



NSAI

Certificate of Registration of Quality Management System to I.S. EN ISO 9001:2008

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 research and development and in-vitro diagnostic products.

Approved by:
Kevin D. Mullaney
Chief Executive Officer

Approved by:
Lisa Greenleaf
Operations Manager

Registration Number: 19.2126
Certification Granted: Aug 14, 1996
Effective Date: Jun 02, 2015
Expiry Date: Jul 12, 2018



National Standards Authority of Ireland, 402 Amherst Street, Nashua, New Hampshire, NH 03063, USA T +1 603 882 4412