

Compatibility List - FluentControl 3.3

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
SPARKControl Magellan	3.0 [†]
Sample Tracking	1.4 SP1
User Management System (UMS)	1.1.13
Licensing	1.1.0
Hydrowasher Driver	4.3.0 [‡]
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
TeMotion II	3.1.2
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
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

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

[†] FluentControl 3.3 is not compatible with version 3.1 of SPARKControl Magellan.

[‡] FluentControl 3.3 is also compatible with Hydrowasher Driver 4.3 SP2(*release pending*)

Balance	MT-SICS Level 0 and 1
Keyence	1.15
UniBootloader	1.0.0.10420
UniBootloader-II	1.3.1
Te-MagS	1.31