

OmniTrust v1.8

Software Update Notification

Minor Release

Date: 16th March 2026

Introduction

OmniTrust consists of three installed applications:

- OmniAccess – system administration, user authentication and authorization configuration
- OmniTrail – audit trail viewer
- Regulated Environment - Malvern Panalytical software service

This SUN outlines the updates and changes applied to these components.

For detailed instructions on how to configure and use OmniTrust and for a list of compatible Malvern Panalytical applications, please refer to the OmniTrust User Guide.

Change summary

Change type	
Electronic record schema	No
OmniTrust configuration file schema	Yes
Audit trail behavior	No
Security control	No

Contents

Introduction.....	1
Change summary.....	1
New features.....	3
Fixed issues.....	3
Known issues.....	3
Installation instructions.....	4
System requirements.....	4
Before installation.....	4
Installation.....	4
After installation.....	5
Appendices.....	5

New features

New features

Active Directory Groups *

Active Directory Groups enables seamless alignment of organizational AD user groups with OmniAccess, eliminating the need to maintain duplicate user entries across multiple systems. Permissions can now be assigned or revoked at the group level, significantly reducing administrative overhead compared to managing users individually.

To enable this feature in OmniAccess, navigate to the 'Options' menu and select 'Active Directory Groups Setting'. 'Groups' will now become available in the 'Authorization Configuration' screen.

Entra ID *

Microsoft Entra ID (formerly Azure Active Directory) is designed to provide cloud-based identity and access management without relying on traditional domain-based Active Directory. It provides secure sign-in, single sign-on, multifactor authentication, and access control for OmniAccess.

To enable this feature in OmniAccess, navigate to the 'Authentication' menu; Entra ID is now available as a source.

Better user search

Domain user search now supports partial matching, making it easier and more flexible to find the right users without needing exact names.

* Note: At time of publishing the Aeris Instrument Suite does not support AD groups and Entra ID

Fixed issues

Fixes

Bug fix - Fixed an issue where roles that contained the same text caused issues adding users

Update - First time installer wizard for OmniAccess now has additional guidance for optimal customer environment setup

Security update - Upgraded installer to fix “.be TEMP folder vulnerable to DLL redirection attacks that allow the attacker to escalate privileges“

Known issues

Issue	Work Around / Details	Comment
Sentinel LDK sleep error - In a few instances when logged into OmniAccess or OmniTrail, if your workstation goes into sleep state an error message (in various forms) may appear on the screen	<ul style="list-style-type: none">• Turn off sleep/hibernate on the workstation• If that is not possible and you see the message, click OK on the message and resume activity.• Or click OK then restart the application (work may not be saved)	Issue in progress with Thales (3 rd party provider)

Installation instructions

System requirements

This software is tested to run correctly on:

- Microsoft Windows 11 24H2 64-bit.

Minimum hardware requirements for the Microsoft Windows operating system:

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

Before installation

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

Installation

The installer contains the License Manager, OmniAccess, OmniTrail and Regulated Environment installers. A setup wizard will take you through the process of installing each in the correct order.

- Navigate to the contents of "OmniTrust 1.8.iso" and double-click on "setup.exe"
- Follow the numbered sequence for installation (License Manager MUST be installed before any of the other packages)



Note: This latest version of OmniTrust will automatically update any existing versions.

After installation

- **OmniAccess** - Start OmniAccess. You will be prompted to activate the license for the software. Please follow the instructions on the screen and use the license activation key as provided by Malvern Panalytical. If you are upgrading from a previous version of OmniAccess, license activation is not required.
- **OmniTrail** - Start OmniTrail. You will be prompted to activate the license for the software. Please follow the instructions on the 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.
- **Regulated Environment (System audit trail server)**
 - Activate the provided license activation key for 'System Audit Trail Server' on the server on which it has been installed.
 - Configure the System Audit Trail Server and clients as detailed in help documentation.

