

Release Notes

Magellan™ V7.5 SP1

1.	PURPOSE	2
2.	VERSION INFORMATION	2
3.	SYSTEM REQUIREMENTS	2
4.	RELEASE NOTES	4
4.1.	General Features	4
4.1.1.	UMS Update	4
4.1.2.	IVDR Compliance	4
4.2.	General Information	4
4.2.1.	Cybersecurity Information	4
4.2.2.	Archive Files with Magellan Tracker	5
4.2.3.	Installation on Windows 10/Windows 11	5
4.3.	Known Anomalies	5
4.3.1.	UMS with Active Directory	5
4.3.2.	Registration under Windows 10/Windows 11	5
4.3.3.	Registration under Windows 10/Windows 11 with Office 365	6
4.3.4.	Compatibility with Firmware F50 (FW-Package 2.01) under Windows 10 / Windows 11	6
4.3.5.	Auto Archive	6
4.3.6.	Upgrade of User Management (Magellan Tracker)	6
4.3.7.	Uninstall Magellan	6
4.3.8.	Kinetic Duration	6
4.3.9.	Automatic Print Limitation	7
4.3.10.	Absorbance dual-wavelength measurement with identical measurement and reference values	7
4.3.11.	Magellan Window vs. Wizard	7
4.3.12.	Copy to Excel and Copy to ASCII	7

1. Purpose

This document describes Release Notes for the following released software package and related hardware and firmware.

2. Version Information

Module	Version
Magellan	7.5 SP1 (Setup 7.5.1.2)
.NET Framework	3.5 4.7.2
User Management System (UMS)	1.1.15
Infinite F50	Series
Firmware F50 (FW-Package 2.01)	MAI (V2.01)
Firmware F50 (FW-Package 3.33)	MAI (V3.33)
Firmware F50 ROBOTIC (FW-Package 3.34)	MAI (V3.34)
Sunrise	Series
Firmware Sunrise	V3.51

3. System Requirements

	Supported	Recommended
PC	Windows compatible PC with a Pentium compatible processor running at 1 GHz (Dual Core)	2 GHz (Dual Core)
Operating System	Windows 10 (32-bit)	
	Windows 10 (64-bit) Editions: Pro Windows® 11 Enterprise LTSC 24H2*	Windows 10 (64-bit) Editions: Pro
Memory	Windows 10 (32-bit): 1 GB RAM	2 GB RAM
	Windows 10 (64-bit): 2 GB RAM	4 GB RAM
	Windows 11 Enterprise LTSC 24H2 (64-bit): 4 GB RAM	4 GB RAM
Free Hard Disk Space	3 GB	5 GB
Monitor	Super VGA Graphics	
Resolution	1024 x 600 and higher	1920 x 1080
Color Depth	256	
Mouse	Microsoft mouse or compatible pointing device	

	Supported	Recommended
Communication	1 x USB 2.0	2 x USB 2.0 1 x RS232 (Serial)
Devices	Windows 10/Windows 11: DirectX 9 graphics device with driver version WDDM 1.0 onwards.	
.NET	.NET Framework 3.5 (in Windows 10/Windows 11 the user will be prompted to install .NET Framework 3.5 if it is not already present) .NET Framework 4.7.2	
Windows Installer	3.1 If this version is not present, the install/upgrade program will install it.	
Microsoft Excel	2007, 2010, 2013, 2016, 2019 (32-bit) Microsoft Excel 365 (32-bit) Only 32-bit editions supported! Starter editions NOT supported!	2010 (32-bit) 2019 (32-bit)

See 30267405 NOTICE MAGELLAN V7.5 SP1 WIN11 COMP

4. Release Notes

Magellan V7.5 SP1 has been developed and documented to meet the In Vitro Diagnostic Medical Devices Regulation (IVDR, Regulation (EU) 2017/746), and is a dedicated software version for Tecan absorbance readers like Infinite F50 and Sunrise.

The new version of Magellan introduces a Cybersecurity service pack for Magellan 7.5 as well as an update UMS component.

4.1. General Features

4.1.1. UMS Update

The UMS component has been updated to V1.1.15.

4.1.2. IVDR Compliance

Magellan V7.5 SP1 is developed to meet the requirements in conjunctions with IVDR, Regulation (EU) 2017/746.

4.2. General Information

4.2.1. Cybersecurity Information



CAUTION

IT IS THE USER'S RESPONSIBILITY TO TAKE THE NECESSARY PRECAUTIONS TO PROTECT THE SYSTEM FROM CYBERSECURITY THREATS.

