

Declaration of Conformity



Sartorius Biohit Liquid Handling Oy hereby certifies that:

Picus® Nxt Electronic Pipette 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. 02/14 dated 04/07/2014 and Competent Body Technical Report/Certificate No. 281023-1 dated 04/09/2015 and Report No.281012-1, dated 08/09/2015 issued by: SGS Fimko Ltd, Särkiniementie 3, FI-00211 Helsinki, Finland, Tel: +358 9 6963 361

Picus® Electronic Pipette 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, Perkkaantie 11, FIN-02601 Espoo, Finland, Tel: +358 424 545 41

Midi Plus Pipetting Controller conforms to the protection requirements of Council Directive EMC Directive 2004/108/EC, relating to Electromagnetic Compatibility, by the Competent Body Technical Report/Certificate No. 02-00924 dated 28/06/2013 issued by: Eleforss Oy, Wahreninkatu 11B, 30100 Forssa.

The following harmonized standards have been applied:

CISPR11:2015

EN 61000-4-2:2009

EN 61000-4-3:2006+AMD2:2010

EN 61000-4-4:2012

EN 61000-4-5:2014

EN 61000-4-6:2014

EN 61000-4-8:2010

EN 61000-4-11:2004

2011/65/EU for RoHS Directive.

This declaration is supported by following EC quality system approval certificates:

ISO 9001:2015 Certificate No. FI17/5101, issued by SGS, Finland, valid 14th December 2020

ISO 13485:2016 Certificate No. FI17/5103, issued by SGS, Finland, valid 14th December 2020

ISO 17025:2005 Certificate No. K041/A10/2016 issued by FINAS, valid 20th March 2024

ISO 14001:2015 Certificate No. FI17/5102, issued by SGS, Finland, valid 14th December 2020

Helsinki, 10th August 2020

Seppo Riikonen

Manager, Quality & Process Development
Sartorius Biohit Liquid Handling Oy