



Personal Molecular Imager™ (PMI™) System

Dedicated to Imaging of Radiolabeled Samples

Product Feature	Benefit
Storage phosphor technology for imaging radioisotopes	<ul style="list-style-type: none"> ▪ The best choice of label for the widest dynamic range and mass spectrometry compatibility
Designed specifically for detection of radiolabeled samples	<ul style="list-style-type: none"> ▪ Best fit for radioisotopes ▪ Safe and controlled laboratory practices
Direct laser excitation of fluorophores and storage phosphors with patented* fiber-optic scanning	<ul style="list-style-type: none"> ▪ Minimizes optical distortion for accurate image data
Self-alignment of external lasers before each scan	<ul style="list-style-type: none"> ▪ Reliable, robust performance; reproducible data
Sample imaging from top by direct contact with scanning beam	<ul style="list-style-type: none"> ▪ Accommodates wide range of samples, including thick agarose gels, microplates, various polyacrylamide gels, blots, and TLC plates
Retractable moisture-sealed trays	<ul style="list-style-type: none"> ▪ Allows imaging of samples of different formats, and facilitates safe and efficient processing of samples
Small footprint (57 x 68 cm)	<ul style="list-style-type: none"> ▪ Optimizes use of limited benchtop
All BaFBr:Eu phosphor screens up to 35 x 43 cm accommodated**	<ul style="list-style-type: none"> ▪ Economical utilization of phosphor screens, new and previously owned
Image acquisition and analysis with Quantity One® 1-D analysis software (included)	<ul style="list-style-type: none"> ▪ Application-directed user interface ▪ Mac and PC compatible ▪ Seamless integration with PDQuest™ and Proteomweaver™ 2-D analysis software and EXQuest™ spot cutter ▪ Options for CFR 21 Part 11



* US patents 4,812,660; 4,822,520; 4,830,875; and 5,266,803.

** Including screens from other manufacturers.



Personal Molecular Imager (PMI) System

Ordering Information

Catalog #	Description
170-9400	Personal Molecular Imager (PMI) System , PC or Mac, 110/240 V, includes Quantity One software, sample tray, USB cable, instructions

Sample Handling Accessories

170-7811	Sample Tray
170-7812	Multi-Sample Tray I , for small aluminum-mounted screens and microplates

Storage Phosphor Screens and Accessories

170-7845	Imaging Screen-K (Kodak)/Tritium , 20 x 25 cm
170-7843	Imaging Screen-K (Kodak) , 20 x 25 cm
170-7841	Imaging Screen-K (Kodak) , 35 x 43 cm
170-7861	Exposure Cassette-K , for 20 x 25 cm Kodak screen
170-7862	Exposure Cassette-K , for 35 x 43 cm Kodak screen
170-7809	Eraser Screen-K , 110/120 V
170-7806	Eraser Screen-K , 220/240 V

Miscellaneous Accessories

170-7869	Replacement Bulb for Eraser Screen-K
161-0722	Bio-Rad Cleaning Concentrate , 50X, 1 kg

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
165-9770	Proteomweaver 4.0 Professional 2-D Analysis Software
165-9780	Proteomweaver 4.0 Enterprise 2-D Analysis Software
165-9798	Proteomweaver 4.0 Enterprise Pro 2-D Analysis Software

* Included with imaging system.

Kodak is a trademark of Eastman Kodak Co. Mac is a trademark of Apple Computer, Inc.

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 905 712 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