

**Compatibility list – vControl 1.2, Build: 3.9.4.1178**

<b>Operating system</b>	<b>Version</b>
Microsoft Windows	Windows 10 Enterprise LTSC 2021

<b>Tecan software packages*</b>	<b>Version</b>
IoT Client	3.1.0.326
MAPlinx Setup	6.3
MAPLinxBASE	Eagle RC2
3DSim	4.0.0.230
Magellan Pro	7.5 (Standard & Tracker) 7.5 SP1 (Standard & Tracker)
Sample Tracking	1.5
Licensing Client	1.2.0
Hydrowasher Driver	4.3.2
SiLa Client	1.4.2
SiLA2 Client	2.3
Optimized Log Extractor G (OLEG)	2.1.1.37327
Tecan IAM	1.1

<b>Other software packages</b>	<b>Version</b>
Microsoft Office	Excel 2016
Microsoft .NET Framework	4.8
Arena SDK (Arena Library only)	1.0.27.9

<b>Tecan instrument modules</b>	<b>Firmware version</b>
TeMotion I	1.10.0.11416
TeMotion II	5.0.0-0-g99f8cf5a
UniBootloader II	1.3.1.0
ARP/ARC	2.1.1.0
LoadingID Application	1.10.0.11408
LoadingID Bootloader	1.0.0.10420
TeGIO	1.3.0.10860
TeGIO-II	1.1.0.0
TeControl Application	1.3.0.11110
TeControl Bootloader	1.3.0.11110
CLLD Main/Support	1.0.0.10483
PMP	1.0.0.10470

<b>Other instrument modules</b>	<b>Firmware version</b>
Inheco MTC/STC	2.83

Other instrument modules	Firmware version
Inheco Black Slot Module	2.83
Inheco Yellow Slot Module	2.87
Inheco Blue Slot Module	2.53
Inheco Teleshake95 AC Application	MTC_SlotTS2.87_04/21
Inheco Teleshake95 AC	Teleshake_AC_FW43
Inheco Teleshake95 AC Bootstrapper	No_BOOT_Attention!
Inheco Incubator Shaker	IncShake_C_V3.50_04/2012
Inheco Incubator Shaker Bootstrapper	BOOT_C_V3.12_05/2007
Inheco ODTC Build	7783.22980
Inheco ODTC Device	ODTC_277_BOOT_001_HW_00003
Balance	MT-SICS Level 0 and 1

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