

RELEASE NOTES

OmniAccess

Release: Version 1.7a, June 2025.

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

Thank you for using OmniAccess.

1 Introduction

OmniAccess is one of the modules of the OmniTrust software solution that is used for administration of the system and configuring user authentication and authorization.

2 Differences between versions

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
- No longer asked to load unsaved changes if changed settings are identical to current ones
- 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 OmniTrust is reported. In OmniAccess 1.5 version 1.4 was reported.
- When using OmniTrust in Mastersizer 3000, OmniAccess now shows the Synchronization tab and the Signing and Reason tab.

Differences between version 1.5 and 1.4

New features:

- None

Fixed issues:

- Updated help documentation.

Differences between versions 1.4 and 1.3

New features:

- When using Central Configuration, upon the deployment of a configuration, a notification is shown informing the user that any running application should be restarted.
- Signing and Reason configuration now allows a set of rules to be associated to a set of roles.
- The OmniAccess Database has been removed.
- It is now possible to duplicate a role with its permissions.
- When enabling Central Configuration, the cache usage is automatically enabled with a default

3 day timeout.

Fixed issues:

- The Synchronization tab is only shown when there is a valid license.
- When using Central Configuration, the roles that are now shown are from the configuration that is centrally deployed.
- When using Central Configuration, disabling the Workstation option on one client and deploying it, will disable this option on other clients. No need to do it manually anymore.
- Some user interface improvements.

Differences between versions 1.3 and 1.2

New features:

- A cache expiry timeout was added to the synchronization configuration when configuring a client.
- The signing and reason configuration only shows the signing events that are applicable to the installed applications.
- After deploying a configuration, a dialog is shown advising to restart the application so that the changes made are applied.

Fixed issues:

- The ApplicationInformation configuration file now supports not having any authorizable functions.
- The OmniAccess wizard does not fail anymore when the logged in user does not have a full name.
- OmniAccess now supports multiple Application Information files for the same application but different versions. The Permissions page will show the applicable version(s) for each permission.
- Some user interface improvements.

Differences between versions 1.2 and 1.1

New features:

- Import and export of all configuration types: authentication, synchronization, signing and reason and authorizations.
- Support for the use of default roles.
- Several user interface improvements.

Fixed issues:

- None

Differences between versions 1.1 and 1.0

New features:

- Graphical user interface to edit authentication, synchronization, and signing & reason configurations.
- Wizard screen to setup initial permissions and options.
- Authorizations report.
- Manual audit entry.

Fixed issues:

- None

3 Remarks and known issues

External files

- When you move configuration files from one computer to another, some compatibility issues may arise. Please check the Help "Configuration > Software overview > Backup existing configurations" for further details.

Single Sign-on

- When Single Sign-on is enabled, Windows Domain or Windows Local will be used. If there is another authentication source enabled, it will only be used for Signing purposes.

Server name in server synchronization configuration

- Currently a server name is not supported in the server synchronization configuration. An IP address must always be provided.

INSTALLATION NOTES

1 Introduction

OmniAccess is a module of the OmniTrust software solution that enables access configuration 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-bits.

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 OmniAccess

The installation procedure, described below, will also upgrade existing versions of OmniAccess 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 or higher, bundled with the DVD.
- Open the OmniAccess 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 OmniAccess. 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 OmniAccess, license activation is not required.

- The first time OmniAccess starts, a wizard screen will help you setup the initial permissions and settings. Check the 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 (x86)\Malvern Panalytical\OmniAccess
Working folder	%AppData%\Malvern Panalytical\OmniAccess