad has a full range of imaging systems.

For more information, visit us on the Web at

www.bio-rad.com/imaging/

Cy and ECL are trademarks of GE Healthcare. Mac is a trademark of Apple Computer, Inc. Pro-Q, SYBR, and SYPRO are trademarks of Molecular Probes, Inc.

**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