



# Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2003 & EN ISO 13485:2012

This is to certify that:

**Bio-Rad Laboratories Limited**

Bio-Rad House  
Maxted Road  
Hemel Hempstead  
HP2 7DX  
United Kingdom

Holds Certificate Number:

MD 636076

and operates a Quality Management System which complies with the requirements of ISO 13485:2003 & EN ISO 13485:2012 for the following scope:

Sales, customer technical support, after sales service and installation, warehouse and distribution of in vitro diagnostic products.

For and on behalf of BSI:

Frank Lee, EMEA Compliance & Risk Director

Original Registration Date: 02/09/2015

Latest Revision Date: 02/09/2015

Effective Date: 02/09/2015

Expiry Date: 01/09/2018

Page: 1 of 3



...making excellence a habit.™

Certificate No: MD 636076

Location	Registered Activities
Bio-Rad Laboratories Limited Bio-Rad House Maxted Road Hemel Hempstead HP2 7DX United Kingdom	Sales, customer technical support, after sales service and installation, warehouse and distribution of in vitro diagnostic products.
DiaMed (GB) Limited Unit 2 Whitehill Business Centre Dalkeith Midlothian EH22 2QB United Kingdom	Sales, customer technical support, after sales service and distribution of in vitro diagnostic products.
Bio-Rad Laboratories BV Winninglaan 3 Temse 9140 Belgium	Sales, customer technical support and after sales service of in vitro diagnostic products.
Bio-Rad Laboratories AB Solna Strandväg 3 17154 Solna Sweden	Sales, customer technical support, after sales service and distribution of in vitro diagnostic products.
Bio-Rad Laboratories Torgbygget Nydalsveien 33 N-0421 Oslo Norway	Sales and after sales service of in vitro diagnostic products.
Bio-Rad Laboratories OY Valimotie 1A FIN-00380 Helsinki Finland	Sales and after sales service of in vitro diagnostic products.
Bio-Rad Laboratories Symbion Science Park Fruebjergvej 3 2100 Kobenhavn O Denmark	Sales of in vitro diagnostic products.
DiaMed Benelux, Winninglaan 3 9140 Temse Belgium	Sales, customer technical support and after sales service of in vitro diagnostic products.

Original Registration Date: 02/09/2015

Effective Date: 02/09/2015

Latest Revision Date: 02/09/2015

Expiry Date: 01/09/2018

Page: 2 of 3

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: + 44 845 080 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.

Certificate No: MD 636076

Location

Registered Activities

Bio-Rad Laboratories SA-NV  
Winninglaan 3  
9140 Temse  
Belgium

Sales, customer technical support and after sales service of in vitro diagnostic products.

DiaMed Fennica OY  
Valmotie 1A  
FIN-00380  
Helsinki  
Finland

Sales and after sales service of in vitro diagnostic products.



Original Registration Date: 02/09/2015

Effective Date: 02/09/2015

Latest Revision Date: 02/09/2015

Expiry Date: 01/09/2018

Page: 3 of 3

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: + 44 845 080 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.