



**Image What Others Only Imagine.**  
The New Fluor-S® MAX2™ Multimager,  
for the Best in Sensitivity and Resolution.

**BIO-RAD**

# the essence of multi



## Chemiluminescence

- Detect all horseradish peroxidase (HRP)-activated substrates including AmpLight™, ECL™, ECL Plus™, SuperSignal™, Ultra, SuperSignal NA, SuperSignal Blaze
- Detect all alkaline phosphatase (AP)-activated substrates including CDP-Star®, CSPD®, AMPPD®
- Perform high-sensitivity detection of chemiluminescence including AttoPhos® and ECL Plus

## Fluorescence

- Compatibility with an extensive range of fluorescent dyes including ethidium bromide, SYBR® Green, Radiant® Red, fluorescein (FITC), rhodamine, and Green Fluorescent Protein, and many more
- Perform true multicolor fluorescence imaging with 8 emission filter positions and interchangeable filters
- Produce perfectly coregistered images for overlay production with patented multiimage alignment

## Densitometry and Colorimetric Imaging

- Perform high-quality imaging and analysis of autoradiograms and colorimetric stains
- Detect numerous colorimetric stains, including copper, zinc, and silver stain

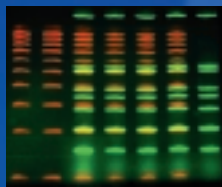


# Imaging

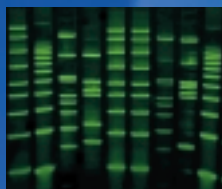
activated substrates,  
SuperSignal®, SuperSignal  
activated substrates, including  
fluorescent substrates,

fluorescent dyes and labels,  
SYBR Gold, SYPRO® dyes,  
Cy™ 2, Cy3, Texas Red®,  
imaging with a choice of  
variable user-accessible filters  
for accurate analysis and  
imaging optics

of 1- and 2-D gels, blots,  
including Coomassie™ Blue,



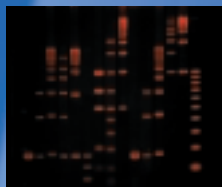
**FITC/Texas Red dual-color  
fluorescence**



**SYBR Green I**



**SYPRO Orange protein stain**



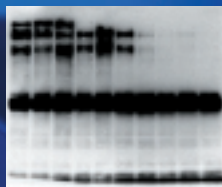
**Ethidium bromide**



**Coomassie Blue**



**CDP-Star array blot**



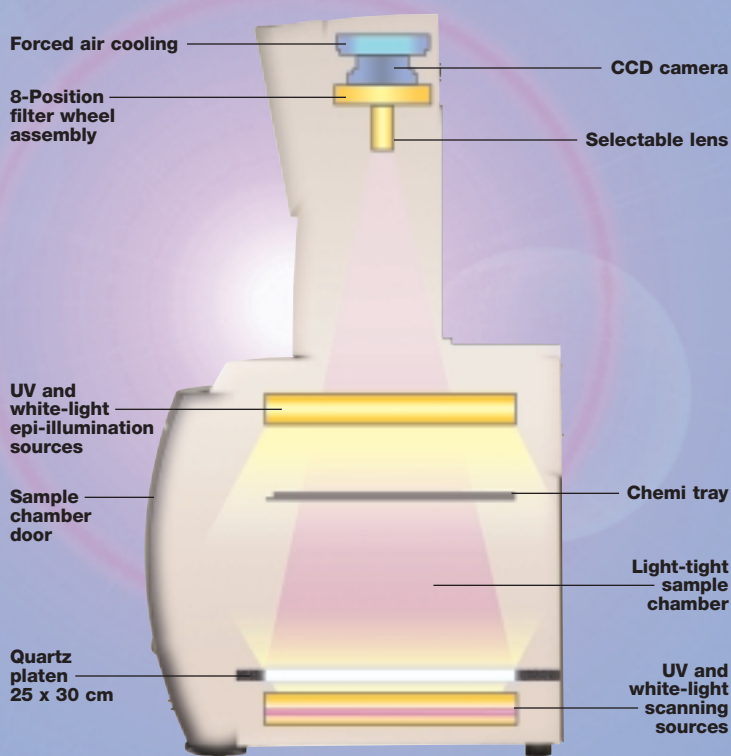
**ECL chemiluminescent  
western blot**

## Advanced Imaging Technology

- Multimode imaging with the ability to select either transmittance illumination to detect transparent samples such as gels and film or epi-illumination to detect opaque samples such as blots
- Rapid scanning transillumination mechanism with higher uniformity than stationary bulb sources
- Broad-bandwidth UV excitation supports the excitation of an extensive range of fluorescent dyes and labels
- Large UV fluorescence Stokes shift (discrimination between the excitation and emission wavelengths) and minimal autofluorescence for improved signal-to-noise ratios and fluorescence sensitivity
- White-light illumination for high-sensitivity detection of colorimetric samples and autoradiographs
- Large, easy-access sample stage (40 x 60 cm) and sample imaging area (25 x 30 cm)
- Flexible optics design compatibility with most f-mount lenses for custom applications
- Advanced dark current correction to remove electronic background noise

