

Malvern Version 5.13 Particle Sizer 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 Ticklt 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.

Lifecycle Documentation:

Requirements Specification
Functional Specification
Design Specification
Design Test Schedules and Reports
Alpha test documentation
Beta test documentation

Malvern Instruments Codes of Practice and Protocols:

Software Development Codes Of Practice ISO 9000:2001 System Design Change Management System Management Procedures Manual

Rights to view this documentation under the terms of a non-disclosure agreement are available as an option in the Malvern Instruments QSpec Validation Elements programme.

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^{RO}
PARTY SOFTWARE USED BY
THEM.

INSTRUMENTS WITH THIS
VERSION OF SOFTWARE
INSTALLED WILL PRODUCE
PRINTED REPORTS WITH THE
CORRESPONDING SOFTWARE
VERSION AND THE
INSTRUMENT SERIAL
NUMBER EMBEDDED IN THE
FOOTER.

