

---

These tables list current specifications for Bio-Rad imaging systems. Refer to the following pages for information about each imaging system. For further information, please call Life Sciences Technical Support at 800-424-6723, opt 2.

## Contents

### Current Products:

Gel Doc™ EZ and Gel Doc™ XR+ Systems with Image Lab™ Software.....	2
ChemiDoc™ XRS+, ChemiDoc™ MP, and ChemiDoc™ Touch Systems with Image Lab Software .....	2
GS-900™ Densitometer with Image Lab Software.....	2
Gel Doc XR+ System with Quantity One® Software .....	3
ChemiDoc XRS+ System with Quantity One Software.....	3

### Previous Products:

GS-800™ Densitometer USB Scanner with Quantity One Software .....	3
Gel Doc XR+ and ChemiDoc XRS+ Systems with Quantity One Software .....	3
Gel Doc™ XR and ChemiDoc™ XRS Systems with Quantity One Software .....	4
PharosFX™ and PharosFX™ Plus USB Scanners with Quantity One Software .....	4
Personal Molecular Imager™ (PMI™) USB Scanner with Quantity One Software.....	4
VersaDoc™ MP System with Quantity One Software .....	4
GS-800 Densitometer SCSI Scanner with Quantity One Software .....	4
PharosFX and PharosFX Plus SCSI Scanners with Quantity One Software .....	4
Personal Molecular Imager (PMI) SCSI Scanner with Quantity One Software .....	4
Computer PCI Slot Requirements .....	5

## Image Lab Software Compatibility: Image Acquisition vs. Image Analysis

See below for the specific operating system and imager serial number cutoff information.

**NOTE:** Image Lab Software, Security Edition (catalog #170-9691, 170-9692, and 170-9693) for 21 CFR part 11 compliance is not compatible with the Macintosh operating system. If you are using this version of the software, you must use a Windows operating system for both image acquisition and analysis.

Imager	Image Acquisition		Image Analysis	
	Windows OS	Mac OS	Windows OS	Mac OS
GS-900 Calibrated Densitometer	•		•	•
Gel Doc EZ System	•	•	•	•
Gel Doc XR+ System	•	•	•	•
ChemiDoc XRS+ System	•	•	•	•
ChemiDoc MP System	•	•	•	•
ChemiDoc Touch System	N/A since Image Lab Touch Software controls image acquisition		•	•

### Current Products:

#### Image Lab Software version 5.2.1:

Gel Doc EZ System

Gel Doc XR+ System (serial numbers starting with 721BR)

ChemiDoc XRS+ System (serial numbers starting with 721BR)

ChemiDoc MP System

ChemiDoc Touch System

Computer	Windows	Mac
Operating system	Win 7 (32- or 64-bit) / Win 8.1 (64-bit)	OS X 10.8 / OS X 10.9
Processor	Intel Core i3 or better	Intel Core i3 or better
Hard disk space	60 GB	60 GB
System memory (RAM)	4 GB	4 GB
Screen resolution	1280 x 1024 or higher; 128 MB video RAM	1280 x 1024 or higher; 128 MB video RAM
USB port	1 free USB 2.0 port	1 free USB 2.0 port

#### Image Lab Software version 5.2.1:

GS-900 Calibrated Densitometer

Computer	Windows	Mac
Operating system	Win 7 (32- or 64-bit) / Win 8.1 (64-bit)	Not compatible
Processor	Intel Core i3 or better	N/A
Hard disk space	60 GB	N/A
System memory (RAM)	4 GB	N/A
Screen resolution	1280 x 1024 or higher; 128 MB video RAM	N/A
USB port	1 free USB 2.0 port	N/A

**Quantity One 1-D Analysis Software version 4.6.9:**

Gel Doc XR+ System (serial numbers starting with 721BR)  
 ChemiDoc XRS+ System (serial numbers starting with 721BR)

**Note:** To run an imager with Quantity One Software, first install instrument drivers from Image Lab Software, then install Quantity One Software.

