

# 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.**

**[Previously certified to BS OHSAS 18001:2007 since 2010-07-05]**

For and on behalf of BSI:

\_\_\_\_\_  
Matt Page, Managing Director Assurance - UK & Ireland

Original Registration Date: 2019-06-07

Latest Revision Date: 2022-06-30

Effective Date: 2022-07-05

Expiry Date: 2025-07-04

Page: 1 of 1



...making excellence a habit.™