




## EC Certificate of Conformity

I, the undersigned, hereby declare that the Product(s)

### **Picus® & Picus® NxT Electronic Pipettes, Charging Stand and Charging Carousel**

conforms to the protection requirements of Council Directive 2004/108/EC, relating to Electromagnetic Compatibility, EN61326 and Safety IEC 61010-1 ed 3.0, by the application of:  
Technical Construction File No. 01/12 dated 26/04/2012 and Competent Body Technical Report/Certificate No. 198349 dated 25/04/2012 and Report No.198349E, dated 31/05/2012 issued by:  
Nemko Oy , Perkkäantie 11, FIN-02601 Espoo, Finland, Tel: +358-424 545 41

This declaration is supported by EC quality system approval certificates:  
ISO 9001 / ISO 13485 Certificate No.108129-2011-AQ-FIN-FINAS issued by DNV on 20. December 2011  
ISO 17025 Certificate No. K041 issued by FINAS on 16. December 2011  
ISO 14001 Certificate No.108133-2011-AE-HEL-FINAS issued by DNV on 20 December 2011

Signed:  Date: 3<sup>rd</sup> June 2014  
Full Name: Jussi Heiniö  
Title: CEO  
Company: Sartorius Biohit Liquid Handling Oy, Laippatie 1, 00880 Helsinki, Finland