

RELEASE NOTES

Malvern Panalytical Software Service (Regulated Environment Service)

Release: Version 1.7a, March 2025.

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

Thank you for using Malvern Panalytical Software service (Regulated Environment Service)

1 Introduction

The System Audit Trail may be stored on a central server in an OmniTrust environment.

2 Differences between versions

2.1 Differences between version 1.7a and 1.7

New features

- None

Fixed issues

- 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:

- None

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 OmniAccess is reported. In OmniAccess 1.5 version 1.4 was reported.
- 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 version 1.5 and 1.4

New features:

- Updated Help documentation of OmniAccess and OmniTrail.

Fixed issues:

- The Malvern Panalytical Software Service window on the central server shows the connected clients.

Differences between version 1.4 and 1.3

New features:

- Supports flexible signing feature.

Fixed issues:

- Synchronization of System Audit Trail events from clients on previous OmniTrust releases.

- Synchronization of System Audit Trail events after a loss and re-establishment of network connection.
- Server name not valid for synchronization configuration.
- User full name is validated on authentication.
- OmniTrust applications are detected upon installation on Centrally Managed systems.

Difference between version 1.3 and 1.2

New features:

- Supports central management (central storage of OmniTrust configuration).

Fixed issues:

- Last system audit trail event sometimes not synced from client to server.
- Unable to use server name in client synchronization configuration.
- System tray UI pops down when a new SAT entry is synchronized.

Differences between version 1.2 and 1.0

New features:

- When there is no OmniTrust license, the system tray pop-up message is improved.
- When a new client connects to a System Audit Trail Server running an older version of OmniTrust, the client will not connect and system tray pop-up message on the client informs the user of the situation.
- The name of the Remote Sync Agent UI executable has changed.

Fixed issues:

- The system tray UI does not present accurate information about the state of remote sync.
- Synchronization does not resume after interruption of network connection.
- System tray icon sometimes disappears.
- Synchronization between client and server can stop in certain situations.

3 Remarks and known issues

Synchronization issues with older versions of OmniTrust software

Previous releases of OmniTrust had issues synchronizing System Audit Trail events when the synchronization client and the synchronization server were from different OmniTrust releases.

It is essential to ensure the synchronization client and synchronization server are on the same OmniTrust release or upgrading the server to the 1.4 or above release solves the issue.

INSTALLATION NOTES

1 Introduction

The System Audit Trail may be stored on a central server in an OmniTrust environment. Installable parts available on the DVD

The following installers should be available on the DVD files

- License Manager 1.3.2
- Regulated Environment 4.3.3

2 Important initial note

This software requires installation of License Manager 1.3.2 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 System Audit Trail Server

The installation procedure described below will also upgrade existing versions of the software.

Upon upgrading System Audit Trail Server 1.7a, the System Audit Trail will automatically convert to the latest version. A conversion event is added to the System Audit Trail, which you should review and approve when you open OmniTrail for the first time after upgrade.

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

- a) Make sure that your computer complies with the system requirements (see Chapter 3).
- b) Ensure any running and/or pending update of Windows or any other software package is concluded before starting the installation.
- c) Log on as user with "Administrator" rights in your local system to obtain appropriate authorization to complete the installation procedure successfully.

Installation

- a) Copy the contents of the DVD to your local hard drive.
- b) Open the Regulated Environment installer from your hard drive using "Run as administrator".
- c) Read and accept the EULA.
- d) Click install and wait for installation to finish.

After installation

- a) Activate the provided license activation key for 'System Audit Trail Server' on the server on which it has been installed.
- b) Configure the System Audit Trail Server and clients as detailed in help documentation.

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\Malvern Panalytical\Regulated Environment
Working folder	%ProgramData%\Malvern Panalytical\Regulated Environment