Appendices

Appendix i: Advised OmniTrust file security settings

#	File type	Extension	Default path %ProgramData%\Malvern Panalytical	Creation	Scope	Minimum required permissions Applicable to Folder and File	Notes
1	*-Application information.config	.config	\Regulated Environment\Application Information	Software installation When an application is installed its corresponding app-info file is created.	System	ReadAndExecute, Write	Application information file with permissions descriptions and application version (this file is not editable in OmniAccess)
2	Authentication.config	.config	\Regulated Environment\Authentication	On Software Install	System	ReadAndExecute, Write	It contains the list of authentication sources that will be used by the Regulated Environment to validate usernames and passwords during login and signature.
3	MalvernPanalyticalUserStore.db3	.db3	\Regulated Environment\Authentication	On Software Install	User	ReadAndExecute, Write	The Malvern Panalytical User Store database file.

4	Authorization files	.config	\Regulated Environment\Authorization	On Software Install	System	ReadAndExecute, Write	Examples: <ul style="list-style-type: none"> • Mastersizer Xplorer- Authorization.config contains the deployed authorization settings (roles, members and permissions) for Mastersizer Xplorer, OmniAccess-Authorization.config for OmniAccess only. • OmniTrail- Authorization.config for OmniTrail only.
5	SigningAndReason.config	.config	\Regulated Environment\Signing	On Software Install	System	ReadAndExecute, Write	File contains the list of possible System Audit trail events and, for each one, settings about whether signing and reason are mandatory or optional.
6	Synchronization.config	.config	\Regulated Environment\System Audit Trail	On Software Install	System	ReadAndExecute, Write	File contains the settings for synchronization of data among Regulated Environment instances on different machines in the network.
7	System_audit_trail.db3	.db3	\Regulated Environment\System Audit Trail	On Software Install	System	ReadAndExecute, Write	The system audit trail database
8	Undeployed .config files	.config	\OmniAccess\Configurations	On Software Install	User	ReadAndExecute, Write (To the folder)	File contains authorization data such as roles, role members, permissions, etc. They are undeployed, but saved so that users can work on them.
9	Log files for OmniAccess	.log	\OmniAccess\Logs	On Software Install	User	ReadAndExecute, Write	Log files can be used for diagnostic and troubleshooting purposes.
10	Log files for OmniTrail	.log	\OmniTrail\Logs	On Software Install	User	ReadAndExecute, Write	Log files can be used for diagnostic and troubleshooting purposes.
11	OmniTrailSettings.json	.json	%USERPROFILE%\AppData\Roaming\Malvern Panalytical\OmniTrail	On Software Install	User	ReadAndExecute, Write	File contains user settings for the OmniTrail application. Preferences such as the number of entries per page visible.

12	OmniAccessSettings.json	.json	\OmniAccess\Configurations	On Software Install	User	ReadAndExecute, Write	File contains user settings for the OmniAccess application. Preferences such as Workstations mode enabled.
13	OmniAccessWorkstations.json	.json	\OmniAccess\Configurations	On Software Install	User	ReadAndExecute, Write	File contains a list of Workstations used in Workstation mode.
14	Filter files	.json	%Public%\Documents\Malvern Panalytical \OmniTrail	By user	User	ReadAndExecute, Write	This is the default folder where filter files are saved and loaded from.



Groewood Road, Malvern,
Worcestershire, WR14 1XZ,
United Kingdom

Tel: +44 1684 892456
Fax: +44 1684 892789

Lelyweg 1,
7602 EA Almelo,
The Netherlands

Tel: +31 546 534 444
Fax: +31 546 534 598

www.malvernpanalytical.com

Disclaimer: Although diligent care has been used to ensure that the information in this material is accurate, nothing herein can be construed to imply any representation or warranty as to the accuracy, correctness or completeness of this information and we shall not be liable for errors contained herein or for damages in connection with the use of this material. Malvern Panalytical reserves the right to change the content in this material at any time without notice. Copyright: © 2025 Malvern Panalytical. This publication or any portion thereof may not be copied or transmitted without our express written permission.