



NSAI

Certificate of Registration of Quality Management System to I.S. EN ISO 13485:2012

The National Standards Authority of Ireland certifies that:

BioTek Instruments, Inc
100 Tigan Street Highland Park
Winooski, VT 05404-0998
USA

has been assessed and deemed to comply with the requirements of the above standard in respect of the scope of operations given below:

The Design, Manufacture, Distribution, Service and Installation of in-vitro diagnostic Microplate Readers, Washers, Dispensers, Stackers, Pipetting Work Stations and Accessories.

Additional sites covered under this multi-site certification are listed on the Annex (File No.MD19.2126)

Approved by:
Kevin D. Mullaney
Chief Executive Officer

Approved by:
Susan Murphy
Risk Management Officer

Registration Number: MD19.2126

Certification Granted: Aug 14, 1996

Effective Date: Jun 02, 2015

Expiry Date: Jul 12, 2018



National Standards Authority of Ireland, 1 Swift Square, Northwood, Santry, Dublin 9, Ireland T +353 1 807 3800



Annex to Certificate number MD19.2126

Scope of Registration :

The Design, Manufacture, Distribution, Service and Installation of in-vitro diagnostic Microplate Readers, Washers, Dispensers, Stackers, Pipetting Work Stations and Accessories.

Activity

Location

Headquarters, Design,
Manufacturing, Warehouse,
Distribution, Service/Repair and
Installation

BioTek Instruments, Inc
100 Tigan Street
Highland Park
PO Box 998
Winooski, VT 05404-0998
USA
File No: MD19.2126

Administration, Distribution,
Warehouse, Service/Repair and
Installation

BioTek Instruments GmbH
Kocherwaldstrasse 34
D-74177 Bad Friedrichshall
Germany
File No: MD19.2126/A

**Verified by:
Operations Manager**