

Computer Specifications

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

| | |
|---|---|
| Gel Doc™ EZ system with Image Lab™ 4 software | 2 |
| Gel Doc XR+ system with Image Lab 4 software | 2 |
| ChemiDoc™ XRS+ (USB or Firewire) system with Image Lab 4 software..... | 2 |
| ChemiDoc MP system with Image Lab 4 software | 3 |
| Gel Doc XR+ and ChemiDoc XRS+ systems with Quantity One® 4.6.9 software | 3 |
| Gel Doc and ChemiDoc systems with Quantity One 4.6.9 or earlier software | 4 |
| Criterion Stain Free™ imager with Image Lab 2 software..... | 4 |
| VersaDoc™ MP 4000 and 5000 cameras with Quantity One 4.6.9 software | 5 |
| VersaDoc 1000, 3000, 4000, and 5000 cameras with Quantity One 4.6.9 software..... | 5 |
| GS-800™ densitometer USB scanner with Quantity One 4.6.9 software..... | 6 |
| GS-800 densitometer SCSI scanner with Quantity One 4.6.9 software | 6 |
| PharosFX™ USB scanner with Quantity One 4.6.9 software..... | 6 |
| FX and FX Pro SCSI scanner with Quantity One 4.6.9 software..... | 7 |
| Personal Molecular Imager™ (PMI™) USB scanner with Quantity One 4.6.9 software..... | 7 |
| Personal Molecular Imager (PMI) SCSI scanner with Quantity One 4.6.9 software..... | 7 |
| Computer PCI slot requirements | 8 |

Gel Doc EZ System with Image Lab 4 Software

| Computer | Windows | Mac |
|---------------------|---|---|
| Operating system | Win XP Pro SP3, Win 7 (32 or 64-bit) | OS 10.6 |
| Processor | Intel 2.0 Ghz; Pentium 4, dual core or better | Core 2 Duo 2.0 GHz or higher |
| Hard disk space | Minimum >20 GB; recommend ≥100 GB | Minimum >20 GB, recommend ≥100 GB |
| System memory (RAM) | Minimum 1 GB; recommend ≥2 GB | Minimum 1 GB, recommend ≥2 GB |
| Screen resolution | 1024 x 768 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 |
| Optional software | Microsoft Office 2003 or 2007 | Microsoft Office 2008 or iWorks |

Gel Doc XR+ System with Image Lab 4 Software

System serial numbers starting with 721BR

Note: To run an imager with Quantity One software, first install instrument drivers from Image Lab 2 software, then install Quantity One software. See 720BR hood chart (page 3) for computer specs.

| Computer | Windows | Mac |
|---------------------|---|---|
| Operating system | Win XP Pro SP3, Win 7 (32 or 64-bit) | OS 10.6 |
| Processor | Intel 2.0 Ghz; Pentium 4, dual core or better | Core 2 Duo 2.0 GHz or higher |
| Hard disk space | Minimum >20 GB; recommend ≥100 GB | Minimum >20 GB, recommend ≥100 GB |
| System memory (RAM) | Minimum 1 GB; recommend ≥2 GB | Minimum 1 GB, recommend ≥2 GB |
| Screen resolution | 1024 x 768 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 |
| Optional software | Microsoft Office 2003 or 2007 | Microsoft Office 2008 or iWorks |

ChemiDoc XRS+ (USB or Firewire) System with Image Lab 4 Software

System serial numbers starting with 721BR

Note: To run an imager with Quantity One software, first install instrument drivers from Image Lab 2 software, then install Quantity One software. See 720BR hood chart (page 3) for computer specs.

