

Octet[®] Instrument Workstation

Pall ForteBio Octet instruments use Microsoft Windows-based computers to control instrument functions and to collect and analyze your experiment data, with Microsoft Office functions generating reports. The communication between the computer and the instrument require the use of USB and serial ports.

For best performance, Pall ForteBio recommends Octet instrument computers meet the following requirements.

Recommended System Requirements

- Windows 7 Professional, 64-bit
- 3.7 GHz Quad Core processor or greater
- 8 GB RAM
- 500 GB hard drive
 - 200 MB available hard disk space required for application files; more to acquire data – a single experiment can range from 1 MB to over 100 MB
- 3 available USB 2.0 ports
 - Octet control features do not support USB 3.0
- 2 available RS-232 serial ports
 - A PCI-Express serial port expansion card is provided with the instrument for adding a second serial port
- Microsoft Office 2013
- Widescreen monitor with 1920x1080 resolution

Notes:

- *Laptops are not recommended for use with data acquisition*
- *A DVD or CD drive is required to install the data acquisition and analysis software*
- *Computer workstations that are not supplied by Pall ForteBio are not covered under the Octet warranty or service contracts.*