

BioFocus[®] 2000 and 3000 Systems Computer and Software Requirements

Hardware Requirements *

| | |
|-----------------------|--|
| Machine Type: | IBM PC (or 100% Compatible) |
| Machine Class: | 80486DX , 80486SX with an 80487SX upgrade, 80386DX with an 80387DX math coprocessor, or 80386SX with an 80387SX math coprocessor |
| Total RAM: | 4,194,304 bytes (commonly referred to as 4 megabytes or 4096 kilobytes)** |
| Disk Drives: | 100 megabyte , or greater 1.44 megabyte, 3.5", internal floppy disk drive, required 1.2 megabyte, 5.25", internal floppy disk drive, optional |
| Video System: | VGA (or 100% compatible) color graphics -or EGA (or 100% compatible) color graphics -or Hercules (or 100% compatible) monochrome graphics |
| Serial Ports: | Two RS-232 ports for asynchronous communications, configured as COM 1 (IRQ4, I/O Port Address 03F8) and COM 2 (IRQ3, I/O Port Address 02F8) |
| Mouse: | Microsoft (or 100% compatible) BUS mouse (IRQ5) (must not use IRQ 0,1,2,3,4,7 or LPT1 IRQ) |

* All hardware documentation, including computer and adapter card settings, must be available.

** This includes 3,801,088 bytes of User RAM (commonly referred to as 3.625 megabytes or 3712 kilobytes) configured as follows: 655,360 bytes of contiguous DOS lower memory, starting at memory address 0000:0000 (commonly referred to as 0.625 megabytes or 640 kilobytes), and 3,145,728 bytes of contiguous extended memory, starting at memory address 10000:0000 (commonly referred to as 3 megabytes or 3072 kilobytes). System RAM comes automatically with the system.

Hardware Options

| | |
|-----------------------|--|
| Parallel Port: | One (1) Centronix (or 100% compatible) for a printer (IRQ7) configured as LPT1 |
| Printer: | Centronix (or 100% compatible) parallel interface. Printer memory should be capable of processing full page graphics at 300 dpi. |

Operating System Requirements

| | |
|----------------------------------|--|
| MS (or PC) DOS: | Version 3.3x, Version 5.x, Version 6.0, 6.1, 6.2 |
| Environment Requirements: | Microsoft Windows 3.1 |
| Environment Option: | Microsoft Excel 4.0 for customizing Bio-Rad Integrator reports |