| Computer | Windows | Mac |
|---|---|---|
| Operating system | Win XP Pro SP3 or Win 7 (32 or 64-bit) | OS 10.6 |
| Processor | Intel 2.0 Ghz; Pentium 4, dual core or better | Core 2 Duo 2.0 GHz or higher |
| Hard disk space | Minimum >20 GB; recommend ≥100 GB | Minimum >20 GB, recommend ≥100 GB |
| System memory (RAM) | Minimum 1 GB; recommend ≥2 GB | Minimum 1 GB, recommend ≥2 GB |
| Screen resolution | 1024 x 768 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 |
| Optional software | Microsoft Office 2003 or 2007 | Microsoft Office 2008 or iWorks |
| Firewire port (only required for ChemiDoc XRS+ system with Firewire camera) | 1 free Firewire port (FW 400; 6 pin) or 1 full size PCI slot for Firewire card Warning: PCI card requires full size PCI slot (see page 8) | 1 free Firewire port (FW 400; 6 pin) |

ChemiDoc MP System with Image Lab 4 Software

| Computer | Windows | Mac |
|---------------------|---|---|
| Operating system | Win XP Pro SP3 or Win 7 (32 or 64-bit) | OS 10.6 |
| Processor | Intel 2.0 Ghz; Pentium 4, dual core or better | Core 2 Duo 2.0 GHz or higher |
| Hard disk space | Minimum >20 GB; recommend \geq 100 GB | Minimum >20 GB, recommend \geq 100 GB |
| System memory (RAM) | Minimum 1 GB; recommend \geq 2 GB | Minimum 1 GB, recommend \geq 2 GB |
| Screen resolution | 1024 x 768 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 |
| Optional software | Microsoft Office 2003 or 2007 | Microsoft Office 2008 or iWorks |

Gel Doc XR+ and ChemiDoc XRS+ Systems with Quantity One 4.6.9 Software

System serial numbers starting with 721BR

Gel Doc XR+ USB camera with Quantity One 4.6.9 software

ChemiDoc XRS+ (USB or Firewire) camera with Quantity One 4.6.9 software

Note: Feedback lens added to all cameras.

Note: To run an imager with Quantity One software, first install instrument drivers from Image Lab 2 or 3 software, then install Quantity One software.

| Computer | Windows | Mac |
|---|---|---|
| Operating system | Win XP Pro SP3 | OS 10.5 |
| Processor | Intel 2.0 Ghz; Pentium 4, dual core or better | Core 2 Duo 2.0 GHz or higher |
| Hard disk space | Minimum >20 GB; recommend \geq 100 GB | Minimum >20 GB, recommend \geq 100 GB |
| System memory (RAM) | Minimum 1 GB; recommend \geq 2 GB | Minimum 1 GB, recommend \geq 2 GB |
| Screen resolution | 1024 x 768 or higher; 128 MB video RAM | 1280 x 1024 or higher; 128 MB video RAM |
| USB port | 2 free USB 2.0 ports; 4 USB ports recommended | 2 free USB 2.0 ports; 4 USB ports recommended |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |
| Optional software | Microsoft Office 2003 or 2007 | Microsoft Office 2008 or iWorks |
| Firewire port (only required for ChemiDoc XRS+ system with Firewire camera) | 1 free Firewire port (FW 400; 6 pin) or 1 full size PCI slot Warning: PCI card requires full size PCI slot (see page 8) | 1 free Firewire port (FW 400; 6 pin) |

System serial numbers starting with 720BR

Gel Doc XR+ USB camera with Quantity One 4.6.9 software

ChemiDoc XRS+ Firewire camera with Quantity One 4.6.9 software

Note: USB hood connection added; later systems included the feedback lens (+).

| Computer | Windows | Mac |
|---|---|---|
| Operating system | Win XP Pro SP3 | OS 10.5 |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | >20 GB |
| System memory (RAM) | >1 GB | >1 GB |
| Cache memory | >2 GB | >2 GB |
| USB ports | 2 free USB 2.0 ports; 4 USB ports recommended | 2 free USB 2.0 ports; 4 USB ports recommended |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |
| Firewire port (only required for ChemiDoc XRS+ system with Firewire camera) | 1 free Firewire port (FW 400; 6 pin) or 1 full size PCI slot Warning: PCI card requires full size PCI slot (see page 8) | 1 free Firewire port (FW 400; 6 pin) |

