

# CERTIFICATE

Number: 2188611

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

## Malvern Panalytical B.V.

Lelyweg 1  
7602 EA Almelo  
The Netherlands

including the implementation meets the requirements of the standard:

# ISO 14001:2015

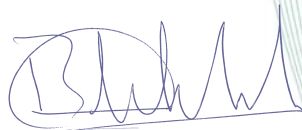
### Scope:

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 2024  
Certificate effective date: 1 December 2021  
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: 2188611

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

## Malvern Panalytical B.V.

Lelyweg 1  
7602 EA Almelo

Certified organization(s) and/or locations:

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

Malvern Panalytical B.V.  
Lelyweg 1  
7602 EA Almelo

Addendum expiry date: 1 December 2024

Addendum effective date: 1 December 2021