

# MALVERN ACCESS CONFIGURATOR: v2.20 (PSW0303-11) SOFTWARE UPDATE NOTIFICATION

## Introduction

This document details the release of software PSW0303-11; version 2.20 of the Malvern Access Configurator (MAC) software.

The MAC is a utility that provides a flexible method of configuring access control for compatible applications from Malvern Panalytical, including the utility itself. The system is based on the assignment of roles to particular Windows users or groups, each of which can be configured to allow certain specific permissions within the compatible Malvern application.

Use of the MAC is required to enable the access control features of the control software for certain Malvern products, including the Mastersizer 3000, OMNISEC, PEAQ-DSC and Kinexus. A guide is available for these products describing how the MAC is used. Please reference to the product software CD-ROM.



### Note:

v2.20 is an English language only release. Any previous localization translations are not active in this release.

## Installation

It is assumed that you have authority to install or update software within your facility. It is also assumed that you have Administrator rights for the system upon which the software is installed, as this is a requirement of the installation process. If you do not have this authority please consult with your I.T. support department before proceeding.

### Minimum system requirements

If the MAC utility is used on a computer controlling a Malvern product, the recommended system requirements will necessarily be those advised for Malvern product itself.

## Installation instructions

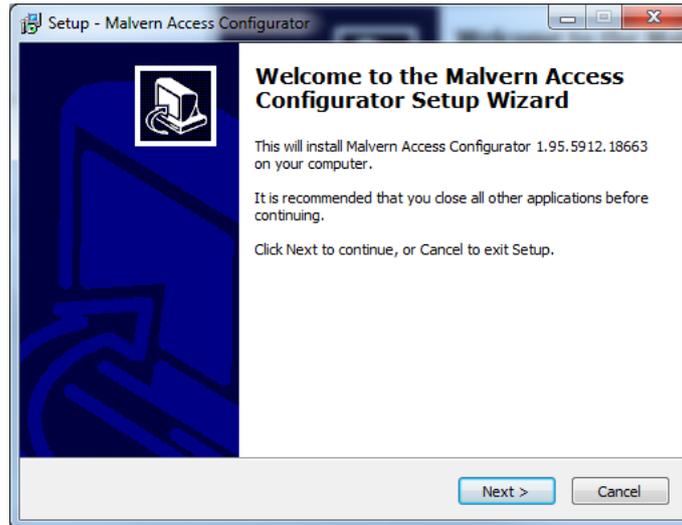
### Prerequisites

The Malvern Access Configurator software requires Windows 7 or Windows 10 to be installed. In addition, the .NET v4.6.1 or later framework must also to be installed.

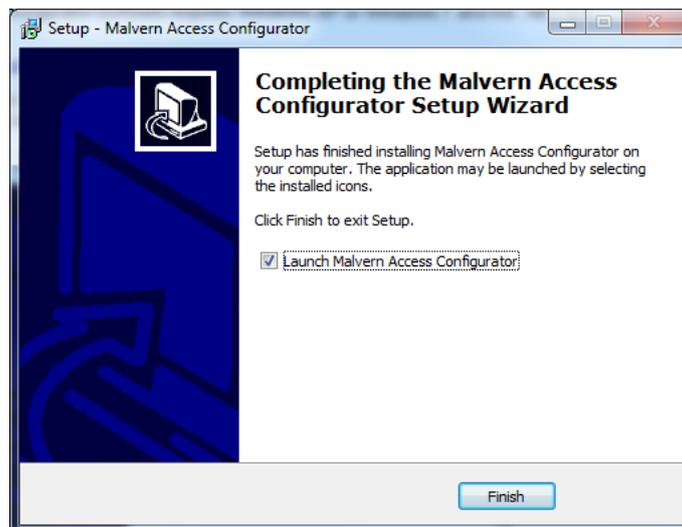
### Installing the software

The software comes as a single executable installer (*setup.exe*). The executable installer will be found on the CD supplied with the Malvern Panalytical product - in the directory called 'Malvern Application Configurator'.

Simply run the installer and follow the onscreen prompts to install the application.



At the end of installation, you will see this window:



If you don't wish to run MAC immediately, remove the tick from the checkbox. Press finish, and the installation is complete.

## Uninstall procedure

The software should be uninstalled using the standard **Add/Remove Programs** feature in the Windows 'Control panel'.

