

ProteOn Manager™ Software Specifications

ProteOn Manager software is supported on Windows operating systems. The software can coordinate ProteOn™ XPR36 protein interaction array system instrument control or can be run, without an instrument connected, for data processing only.

Software Used for Instrument Control

The ProteOn™ XPR36 system includes a controller and display configured to optimally run ProteOn Manager software and the ProteOn XPR36 instrument. **Do not substitute products without approval from Bio-Rad.**

Software Used for Data Processing Only

ProteOn Manager software can be run on a computer not connected to a ProteOn XPR36 instrument, for data processing in a remote location.

System Specifications for Instrument Control

| | |
|----------------------------|--|
| Processor/model | Pentium 4 (3 GHz processor minimum) |
| Operating system | Windows XP Professional service pack (SP2) |
| RAM | 2 GB |
| Hard disk space | 120 GB minimum, serial ATA/150 minimum |
| Display | 19" flat-panel display |
| Communication ports | 6 available USB ports (included as standard) |
| Graphics | Embedded Intel Extreme Graphics 2 |
| Network card #1 | Integrated gigabit Ethernet |
| Network card #2 | Intel PRO/1000 GT desktop adaptor |
| DVD drives | 16x/8x/16x DVD+RW, 12x/4x/16x DVD-RW |
| Power | 115 or 230 VAC (universal manual switch), 200 W, 50 or 60 Hz |
| Software | Microsoft Office Professional (Office 2003) |
| Language | US English |

System Requirements for Data Processing

| | |
|---------------------------|--|
| Desktop | |
| Processor/model | Pentium 4 (3 GHz processor minimum) |
| Operating system | Windows XP Professional SP2 |
| RAM | 2 GB (1 GB minimum) |
| Hard disk space | 80 GB |
| Display* | 1,280 x 1,024 resolution (absolutely required) |
| Communication port | 1 available USB port |
| Laptop | |
| Processor/model | Pentium 4 (3 GHz processor minimum) |
| Operating system | Windows XP Professional SP2 |
| RAM | 2 GB (1 GB minimum) |
| Hard disk space | 80 GB |
| Display* | 1,280 x 1,024 resolution (absolutely required) |
| Communication port | 1 available USB port |

* Widescreen laptops and monitors are not supported at this time due to the aspect ratio. For more information, contact Bio-Rad's Technical Support.

Pentium is a trademark of Intel Corporation. Windows and Windows XP are trademarks of Microsoft Corporation.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 905 712 2771 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
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 **United Kingdom** 020 8328 2000