Gel Doc and ChemiDoc Systems with Quantity One 4.6.9 or Earlier Software

System serial numbers starting with 76S

Gel Doc XR Firewire camera with Quantity One software

ChemiDoc XRS Roper LVDS camera with Quantity One software

Note: Always install drivers before Roper PCI card.

| Computer | Windows | Mac |
|--|--|---|
| Operating system | Win XP Pro SP3, Vista SP2 (32-bit only) | OS 10.4 or 10.5 (requires power PC) |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | >20 GB |
| System memory (RAM) | >1 GB | >1 GB |
| Cache memory | >2 GB | >2 GB |
| USB ports | 1 free USB 1.0 or greater ports for HASP dongle | 1 free USB 1.0 or greater ports for HASP dongle |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |
| Firewire port (Gel Doc XR system with Firewire camera only) | 1 free Firewire port (FW 400; 6 pin) or 1 full size PCI slot Warning: PCI card requires full size PCI slot | 1 free Firewire port (FW 400; 6 pin) |
| Full size PCI slot (for ChemiDoc XRS+ system with LVDS camera) | 1 full size PCI slot for Roper LVDS video card Warning: PCI card requires full size PCI slot (see page 8) | 1 full size PCI slot for Roper LVDS video card Warning: PCI card requires full size PCI slot (see page 8) |

System serial numbers starting with 76S or 75S

Gel Doc EQ Scion camera

ChemiDoc XRS Roper LVDS or Scion camera

Note: Always install drivers before Roper PCI card.

| Computer | Windows | Mac |
|--------------------------|--|--|
| Operating system | Win XP Pro SP3 (32-bit only) | OS 10.4 or 10.5 (ChemiDoc XRS LVDS requires power PC) |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | >20 GB |
| System memory (RAM) | >1 GB | >1 GB |
| Cache memory | >2 GB | >2 GB |
| Serial port | 1 free serial port | 1 free USB port; Keyspan serial to USB converter |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |
| Full size PCI slot | 1 full size PCI slot for Scion or Roper LVDS video card Warning: PCI card requires full size PCI slot (see page 8) | 1 full size PCI slot for Scion or Roper LVDS video card Warning: PCI card requires full size PCI slot (see page 8) |

Criterion Stain Free Imager with Image Lab 2 Software

| Computer | Windows | Mac |
|---------------------|--|--|
| Operating system | Win XP Pro SP3, Vista SP2 (32-bit only) | OS 10.4 or 10.5 |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | >20 GB |
| System memory (RAM) | >1 GB | >1 GB |
| Cache memory | >2 GB | >2 GB |
| USB ports | 1 free USB 2.0 port; 4 USB ports recommended | 1 free USB 2.0 port; 4 USB ports recommended |

VersaDoc MP 4000 and 5000 Cameras with Quantity One 4.6.9 Software

Note: Always install drivers before Roper PCI card.

| Computer | Windows | Mac |
|--------------------------|--|--|
| Operating system | Win XP Pro SP3, Vista SP2 or Win 7 (32-bit only) | OS 10.4 or 10.5 (requires power PC) |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | >20 GB |
| System memory (RAM) | >1 GB | >1 GB |
| Cache memory | >2 GB | >2 GB |
| USB ports | 2 USB 2.0 ports; 4 USB ports recommended | 2 USB 2.0 ports; 4 USB ports recommended |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |
| Full size PCI slot | 1 full size PCI slot for video card Warning: PCI card requires full size PCI slot (see page 8) | 1 full size PCI slot for video card Warning: PCI card requires full size PCI slot (see page 8) |

VersaDoc 1000, 3000, 4000, and 5000 Cameras with Quantity One 4.6.9 Software

Note: Contact technical support about updating the PCI video card firmware for older systems.

