

## Compatibility List - FluentControl 3.8 SP1

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

  

Tecan software packages*	Version
IoT Client	3.1
Magellan Pro	7.5 (Standard & Tracker)
SPARKControl Magellan	3.2 SP1
Sample Tracking	1.5
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	1.5
SiLA2 Client	2.3
Driver Framework	3.3.14 or higher†
Optimized Log Extractor G (OLEG)	2.1.0.36743
MAPlinx	Crocodile
MAPlinx Setup	6.2

  

Other software packages	Version
eGalaxTouch	5.14.0.24318
DisplayLink USB Graphics Software for Windows	12.1
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
Te-GIO-II	1.1.0.0
Te-Shake	1.10

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

† The Driver Framework is compatible with Windows 10 and Windows 11. Some drivers require a software package from a 3rd party manufacturer. For such drivers, consult the documentation of the 3rd party manufacturer to ensure appropriate operating system compatibility.

Te-Stack	1.50
Wash and Refill Center	2.11
Te-VacS	2.10
MIO2	2.10
Fluent Carousel	0.9A
Loading ID Fluent	1.10.0.11408
Balance	MT-SICS Level 0 and 1
Keyence	1.15
UniBootloader	1.0.0.10420
UniBootloader-II	1.3.1
Te-MagS	1.31