

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 No:

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, 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:



Carlos Pitanga, Chief Operating Officer Assurance – Americas

Original Registration Date: 2009-06-12

Latest Revision Date: 2018-07-12

Effective Date: 2018-06-12

Expiry Date: 2021-06-11



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 660-665 Eskdale Road Winnersh Triangle Wokingham Berkshire RG41 5TS 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 Nanhui District Shanghai 201300 China	Production of equipment, information software and reagents for R&D laboratory use in cell biology, biochemistry and physiology.
Molecular Devices (Austria) GmbH Lagerhausstrasse 45 5071 Wals Austria	Design, development, and production 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: 2018-07-12

Effective Date: 2018-06-12

Expiry Date: 2021-06-11

Page: 2 of 2