

Compatibility List - FluentControl 3.7

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
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	3.1.12 or higher
Optimized Log Extractor G (OLEG)	2.1.0.36743
M.A.P.	Bull – RC3

Other software packages	Version
Microsoft Office	Excel 2016
eGalaxTouch	5.14.0.24318
DisplayLink USB Graphics Software for Windows	11.5 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	4.1.1
CxP	1.3.0.11238
cLLD Main / Support	1.0.0.10483
PMP	1.0.0.10470
Te-GIO	1.3.0.10860
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

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

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