



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™ analysis software and EXQuest™ spot cutter ▪ Options for CFR 21 Part 11



* U.S. 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
170-7819	Multi-Sample Tray II, for scanning gels mounted to glass plates

Storage Phosphor Screens and Accessories

170-7845	Imaging Screen-K (Kodak)/Tritium, 20 x 24 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

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

* Included with imaging system.

Kodak is a trademark of Eastman Kodak Company. Mac is a trademark of Apple, Inc.



BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **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 64 9 415 2280 **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 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000