

The following computer hardware and operating system specifications are listed to ensure proper operation of both SoftMax Pro and SoftMax Pro GxP software.

WINDOWS[®] COMPATIBLE COMPUTERS

Recommended	Minimum
<ul style="list-style-type: none"> • Intel[®] Pentium[®] 4, 2+ GHz processor or equivalent <ul style="list-style-type: none"> ○ 2.8+ GHz for FlexStation[®] 3 • Windows[®] XP, Vista[®] (Note 1) • Color monitor at 1024x768 resolution display • 1 GB RAM (Note 2) <ul style="list-style-type: none"> ○ 2 GB RAM for FlexStation 3 or StakMax[™] • 150 MB available hard disk space • CD-ROM drive • Available serial port (COM1-COM9) <p style="text-align: center;">- or -</p> • Keyspan-by-Tripp Lite High-Speed USB Serial Adapter (Note 3) 	<ul style="list-style-type: none"> • Intel Pentium[®] 4, 1.6+ GHz processor or equivalent <ul style="list-style-type: none"> ○ 2.8 GHz for FlexStation 3 • Windows XP, Vista (Note 1) • Color monitor at 1024x768 resolution display • 512 MB RAM (Note 2) <ul style="list-style-type: none"> ○ 1 GB RAM for FlexStation 3 or StakMax • 90 MB available hard disk space • CD-ROM drive • Available serial port (COM1) <p style="text-align: center;">- or -</p> • Keyspan-by-Tripp Lite High-Speed USB Serial Adapter (Note 3)

MAC[®] COMPUTERS

Recommended	Minimum
<ul style="list-style-type: none"> • Intel Core Duo • Mac OS X 10.5.x (Note 1) • Color monitor at 1024x768 resolution display • 1 GB RAM (Note 2) <ul style="list-style-type: none"> ○ 2 GB RAM for FlexStation 3 or StakMax • 150 MB available hard disk space • CD-ROM drive • Available serial port (8-pin DIN) <p style="text-align: center;">- or -</p> • Keyspan-by-Tripp Lite High-Speed USB Serial Adapter (Note 3) 	<ul style="list-style-type: none"> • PowerPC processor (G3 > 600 MHz) • Mac OS X 10.4.x • Color monitor at 800x600 resolution display • 512 MB RAM (Note 2) <ul style="list-style-type: none"> ○ 1 GB RAM for FlexStation 3 or StakMax • 90 MB available hard disk space • CD-ROM drive • Available serial port (8-pin DIN) <p style="text-align: center;">- or -</p> • Keyspan-by-Tripp Lite High-Speed USB Serial Adapter (Note 3)

Note 1: Windows Vista and Mac OS 10.5 compatibility began with SoftMax Pro v5.3.

Note 2: Users integrating with a robotics system, manipulating large data files (>20 plates per file), or reading in FLEX mode are advised to use a computer system that meets or exceeds the Recommended specifications. For users with data files containing multiple plate section of kinetic data, selecting “Suspend Recalculation” from the *Edit* menu may improve system performance.

Note 3: Please visit the manufacturer’s website to download and install the latest software drivers for the USB serial adapter for full OS compatibility.