WHEN INSTALLING VIRUS PROTECTION SOFTWARE OR SECURITY-RELATED UPDATES OF THE OPERATING SYSTEM PLEASE FOLLOW THE RECOMMENDATIONS OF THE LOCAL IT DEPARTMENT. BACK UP MAGELLAN FILES BEFORE STARTING UPDATES. RUN INSTALLATION QUALIFICATION (IQ), OPERATION QUALIFICATION (OQ) AND PERFORMANCE QUALIFICATION (PQ) OF MAGELLAN AFTER EVERY SYSTEM MODIFICATION.

Back up your data:

The importance of backing up files to be able to recover in case of a ransomware attack cannot be overstated. Be sure to create destruction-resistant backups of your critical systems and data. There are many tools and services available for file backup, restoration (of previous file versions) and recovery. Also, be sure to regularly test that the backups are working.

Use Magellan archive functionality or your product of choice to backup Magellan data files.

Manage and control privileged access to data:

To minimize risk of credential compromise and abuse, implement unique local administrator passwords on all systems, separate and protect privileged accounts and reduce broad permissions on file repositories.

Use standard windows operators (not administrators) to operate Magellan.

Use Magellan Tracker functionality to control access and further secure the application.

Operation Qualification

In addition to the installation qualification check, Tecan also recommends that the calculation abilities of Magellan are tested.

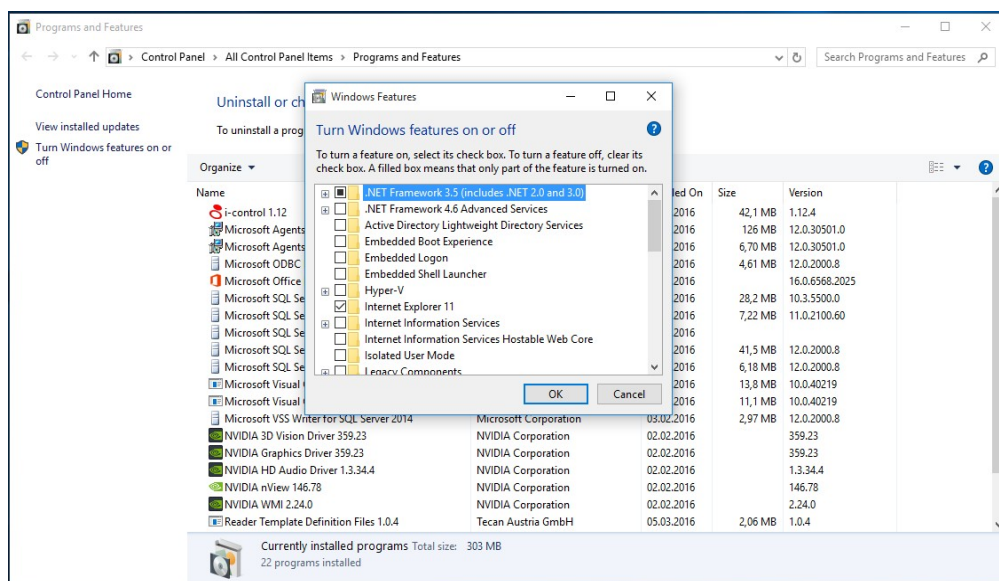
4.2.2. Archive Files with Magellan Tracker

To archive Magellan files (remove files from folder in Magellan default path and transfer them to folder in Archive path) Windows Administrator rights are necessary.

4.2.3. Installation on Windows 10/Windows 11

To install Magellan on a PC with the operating system Windows 10/11, .NET Framework 3.5, has to be installed. This Framework includes the Framework version .NET 2.0 that is required to operate Magellan; later .NET Framework versions (> 4.0) do not include V2.0 anymore.

The framework version V3.5 is not included on the Magellan CD and therefore has to be downloaded or installed from the operating system installation CD. It can be installed via Control Panel > Programs and Features. Open Turn Windows features on or off and select .NET Framework 3.5.



