

# Certificate of Registration

## ENVIRONMENTAL MANAGEMENT SYSTEM - ISO 14001:2015

This is to certify that:

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

Holds Certificate Number:

EMS 684720

and operates an Environmental Management System which complies with the requirements of ISO 14001:2015 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:



Matt Page, Senior Vice President, EMEA Assurance

Original Registration Date: 2007-06-18

Latest Revision Date: 2025-03-27

Effective Date: 2025-06-18

Expiry Date: 2028-06-17



Page: 1 of 1

...making excellence a habit.™