

# Certificate of Registration

INFORMATION SECURITY MANAGEMENT SYSTEM - ISO/IEC 27001:2013

This is to certify that:

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

Holds Certificate Number:

IS 751368

and operates an Information Security Management System which complies with the requirements of ISO/IEC 27001:2013 for the following scope:

**The development, supply and support from the Company's digital solutions department known internally as Frontier Analytics and the associated controls serviced by its internal and external partners. It also relates where appropriate to external risk sources including functions which are outsourced in accordance with the Statement of Applicability v2.0 dated 21/10/2021.**

For and on behalf of BSI:

\_\_\_\_\_  
Matt Page, Managing Director Assurance - UK & Ireland

Original Registration Date: 2022-03-11

Latest Revision Date: 2022-03-11

Effective Date: 2022-03-11

Expiry Date: 2025-03-10

Page: 1 of 2



...making excellence a habit.™

Certificate No: IS 751368

Location	Registered Activities
Malvern Panalytical Ltd Groewood Road Enigma Business Park Malvern WR14 1XZ United Kingdom	The development, supply and support from the Company's digital solutions department known internally as Frontier Analytics and the associated controls serviced by its internal and external partners. It also relates where appropriate to external risk sources including functions which are outsourced in accordance with the Statement of Applicability v2.0 dated 21/10/2021.
Malvern Panalytical Ltd DeskLodge House Redcliffe Way Bristol BS1 6NL United Kingdom	The development, supply and support from the Company's digital solutions department known internally as Frontier Analytics and the associated controls serviced by its internal and external partners. It also relates where appropriate to external risk sources including functions which are outsourced in accordance with the Statement of Applicability v2.0 dated 21/10/2021.

Original Registration Date: 2022-03-11

Latest Revision Date: 2022-03-11

Effective Date: 2022-03-11

Expiry Date: 2025-03-10

Page: 2 of 2

This certificate relates to the information security management system, and not to the products or services of the certified organisation. The certificate reference number, the mark of the certification body and/or the accreditation mark may not be shown on products or stated in documents regarding products or services. Promotion material, advertisements or other documents showing or referring to this certificate, the trademark of the certification body, or the accreditation mark, must comply with the intention of the certificate. The certificate does not of itself confer immunity on the certified organisation from legal obligations.

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.