

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, 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:


Carlos Pitanga, Chief Operating Officer Assurance – Americas

Original Registration Date: 2009-06-12

Latest Revision Date: 2018-12-19

Effective Date: 2018-06-12

Expiry Date: 2021-06-11



Page: 1 of 3

...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.
Molecular Devices 47661 Fremont Boulevard Fremont California 94538 USA	Manufacture of Octet, BLItz and Pioneer instruments and consumables, MicroReactor cassettes and Genedisc calibration tools including field service, technical support, research and development

Original Registration Date: 2009-06-12

Effective Date: 2018-06-12

Latest Revision Date: 2018-12-19

Expiry Date: 2021-06-11

Page: 2 of 3

This certificate remains the property of BSI and shall be returned immediately upon request.

An electronic certificate can be authenticated [online](http://www.bsigroup.com/ClientDirectory). Printed copies can be validated at www.bsigroup.com/ClientDirectory To be read in conjunction with the scope above or the attached appendix.

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.

Certificate No: **FS 534246**

Location

Molecular Devices
No. 88 Shanghai Ke Road
Zhangjiang HI-Tech Park
Shanghai
Shanghai
201210
China

Registered Activities

Manufacture of biosensors and biological experiment analysis instrument and related components.



Original Registration Date: 2009-06-12

Latest Revision Date: 2018-12-19

Effective Date: 2018-06-12

Expiry Date: 2021-06-11

Page: 3 of 3