

RELEASE NOTES

OmniAccess

Release: Version 1.6, August 2023.

© 2023 Malvern Panalytical B.V. All rights reserved.

Thank you for using Malvern Panalytical's 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 System requirements

The requirements with respect to operating systems, applications, computer hardware and other system aspects are described in the OmniAccess installation notes.

3 Differences between versions

3.1 Differences between version 1.6 and 1.5

This release includes bug fixes that were reported from version 1.5:

- 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.

3.2 Differences between version 1.5 and 1.4

Updated help documentation.

3.3 Differences between versions 1.4 and 1.3

OmniAccess 1.4 includes the following 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.

This release includes bug fixes that were reported from version 1.3, here are the most relevant:

- 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.

3.4 Differences between versions 1.3 and 1.2

OmniAccess 1.3 includes the following 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.

This release includes bug fixes that were reported from version 1.2, here are the most relevant:

- 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.

3.5 Differences between versions 1.2 and 1.1

OmniAccess 1.2 includes the following 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.

3.6 Differences between versions 1.1 and 1.0

OmniAccess 1.1 includes the following 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.

4 Remarks and known issues

4.1 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.

4.2 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.

4.3 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.