

SureScan Scan Control SW Revisions and PC/OS Compatibility

ScanControl Revision (Release Date)	HP PC Model	Microsoft Operating System	
		Edition	Version
9.1.15.0 (Sep-2022)	Z2 G9	Windows 11-64 bit Pro	21H2
	Z2 G5	Windows 11-64 bit Pro	23H2
		Windows 10-64 bit Pro	22H2
			20H2
9.1.13.1 (Jan-2020)	Z2 G5	Windows 10-64 bit Pro	20H2
	Z2 G4	Windows 10-64 bit Pro	20H2
			1909
			1903
	Z240	Windows 10-64 bit Pro	1909
9.1.11.13 (Jan-2018)	Z2 G4	Windows 10-64 bit Pro	1903
			1803
	Z240	Windows 10-64 bit Pro	1903
			1703
9.1.11.7 (Sep-2015)	Z240	Windows 10-64 bit Pro	1703
	Z230	Windows 7-64 bit Pro	SP1
9.1.9.1 (Jan-2015)	Z230	Windows 7-64 bit Pro	SP1
9.1.7.1 (Dec-2012)	Z220	Windows 7-64 bit Pro	SP1
9.1.5.1 (Mar-2012)			
9.1.3.1 (Feb-2012)			
9.1.1.1 (Aug-2011)	Z210	Windows 7-64 bit Pro	SP1