

# Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

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

Holds Certificate Number:

FS 684721

and operates a Quality Management System which complies with the requirements of ISO 9001:2015 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:

Andrew Launn, EMEA Systems Certification Director

Original Registration Date: 1996-08-07

Latest Revision Date: 2021-07-07

Effective Date: 2020-09-01

Expiry Date: 2023-08-31

Page: 1 of 2



...making excellence a habit.™

Certificate No: FS 684721

Location

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

Malvern Panalytical Inc  
22 Industrial Drive East  
Northampton  
Massachusetts  
01060  
USA

Registered Activities

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

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



Original Registration Date: 1996-08-07

Latest Revision Date: 2021-07-07

Effective Date: 2020-09-01

Expiry Date: 2023-08-31

Page: 2 of 2

This certificate was issued electronically and remains the property of BSI and is bound by the conditions of contract.  
An electronic certificate can be authenticated [online](#).  
Printed copies can be validated at [www.bsigroup.com/ClientDirectory](http://www.bsigroup.com/ClientDirectory)

Information and Contact: BSI, Kitemark Court, Davy Avenue, Knowlhill, Milton Keynes MK5 8PP. Tel: + 44 345 080 9000  
BSI Assurance UK Limited, registered in England under number 7805321 at 389 Chiswick High Road, London W4 4AL, UK.  
A Member of the BSI Group of Companies.