






## Imaging System Selection Guide

Application	Molecular Imager Systems							
	ChemiDoc™ MP	Gel Doc™ EZ	GS-800™ Densitometer	Gel Doc™ XR+	ChemiDoc™ XRS+	PharosFX™	PharosFX™ Plus	Personal Molecular Imager™
								
<b>Nucleic Acid Detection</b>								
Ethidium bromide stain	5	4	—	4	4	5	5	—
SYBR® Green I stain	5	4	—	4	4	5	5	—
SYBR® Safe stain	5	4	—	4	4	5	5	—
Coumarin stain	5	4	5	—	—	—	—	—
Fast Blast™ DNA stain	4	4	5	4	4	—	—	—
<b>Protein Detection, 1-D Gels</b>								
<b>Protein Detection, 2-D Gels</b>								
<b>Blot Detection</b>								
<b>Micro- and Macroarray** Detection</b>								
<b>Colony Counting</b>								
<b>Isotopic Detection</b>								

— Not recommended; 1–5, recommendation level (5 = highest).

\* With spot features ≥400 µm



**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

---