

MALVERN ACCESS CONFIGURATOR: v2.21 (PSW0303-12) SOFTWARE UPDATE NOTIFICATION

Introduction

This document details the release of software PSW0303-12; version 2.21 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 Panalytical products, including the Mastersizer 3000, OMNISEC and PEAQ-DSC. A guide is available for these products describing how the MAC is used.



Note:

v2.21 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 Panalytical product, the recommended system requirements will be those advised for the Malvern Panalytical product itself.

Installation instructions

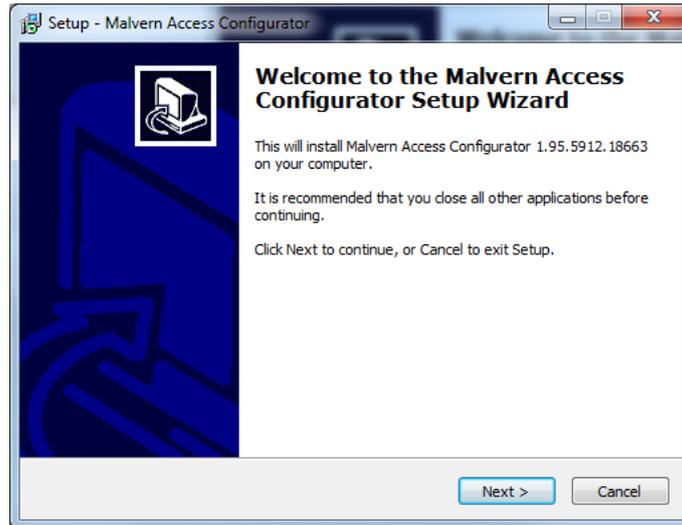
Prerequisites

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

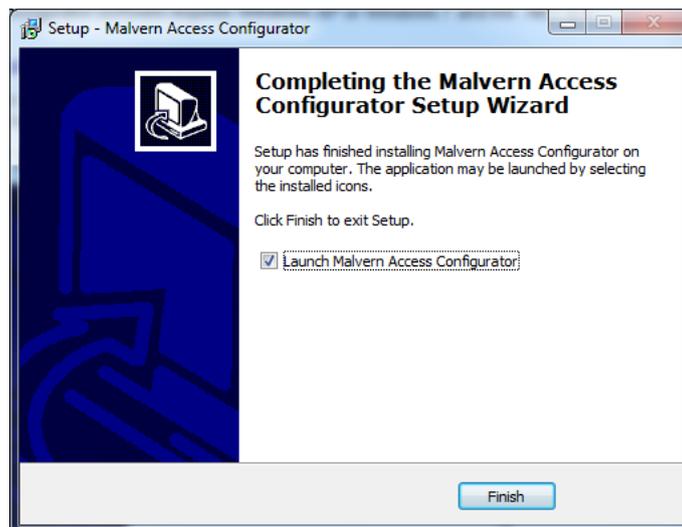
Installing the software

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

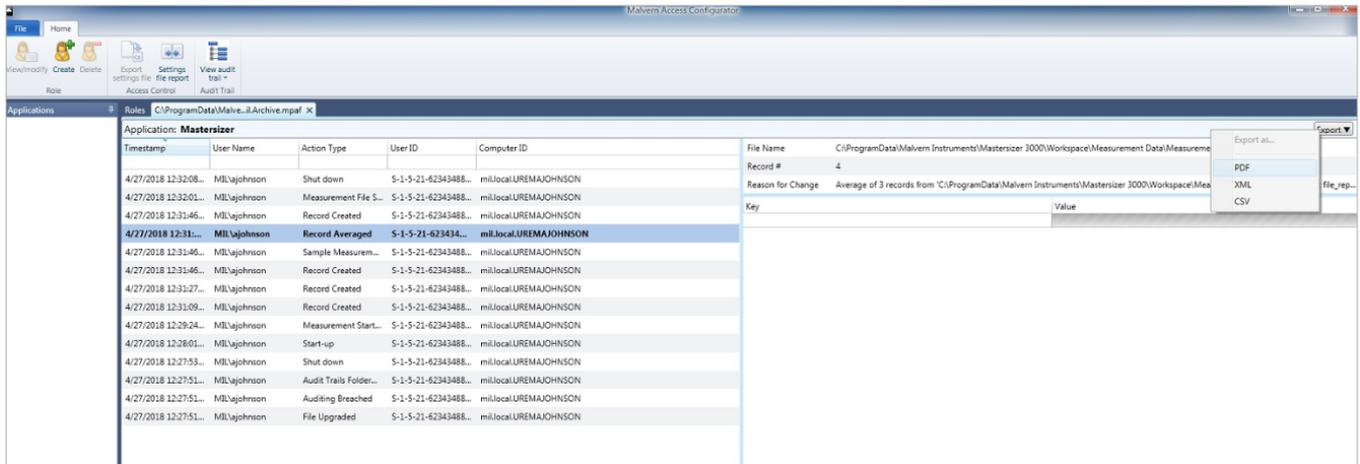
MAC v2.00 or higher is not compatible with permission files generated with versions of the application prior to v1.50.

