



BioTek Instruments, Inc.
Highland Park, P.O. Box 998
Winooski, Vermont 05404-0998, USA

Tel: 888-451-5171
Fax: 802-655-7941
Outside the USA: 802-655-4740
Email: customercare@biotek.com
www.biotek.com



CERTIFICATE OF SOFTWARE VALIDATION

As a responsible official of BioTek Instruments, Inc., I hereby certify that all versions of Gen5 software have been fully validated to operate per written product specifications. This validation was conducted utilizing several controlled test protocols, which challenged the data transfer, data reduction and instrument interface functionality. The integrity of the Microsoft[®] Excel spreadsheet data transfer and operation was also validated. Additionally, the Electronic Signature aspects of the software (Gen5 Secure) were fully validated per US FDA's 21 CFR Part 11 software requirements. All software validation activities are certified to meet the US FDA and ISO 9001/ISO 13485 requirements.

The software validation documentation, Master Software Copy, and Source Code are available for review by authorized agents of a customer, US FDA, Health Canada, the European Community or other qualified governmental bureaus at:

BioTek Instruments, Inc.
100 Tigan Street
Highland Park
P.O. Box 998
Winooski, Vermont 05404-0998
USA

Field updates of software are handled based on the type of update. That is, problem fixes are distributed free-of-charge to the Sales Order site, while enhancements are shipped on a pro-rated basis depending on age of product.

I hereby certify that the above information is true and accurate:

Michael N. Sevigny
Director Quality Assurance and Regulatory Affairs
sevignym@biotek.com
802-861-8618

5 August 2015

