



# Molecular Imager<sup>®</sup> PharosFX<sup>™</sup> and PharosFX<sup>™</sup> 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"> <li>Minimizes optical distortion for accurate image data</li> </ul>
Sample imaging from top by direct contact with scanning beam	<ul style="list-style-type: none"> <li>Accommodates wide range of samples, including thick agarose gels, microplates, various polyacrylamide gels, blots, and TLC plates</li> </ul>
Self-alignment of external lasers before each scan	<ul style="list-style-type: none"> <li>Reliable, robust performance; reproducible data</li> </ul>
User-configured custom emission filters in scanner hardware and software	<ul style="list-style-type: none"> <li>Unlimited choice of fluorophores in visible spectrum</li> </ul>
Transillumination screen	<ul style="list-style-type: none"> <li>Performs gel documentation for colorimetric stains</li> </ul>
Retractable moisture-sealed trays	<ul style="list-style-type: none"> <li>Allows imaging of samples of different formats, and facilitates safe and efficient processing of samples</li> </ul>
Small footprint (57 x 68 cm)	<ul style="list-style-type: none"> <li>Optimizes use of limited benchtop</li> </ul>
Image acquisition and analysis with Quantity One <sup>®</sup> 1-D analysis software (included)	<ul style="list-style-type: none"> <li>Application-directed user interface</li> <li>Mac and PC compatible</li> <li>Seamless integration with PDQuest<sup>™</sup> analysis software and EXQuest<sup>™</sup> spot cutter</li> <li>Options for CFR 21 Part 11</li> </ul>
<b>PharosFX Plus</b> All BaFBr:Eu phosphor screens up to 35 x 43 cm accommodated**	<ul style="list-style-type: none"> <li>Economical utilization of phosphor screens, new and previously owned</li> </ul>



\* U.S. 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

- 170-7893 **635 nm External Laser Upgrade**, 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<sup>®</sup> 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
- 170-7814 **Microplate Adaptor**, for multi-sample tray I

Catalog # Description

### 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 **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**

\* Included with imaging system.

Alexa Fluor, Pro-Q, SYPRO, and Texas Red are trademarks of Invitrogen Corporation. SYBR is a trademark of Life Technologies Corporation. AttoPhos is a trademark of Promega Corporation. Cy and ECL Plus are trademarks of GE Healthcare group companies. Kodak is a trademark of Eastman Kodak Company. Mac is a trademark of Apple, Inc.

Bio-Rad Laboratories, Inc. is licensed by Invitrogen Corporation to sell SYPRO products for research use only, under U.S. patent 5,616,502.

Bio-Rad Laboratories, Inc. is licensed by Life Technologies Corporation to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://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