



Molecular Imager[®] PharosFX[™] and PharosFX[™] Plus Systems

Efficiency and Flexibility for Accurate Molecular Imaging

Product Feature	Benefit
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
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
Self-alignment of external lasers before each scan	<ul style="list-style-type: none"> Reliable, robust performance; reproducible data
User-configured custom emission filters in scanner hardware and software	<ul style="list-style-type: none"> Unlimited choice of fluorophores in visible spectrum
Transillumination screen	<ul style="list-style-type: none"> Performs gel documentation for colorimetric stains
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
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
PharosFX Plus 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



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

** Including screens from other manufacturers.



Molecular Imager **PharosFX** and **PharosFX Plus** Systems

Ordering Information

Catalog # Description

Systems

170-9450 **Molecular Imager PharosFX System**, PC or Mac, 100/240 V, includes Quantity One software, sample tray set, fluorescence filters (170-7866, 170-7896), USB2 cable, instructions

170-9460 **Molecular Imager PharosFX Plus System**, PC or Mac, 110/240 V, includes Quantity One software, sample tray set, fluorescence (170-7866, 170-7896) and phosphor imaging filters, USB2 cable, instructions

Lasers

170-7890 **External Laser**, 488 nm, includes 170-9459 filter

170-7893 **635 nm External Laser Upgrade**, for 170-7890, includes 170-7865 filter

170-7892 **External Lasers**, 488 nm and 635 nm, includes 170-9459 and 170-7865 filters

Filters

170-9459 **Filter 530 nm BP**, for ECL Plus, AttoPhos, SYBR Green I, Alexa Fluor 488, FITC, Cy2, and Pro-Q Emerald dyes

170-7863 **Filter 555 nm LP**, for Texas Red dye

170-7866 **Filter 605 nm BP**, for Flamingo™, ethidium bromide, SYPRO Red, SYPRO Ruby, Alexa Fluor 532 and 546, and Cy3 dyes

170-7896 **Filter 640 nm BP**, for Texas Red dye

170-7865 **Filter 695 nm BP**, for Cy5 and Alexa Fluor 635 dyes

170-7867 **Blank Filter Holder**

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

170-7813 **Sample Holders**, for gels

170-7814 **Microplate Adaptor**, for multi-sample tray I

Catalog # Description

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

170-7869 **Replacement Bulb for Eraser Screen-K**

Miscellaneous Accessories

161-0722 **Bio-Rad Cleaning Concentrate**, 50x, 1 kg

165-3414 **Gel Clip**, holds any gel size

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**

165-9775 **Proteomweaver 4.0 Professional 2-D Analysis Software**, with multifuorescence analysis (MFA)

165-9790 **Proteomweaver 4.0 Enterprise 2-D Analysis Software**, with multifuorescence analysis (MFA)

165-9798 **Proteomweaver 4.0 Enterprise Pro 2-D Analysis Software**

* Included with imaging system.

Alexa Fluor, Pro-Q, SYBR, SYPRO, and Texas Red are trademarks of Molecular Probes, Inc. AttoPhos is a trademark of Promega Corporation. Cy and ECL Plus are trademarks of GE Healthcare. 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