

# RELEASE NOTES

## OmniTrust

Release: Version 1.7a, June 2025.

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

Thank you for purchasing Malvern Panalytical's OmniTrust.

### 1 Introduction

OmniTrust is Malvern Panalytical's instrument-agnostic solution for regulatory compliance and data integrity.

Version 1.7a consists of the following software modules:

- OmniAccess 1.7a - 1.7.8
- OmniTrail 1.7a - 1.7.8
- Malvern Panalytical Software Service (Regulated Environment Service) 4.3.4

The OmniTrust software modules require installation of License Manager 1.3.2 or higher.

### 2 System requirements

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

### 3 Differences between versions

#### 3.1 Differences between versions 1.7a and 1.7

OmniTrail has been updated to version 1.7a.

OmniAccess has been updated to version 1.7a.

Malvern Panalytical Software Service (Regulated environment Service) has been updated to version 1.7a.

See the release notes of the updated versions for the changes.

## INSTALLATION NOTES

The installation notes can be found in the Release and Installation Notes of the software modules mentioned in this document.