

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, Managing Director Assurance - UK & Ireland

Original Registration Date: 2007-06-18

Latest Revision Date: 2022-06-10

Effective Date: 2022-06-18

Expiry Date: 2025-06-17

Page: 1 of 1



...making excellence a habit.™