## Compatibility

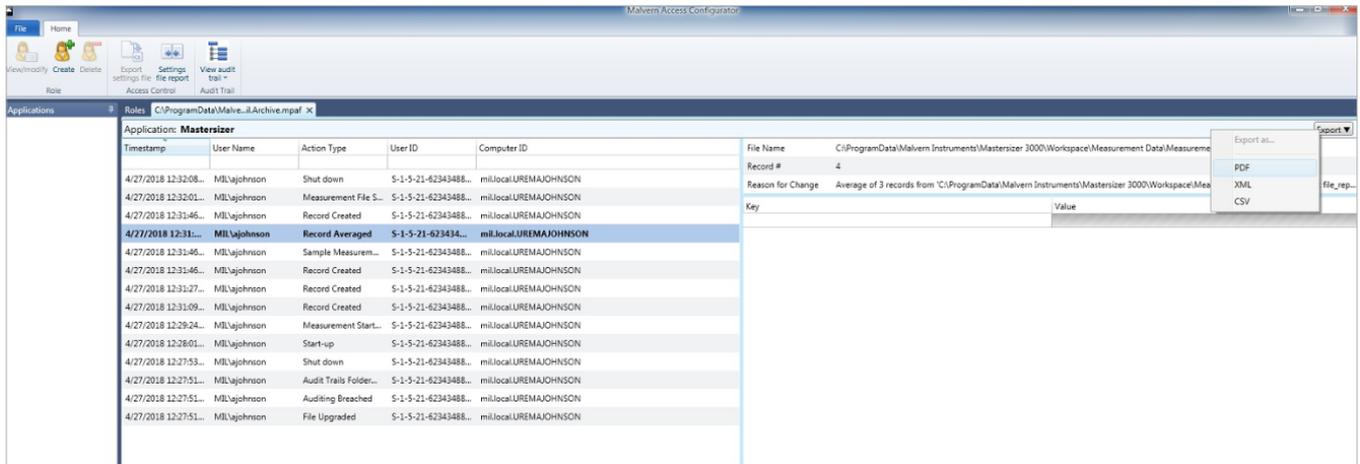
MAC v2.00 or higher is not compatible with permission files generated with versions of the application prior to v1.50.

## Main new feature description

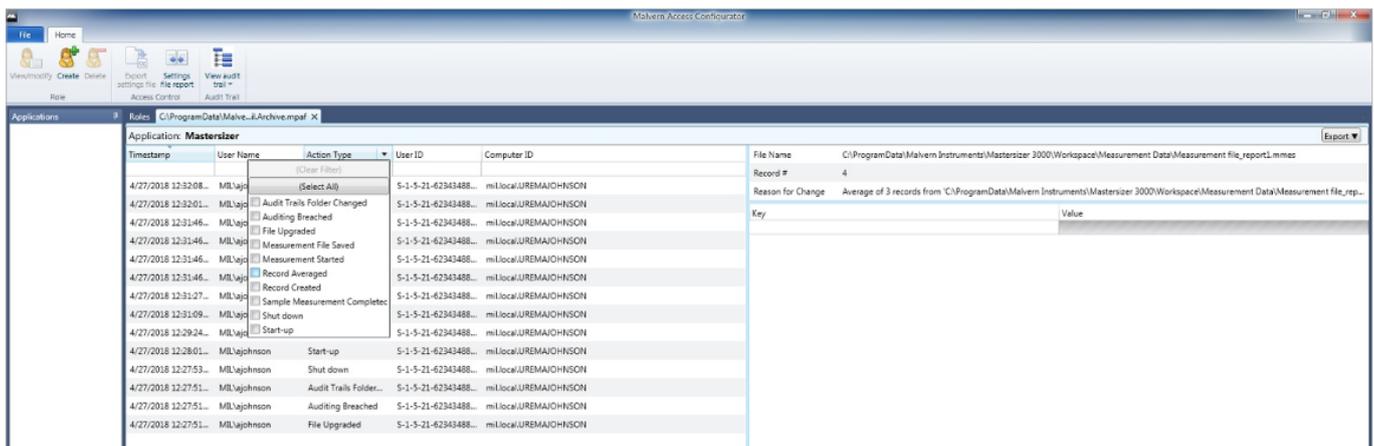
### Searching and exporting the audit history

The audit trail viewer within the MAC application has been updated to enable easier review of audit entries.

When an Audit Trail is opened, the detail for each entry can be seen on the righthand side of the screen, enabling easier review of audit information:



Once an Audit Trail has opened, it is also possible to filter the audit entries based on Date/Time, user name, action type and User ID:



To Export the entire audit trail, or the list of entries obtained with filters applied, click on the export button dropdown in the top-right corner of the Audit Trail viewer ( **Export ▼** ). The audit trail can be exported in PDF, XML or CSV formats. By default, the file will be saved into the ...*Malvern Instruments\Mastersizer 3000\Audit Exports\* folder.

Example of the exported files are shown below.

### PDF export



### Audit Trail Export - Mastersizer 3000

**Date range**  
29/03/2018 09:50:35 to 29/03/2018 12:27:34

**Filtered Audit Trail**

Timestamp	User Name	User ID	Computer ID	Action Type
29/03/2018 10:22:48	MIL\mutbeem	S-1-5-21-623434881-1009198268-2996851928-1573	mil.local.UKREMRNUTB EEM	Start-up
29/03/2018 10:22:37	MIL\MallInstall	S-1-5-21-623434881-1009198268-2996851928-21636	mil.local.UKREMRNUTB EEM	Shut down
29/03/2018 09:50:35	MIL\MallInstall	S-1-5-21-623434881-1009198268-2996851928-21636	mil.local.UKREMRNUTB EEM	Auditing Enabled

### XML export

### Audit Trail Export

**Audit Trail export details**

Time of export	2018-04-30 11:14:00
Exported by User	MIL\AHorton
Application Version	Mastersizer Software v3.70.1803.270 Alpha (non-validated)

**Date-Range**

From: 11/04/2018 10:24:57  
To: 30/04/2018 11:13:58

**Filtered Audit Trail**

The filtered Audit Trail. If no table is displayed then the filters applied returned no results.

