



Agilent Seahorse Software Specifications

Instrument Control and Data Acquisition Software

The Agilent Seahorse XFe and Agilent Seahorse XF Analyzer Controller (computer) enables instrument operation and data acquisition. The XFe Controller specifications listed in [Table 1](#) are factory default, and must never be modified. Unauthorized modifications to the XFe Controller can result in unstable performance and void the instrument warranty.

Table 1 XFe Controller specifications

| Instrument type | Specifications | Software name |
|--|--|--------------------------------|
| Seahorse XFe96 Analyzer Seahorse XFe24 Analyzer | Operating system: Windows 7, 32-bit Processor: Intel Core i3-3220 (latest) System memory: 4 GB RAM | Seahorse Wave Controller |
| 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 | Seahorse XF96 Seahorse XF24 |

Data Analysis Software: Wave Desktop 2.4

Wave Desktop is the assay design, data analysis, and file management software for the Agilent Seahorse XF Analyzers. Wave Desktop 2.4 supports:

- Analysis of data files from all Seahorse XF Analyzers (XFe96, XFe24, XFp, XF96, and XF24).
- Create and customize assay templates for the Seahorse XFe24, XFe96, and XFp Analyzers.
- Export data to (1) Microsoft Excel, (2) GraphPad Prism, or (3) Seahorse XF Report Generators.
- Supports Microsoft Excel (32 and 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): 2 GB minimum* Screen resolution: 1280 x 800 minimum Supported Excel versions: 2010, 2013, and 2016 |





| Computer | Specifications |
|---|--|
| Macintosh PC (requires use of a virtual machine) | Operating system: Mac OS w. Parallels 9 (or higher) and Windows 7, 10 Processor: Intel Core i3 (or better) Hard disk space: 175 GB System memory (RAM): 2 GB minimum* Screen resolution: 1280 x 800 minimum Supported Excel versions: 2011 and 2016 |

Optional Data Analysis Software: XF Reader Software

(For XF96 and 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 XP (Service Pack 3), Vista, or 7 (32 and 64-bit) Processor: Intel Core 2 Duo (2.2 GHz) Hard disk space: 175 GB System memory (RAM): 2 GB minimum Screen resolution: 1024 x 768 minimum Supported Excel versions: 2003, 2007, and 2010 |
| Incompatible elements | Operating systems: Windows 8, 8.1, 10; All Mac OS Microsoft Excel: 2013, 2016, Office 365, Excel for Mac, Excel file types *.xlsx and *.xlsm |

Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

© Agilent Technologies, Inc. 2017
Printed in USA
Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19808-1610 USA

For Research Use Only
Not for use in diagnostic procedures.



S7891-10000