

Agilent Seahorse Software Specifications



Instrument control & data acquisition software:

All Agilent Seahorse XFe, XFp, XF HS Mini, and XF Analyzer Controller (computer) specifications are listed below. The *XFe Analyzer Controller* has 2 factory default configurations – Windows 7, and Windows 10 on newer models.

Important – The factory default configuration must never be modified. Unauthorized modifications can result in unstable performance and void the instrument warranty. Contact Seahorse Technical Support if you have any questions about your XFe Analyzer Controller model.

| Instrument Type | Specifications | Software Name |
|--|--|--------------------------------|
| Seahorse XFe96 Analyzer Seahorse XFe24 Analyzer | Operating System: Windows 10, 64-bit Processor: Intel Core i5 System Memory: 16 GB RAM Latest Software: Wave Controller 2.6 | Seahorse Wave Controller |
| Seahorse XFe96 Analyzer Seahorse XFe24 Analyzer | Operating System: Windows 7, 32-bit Processor: Intel Core i3 System Memory: 4 GB RAM Latest Software: Wave Controller 2.4 | Seahorse Wave Controller |
| Seahorse XF HS Mini | Operating System: Windows 10 Embedded | Seahorse XF HS Mini |
| Seahorse XFp Analyzer | Operating System: Windows 7 Embedded | Seahorse XFp Analyzer |
| Seahorse XF96 Analyzer Seahorse XF24 Analyzer | Operating System: Windows Vista or 7, 32-bit Processor: Intel Celeron or Intel Core 2 Duo (2.2 GHz) System Memory: 2 GB RAM (minimum) | Seahorse XF96 Seahorse XF24 |



Data analysis software: Wave Desktop 2.6

Wave Desktop is the assay design & data analysis software for all Seahorse Analyzers and supports:

- Analysis of data files from all Seahorse Analyzers (XFe96, XFe24, XFp, XF96 & XF24).
- Create and customize assay templates for XFe96, XFe24 & XFp Analyzers.
- Export data to (1) Microsoft Excel, (2) GraphPad Prism, or (3) XF Report Generators.
- Supports Microsoft Excel (32 & 64-bit) for both Windows and Macintosh PCs.

| Computer | Specifications |
|--|--|
| Windows PC | Operating System: Windows 7, 8.1, and 10 Processor: Intel Core i3 (or better) Hard Disk Space: 175 GB System Memory (RAM): 4 GB (minimum*) Screen Resolution: 1280 x 800 (minimum) Supported Excel Versions: 2010, 2013 & 2016 |
| Macintosh PC <i>(requires use of a virtual machine)</i> | Operating System: Mac OSx 10.11 or higher Virtual Machine: Parallels 12 & Windows 7, 8.1 and 10 Processor: Intel Core i3 (or better) Hard Disk Space: 175 GB System Memory (RAM): 4 GB (minimum*) Screen Resolution: 1280 x 800 (minimum) Supported Excel Versions: 2011 & 2016 |

**For the optimal experience, 8 GB (or higher) of System Memory (RAM) is recommended.*



Optional Data Analysis Software: XF Reader Software

(for XF96 & XF24 Analyzers only)

- Supports analysis of data acquired by Agilent Seahorse XF96 and XF24 Analyzers *only*.
- Supports Microsoft Excel 32-bit *only*.

| Computer | Specifications |
|-----------------------|---|
| Windows PC (only) | Operating System: Windows 7, 8.1, and 10 Processor: Intel Core i3 (or better) Hard Disk Space: 175 GB System Memory (RAM): 4 GB (minimum*) Screen Resolution: 1280 x 800 (minimum) Supported Excel Versions: 2010, 2013 & 2016 |
| Incompatible Elements | Operating Systems: Windows 8, 8.1, 10; All Mac OS Virtual Machine: 2013, 2016, Office 365, Excel for Mac, Excel file types .xlsx & .xlsm |

www.agilent.com/chem/discoverxf

For Research Use Only.
Not for use in diagnostic procedures.

DE: 5757060185

This information is subject to change without notice.

© Agilent Technologies, Inc. 2020
 Printed in the USA, January 10, 2020
 57891-10000 Rev B