



NSAI

Certificate of Registration of Quality Management System to I.S. EN ISO 13485:2016

The National Standards Authority of Ireland certifies that:

BioTek Instruments, Inc
100 Tigan Street
Highland Park
PO Box 998
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, Automated Microscopes, Washers, Dispensers, Stackers, Incubators, Pipetting Work Stations and Accessories.

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

Approved by:
Geraldine Larkin
Chief Executive Officer

Approved by:
Caroline Dore Geraghty
Director of Medical Devices /
Head of Notified Body

Registration Number: MD19.2126
Certification Granted: August 14, 1996
Effective Date: July 13, 2021
Expiry Date: July 12, 2022





NSAI

Annex to Certificate Number: MD19.2126

Scope of Registration:

The Design, Manufacture, Distribution, Service and Installation of In-Vitro Diagnostic Microplate Readers, Automated Microscopes, Washers, Dispensers, Stackers, Incubators, Pipetting Work Stations and Accessories.

Activity

Location

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

BioTek Instruments, Inc
100 Tigan Street
Highland Park
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**