Timestamp	User Name	User ID	Computer ID	Action Type
30/04/2018 11:13:49	MIL\AHorton	S-1-5-21-623434881-1009198268-2996851928-29753	mil.local.UKMAL1T3399	Start-up
30/04/2018 11:04:55	MIL\AHorton	S-1-5-21-623434881-1009198268-2996851928-29753	mil.local.UKMAL1T3399	Start-up
+ 30/04/2018 10:22:27	MIL\AHorton	S-1-5-21-623434881-1009198268-2996851928-29753	mil.local.UKMAL1T3399	Record Created
+ 30/04/2018 10:22:26	MIL\AHorton	S-1-5-21-623434881-1009198268-2996851928-29753	mil.local.UKMAL1T3399	Record Averaged
+ 30/04/2018 10:22:26	MIL\AHorton	S-1-5-21-623434881-1009198268-2996851928-29753	mil.local.UKMAL1T3399	Sample Measurement Completed
+ 30/04/2018 10:22:26	MIL\AHorton	S-1-5-21-623434881-1009198268-2996851928-29753	mil.local.UKMAL1T3399	Record Created
+ 30/04/2018 10:22:03	MIL\AHorton	S-1-5-21-623434881-1009198268-2996851928-29753	mil.local.UKMAL1T3399	Record Created
+ 30/04/2018 10:21:40	MIL\AHorton	S-1-5-21-623434881-1009198268-2996851928-29753	mil.local.UKMAL1T3399	Record Created
+ 30/04/2018 10:21:17	MIL\AHorton	S-1-5-21-623434881-1009198268-2996851928-29753	mil.local.UKMAL1T3399	Record Created

In the XML export, click on the  next to the action to see any expanded mode properties.

### CSV export

	A	B	C	D	E	F	G	H	I
1	Audit Trail	Timestamp	User Name	User ID	Computer ID	Action Type		Filters	Action Type
2		13/06/2018 11:11	MIL\AHorton	S-1-5-21-623434	mil.local.UKM	Login Successful			Auditing Breached
3		13/06/2018 10:36	MIL\AHorton	S-1-5-21-623434	mil.local.UKM	Login Successful			Insufficient Records Created
4		12/06/2018 15:02	MIL\AHorton	S-1-5-21-623434	mil.local.UKM	Login Successful			Measurement Started
5		12/06/2018 11:39	MIL\AHorton	S-1-5-21-623434	mil.local.UKM	Login Successful			Record Averaged
6		12/06/2018 11:38	MIL\AHorton	S-1-5-21-623434	mil.local.UKM	Login Successful			Login Successful

## New features list

This version of the MAC software includes the following new features.

Reference	Issue
84725	MAC must be able to open and view audit trails produced by Mastersizer 3000 v3.70 or higher
84731	Provide the ability to export and print audit trails
84733	Provide security permissions for audit trail export functions
84734	Improving the performance of caching LDAP / Active Directory queries
12828	Export destinations for security permission files should be remembered by application context.
84811	Version control should be provided for exported permissions files
73006	Provide means of comparing the security configurations for a given product
82270	Allow MAC to view OMNISEC per-item audit trails

## Fixed issues list

The main issues fixed in this release of the MAC software are confirmed in the table below.

Reference	Issue
80692	Administrator check fails on some customer networks
86927	'Valid From' fields can be set to a time in the past with user and permission configuration dialogues
83093	Translation errors cause an incorrect space to be inserted in xml file names (**. xml or **.xml)
76298	Multiple windows can be resized such that they are too small to see any information
76297	Role columns cannot be resized after a user opens the Role Detail window

## Known issues list

The issues found in this release but not rectified are listed below.

Reference	Issue
92511	Unable to install previous software versions without first manually uninstalling v2.20
91998	Valid from/to dates are not saved in the Role settings the first time a value is entered
91968	Valid to/from dates remain in the date window when a change is cancelled
91967	Timed permissions do not activate unless the software is restarted

As mentioned earlier in this document, this release is an English language only version. Some translations had been reported as not being correct. It is intended that an update to the translations will occur in a later release.

Also, some help files have not been fully updated to incorporate recent release features, these will also be updated in a future MAC release.



**Malvern  
Panalytical**

## MALVERN PANALYTICAL

Malvern Panalytical Ltd.  
Groewood Road, Malvern,  
Worcestershire, WR14 1XZ,  
United Kingdom

Malvern Panalytical B.V.  
Lelyweg 1, 7602 EA Almelo,  
Netherlands

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

Tel: +33 546 534 444  
Fax: +33 54 534 598

[info@malvernpanalytical.com](mailto:info@malvernpanalytical.com)  
[www.malvernpanalytical.com](http://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: © 2018 Malvern Panalytical. This publication or any portion thereof may not be copied or transmitted without our express written permission.

