

Certificate of Registration

OCCUPATIONAL HEALTH & SAFETY MANAGEMENT SYSTEM - ISO 45001:2018

This is to certify that:

Malvern Panalytical Ltd
Groewood Road
Enigma Business Park
Malvern
WR14 1XZ
United Kingdom

Holds Certificate Number:

OHS 684719

and operates an Occupational Health and Safety Management System which complies with the requirements of ISO 45001:2018 for the following scope:

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

For and on behalf of BSI:

Andrew Launn, EMEA Systems Certification Director

Original Registration Date: 2010-07-05

Effective Date: 2019-07-05

Latest Revision Date: 2020-04-20

Expiry Date: 2022-07-04

Page: 1 of 1



003

...making excellence a habit.™