

# 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: 2021-11-03

Effective Date: 2020-09-01

Expiry Date: 2023-08-31

Page: 1 of 3



...making excellence a habit.™

Certificate No: FS 684721

Location	Registered Activities
Malvern Panalytical Ltd Groewood Road Enigma Business Park Malvern WR14 1XZ United Kingdom	The design, development, manufacture, sale and support of material analysis and characterization systems and software with associated products and services.
Malvern Panalytical Inc 22 Industrial Drive East Northampton Massachusetts 01060 USA	The design, development, manufacture, sale and support of material analysis and characterization systems and software with associated products and services.
Malvern Panalytical, Inc. Canada 4921 Place Olivia St-Laurent Québec H4R 2VR Canada	Sales and service office.
Malvern Panalytical, Inc. Mexico Malvern Panalytical, a division of Spectris Mexico S de RL de CV Av. Pedro Ramirez Vazquez 200-13 Nivel 1, Parque Corporativo Valle Oriente, Col. Valle Oriente San Pedro Garza Garcia Nuevo León 66269 Mexico	Administrative work (Finance, Logistics, Marketing, Inside Sales, HR, Service Backoffice), Applications/ Laboratory, Repair.
Malvern Panalytical, Inc. 2400 Computer Drive Westborough Massachusetts 01581 USA	Customer Service & Support, Apps Lab, Repair Depot, Customer Training and Warehouse.

Original Registration Date: 1996-08-07

Latest Revision Date: 2021-11-03

Effective Date: 2020-09-01

Expiry Date: 2023-08-31

Page: 2 of 3

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)

Certificate No: FS 684721

Location

Malvern Panalytical, Inc. Sugar land  
12300 Dairy Ashford Rd.  
Suite 400  
Sugar Land  
Texas  
77478  
USA

Registered Activities

Proof-of-principle analysis of customer samples in support of sales



Original Registration Date: 1996-08-07

Latest Revision Date: 2021-11-03

Effective Date: 2020-09-01

Expiry Date: 2023-08-31

Page: 3 of 3

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.