



Agilent Technologies

Compatibility of Chemstation software with Windows Operating System

This document is believed to be accurate and up-to-date. However, Agilent Technologies, Inc. cannot assume responsibility for the use of this material.
The information contained herein is intended for use by informed individuals who can and must determine its fitness for their purpose.



Title :Compatibility of Chemstation software with Windows Operating System

	A.03.0x	A.04.0x	A.05.0x	A.06.0x	A.07.0x	A.08.0x	A.09.0x	A.10.0x	B.01.01	B.01.03	B.01.03 SR1	B.02.01	B.02.01 SR 1
Win 3.1 / WFW 3.11	Yes	Yes	No	No	No	No	No	No	No	No	No	No	No
Win 95	No	Yes	Yes	Yes	No	No	No	No	No	No	No	No	No
Win NT 4.0	No	No	Yes	Yes	Yes	Yes	Yes	No	No	No	No	No	No
SP3	No	No	Yes	Yes	No	No	No	No	No	No	No	No	No
SP4	No	No	No	Yes	Yes	No	No	No	No	No	No	No	No
SP5	No	No	No	No	No	No	No	No	No	No	No	No	No
SP6a	No	No	No	No	Yes	Yes	Yes	No	No	No	No	No	No
Win 2000 Pro	No	No	No	No	No	No	Yes	Yes	Yes	Yes	Yes	Yes	Yes
SP2	No	No	No	No	No	No	Yes	Yes	No	No	No	No	No
SP3	No	No	No	No	No	No	A.09.03 Only	Yes	No	No	No	No	No
SP4	No	No	No	No	No	No	A.09.03 Only	Yes	Yes	Yes	Yes	Yes	Yes
	A.03.0x	A.04.0x	A.05.0x	A.06.0x	A.07.0x	A.08.0x	A.09.0x	A.10.0x	B.01.01	B.01.03	B.01.03 SR1	B.02.01	B.02.01
Win XP Pro SP1A	No	No	No	No	No	No	No	Yes	Yes	No	No	No	No

This document is believed to be accurate and up-to-date. However, Agilent Technologies, Inc. cannot assume responsibility for the use of this material.
The information contained herein is intended for use by informed individuals who can and must determine its fitness for their purpose.

Win XP Pro SP2	No	No	No	No	No	No	No	No	Yes	Yes	Yes	Yes	Yes
-------------------	----	----	----	----	----	----	----	----	-----	-----	-----	-----	-----