

## Certificate of Conformity

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

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.

This declaration is supported by following EC quality system approval certificates: ISO 9001:2015 Certificate No. FI17/5101, issued by SGS, Finland, valid 14<sup>th</sup> December 2020 ISO 13485:2016 Certificate No. FI17/5103, issued by SGS, Finland, valid 14<sup>th</sup> December 2020 ISO 17025:2017 Certificate No. K041 issued by FINAS, valid 20<sup>th</sup> March 2024 ISO 14001:2015 Certificate No. FI17/5102, issued by SGS, Finland, valid 14<sup>th</sup> 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 17025:2017 (Cert. No. KO41)

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 10<sup>th</sup> August 2020

Full Name: Seppo Riikonen

Title: Manager, Quality & Process Development

Company: Sartorius Biohit Liquid Handling Oy, Laippatie 1, 00880 Helsinki, Finland