

Agilent Seahorse Software Specifications



Instrument control & data acquisition software:

All Agilent Seahorse XF Pro, 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	Latest Software Version
Seahorse XF Pro Analyzer	Operating System: Windows 10 Enterprise LTSC (Version 1809) 64-bit Processor: Intel Core i5 System Memory: 16 GB RAM	XF Pro Controller 10.0.1.72
Seahorse XFe96 Analyzer	Operating System: Windows 10 Enterprise LTSC (Version 1809) 64-bit Processor: Intel Core i5 System Memory: 16 GB RAM	Wave Controller 2.6.3
Seahorse XFe24 Analyzer	Operating System: Windows 10 Enterprise LTSC (Version 1809) 64-bit Processor: Intel Core i5 System Memory: 16 GB RAM	Wave Controller 2.6.1
Seahorse XFe96 Analyzer	Operating System: Windows 7, 32-bit Processor: Intel Core i3 System Memory: 4 GB RAM	Wave Controller 2.4.3
Seahorse XFe24 Analyzer	Operating System: Windows 7, 32-bit Processor: Intel Core i3 System Memory: 4 GB RAM	Wave Controller 2.4.2
Seahorse XF HS Mini	Operating System: Windows 10 Embedded (Version 1809)	XF HS Mini Controller software 3.0.0.41
Seahorse XFp Analyzer	Operating System: Windows 7 Embedded Operating System: Windows 10 Embedded	XFp 1.1.1.3 XFp 2.0.0.72
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)	XF96 1.4.2 software XF24 1.8.1 software



Data analysis software: Agilent Seahorse Analytics

Agilent Seahorse Analytics is a web-based software application for Seahorse data analysis, interpretation, and collaboration and supports:

- Analysis of data generated by all Agilent Seahorse XF Analyzers.
- Free XF data file storage up to 10 GB.
- Greater control and flexibility for data export to Microsoft Excel and GraphPad Prism.

Specification	Detail
Internet Browser Compatibility	Google Chrome Safari Mozilla Firefox
	Microsoft Edge (Internet Explorer is not recommended)
Export Options	Microsoft Excel (.xlsx)GraphPad Prism (.pzfx)Image (.jpg and .png)

Other data analysis software: Wave Pro 10.0.1

Wave Pro supports:

- Analysis of data generated by the Seahorse XF Pro and XF HS Mini Analyzers
- Create and customize assay templates for Seahorse XF Pro and XF HS Mini Analyzers
- Export data for (1) Microsoft Excel, (2) GraphPad Prism, or (3) XF Report Generators

Computer	Specifications	
Windows PC	Operating system: 10 Processor: Intel Core i3 (or better) Hard disk space: 175 GB	System memory (RAM): 4 GB (minimum*) Screen resolution: 1280 × 800 (minimum) Supported Excel versions: 2010, 2013, and 2016

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



Wave Desktop 2.6

Wave Desktop supports:

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

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

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

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.



