

Compatibility List - FluentControl 3.5

Operating system	Version
Microsoft Windows	Windows 10 Enterprise LTSC 2019 Windows 10 Enterprise LTSC 2021

Tecan software packages*	Version
IoT Client	2.7.5
Magellan Pro	7.5 (Standard & Tracker)
SPARKControl Magellan	3.2
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	contact Tecan
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	11.1 M1
IDS uEye	4.91
Arena SDK (Arena Library only)	1.0.27.9

Tecan instrument modules	Firmware version
TeControl2 / TeControl2 Bootloader	1.3.0.11110 / 1.3.0.11110
TeMotion	1.10.0.11416
TeMotion II	4.0.1
CxP	1.3.0.11238
cLLD Main / Support	1.0.0.10483
PMP	1.0.0.10470
TeGIO	1.3.0.10860
TeGIO-II	1.1.0.0
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

* 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