



RELEASE NOTES

OmniAccess

Release: Version 1.4, November 2021.

© 2021 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 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.

3.2 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

© 2021 Malvern Panalytical



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.

3.3 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.

3.4 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.