

CERTIFICATE

Number: 75160

The management system of the organization(s) and locations mentioned on the addendum belonging to:

Malvern Panalytical B.V.

Lelyweg 1
7602 EA Almelo
Netherlands

including the implementation meets the requirements of the standard:

ISO 9001:2015

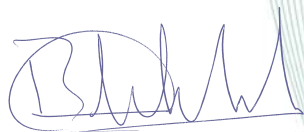
Scope:

The design, development, manufacture, sell, supply and support of professional material analysis and material characterization systems with associated products, software and services.

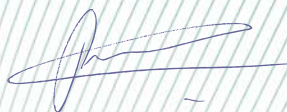
Certificate expiry date: 1 December 2027
Certificate effective date: 1 December 2024
Certified since*: 1 August 1991

This certificate is valid for the organization(s) and/or locations mentioned on the addendum.

DEKRA Certification B.V.



B.T.M. Holtus
Managing Director



R.C. Verhagen
Certification Manager

© Integral publication of this certificate and adjoining reports is allowed.

* Against this certifiable standard / possibly by another certification body.



ADDENDUM

To certificate: 75160

The management system of the organization(s) and/or location(s) of:

Malvern Panalytical B.V.

Lelyweg 1
7602 EA Almelo
Netherlands

Certified organization(s) and/or locations:

Malvern Panalytical B.V.
De Schakel 18
5651 GH Eindhoven
Netherlands

Malvern Panalytical B.V.
Lelyweg 1
7602 EA Almelo
Netherlands

Addendum expiry date: 1 December 2027
Addendum effective date: 1 December 2024

