

Mastersizer

3000

Software Suite PSS0223-02

Certificate of Conformance with Validation Procedures

We hereby certify that the above-referenced software has been developed, tested and released in conformance with Malvern Instruments ISO9001: 2008 and TickIT compliance procedures.

These procedures, which provide a fundamental basis for the end user's Validation of the software, include the production of the following documentation tracing the conformance of the software to design requirements in accordance with GAMP guidelines.

Malvern Instruments Codes of Practice and Protocols:
Software Development Codes Of Practice
ISO 9001:2008 System
Business Management System

This documentation is open to inspection as part of a paid audit.




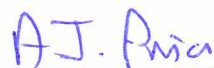

The following items are available to help Quality Assurance personnel to complete the validation of Mastersizer systems:

- **Installation Qualification/ Operational Qualification (IQ/OQ) procedures**
- **21 CFR Part 11 Feature Key** providing users with a means to achieve technical compliance to the requirements of 21 CFR Part 11
- **21 CFR Part 11 Guide** offering guidance as to the features of the software which can support operation within a 21CFR11 compliant environment.

Please Note:

The Scope of the assurance offered by this Certificate is limited to Malvern Documentary Procedures only.

The issue of this certificate does not relieve users of their ultimate responsibilities for validation of 3rd Party Software used by them.

 Malvern  001  ISO 9001:2008 The TickIT Guide Issue 5.5	Date: 16 March 2012
	Andy Prior, Project Manager 
	Simon Wihl, Design Authority 
	Steve Anderson, Technical Authority 