

Compatibility List - FluentControl 3.7 Service Pack 1

Operating system	Version
Microsoft Windows	Windows 10 Enterprise (IoT) LTSC 2021 Windows 11 Enterprise (IoT) LTSC 2024

Tecan software packages*	Version
IoT Client	3.0
Magellan Pro	7.5 (Standard & Tracker)
SPARKControl Magellan	3.2 SP1
Sample Tracking	1.4 SP2 [†]
User Management System (UMS)	1.1.13
Identity Access Management (IAM)	1.1
Licensing Client	1.2.0
Hydrowasher Driver	4.3.2
InhecoMTC Driver	2.0
SiLA Client	1.4.2
SiLA2 Client	2.3
Driver Framework	3.2.30 or higher [‡]
Optimized Log Extractor G (OLEG)	2.1.0.36743
MAPlinx	Bull SP1
MAPlinx Setup	6.0 [§]

Other software packages	Version
Microsoft Office	Excel 2016
eGalaxTouch	5.14.0.24318
DisplayLink USB Graphics Software for Windows	11.6 M1
IDS uEye	4.91
Arena SDK (Arena Library only)	1.0.27.9

Tecan instrument modules	Firmware version
Te-Control II / Te-Control II Bootloader	1.3.0.11110 / 1.3.0.11110
Te-Motion	1.10.0.11416
Te-Motion II	4.2.1
CxP	1.3.0.11238
cLLD Main / Support	1.0.0.10483
PMP	1.0.0.10470
Te-GIO	1.3.0.10860

* Components/drivers, that are part of FluentControl, are not explicitly listed.

[†] Sample Tracking 1.4 SP2 is not compatible with IAM. Please contact Tecan if Sample Tracking and User Administration is needed on the system.

[‡] Some drivers require a software package from a 3rd party manufacturer. For such drivers, consult the documentation of the 3rd party manufacturer to ensure Windows version compatibility.

[§] MAPlinx Setup 6.0 is not compatible with Windows 11. For Resolvex i300 systems, Windows 10 must be used.

Te-GIO-II	1.1.0.0
Te-Shake	1.10
Te-Stack	1.50
Wash and Refill Center	2.11
Te-VacS	2.10
MIO2	2.10
Fluent Carousel	0.9A
Loading ID Fluent	1.10.0.11.408
Balance	MT-SICS Level 0 and 1
Keyence	1.15
UniBootloader	1.0.0.10420
UniBootloader-II	1.3.1
Te-MagS	1.31