

Compatibility List - FluentControl 3.4 SP1

Operating system	Version
Microsoft Windows	Windows 10 Enterprise LTSC 2019 Windows 10 Enterprise LTSC 2021
Tecan software packages*	Version
IoT Client	2.6
Magellan Pro	7.5 (Standard & Tracker)
SPARKControl Magellan	3.2
Sample Tracking	1.4 SP2
User Management System (UMS)	1.1.13
Licensing	1.1.0
Hydrowasher Driver	4.3.2
InhecoMTC Driver	2.0
SiLA Driver	1.4.2
SiLA2 Driver	2.0
Driver Framework	3.1.12 or higher
Other software packages	Version
Microsoft Office	Excel 2016
eGalaxTouch	5.14.0.15315
DisplayLink USB Graphics Software for Windows	10.3 M0
IDS uEye	4.91
Arena SDK (Arena Library only)	1.0.27.9
Tecan instrument modules	Firmware version
TeControl / TeControl Bootloader	1.3.0.11110 / 1.3.0.11110
TeMotion	1.10.0.11416
TeMotion II	3.3.0
CxP	1.3.0.11238
cLLD Main / Support	1.0.0.10483
PMP	1.0.0.10470
TeGIO	1.3.0.10860
TeGIO-II	1.2
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
Balance	MT-SICS Level 0 and 1

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

Keyence	1.15
UniBootloader	1.0.0.10420
UniBootloader-II	1.3.1
Te-MagS	1.31