

MALVERN ACCESS CONFIGURATOR: v2.10 software

CERTIFICATE OF CONFORMANCE

with Validation Procedures

-  PARTICLE SIZE
-  MOLECULAR SIZE
-  MOLECULAR WEIGHT
-  MOLECULAR STRUCTURE
-  RHEOLOGICAL PROPERTIES
-  LABEL FREE BINDING
-  MICROCALORIMETRY

Malvern Instruments Ltd

Enigma Business Park
Grovewood Road
Malvern.
Worcestershire. UK. WR14 1XZ

Tel: +44 (0)1684 892456
Fax: +44 (0)1684 892789

Malvern Instruments Worldwide

Sales and service centres in
over 65 countries.

For details visit
www.malvern.com/contact

© Malvern Instruments Ltd 2017

more information at
www.malvern.com

We hereby certify that the above referenced software has been developed, tested and released in conformance with Malvern Instruments' ISO9001:2008 and TickIT Plus compliance procedures. These are maintained as part of Malvern's:

- Software Development Process
- ISO 9001:2008 Business Management System
- ISO 14001:2004 Environment Management System
- OHSAS 18001:2007 H&S Management System

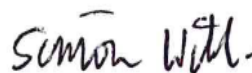
These procedures define Malvern's Design Qualification (DQ) process, providing a fundamental basis for an end user's validation of the software. This includes the production of documentation tracing the conformance of the software to design requirements in accordance with GAMP 5 guidelines. This documentation is open to inspection as part of a paid audit.

The Malvern Access Controller is designed to be installed with the application software provided with Malvern's Mastersizer 3000, Kinexus, OMNISEC and PEAQ-DSC instruments. The following items are available to help users to complete the validation of these instruments:

- Installation Qualification and Operational Qualification procedures, covering both hardware and software qualification in line with USP<1058> requirements;
- A 21 CFR Part 11 Feature Key providing users with access to software features designed to aid technical compliance to the requirements of 21 CFR Part 11;
- A 21 CFR Part 11 Guide and Gap Analysis offering guidance regarding how Malvern software can be configured within a 21CFR Part 11 compliant environment;
- A Malvern Access Configurator manual, providing advice on configuring and deploying a security system;
- A 21 CFR Part 11 Test Schedules providing a series of tests which may be used to validate a software's 21CFR Part 11 features.



Steve Anderson
Chief Engineer



Simon Wihl
Software Manager



Paul Kippax
Product Manager

Date: July 27, 2017

Please Note: This certificate relates to Malvern Instrument's business management systems only. It does not cover any validation procedures implemented by the end user of the software.