Computer	Windows	Mac
Operating system	Win XP Pro SP3 / Win 7 (32-bit)	OS X 10.4 / OS X 10.5
Processor	Intel Core i3 or equivalent	PowerPC
Hard disk space	60 GB	60 GB
System memory (RAM)	4 GB	4 GB
Screen resolution	1280 x 1024 or higher; 128 MB video RAM	1280 x 1024 or higher; 128 MB video RAM
USB port	1 free USB 2.0 port	1 free USB 2.0 port
Quantity One HASP dongle	1 free USB 1.0 or better port	1 free USB 1.0 or better port

**Previous Products:****Quantity One 1-D Analysis Software version 4.6.9:**

GS-800 Densitometer USB Scanner

Computer	Windows	Mac
Operating system	Win XP Pro SP3	Not compatible
Processor	Intel 2.0 GHz; Pentium 4, Dual-Core, or equivalent	N/A
Hard disk space	Minimum >20 GB	N/A
System memory (RAM)	Minimum >1 GB	N/A
Cache memory	Minimum >2 GB	N/A
Screen resolution	1280 x 1024 or higher; 128 MB video RAM	N/A
USB port	1 free USB 2.0 port	N/A
Quantity One HASP dongle	1 free USB 1.0 or better port	N/A

**Quantity One 1-D Analysis Software version 4.6.9:**

Gel Doc XR+ System (serial numbers starting with 720BR)  
 ChemiDoc XRS+ (FireWire) System

Computer	Windows	Mac
Operating system	Win XP Pro SP3	OS X 10.4 / OS X 10.5
Processor	Intel 2.0 GHz; Pentium 4, Dual-Core, or equivalent	PowerPC
Hard disk space	Minimum >20 GB	Minimum >20 GB
System memory (RAM)	Minimum 1 GB	Minimum 1 GB
Screen resolution	1280 x 1024 or higher; 128 MB video RAM	1280 x 1024 or higher; 128 MB video RAM
USB port	1 free USB 2.0 port	1 free USB 2.0 port
Quantity One HASP dongle	1 free USB 1.0 or better port	1 free USB 1.0 or better port
FireWire port (only required for ChemiDoc XRS+ FireWire camera)	1 free FireWire port (FW 400; 6 pin) or 1 full size PCI slot	1 free FireWire port (FW 400; 6 pin)

### Quantity One 1-D Analysis Software version 4.6.9:

Gel Doc XR (FireWire) System (serial numbers starting with 76S)  
 ChemiDoc XRS (Roper LVDS camera) System (serial numbers starting with 76S)

**Note on ChemiDoc XRS (Roper LVDS camera):** Always install drivers before Roper PCI card.

Computer	Windows	Mac
Operating system	Win XP Pro SP3	OS X 10.4 / OS X 10.5
Processor	Intel 2.0 GHz; Pentium 4, Dual-Core, or equivalent	PowerPC
Hard disk space	Minimum >20 GB	Minimum >20 GB
System memory (RAM)	Minimum >1 GB	Minimum >1 GB
Cache memory	Minimum >2 GB	Minimum >2 GB
Screen resolution	1280 x 1024 or higher; 128 MB video RAM	1280 x 1024 or higher; 128 MB video RAM
Serial or USB port	1 free serial port	1 free USB port; Keyspan serial to USB convertor
Quantity One HASP dongle	1 free USB 1.0 or better port	1 free USB 1.0 or better port
FireWire port (only required or GelDoc XR FireWire camera)	1 free FireWire port (FW 400; 6 pin) or 1 full size PCI slot	1 free FireWire port (FW 400; 6 pin)
Full size PCI slot	1 full size PCI slot for video card	1 full size PCI slot for video card

### Quantity One 1-D Analysis Software version 4.6.9:

PharosFX and PharosFX Plus USB Scanners\*  
 Personal Molecular Imager (PMI) USB Scanner

Computer	Windows	Mac
Operating system	Win XP Pro SP3 / Win 7 (32- or 64-bit)	OS X 10.4 / OS X 10.5
Processor	Intel Core i3 or equivalent	PowerPC
Hard disk space	60 GB	60 GB
System memory (RAM)	4 GB	4 GB
Screen resolution	1280 x 1024 or higher; 128 MB video RAM	1280 x 1024 or higher; 128 MB video RAM
USB port	1 free USB 2.0 port	1 free USB 2.0 port
Quantity One HASP dongle	1 free USB 1.0 or better	1 free USB 1.0 or better port

\*PharosFX and PharosFX Plus USB Scanners can use Windows 7 64-bit only with updated HASP drivers. Contact Technical Support at 800-424-6723, opt 2.

