

# RELEASE NOTES

## OmniTrail

Release: Version 1.7a, June 2025.

© 2025 Malvern Panalytical B.V. All rights reserved.

Thank you for using OmniTrail.

### 1 Introduction

OmniTrail is an OmniTrust software module to view and audit system audit trail events in an OmniTrust environment.

OmniTrail access and authorizations are configured through the OmniTrust module OmniAccess.

### 2 Differences between versions

#### Differences between version 1.7a and 1.7

New features

- None

Fixed issues

- Solved an issue that sometimes (depending on the combination of licenses) prevented a user to see all relevant audit trail entries
- Merged the release and installation notes into one PDF document

#### Differences between version 1.7 and 1.6

New features

- Added support for SuperQ
- Updated version of license manager
- Security updates to installer

Fixed issues

- Special characters used in filenames no longer get displayed as %xx in the system audit trail. E.g., # was displayed as %23.

#### Differences between version 1.6 and 1.5

New features

- None

Fixed issues

- When writing to the system audit trail, the correct version (1.6) of OmniTrail is reported. In OmniTrail 1.5 version 1.4 was reported.

#### Differences between versions 1.5 and 1.4

New features

- None

Fixed issues

- Updated help documentation.

#### Differences between versions 1.4 and 1.3

New features

- Support showing the new Signing and Reason configuration.
- When using Central Management, upon deploying a configuration, a notification is shown

telling the users to restart any running applications.

Fixed issues

- Minor UI issues.

### **Differences between versions 1.3 and 1.2**

New features

- It is now possible to see information regarding the ElectronicRecordNotFound event.
- A conversion event review can only be done by users with permissions to do so.
- Conversion event details can now be found on the additional data.
- The user can now view the configurations that were used on each entry.

Fixed issues

- The custom filter of the timestamp columns does not have any character restrictions, allowing the user to input different formats.
- Installer message fixed upon uninstalling OmniTrail.
- The origin of the configuration files can now be found on its own column.

### **Differences between versions 1.2 and 1.1**

New features

- There is a new Special Filter 'Unapproved records' which shows electronic records that were created but not approved.
- Three new filters are installed to the default filter folder of OmniTrail. These are helpful to detect multiple aborts of consecutive records by the same user.
- Two different authorization configuration files can now be compared, and the differences viewed and printed. There are two comparison views available (Table and Tree view).
- The details of a single OmniTrail entry can now be printed by clicking a button in the details panel.
- A manual audit entry can now be created by a user, allowing the input of any text desired. This is then stored as an event in the System Audit Trail Database.
- New feature implemented to allow a user to review an event of conversion of the database and print a report. This happens when there was an update in the System Audit Trail Database.
- New column 'Electronic record type' shows the type of the electronic record associated with the entry.
- New column 'Electronic record location' shows the location of the electronic record when the event was registered in the System Audit.

Fixed issues:

- None

### **Differences between versions 1.1 and 1.0**

New features

- The event time stamp column now has a thicker right border to be better highlighted since it does not move when scrolling to the right.
- The time stamps can now be shown in UTC.
- When working in UTC mode, this will be displayed as UTC instead of +00:00.
- There is a new Special Filter 'Invalid hashes' which shows entries have been tempered with.
- A new live mode now allows the user to always see new entries as they come, refreshing OmniTrail automatically.
- New column 'User Identifier' shows the user identification with more detail (Active Directory, Local, etc.).
- Filters can now have a name associated.
- Configuration files can now also be downloaded from the details view.
- On the filter description, the 'Multiple Values Selected' when filtering multiple values is now clickable, showing in detail what values were filtered.
- There is a new Special Filter 'Missing Entries' which shows entries that are missing from the

database.

#### Fixed issues

- In the details panel, the word "Reason" would sometimes appear twice.
- The hamburger menu now has a tooltip.
- The entries report and reviewed entries report time stamp are now uniformized.
- A loss of connection from OmniTrail to the Regulated Environment Service no longer causes an application crash. A clear message is now shown to the user when this happens.
- Launching OmniTrail from the desktop icon now brings it to the front of all windows correctly.
- When having multiple windows, the tasks list now shows all opened windows correctly.

### **3 Remarks and known issues**

#### **Live updates**

- It is not advised to leave Live Updates turned on if not for observing the incoming events. The reason for this is that, in extremely rare situations, an application error may occur after 1 or 2 days without turning the Live Updates off.

# INSTALLATION NOTES

## 1. Introduction

OmniTrail is an OmniTrust software module to view and audit system audit trail events in an OmniTrust environment.

## 2. Important initial note

This software requires installation of License Manager 1.3.2 or higher and Regulated Environment 4.3.3 or higher.

## 3. System requirements

This software is tested to run correctly on:

- Microsoft Windows 10 1909 64-bit or above.
- Microsoft Windows 11 23H2 64-bit.

To run this software a PC configuration that meets the minimum hardware requirements for the Microsoft Windows operating system will be sufficient.

For Microsoft Windows 10 see

<http://support.microsoft.com/en-us/help/4028142/windows-10-system-requirements>.

For Microsoft Windows 11 see

<https://support.microsoft.com/en-us/windows/windows-11-system-requirements-86c11283-ea52-4782-9efd-7674389a7ba3>.

## 4. Upgrading OmniTrail

The installation procedure described below, will upgrade existing versions of OmniTrail to version 1.7a.

## 5. Installation procedure

It is important that you follow the installation sequence as given in the three sections below. If you do not follow this sequence the software may not work correctly.

### Before installation

- Make sure that your computer complies with the system requirements (see paragraph 4).
- Log on as user with "Administrator" rights in your local system to obtain appropriate authorization to complete the installation procedure successfully.

### Installation

- Install License Manager 1.3.2 or higher, bundled with the DVD
- Install Regulated Environment 4.3.3, bundled with the DVD
- Open the OmniTrail installer
- Read and accept the EULA
- Click install and wait for installation to finish
- No other user input is required to finish the installation

### After installation

- Remove the DVD and store it in a safe place.
- Start OmniTrail. You will be prompted to activate the license for the software. Please follow the instructions on your screen and use the license activation key as provided by Malvern Panalytical.  
If you are upgrading from a previous version of OmniTrail, license activation is not required.
- A System Audit Trail Conversion event will be created. This system event contains information

about the upgrade of the Database schema that occurred during the upgrade process. When opening OmniTrail for the first time after the upgrade, if you have permissions, you will have the opportunity to review this event that contains useful information about the conversion.

## 6. Installation remarks

Please read the following carefully because the remarks made might apply to your situation.

### Installation folders

The default folders are:

Installation folder	C:\Program Files (x86)\Malvern Panalytical\OmniTrail
Working folder	%AppData%\Malvern Panalytical\OmniTrail