

Compatibility list – Freedom EVOware 2.9

Operating system	Version
Microsoft Windows	Windows 10 Enterprise LTSC 21H2 Windows 11 IOT Enterprise LTSC 2024

Freedom EVOware add-ons	Version
Sample Oriented	1.3
Sample Tracking	1.5
Normalization Wizard	2.2
Hit-Picking Wizard	1.0 SP3
TouchTools Suite	3.3 SP3
Te-Chrom	1.1
Buffer Creation Wizard	1.2
QC Kit	1.3
Weigh Station Wizard	1.0 SP2
PCR Wizard	1.1 SP1

Tecan software packages	Version
Instrument Software	8.1
Magellan Pro	7.5
I-control	2.0
EVOware Reader Driver	1.12.2
HydroControl	4.3.1
EVOware Hydro Washer Driver	4.3.3
Magellan SPARK Control	3.2 SP1

Tecan Common Components	Version
Communication	2.3.0-14
Logging	2.3.3-01
PosID-3	2.3.2-01
PMP	2.3.1-11
UMS	1.1.14
IOT Client	2.8 SP1

Other Tecan software	Version
WinWash Plus	2.22
RdrOle	4.51
RdrOle XFluor4 Safire2 add-on	4.62

Other software packages	Version
Microsoft Outlook	16.0 (Outlook 2016, 2019)
Microsoft Office	Excel 2016

Tecan instrument modules	Firmware version
DC Servo 2	1.21

LiHa	1.50,1.51,1.60,1.70*, 1.72, 1.80
MCA384	1.01, 1.10, 1.11
MCA96	1.20, 1.21
MIO	1.41
MIO2	2.02, 2.10
MPO	2.00, 3.20
PMP	1.02
PnP	1.08
PnP2	2.00
PosID-2	2.51
PosID-3	1.21
RoMa (DC Servo 1)	1.13
RoMa (DC Servo 2)	2.20, 2.21
SAFY	1.30
SPO	3.20
Te-Link	1.00
Te-MagS	1.30
Te-MO 3/5	2.02
Te-MO 3/3	2.02
Te-MO WRC	2.11
Te-PS	1.00
Te-Servo	1.00
Te-Shake	1.10
Te-Stack	1.40, 1.50
Te-VacS	2.10
TeCU	1.40
XP2000	1.10
XMP	1.0.1
ZAAP	1.1, 1.2
Readers & Washers	
HydroSpeed	1.70 and 1.71
HydroFlex	4.20

External Tecan devices	Firmware version
Carousel/Carousel HS	3.00
Carousel LPT220 EVO	7.2.0.0
Carousel LPT280 EVO	8.0.0.1
Washer PW384 (with PW384GP driver)	2.06
Washer PW384 (via WinWash Plus)	See section Other Tecan software
Washer 96PW	3.0
Washer Columbus	4.27
Sunrise (via RdrOle)	See section Other Tecan software
SpectraFluor (via RdrOle)	See section Other Tecan software
SpectraFluor Plus (via RdrOle)	See section Other Tecan software
Genios (via RdrOle)	See section Other Tecan software
Genios Pro (via RdrOle)	See section Other Tecan software
Safire (via RdrOle)	See section Other Tecan software
Ultra (via RdrOle)	See section Other Tecan software

3 rd party devices	Firmware version
Barcode scanner MS954	-
Barcode scanner LS1220	-
Centrifuge Hettich Rotanta 46 RS	4.070
Centrifuge Hettich Rotanta 460 RSC	1.18
Balance Mettler AG285	MT-SICS 0
Balance Mettler SAG 285/01	MT-SICS 0
Waterbath Lauda	2.2