

# Compliance Declaration: Malvern Panalytical Customer Validation Documents (CVDs)

Enigma Business Park  
Grovewood Road  
Malvern  
WR14 1XZ  
United Kingdom

Malvern Panalytical  
Lelyweg 1 (7602 EA)  
P.O. Box 13  
7600 AA Almelo  
Netherlands

[info@malvernpanalytical.com](mailto:info@malvernpanalytical.com)  
[www.malvernpanalytical.com](http://www.malvernpanalytical.com)

## Customer Validation Documents

Customer Validation Documents (CVD) are documents used by customers to validate that Malvern Panalytical instruments, accessories and software are installed and operating correctly – they include: Installation Qualification (IQ), Operational Qualification (OQ) and Regulated Environment Compliance Test document (CTD).

We hereby certify that Malvern Panalytical instrument CVDs have been developed, tested and released to be compliant with the following procedures:

- ISO9001:2015 certification

These procedures ensure that CVDs are version controlled, tested and approved using a recorded and auditable method by qualified company representatives applicable to the product.

Digitally signed approval confirmation material for specific CVDs, and further details on these processes, is open to inspection for audit by a customer on request.

Signed:



Kristel Zonder, Global Quality Director

21 August 2025