

Certificate of Conformity



Sartorius Biohit Liquid Handling Oy hereby confirm that the product line(s):

Picus & NxT Electronic pipettes+ Charging Stands
eLINE Electronic pipettes + Charging Stands
Biohit Midi Plus
Tacta Mechanical pipettes
mLINE Mechanical pipettes
Proline Plus Mechanical pipettes
Proline Mechanical pipettes

Are manufactured in:
Sartorius Biohit Liquid Handling Oy
Tietokatu 4, 87400 Kajaani, FINLAND

Country of origin is Finland.

All product lines are designed and manufactured according to the Essential Requirements listed in Annex I of EC Directive 98/79/EC (In Vitro Medical Devices).

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 2020

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

Furthermore the Calibration Laboratory for Volumetric Instruments at Sartorius Biohit has been officially accredited by the Finnish Accreditation Service (FINAS) according to the **ISO/IEC 17025** (Cert. No. K041/2012).

All Sartorius Biohit products are individually tested according to the ISO 8655 and all delivered units have serial numbered QC-certificate included. The QC-certificate indicates that the unit has passed the given specifications.

Signed: 

Helsinki 16th January 2019

Full Name: **Seppo Riikonen**
Title: **Director, Quality & Process Development**
Company: **Sartorius Biohit Liquid Handling Oy, Laippatie 1, 00880 Helsinki, Finland**