

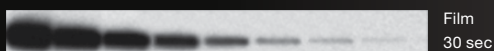
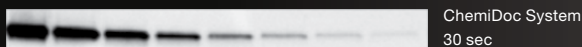


# ChemiDoc™ Imaging System

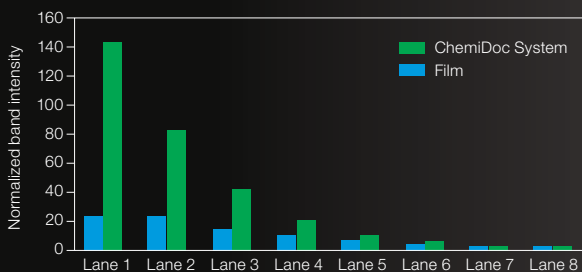
## The best of chemiluminescence.

Gain a larger dynamic range and the benefits of digital quantification while keeping the sensitivity of film.

### The sensitivity of film



### A larger dynamic range

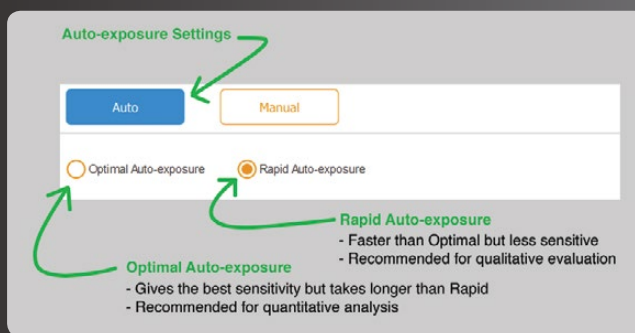


## Not your traditional software.

Image Lab™ Touch Software:

- Teaches as you tap and explore
- Picks optimal light source and filters
- Permits manual entry and automated calculation of exposure times

### ChemiDoc System Help Screen



## A path to fluorescence.

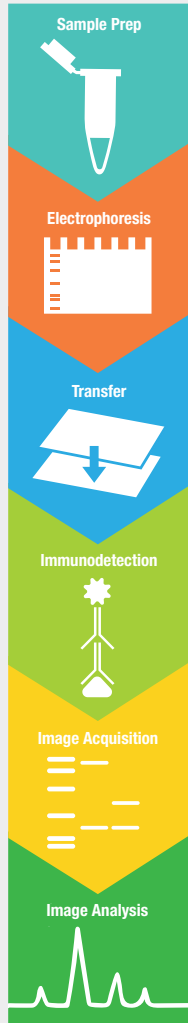
A simple on-site hardware update provides:

- Instant access to RGB and NIR fluorescent western blotting
- The ability to detect three different proteins on a single blot
- Compatibility with a wider range of fluorophores than traditional fluorescence scanners



Place business card here

## 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

