

Certificate of Conformance – Software Products

This Certificate of Conformance certifies that all software products developed and released by Malvern Panalytical conform to the specifications and quality standards defined by the company and applicable regulatory requirements.

Malvern Panalytical follows a ISO 9001:2015 certified quality management system and ensures that each software product has been:

- Developed in accordance with documented design and development procedures.
- Verified and validated through structured testing protocols.
- Released under controlled conditions by authorized personnel.
- Version controlled and is traceable.

This certificate confirms that the software product delivered to the customer meets the intended functional and quality requirements at the time of release.

Signed,



Kristel Zonder
Global Quality Director