

The following minimum & recommended computer hardware/OS specifications are listed to ensure proper operation of both SoftMax® Pro Standard & SoftMax Pro Enterprise edition software.

WINDOWS COMPATIBLE COMPUTERS

Recommended	Minimum
<ul style="list-style-type: none"> • Intel Pentium® IV, 1.8+ GHz processor • Microsoft Windows® 2000 or XP • Color monitor @ 1024x768 resolution display • 256 MB RAM • 150MB of available hard disk space • CD-ROM drive • Available serial port (COM1 – COM9) 	<ul style="list-style-type: none"> • Pentium® II, 300+MHz processor • Microsoft Windows® 2000 or XP • Color monitor @ 800x600 resolution display • 128 MB RAM • 75MB of available hard disk space • CD-ROM drive • Available serial port (COM1)

NOTE: If you will be creating large files (>20 plates per file) or if you will be using any robotic applications, Molecular Devices recommends that you increase the your computer's RAM to 256 MB or greater.

NOTE: For best system performance, it is recommended that users with data files containing multiple plate sections of kinetic data disable Recalculation by selecting "Suspend Recalculation" from the Edit menu.

APPLE MACINTOSH COMPUTERS

Recommended	Minimum
<ul style="list-style-type: none"> • PowerPC processor (G4/G5 at any speed) • Macintosh OS 9.2.2; OS X 10.3 in Classic mode • Color monitor @ 1024x768 resolution display • 256 MB RAM • 150 MB of available hard disk space • CD-ROM drive • Available serial port (8-pin DIN) • USB-to-Serial Adapter and adapter software 	<ul style="list-style-type: none"> • PowerPC processor (G3 > 300MHz) • Macintosh OS 9.2; OS X 10.2.1 in Classic Mode • Color monitor @ 800x600 resolution display • 128 MB RAM • 75 MB of available hard disk space • CD-ROM drive • Available serial port (8-pin DIN) • USB-to-Serial Adapter and adapter software

NOTE: Mac OS X customers can operate and use SoftMax Pro v4.x software in Classic Mode.

NOTE: If you will be creating large files (>20 plates per file) or if you will be using any robotic applications, Molecular Devices recommends that you increase the your computer's RAM to 256 MB or greater.

NOTE: While **Enterprise Administrator 1.1** is a Windows-only application, the proprietary Administrator Data File (edb) file generated can be used by Macintosh users of SoftMax Pro Enterprise Edition either locally or from a network volume.