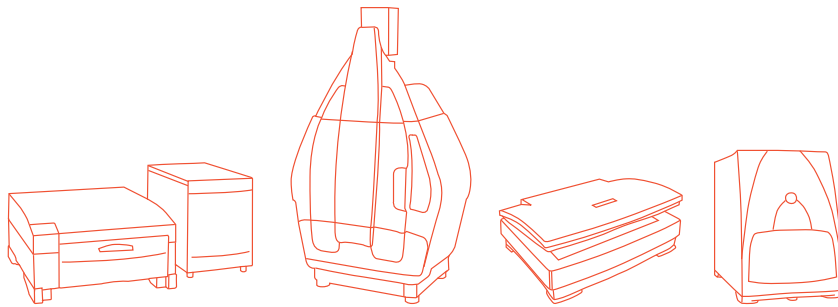




## Bio-Rad Imaging Systems Family







Unlike other imagers, Bio-Rad imaging systems were developed by scientists for scientists. Bio-Rad understands your challenges and knows how to overcome them. That's why Bio-Rad imaging systems have consistently won the prestigious Life Science Industry Awards for most outstanding image analysis system. Life scientists choose the most reliable and best performing products to receive these awards in categories spanning all of the life science disciplines.

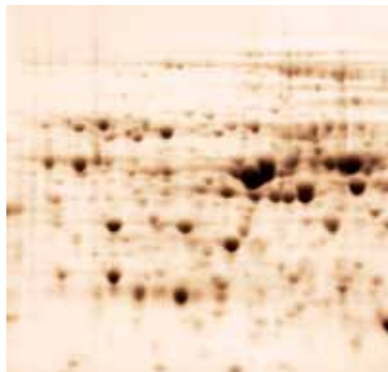


Bio-Rad Laboratories, Inc.

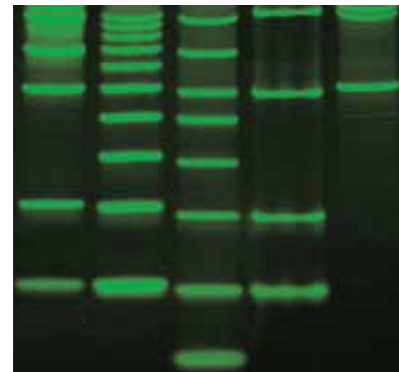
2006, 2007, and 2008 Award Winner  
Best Image Analysis Systems

2010 Award Winner  
Top Ten Innovations

COMPACT



EFFICIENT



DNA Gels



Protein Gels



3-Color Fluorescence



2-D Protein Gels



X-Ray Film



Colony Counting



2-D DIGE



Phosphor Imaging



Chemiluminescent Blots



Colorimetric Blot Detection



Stain-free imaging

### Gel Doc™ EZ System



**Imaging and analysis of protein and DNA gels in seconds with the push of just one button**

#### About the system

- A simple and compact system that is customizable, easy to use, and that can fit easily on a laboratory bench
- Offers remarkable flexibility with its four tray options for different applications
- Provides expedited workflow and increased free time. A 2-hour Coomassie staining protocol can be condensed to 5 minutes using the Stain-Free tray on the Gel Doc EZ system
- Provides high-quality, publication-ready images without any manipulation or user-introduced errors
- Possesses the sophisticated and intuitive Image Lab™ software, which provides images and analyzed results that are highly reproducible, within a matter of seconds

### Gel Doc™ XR+ System

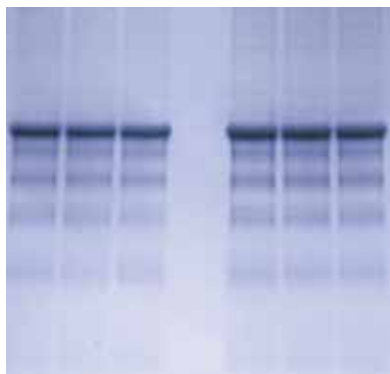


**Affordable system for routine imaging of DNA and protein gels**

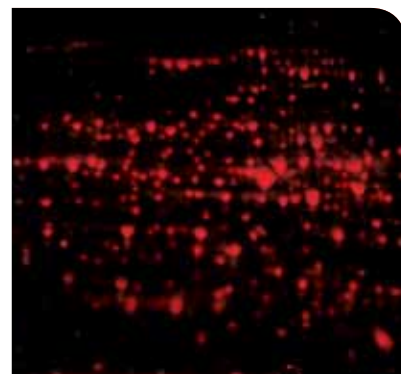
#### About the system

- Image Lab software provides automated and reproducible results from image acquisition to analysis
- Autofocus at any zoom level ensures high-quality images that are highly reproducible
- Trans-blue light plate allows on-plate band excision using safe DNA staining (optional)
- Can be upgraded to the ChemiDoc™ XRS+ system
- Installation qualification/operational qualification (IQ/OQ) protocols (optional)
- Quantity One® 1-D analysis software 21 CFR Part 11 module (optional)

