

## Compatibility list – vControl 1.1 SP1, Build: 3.8.12.735

| Operating system  | Version                         |
|-------------------|---------------------------------|
| Microsoft Windows | Windows 10 Enterprise LTSC 2021 |

| Tecan software packages*         | Version                  |
|----------------------------------|--------------------------|
| IoT Client                       | 3.1.0.326                |
| MAPlinx Setup                    | 6.1 SP1                  |
| MAPlinxBase                      | Cobra SP2                |
| 3DSim                            | 3.13.0.137               |
| Magellan Pro                     | 7.5 (Standard & Tracker) |
| Sample Tracking                  | 1.4 SP2 <sup>+</sup>     |
| Licensing Client                 | 1.2.0                    |
| Hydrowasher Driver               | 4.3.2                    |
| SiLa Client                      | 1.4.2                    |
| SiLA2 Client                     | 2.3                      |
| Optimized Log Extractor G (OLEG) | 2.1.0.36743              |
| User Management System (UMS)     | 1.1.13                   |
| Tecan IAM                        | 1.1                      |

| Other software packages        | Version    |
|--------------------------------|------------|
| Microsoft Office               | Excel 2016 |
| Microsoft .NET Framework       | 4.8        |
| Arena SDK (Arena Library only) | 1.0.27.9   |

| Tecan instrument modules | Firmware version   |
|--------------------------|--------------------|
| TeMotion I               | 1.10.0.11416       |
| TeMotion II              | 4.2.1.0            |
| UniBootloader II         | 1.3.1.0            |
| ARP/ARC                  | 2.1.1-0-g7fe9e5cd  |
| LoadingID Application    | 1.10               |
| LoadingID Bootloader     | 1.0.0.10420        |
| TeGIO                    | 1.3.0.10860        |
| TeGIO-II                 | V1.1.0-0-g8e1cb355 |
| TeControl Application    | 1.3.0.11110        |
| TeControl Bootloader     | 1.3.0.11110        |
| MTC/STC                  | 2.83               |
| Black Slot Module        | 2.83               |
| Yellow Slot Module       | 2.87               |

| <b>Tecan instrument modules</b> | <b>Firmware version</b>  |
|---------------------------------|--------------------------|
| Blue Slot Module                | 2.53                     |
| Teleshake95AC Application       | MTC_SlotTS2.87_04/21     |
| Teleshake95 AC                  | Teleshake_AC_FW43        |
| Teleshake95AC Bootstrapper      | No_BOOT_Attention!       |
| Inheco Incubator                | IncShake_C_V3.50_04/2012 |
| Inheco ODTC                     | 277                      |
| CLLD Main/Support               | 1.0.0.10483              |
| PMP                             | 1.0.0.10470              |
| HEPA Filter                     | 3.1.8.756                |
| UV Light                        | 3.0.20.756               |
| Interior Light                  | 2.0.19.756               |
| RGB Light                       | 3.0.19.756               |
| Cooling Fans                    | 3.0.20.756               |
| WXS Balance                     | MT-SICS Level 0 and 1    |

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

\*Sample Tracking 1.4 SP2 is not compatible with IAM. Please contact Tecan if Sample Tracking and User Administration is needed on the system.