Main new feature description

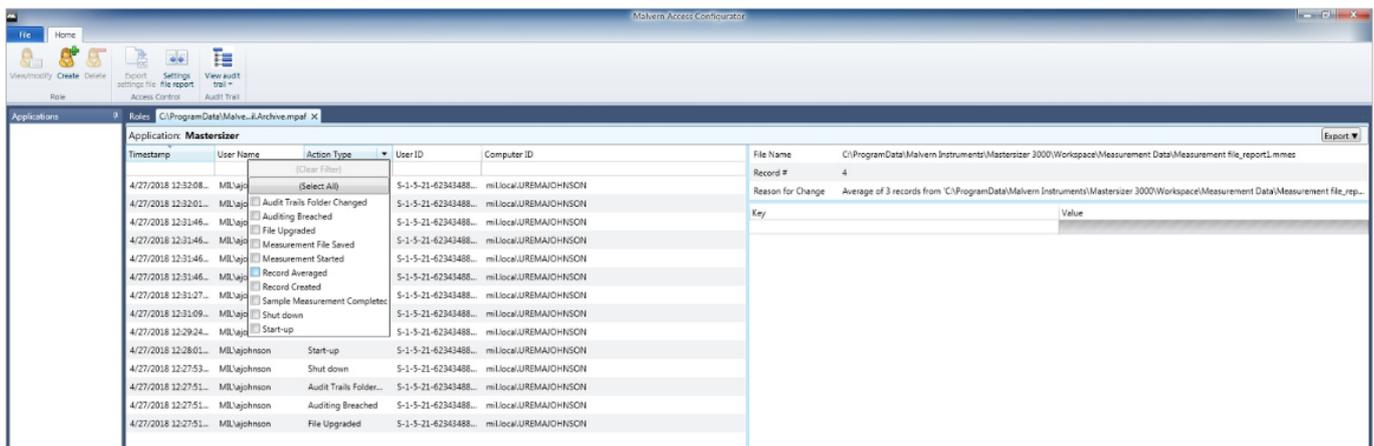
Searching and exporting the audit history

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 Timestamp, 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 no new features.

Fixed issues list

The no issues were fixed in this release of the MAC software.

Security Advisories

The following section details any security updates that have been addressed in this release, including fixes for identified vulnerabilities.



Note:

We always recommend updating to the latest software version which will provide you with new features, bugfixes and most importantly, security updates.



Note:

Other products may also be affected by any issue described here. We recommend you regularly check the Software Updates Notifications (SUNs) for all of your Malvern Panalytical products, and register on our website to receive updates.

Reference	Description	Recommendation
147839	Version 2.20 and earlier of the Malvern Access Configurator (MAC) contains a vulnerability in the audit trail viewing feature which could allow an attacker to craft a malicious audit trail file. If a user is tricked into loading a malicious audit trail file, this could result in arbitrary code execution. Version 2.21 introduces a fix to completely mitigate this vulnerability.	Upgrade to version 2.21 of the MAC. Never open files from an untrusted source, even if they appear to be non-executable.



**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
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.

