nga global assurance

This is to certify that the Quality Management System of:

Bio-Rad Laboratories (Singapore) Pte. Ltd.

1 Kaki Bukit View #03-01 Techview 415941 Republic of Singapore

applicable to:

Manufacture of analytical instruments and chemicals

has been assessed and approved by National Quality Assurance, U.S.A., against the provisions of:

ISO 9001: 2008

KM Bun

For and on behalf of NQA, USA



Certificate Number: 12400 EAC Code: 19 Certified Since: October 12, 2007 Valid Until: September 14, 2018 Reissued: November 11, 2015 Cycle Issued: November 9, 2015

Page 1 of 1

This approval is subject to the company maintaining its system to the required standard, which will be monitored by NQA, USA, 289 Great Road, Suite 105, Acton, MA 01720, an accredited organization under the ANSI-ASQ National Accreditation Board.

nqa global assurance

This is to certify that the Quality Management System of:

Bio-Rad Laboratories (Singapore) Pte. Ltd.

1 Kaki Bukit View #03-01 Techview 415941 Republic of Singapore

applicable to:

Manufacture of analytical instruments and chemicals

has been assessed and approved by National Quality Assurance, U.S.A., against the provisions of:

ISO 13485:2003

KM Bun

For and on behalf of NQA, USA



Certificate Number: 13217 EAC Code: 19 Certified Since: December 23, 2009 Valid Until: November 8, 2018 Reissued: November 11, 2015 Cycle Issued: November 9, 2015

Page 1 of 1

This approval is subject to the company maintaining its system to the required standard, which will be monitored by NQA, USA, 289 Great Road, Suite 105, Acton, MA 01720, an accredited organization under the ANSI-ASQ National Accreditation Board.