Note: Always install drivers before Roper PCI video card.

| Computer | Windows | Mac |
|--------------------------|--|--|
| Operating system | Win XP Pro SP3, Vista SP2 (32-bit only) | OS 10.4 or 10.5 (requires power PC) |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | >20 GB |
| System memory (RAM) | >1 GB | >1 GB |
| Cache memory | >2 GB | >2 GB |
| USB or serial ports | 1 USB 2.0 ports; 4 USB ports recommended Note: Older hoods require serial port | 2 USB 2.0 ports; 4 USB ports recommended Note: Older hoods require serial to USB converter |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |
| Full size PCI slot | 1 full size PCI slot for video card Warning: PCI card requires full size PCI slot (see page 8) | 1 full size PCI slot for video card Warning: PCI card requires full size PCI slot (see page 8) |

GS-800 Densitometer USB Scanner with Quantity One 4.6.9 Software

| Computer | Windows | Mac |
|--------------------------|---|-------------------------------------|
| Operating system | Win XP Pro SP3, or Vista SP2 (32-bit only) | Not compatible |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | N/A |
| System memory (RAM) | >1 GB | N/A |
| Cache memory | >2 GB | N/A |
| USB ports | 2 free USB 2.0 ports; 4 USB ports recommended | N/A |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |

GS-800 Densitometer SCSI Scanner with Quantity One 4.6.9 Software

Note: Adaptec ASPI layer required for WinXP computers. Call technical support for details.

| Computer | Windows | Mac |
|--------------------------|---|---|
| Operating system | Win XP Pro SP3 (32-bit only) | Mac OS 10.4 or 10.5 |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | >20 GB |
| System memory (RAM) | >1 GB | >1 GB |
| Cache memory | >2 GB | >2 GB |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |
| Full size PCI slot | 1 full size PCI slot for SCSI card Warning: PCI card requires full size PCI slot (see page 8) | 1 full size PCI slot for SCSI card Warning: PCI card requires full size PCI slot (see page 8) |

PharosFX USB Scanner with Quantity One 4.6.9 Software

| Computer | Windows | Mac |
|--------------------------|---|---|
| Operating system | Win XP Pro SP3, or Vista SP2 | Mac OS 10.4 or 10.5 |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | >20 GB |
| System memory (RAM) | >1 GB | >1 GB |
| Cache memory | >2 GB | >2 GB |
| USB ports | 2 free USB 2.0 ports; 4 USB ports recommended | 2 free USB 2.0 ports; 4 USB ports recommended |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |

FX and FX Pro SCSI Scanner with Quantity One 4.6.9 Software

Note: Adaptec ASPI layer required for WinXP computers. Call technical support for details.

| Computer | Windows | Mac |
|--------------------------|---|---|
| Operating system | Win XP Pro SP3; Win 7 (32-bit only) | Mac OS 10.4 or 10.5 |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | >20 GB |
| System memory (RAM) | >1 GB | >1 GB |
| Cache memory | >2 GB | >2 GB |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |
| Full size PCI slot | 1 full size PCI slot for SCSI card Warning: PCI card requires full size PCI slot (see page 8) | 1 full size PCI slot for SCSI card Warning: PCI card requires full size PCI slot (see page 8) |

Personal Molecular Imager (PMI) USB Scanner with Quantity One 4.6.9 Software

| Computer | Windows | Mac |
|--------------------------|---|---|
| Operating system | Win XP Pro SP3, or Vista SP2; Win 7 (32-bit only) | Mac OS 10.4 or 10.5 |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | >20 GB |
| System memory (RAM) | >1 GB | >1 GB |
| Cache memory | >2 GB | >2 GB |
| USB ports | 2 free USB 2.0 ports; 4 USB ports recommended | 2 free USB 2.0 ports; 4 USB ports recommended |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |

Personal Molecular Imager (PMI) SCSI Scanner with Quantity One 4.6.9 Software

Note: Adaptec ASPI layer required for WinXP computers. Call technical support for details.

