

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:



Matt Page, Senior Vice President, EMEA Assurance

Original Registration Date: 1996-08-07

Latest Revision Date: 2026-07-30

Effective Date: 2026-09-01

Expiry Date: 2029-08-31



Page: 1 of 2

...making excellence a habit.™

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

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 Seventh and Eighth Floors, The Acre, 90 Long Acre, London, WC2E 9RA, UK.

A Member of the BSI Group of Companies.

Certificate No: FS 684721

Location	Registered Activities
Malvern Panalytical Ltd Groveswood 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 and support of material analysis and characterisation systems and software, with associated products and services.
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	The sale and support of material analysis and characterisation systems and software, with associated products and services.
Malvern Panalytical, Inc. 2400 Computer Drive Westborough Massachusetts 01581 USA	The sale and support of material analysis and characterisation systems and software, with associated products and services.
Malvern Panalytical, Inc. Sugar land 12300 Dairy Ashford Rd. Suite 400 Sugar Land Texas 77478 USA	The sale and support of material analysis and characterisation systems and software, with associated products and services.

Original Registration Date: 1996-08-07

Latest Revision Date: 2026-07-30

Effective Date: 2026-09-01

Expiry Date: 2029-08-31

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

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 Seventh and Eighth Floors, The Acre, 90 Long Acre, London, WC2E 9RA, UK.

A Member of the BSI Group of Companies.