

Agilent 2100 Expert Software



Compatibility matrix for supported operating systems

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

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

	B.02.11	B.02.10	B.02.09	B.02.08
Windows XP	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. Expected to be compatible.	Compatible	Compatible
Windows 10 Build 1511	Not tested. Expected to be compatible.	Compatible	Compatible	Not supported.
Windows 10 Version 1607	Not tested. Expected to be compatible.	Compatible	Compatible	Not supported.
Windows 10 Version 1709	Not tested. Expected to be compatible.	Compatible	Not tested. Expected to be compatible.	Not supported.
Windows 10 Version 1803	Not tested. Expected to be compatible.	Compatible	Not tested. Expected to be compatible.	Not supported.
Windows 10 Version 1809	Compatible	Compatible	Not tested. Expected to be compatible.	Not supported.
Windows 10 Version 1903	Not tested. Expected to be compatible.	Not tested. Expected to be compatible.	Not tested. Expected to be compatible.	Not supported.
Windows 10 Version 1909	Not tested. Expected to be compatible.	Not tested. Expected to be compatible.	Not tested. Expected to be compatible.	Not supported.
Windows 10 Version 2004	Compatible	Not tested. Expected to be compatible.	Not tested. Expected to be compatible.	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.

		B.02.11	B.02.10	B.02.09	B.02.08
End of Support	G2938A (B.02.06 required)	Not compatible	Not compatible	Not compatible	Not compatible
Support Phase	G2938B	Compatible	Compatible	Compatible	Compatible
	G2938C	Compatible	Compatible	Compatible	Compatible
	G2939A	Compatible	Compatible	Compatible	Compatible
Active Phase	G2939B	Compatible	Compatible	Compatible	Not compatible