| Computer | Windows | Mac |
|--------------------------|---|---|
| Operating system | Win XP Pro SP3; Win 7 (32-bit only) | Mac OS 10.4 or 10.5 |
| Processor | 32-bit; 64-bit might not work | N/A |
| Hard disk space | >20 GB | >20 GB |
| System memory (RAM) | >1 GB | >1 GB |
| Cache memory | >2 GB | >2 GB |
| Quantity One HASP dongle | 1 USB 1.0 or better or parallel port | 1 USB 1.0 or better or Mac ADB port |
| Full size PCI slot | 1 full size PCI slot for SCSI card Warning: PCI card requires full size PCI slot (see page 8) | 1 full size PCI slot for SCSI card Warning: PCI card requires full size PCI slot (see page 8) |

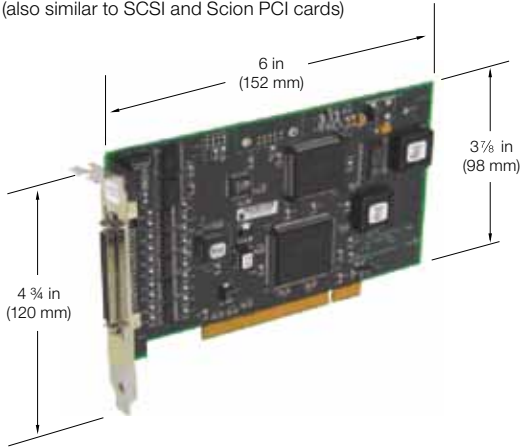
Computer PCI Slot Requirements

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

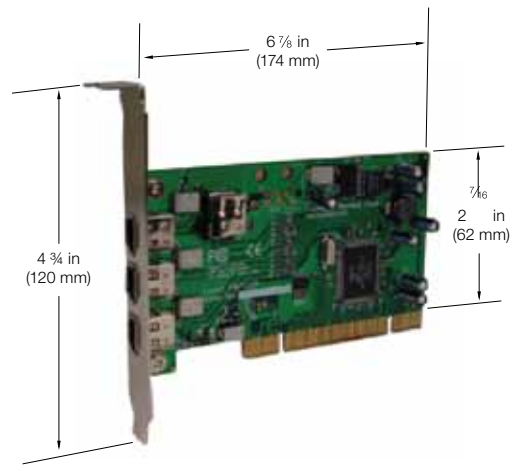
Warning: All PCI cards require full size PCI slots.

VersaDoc 4000 LVDS PCI Card

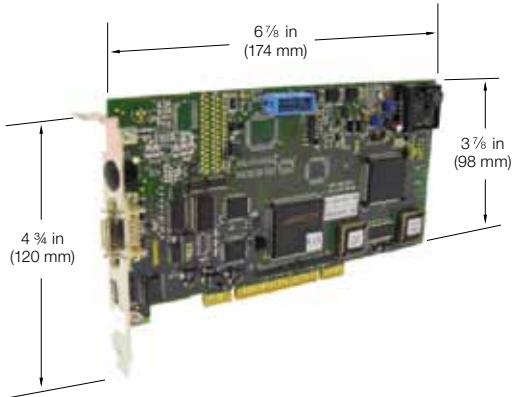
(also similar to SCSI and Scion PCI cards)



ChemiDoc XRS or Gel Doc XR Firewire PCI Card



VersaDoc 5000 LVDS PCI Card



Adaptec is a trademark of PMC-Sierra, Inc. Firewire and Mac are trademarks of Apple, Inc. HASP is a trademark of Aladdin Knowledge Systems LTD Corporation. Intel Core and Pentium are trademarks of Intel Corporation. Keyspan is a trademark of Innosys, Inc. Microsoft Office, Windows, and Windows Vista are trademarks of Microsoft Corporation.



BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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 31 3689 6600
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** 30 210 777 4396 **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 **Malaysia** 60 3 2117 5260 **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 3170 **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** 66 2 6518311 **United Kingdom** 020 8328 2000