

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016

This is to certify that:

Meso Scale Diagnostics, LLC
1601 Research Blvd.
Rockville
Maryland
20850
USA

Holds Certificate Number:

FM 579698

and operates a Quality Management System which complies with the requirements of ISO 13485:2016 for the following scope:

Design and development of in-vitro diagnostic products including instruments, plates, reagents and kits.

For and on behalf of BSI:

Graeme Tunbridge, Senior Vice President Medical Devices

Original Registration Date: 2012-05-03

Latest Revision Date: 2024-02-21

Effective Date: 2024-05-03

Expiry Date: 2027-05-02

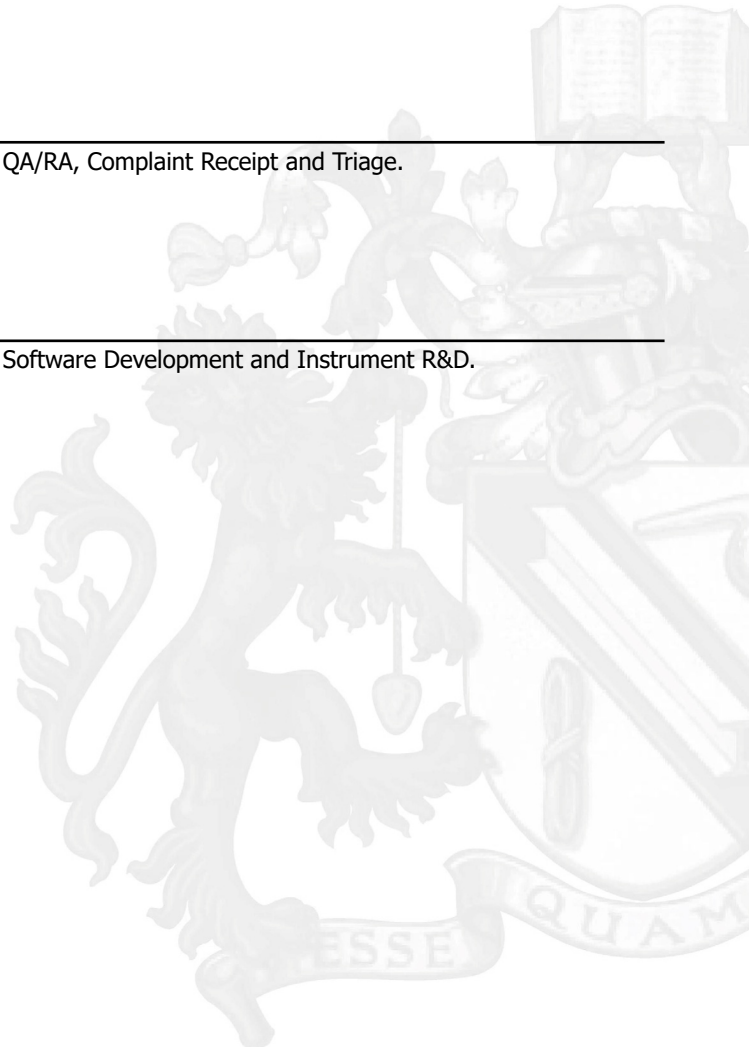
Page: 1 of 2



...making excellence a habit.™

Certificate No: FM 579698

Location	Registered Activities
Meso Scale Diagnostics, LLC 1601 Research Blvd. Rockville Maryland 20850 USA	Head Quarter Activities including- Finance, IT, Compliant Receipt and Triage, Legal, Human Resources, Project Management Office, Technical Support and Infrastructure Support.
Meso Scale Diagnostics, LLC 16020 Industrial Drive Gaithersburg Maryland 20877 USA	Consumables/ Assay R&D, Document and Record Controls.
Meso Scale Diagnostcs, LLC 1701 Research Blvd. Rockville Maryland 20850 USA	QA/RA, Complaint Receipt and Triage.
Meso Scale Diagnostics LLC 9120 Gaither Road Gaithersburg Maryland 20878 USA	Software Development and Instrument R&D.



Original Registration Date: 2012-05-03

Effective Date: 2024-05-03

Latest Revision Date: 2024-02-21

Expiry Date: 2027-05-02

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

Issuing Body: BSI Group The Netherlands B.V., John M. Keynesplein 9, 1066 EP Amsterdam, The Netherlands

BSI Group The Netherlands B.V. is registered in The Netherlands under number 33264284 | A Member of the BSI Group Holdings B.V.

Contact Office: 12950 Worldgate Drive, Suite 800, Herndon, VA 20170-6007 USA.