

# Agilent 2100 Expert Software



## Compatibility matrix for supported operating systems

NOTE: With Windows 10 and Windows 11, this application requires the feature “.NET Framework 3.5” activated.

NOTE: We do not expect differences between “Enterprise”, “Education”, “Professional” or “Home” versions.

	B.02.12	B.02.11	B.02.10	B.02.09	B.02.08 (End of support)
Windows XP	Not tested. Not recommended.	Not tested. Not recommended.	Not tested. Not recommended.	Not tested. Not recommended.	Not tested. Expected to be compatible.
Windows 7 Service Pack 1	Not tested. Not recommended.	Not tested. Not recommended.	Not tested. Expected to be compatible.	Compatible	Compatible
Windows 10 Build 1511	Not tested. Not recommended.	Not tested. Expected to be compatible.	Compatible	Compatible	Not supported.
Windows 10 Version 1607	Not tested. Not recommended.	Not tested. Expected to be compatible.	Compatible	Compatible	Not supported.
Windows 10 Version 1709	Not tested. Not recommended.	Not tested. Expected to be compatible.	Compatible	Not tested. Expected to be compatible.	Not supported.
Windows 10 Version 1803	Not tested. Not recommended.	Not tested. Expected to be compatible.	Compatible	Not tested. Expected to be compatible.	Not supported.
Windows 10 Version 1809	Not tested. Not recommended.	Compatible	Compatible	Not tested. Expected to be compatible.	Not supported.
Windows 10 Version 1903	Not tested. Not recommended.	Not tested. Expected to be compatible.	Not tested. Expected to be compatible.	Not tested. Expected to be compatible.	Not supported.
Windows 10 Version 1909	Compatible.	Not tested. Expected to be compatible.	Not tested. Expected to be compatible.	Not tested. Expected to be compatible.	Not supported.
Windows 10 Version 2004	Not tested. Expected to be compatible.	Compatible	Not tested. Expected to be compatible.	Not tested. Expected to be compatible.	Not supported.
Windows 10 Version 21H2	Compatible	Compatible	Not tested. Expected to be compatible.	Not tested. Expected to be compatible	Not supported.
Windows 10 Version 22H2	Compatible	Not tested. Expected to be compatible.	Not tested. Expected to be compatible.	Not tested. Expected to be compatible	Not supported.

# Agilent 2100 Expert Software



## Compatibility matrix for supported operating systems

NOTE: With Windows 10 and Windows 11, this application requires the feature “.NET Framework 3.5” activated.

NOTE: We do not expect differences between “Enterprise”, “Education”, “Professional” or “Home” versions.

	B.02.12	B.02.11	B.02.10	B.02.09	B.02.08 (End of support)
Windows 11 Version 21H2	Compatible	Not supported.	Not supported.	Not supported.	Not supported.
Windows 11 Version 22H2	Compatible	Not supported.	Not supported.	Not supported.	Not supported.
Windows 11 Version 23H2	Not tested. Expected to be compatible	Not supported.	Not supported.	Not supported.	Not supported.
Windows 11 Version 24H2	Compatible	Not supported.	Not supported.	Not supported.	Not supported.

# Agilent 2100 Expert Software



## Compatibility matrix for supported hardware

NOTE: Max. of two instruments per PC are supported. 1x RS-232 and 1x USB/Serial connection. 2x USB/Serial connection for 2100 Bioanalyzer bundle laptop HP ProBook 850 G8 and higher

		B.02.12	B.02.11	B.02.10	B.02.09	B.02.08 (End of support)
End of support	G2938A (B.02.06 required)	Not compatible	Not compatible	Not compatible	Not compatible	Not compatible
End of guaranteed support/ support on best-effort basis	G2938B	Compatible	Compatible	Compatible	Compatible	Compatible
	G2938C	Compatible	Compatible	Compatible	Compatible	Compatible
	G2939A	Compatible	Compatible	Compatible	Compatible	Compatible
Guaranteed support	G2939B	Compatible	Compatible	Compatible	Compatible	Not compatible