



# ChemiDoc™ MP Imaging System

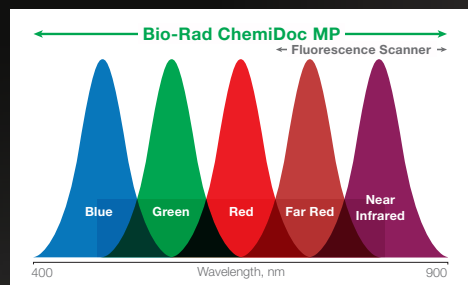
The best of chemiluminescence and fluorescence in one imager.

Leave behind traditional film and laser scanners and get the superior quantitation of digital, the multiplexing of fluorescence, and the brightest new fluorophores all in a single imaging system.



## More flexibility in dye selection.

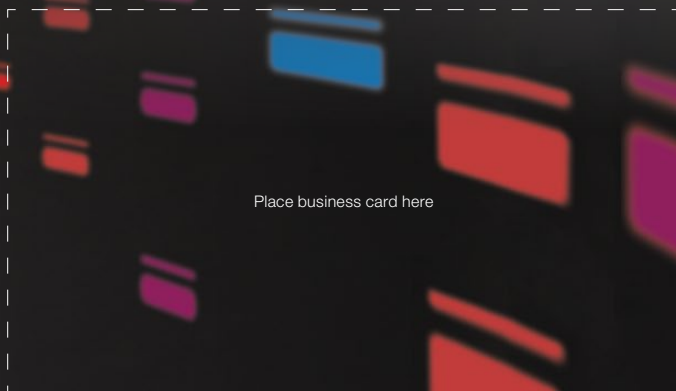
Detect up to three different proteins in a single experiment with a wide range of supported fluorophore options.



## Not your traditional software.

Image Lab™ Touch Software:

- Teaches as you tap and explore
- Picks optimal light source and filters
- Suggests exposure times



## More than an imaging system. A fully supported western blotting workflow.



### Sample Prep:

SureBeads™ Protein A/G–Conjugated Beads

*Easy, fast, and cost-effective for better immunoprecipitation.*

### Electrophoresis:

Stain-Free Gels

*Rapid protein separation and normalization without the hassles of staining.*

### Transfer:

Trans-Blot® Turbo™ Transfer System

*Get to the finish line faster with transfers as short as 3 minutes.*

### Immunodetection:

PrecisionAb™ Validated Western Blotting Antibodies

*Sensitivity, specificity, reproducibility — antibodies you can rely on.*

### Image Acquisition:

ChemiDoc MP Imaging System

*Fluorescence and chemiluminescence detection without compromises.*

### Image Analysis:

Image Lab Software Family

*Easy to use, automated, and powerful.*

Visit [bio-rad.com/WesternWorkflow](http://bio-rad.com/WesternWorkflow) to learn more.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00  
**France** 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 972 03 963 6050  
**Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666 **New Zealand** 64 9 415 2280  
**Norway** 47 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 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026674 55 05 **Taiwan** 886 2 2578 7189  
**Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