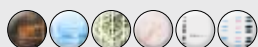
**AUTOMATED**



**POWERFUL**



## ChemiDoc™ XRS+ System



**Affordable system for chemiluminescent western blotting and routine gel imaging**

### About the system

- System delivers both high sensitivity and excellent resolution
- Image Lab software provides automated and reproducible results from image acquisition to analysis
- Autofocus at any zoom level ensures high-quality images that are highly reproducible
- Trans-blue light plate allows on-plate band excision using safe DNA staining (optional)
- IQ/OQ (optional)
- Quantity One 1-D analysis software module is available for compliance with U.S. FDA 21 CFR Part 11 requirements (optional)

## ChemiDoc™ MP System



**Designed for chemiluminescence and multiplex fluorescence detection, this imager can also be used for routine gel imaging**

### About the system

- Our most sensitive charge-coupled device (CCD)-based imaging system
- Image Lab software provides automated and reproducible results from image acquisition to analysis
- Autofocus at any zoom level ensures high-quality images that are highly reproducible
- 3-color multiplex capability
- Automated, preprogrammable filter and light operation determined by the application selected
- Trans-blue light plate allows on-plate band excision using safe DNA staining (optional)
- Total protein normalization with stain-free technology

## GS-800™

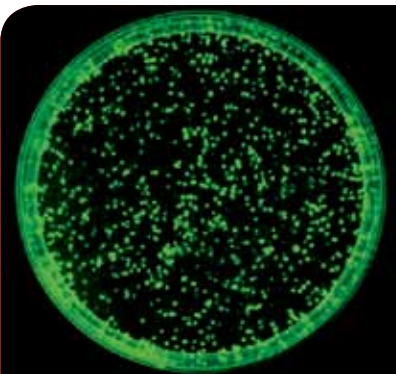
## Calibrated Densitometer



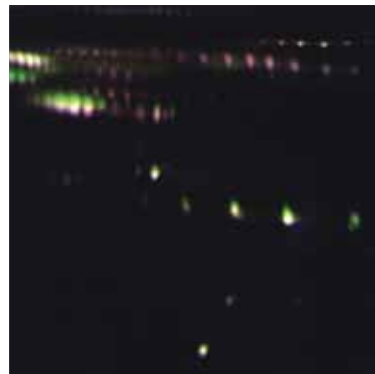
**Affordable, high-resolution scanner for stained gels and blots**

### About the system

- Self-calibrating with internal standards for run-to-run consistency
- IQ/OQ (optional)
- Quantity One 1-D analysis software 21 CFR Part 11 module (optional)



REPRODUCIBLE



FAST



## PharosFX™ Plus System



**Laser-scanning system allows high-resolution 3-color imaging and phosphor imaging**

### About the system

- A high-resolution laser scanning system for fluorescent and radioisotopic applications
- 3-color multiplex capability
- Quantity One 1-D analysis software 21 CFR Part 11 module (optional)
- Fully upgradable from single laser to 3-laser capability

## Personal Molecular Imager™ (PMI™) System



**Basic phosphor imaging system for radioisotopic applications**

### About the system

- A high-resolution laser scanning system for radioisotopic applications
- Used with storage phosphor screens for  $^{32}\text{P}$ ,  $^{33}\text{P}$ ,  $^{35}\text{S}$ ,  $^{14}\text{C}$ , and  $^3\text{H}$
- Quantity One 1-D analysis software 21 CFR Part 11 module (optional)

# Imaging System Packages

For most researchers, no single imager addresses every experimental need. Therefore, Bio-Rad has created unified image acquisition and analysis software to allow one computer to operate a combination of instruments.

These are our most often requested imaging system combinations.

## Gel, Blot, and Isotopic Imaging Package



▪ ChemiDoc XRS+ system

▪ Personal Molecular Imager (PMI) system

## Multiplex Blot, Gel, and Isotopic Imaging Package



▪ ChemiDoc MP system

▪ Personal Molecular Imager (PMI) system

## 2-D Imaging Package



▪ PharosFX Plus system with blue and red lasers

▪ PDQuest™ Advanced 2-D analysis software

▪ EXQuest™ spot cutter

Life Science Industry Awards is a service mark of The Scientist.  
Coomassie is a trademark of BASF Aktiengesellschaft.



**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