



Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2008

This is to certify that:

Molecular Devices
1311 Orleans Drive
Sunnyvale
California
94089
USA

Holds Certificate No:

FS 534246

and operates a Quality Management System which complies with the requirements of ISO 9001:2008 for the following scope:

Instrumentation, Software and Reagents for R&D Laboratory Use in Cell Biology, Biochemistry, Physiology.

For and on behalf of BSI:

Reg Blake, VP Regulatory Affairs, BSI Group America Inc.

Original Registration Date: 06/12/2009

Latest Revision Date: 06/08/2015

Effective Date: 06/12/2015

Expiry Date: 06/11/2018

Page: 1 of 2



...making excellence a habit.™

Certificate No: **FS 534246**

Location	Registered Activities
Molecular Devices 1311 Orleans Drive Sunnyvale California 94089 USA	Instrumentation, Software and Reagents for R&D Laboratory Use in Cell Biology, Biochemistry, Physiology.
Molecular Devices 402 Boot Road Downingtown Pennsylvania 19335 USA	Software for R&D Laboratory Use in Cell Biology, Biochemistry, Physiology.
Molecular Devices 660-665 Eskdale Road Winnersh Triangle Wokingham Berkshire RG41 5TS United Kingdom	Service and Distribution of Instrumentation, 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	Instrumentation, Software and Reagents for R&D Laboratory Use in Cell Biology, Biochemistry, Physiology.
Molecular Devices (Austria) GmbH Lagerhausstrasse 45 5071 Wals Austria	Instrumentation, Software and Reagents for R&D Laboratory Use in Cell Biology, Biochemistry, Physiology.

Original Registration Date: 06/12/2009

Latest Revision Date: 06/08/2015

Effective Date: 06/12/2015

Expiry Date: 06/11/2018

Page: 2 of 2

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: + 44 845 080 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.