### Quantity One 1-D Analysis Software version 4.6.9:

VersaDoc MP (4000 and 5000 cameras) System  
 GS-800 Densitometer SCSI Scanner  
 PharosFX and PharosFX Plus SCSI Scanners  
 Personal Molecular Imager (PMI) SCSI Scanner

**Note on GS-800 SCSI, FX and FX Plus SCSI, PMI SCSI:** Adaptec ASPI layer required for Win XP computers. Call technical support for details.

Computer	Windows	Mac
Operating system	Win XP Pro SP3	OS X 10.4 / OS X 10.5
Processor	Intel 2.0 GHz; Pentium 4, Dual-Core, or equivalent	PowerPC
Hard disk space	Minimum >20 GB	Minimum >20 GB
System memory (RAM)	Minimum >1 GB	Minimum >1 GB
Cache memory	Minimum >2 GB	Minimum >2 GB
Screen resolution	1280 x 1024 or higher; 128 MB video RAM	1280 x 1024 or higher; 128 MB video RAM
USB port	1 free serial port	1 free USB port
Quantity One HASP dongle	1 free USB 1.0 or better port	1 free USB 1.0 or better port
Full size PCI slot	1 full size PCI slot for video card	1 full size PCI slot for video card

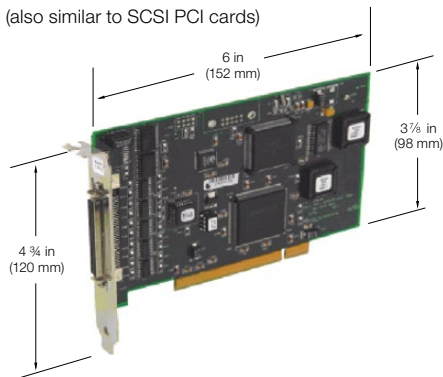
## Computer PCI Slot Requirements

The computer specifications are similar for all FireWire, Roper, and SCSI PCI cards, but only recent models are pictured in this document.

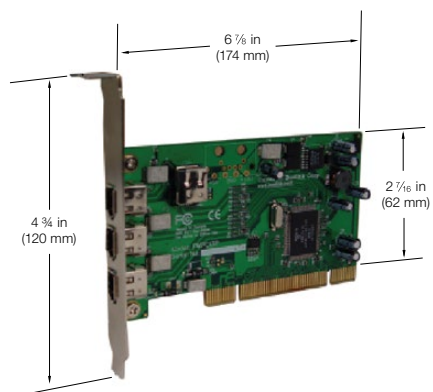
**Warning:** All PCI cards require full size PCI slots.

### VersaDoc 4000 LVDS PCI Card

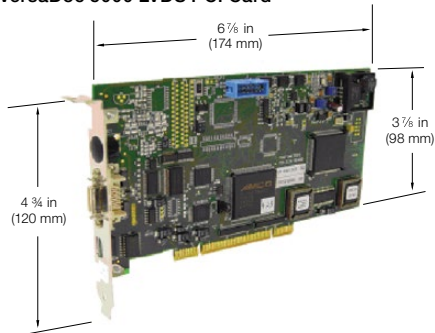
(also similar to SCSI PCI cards)



### ChemiDoc XRS or Gel Doc XR FireWire PCI Card



### VersaDoc 5000 LVDS PCI Card



Adaptec is a trademark of PMC-Sierra, Inc. FireWire, Mac, Macintosh, and OS X are trademarks of Apple Inc. HASP is a trademark of Aladdin Knowledge Systems, Ltd. Intel, Intel Core, and Pentium are trademarks of Intel Corporation. Windows is a trademark of Microsoft Corporation. PowerPC is a trademark of International Business Machines Corporation.



**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 43 1 877 89 01 Belgium 03 710 53 00 Brazil 55 11 3065 7550  
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 49 89 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 81 3 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 (0) 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 026674 55 05  
Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000