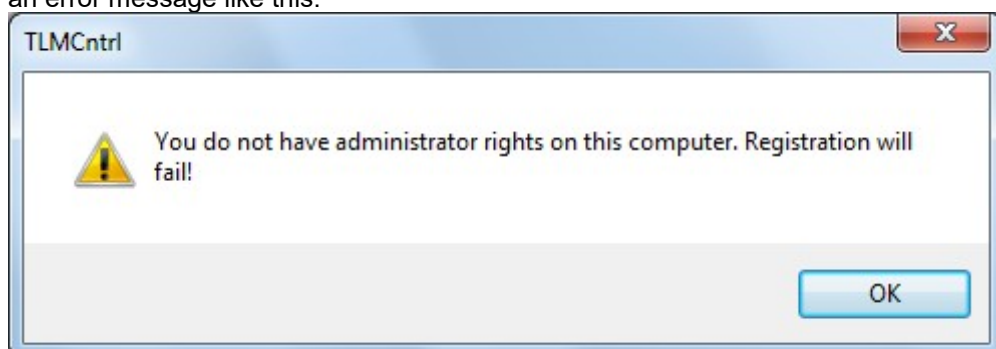
4.3. Known Anomalies

4.3.1. UMS with Active Directory

If UMS is used together with Active Directory, the logins of unknown users and logins with wrong passwords are not logged in the UMS Audit Trail nor in the Magellan System Audit Trail.

4.3.2. Registration under Windows 10/Windows 11

When registering Magellan under Windows 10/11, Tecan recommends starting Magellan with the option "Run as Administrator", otherwise the registration process might lead to an error message like this:



Nevertheless, the registration is still valid. For confirmation of the successful registration just view the About Dialog and check that the License is correct.

4.3.3. Registration under Windows 10/Windows 11 with Office 365

When using a PC with Office 365, the proposed solution in Chapter 4.3.2. prevents a TLM error but leads to the following subsequent error: Outlook, used for sending the licensing request to Tecan within the registration procedure, cannot be started automatically.

Workaround: save license request, attach it to an e-mail manually and send it to Tecan.

4.3.4. Compatibility with Firmware F50 (FW-Package 2.01) under Windows 10 / Windows 11

When working with Infinite F50 and firmware package 2.01 under Windows 10/11, PC-related USB connection problems may appear. Contact Tecan for further support.

4.3.5. Auto Archive

If the auto archive storage path is defined on a network directory and the archiving is triggered when the connection is temporarily unavailable, the System Audit Trail will not give any indication of that error in the log file. In such a case, the Microsoft operating system buffers the data to be written and writes the data as soon as the connection is stable again. In such a case, Tecan advises to make sure to check if the archiving was finished successful by the operating system.

4.3.6. Upgrade of User Management (Magellan Tracker)

If the default roles are modified, only the index of the corresponding right is displayed in the AuditTrail. If new roles are created the description of the right is also displayed in the AuditTrail.

Upgrade of Unused Users from Older Magellan Version:

If a user has never logged into the previous Magellan installation, the user is migrated to Magellan Tracker 7.x, but at first login the password does not have to be reset.

Upgrade of Groups, Roles and Option Settings from Older Magellan Version

Default groups or roles are reset to their default rights. User defined groups/roles are migrated with the name but without the defined rights. Option settings are reset to default values. To avoid losing that information, write down the changes already made to the groups/roles and option settings of your current Magellan version before the migration and reset those values after the upgrade to Magellan Tracker 7.x.

4.3.7. Uninstall Magellan

During the uninstall procedure of Magellan two dialog boxes for the Tracker version (one dialog box for the Standard version) are displayed. These dialog boxes handle the removal of User Management Administration and Log-files. They might appear behind the main de-installation wizard window. Please move those windows aside to get access to those dialogs. Our recommendation is to move the dialog to the side of the screen at the start of de-installation.

Workaround:

Move dialog to the side of the screen at the beginning of the uninstall procedure.

4.3.8. Kinetic Duration

For measurements in combination with a defined kinetic duration instead of a defined number of cycles, minimal variations in the measurement processing time may result in a different total number of cycles for each measurement label.

4.3.9. Automatic Print Limitation

Before starting an automatic print, Magellan checks the availability of the standard printer. Because of limitations of the built-in Windows function (name of the printer is limited to 32 characters) to perform this check, the check fails when the name of the standard printer is longer than 32 characters and no automatic print is performed.

Workaround:

Change the standard printer name to less than 32 characters.

4.3.10. Absorbance dual-wavelength measurement with identical measurement and reference values

When an absorbance dual-wavelength measurement is performed and the difference value (measurement value minus reference value) is exactly zero, Magellan might crash when further division is done with this zero value.

Workaround:

Unfortunately, no workaround is currently available. The probability for this is considered to be very low.

4.3.11. Magellan Window vs. Wizard

It might be possible that in rare cases the wizard is not displayed correctly in front of the main window (grey background). In such cases the wizard has moved behind the main window.

Workaround:

Click on the grey area of the main window to move the wizard to the front.

4.3.12. Copy to Excel and Copy to ASCII

If only individual wells are measured (contain data) copying to Excel does not work if the upper left well does not contain data.

Workaround:

Measure upper left well or entire plate if you want to use Copy to Excel/ASCII or use the Export options.