

SoftMax Pro Software, Version 7.0.1

Computer System Requirements

SoftMax® Pro Software can be installed on a computer with the following system specifications.

Table 1-1: Minimum and Recommended Requirements for SoftMax Pro Software

Item	Minimum	Recommended
Processor	Single-core, 2 GHz or faster	Quad core or faster
Operating system (SoftMax Pro Software requires the Professional Edition of the operating systems listed.)	Windows 7 (SP1 or newer), 32-bit or 64-bit (x86 or x64) and .NET Framework 4.6 (.NET Framework 4.6 is installed automatically by the SoftMax Pro Software installer if required.)	Windows 7 (SP1 or newer), 32-bit or 64-bit (x86 or x64), Windows 8, 32-bit or 64-bit (x86 or x64), or Windows 10, 32-bit or 64-bit (x86 or x64) The SoftMax Pro MiniMax Imaging Edition must be installed on a 64-bit operating system.
Data connection	RS-232 serial port or USB 2.0 port Network port (depending on the instrument)	USB 2.0 ports For instruments that require an RS-232 serial port, you can use a USB 2.0 port with a Keyspan USB-to-serial adapter. Keyspan USB-to-serial adapters have been field tested and given approval by Molecular Devices. Network port
Memory	2 GB RAM	4 GB RAM If running on a virtual machine, Molecular Devices recommends a minimum of 6 GB RAM. For automation, Molecular Devices recommends a minimum of 8 GB RAM.
Hard drive	500 MB of available space	1 GB of available space, or more
Graphics display	Graphics display adapter 1024 x 768 or higher-resolution display	32-bit graphics display with 256 MB video RAM 1280 x 1024 or higher-resolution display
Software activation	Internet connection or external USB drive	Internet connection or external USB drive

To prevent data loss, turn off all sleep and hibernation settings for the hard disk, the CPU, and the USB ports. Also, disable automatic Windows Updates. You can update Windows manually when the instrument is not being used by the software. You can set these options in Windows Control Panel.

The SpectraMax MiniMax 300 Imaging Cytometer is shipped with a computer that meets greater minimum computer system specifications than those required for the standard installation of the SoftMax Pro Software. The supplied computer must be used as the acquisition computer connected to the instrument.



Note: Installing and using the SoftMax Pro Software on the Windows XP operating system is no longer supported. The software is neither tested nor validated on Windows XP.

