

# Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

Molecular Devices  
3860 N. First Street  
San Jose  
California  
95134  
USA

Holds Certificate Number:

FS 534246

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

**Design, development, production, installation, distribution and servicing of equipment, information software and reagents for R&D laboratory use in cell biology, biochemistry, and physiology.**

For and on behalf of BSI:

\_\_\_\_\_  
Matt Page, Senior Vice President, EMEA Assurance

Original Registration Date: 2009-06-12

Effective Date: 2024-06-12

Latest Revision Date: 2026-04-27

Expiry Date: 2027-06-11

Page: 1 of 2



...making excellence a habit.™

Certificate No: FS 534246

Location	Registered Activities
Molecular Devices 3860 N. First Street San Jose California 95134 USA	Design, development, production, installation, and servicing of equipment, information software and reagents for R&D laboratory use in cell biology, biochemistry, and physiology.
Molecular Devices 402 Boot Road Downingtown Pennsylvania 19335 USA	Design, development of information software for R&D laboratory use in cell biology, biochemistry, and physiology.
Molecular Devices 1180 Eskdale Rd Winnersh Triangle Wokingham Berkshire RG41 5TU United Kingdom	Service and distribution of instrumentation, information software and reagents for R&D laboratory use in cell biology, biochemistry, and physiology.
Molecular Devices 555 Yuan Xi Road Nanhui Industrial Zone Pudong New Area Shanghai 201300 China	Design, development and production of equipment for R&D laboratory use in cell biology, biochemistry, and physiology.
Molecular Devices (Austria) GmbH Urstein Süd 17 Puch bei Hallein Salzburg 5412 Austria	Design, development, production and servicing of equipment, information software, and reagents for R&D laboratory use in cell biology, biochemistry, and physiology.

Original Registration Date: 2009-06-12

Latest Revision Date: 2026-04-27

Effective Date: 2024-06-12

Expiry Date: 2027-06-11

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

A Member of the BSI Group of Companies.