

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

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

Holds Certificate No:

FM 579699

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

The design, development, manufacture, distribution and service of instruments, plates, reagents, assays and kits for research purposes.

For and on behalf of BSI:

Carlos Pitanga, Managing Director Assurance, Americas

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 3



...making excellence a habit.™

Certificate No: **FM 579699**

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 9161 Industrial Ct Gaithersburg Maryland 20877 USA	Plate manufacturing.
Meso Scale Diagnostics, LLC 9105-9125 Arbuckle Drive Gaithersburg Maryland 20877 USA	Warehouse activities and Machine Shop.
Meso Scale Diagnostics, LLC 16020 Industrial Drive Gaithersburg Maryland 20877 USA	Shipping/ Receiving, Consumables/ Assay R&D, Document and Record Controls, Assay Service Group (assay design and development and testing service for research purposes only).
Meso Scale Diagnostcs, LLC 1701 Research Blvd. Rockville Maryland 20850 USA	Consumable Quality Control, Shipping Receiving, Kit Assembly, Reagent Manufacturing (Bulk, Diluent and Plate Manufacturing), Purchasing/Supplier Management, QA/RA, Order Entry, Warehouse, Complaint Receipt and Triage.
Meso Scale Diagnostics LLC 9120 Gaither Road Gaithersburg Maryland 20878 USA	Instrument Service - refurbishment of instruments/parts, Quality Control for Instruments, Instrument manufacturing, Warehouse, Software Development, and Instrument R&D and Machine Shop.

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 3

Certificate No: **FM 579699**

Location

Registered Activities

Meso Scale Diagnostics, LLC
9125 Industrial Court
Gaithersburg
Maryland
20877
USA

Manufacturing and R&D of plate components.

Meso Scale Diagnostics, LLC
9346 Gaither Road
Gaithersburg
Maryland
20877
USA

Automation.



Original Registration Date: 2012-05-03

Effective Date: 2024-05-03

Latest Revision Date: 2024-02-21

Expiry Date: 2027-05-02