## Powerful, Easy-to-Use Software

- Turnkey application-driven acquisition, requiring no user knowledge of sample excitation or collection settings
- Automated control of scanning system through a flexible decoupled computer interface
- Intelligent multiplatform software that operates in Macintosh® and Windows® 98, 2000, and NT environments
- Choice of application-specific programs for fingerprinting, 2-D, and sequence analysis



## Specifications

Cooling system	Thermoelectric
Cooling temperature	-25°C
Pixel size	9 x 9 µm
Pixel density	1,536 x 1,024 (H x W)
Dynamic range	3.4 orders
Data density	12 bit (4,096 gray levels)
Illumination modes	Scanning and epi-illumination
Excitation wavelengths	Broad-bandwidth UV and white light
Standard emission filters	520 long pass, 530DF30, 610 long pass
Maximum sample size	40 x 60 cm
Maximum imaging area	25 x 30 cm
Software compatibility	Windows 98, 2000, NT, and Macintosh
Dimensions	68 x 54 x 109 cm (W x D x H)
Weight	48 kg

## Ordering Information

Catalog #	Description
170-7740	Fluor-S MAX2 Multimager System, PC, 100/240 V, includes 50 mm and 28–80 mm zoom lens, sample tray, PC SCSI cable, cleaning kit, Quantity One® software, instructions
170-7741	Fluor-S MAX2 Multimager System, Mac®, 100/240 V, includes 50 mm and 28–80 mm zoom lens, sample tray, Mac SCSI cable, cleaning kit, Quantity One software, instructions

## Fluorescent Reagents and Kits

### AmpLight Fluorescent and Chemiluminescent Detection Kits

These kits employ novel HRP-based chemistries which have been optimized for highly sensitive western detection using digital imaging systems.

### SYPRO Orange Fluorescent Protein Stain

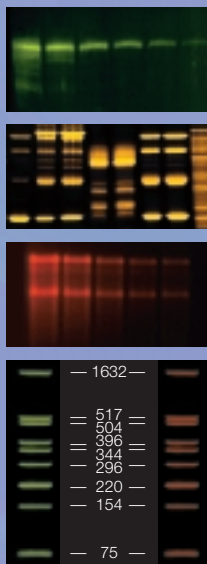
Highly sensitive and easy-to-use stain for visualizing 1-D and 2-D SDS-PAGE and native protein gels.

### Radiant Red Fluorescent RNA Stain

A sensitive fluorescent stain for visualizing RNA following electrophoresis in denaturing agarose gels.

### 2-Color Fluorescent DNA Standards

Our DNA standards allow the ultimate in fragment sizing accuracy. Standards are pre-labeled with either fluorescein or Texas Red and are available in a range of sizes, including convenient molecular weight rulers. Complementary sample end-labeling kits are also available.



AMPPD, CDP-Star, and CSPD are trademarks of Tropix, Inc. AttoPhos is a trademark of JBL Scientific, Inc. Coomassie is a trademark of Imperial Chemical Industries PLC. Cy, ECL, and ECL Plus are trademarks of Amersham International PLC. Macintosh and Mac are trademarks of Apple Computer, Inc. SuperSignal, SuperSignal Ultra, SuperSignal Blaze, and SuperSignal NA are trademarks of Pierce Chemical Co. SYBR, SYPRO, and Texas Red are trademarks of Molecular Probes, Inc. Windows, Windows 98, 2000, and NT are trademarks of Microsoft, Inc.

## Integrated Imaging Systems and Software



**Fluor-S MAX2 Multimager System**



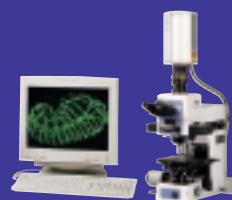
**Molecular Imager® FX Pro Plus™ System**



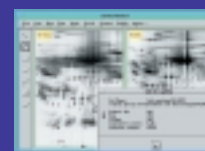
**GS-800™ Imaging Densitometer**



**Gel Documentation Systems**



**Confocal Imaging Systems**



**Specialized RFLP, 2-D, and Sequencing Software**

**BIO-RAD**

**Bio-Rad  
Laboratories**

Life Science  
Group

**Web Site** [www.bio-rad.com](http://www.bio-rad.com) **U.S.** (800) 4BIORAD **Australia** 02 9914 2800 **Austria** (01)-877 89 01 **Belgium** 09-385 55 11 **Canada** (905) 712-2771  
**China** 86-10-62051850/51 **Denmark** 45 39 17 99 47 **Finland** 358 (0)9 804 2200 **France** 01 43 90 46 90 **Germany** 089 318 84-0  
**Hong Kong** 852-2789-3300 **India** (91-11) 461-0103 **Israel** 03 951 4127 **Italy** 39-02-216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460  
**Latin America** 305-894-5950 **Mexico** 514-2210 **The Netherlands** 0318-540666 **New Zealand** 64-9-4152280 **Norway** 22-74-18-70 **Russia** 7 095 979 98 00  
**Singapore** 65-2729877 **Spain** 34-91-661-7085 **Sweden** 46 (0)8-55 51 27 00 **Switzerland** 01-809 55 55 **United Kingdom** 0800-181134