

PEAQ-ITC Analysis SOFTWARE: v1.30

CERTIFICATE OF CONFORMANCE




We hereby certify that the above referenced software has been developed, tested and released in conformance with Malvern Panalytical's ISO9001:2015 compliance procedure. These are maintained as part of Malvern Panalytical's:

- Software Development Process
- ISO 9001:2015 Business Management System

These procedures provide a fundamental basis for an end user's validation of the software, and include the production of documentation tracing the conformance of the software to design requirements in accordance with GAMP guidelines. 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 their PEAQ-ITC systems:

- Installation Qualification and Operational Qualification procedures.

Approvals		
Role	Name	Signature
Acceptance	Ronan O'Brien	
Approval	Steve Anderson	
Release	James Wraight	
Release Date:	06/04/2020	

Please Note:

The Scope of the assurance offered by this Certificate is limited to Malvern Panalytical 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.