

## Malvern Version 5.22 Software for the Mastersizer 2000

### 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 ISO9000: 2001 and TickIT Compliance Procedures.

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

<p><b>Lifecycle Documentation:</b></p> <p>Requirements Specification Functional Specification Design Specification Design Test Schedules and Reports Alpha test documentation Beta test documentation</p>	<p><b>Malvern Instruments Codes of Practice and Protocols:</b></p> <p>Software Development Codes Of Practice ISO 9000:2001 System Design Change Management System Management Procedures Manual</p>
---	--




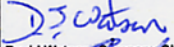



Rights to view this documentation are available as an option in the Malvern Instruments QSpec Validation Elements programme to assist users and regulatory auditors.

The QSpec Validation Elements programme includes the following **separately** available items to help Quality Assurance personnel to complete the full validation of their Mastersizer 2000 systems:

- Installation Qualification/ Operational Qualification (IQ/OQ )
- IQ/OQ Preview and Validation Planning Folder
- Software Update Verification
- Escrow Licenses
- QSpec Elements Rights-to-View Certification and Registration  
*Includes the provision of a Logbook which acts as a validation Workbook for individual Mastersizer installations*
- User Requirements Specification Matrix  
*This is a detailed specification of the Mastersizer 2000 describing the system and its capabilities and detailing pass/fail criteria which can form the basis of a User Performance Qualification.*
- 21 CFR Part 11 Feature Keys  
*These provide users with the technical means to achieve 21CFR Part 11 compliance.*

**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 3<sup>RD</sup> PARTY SOFTWARE USED BY THEM.

   ISO 9001:2000 The TickIT Guide Issue 5	Date: 17 <sup>th</sup> September 2004  David Watson, Company Chief Engineer
	 Steve Anderson, Engineering Manager
	 Andy Prier, Software Manager
	 Rod Jones, QSpec Validation Elements Program Manager