



# 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:  
Geraldine Larkin  
Chief Executive Officer

Approved by:  
Susan Murphy  
European Medical Device  
Operations Manager

---

Registration Number: MD19.2126  
Certification Granted: Aug 14, 1996  
Effective Date: Jul 13, 2018  
Expiry Date: Mar 31, 2019





## **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  
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**