



Certificate FI17/5103

The management system of

# Sartorius Biohit Liquid Handling Oy

Laippatie 1  
00880 Helsinki  
Finland

has been assessed and certified as meeting the requirements of



## ISO 13485:2016 EN ISO 13485:2016

For the following activities

**Design, development, manufacture, marketing, sales,  
training and service of liquid handling products, their  
disposables for research, health care and industrial  
laboratories**

This certificate is valid from 15<sup>th</sup> of December 2020  
until 14<sup>th</sup> of December 2023  
Issue 4. Certified since June 1997

Certified activities performed by additional sites listed on subsequent pages.

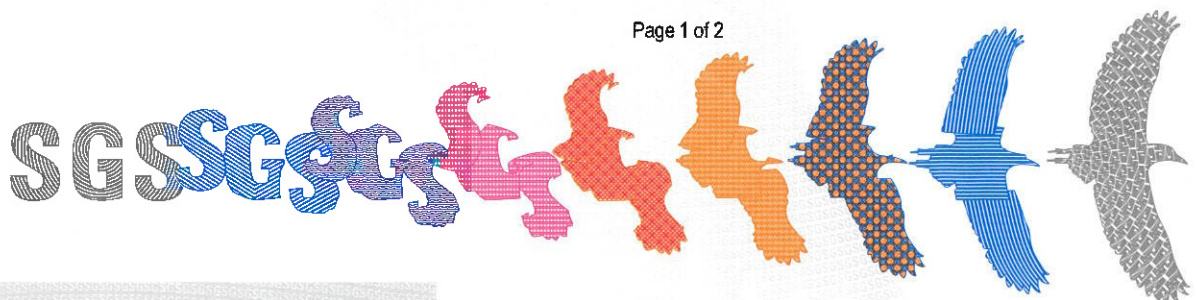
Authorised by

Harri Kaartinen  
Manager

SGS Fimko Ltd Certification and Business Enhancement  
Takomotie 8, FI-00380 Helsinki, Finland  
t +358 (0)9 6963 61 - f +358 (0)9 692 5474 www.sgs.fi



Page 1 of 2



This document is issued by the Company subject to its General Conditions of Certification Services accessible at [www.sgs.com/terms\\_and\\_conditions.htm](http://www.sgs.com/terms_and_conditions.htm). Attention is drawn to the limitations of liability, indemnification and jurisdictional issues established therein. The authenticity of this document may be verified at <http://www.sgs.com/en/certified-clients-and-products/certified-client-directory>. Any unauthorized alteration, forgery or falsification of the content or appearance of this document is unlawful and offenders may be prosecuted to the fullest extent of the law.

# Sartorius Biohit Liquid Handling Oy

## ISO 13485:2016 EN ISO 13485:2016



Issue 4

**Certified activities are performed by the sites on the list.**

**Sites:**

- |  |   |
|--|---|
| <p>Laippatie 1<br/>00880 Helsinki<br/>Finland</p>    | <p>Design, development, marketing, training and service of liquid handling products, their disposables for research, health care and industrial laboratories</p>  |
| <p>Tulppatie 1<br/>00880 Helsinki<br/>Finland</p>    | <p>Design and development of molds for research, health care and industrial laboratories and manufacture of liquid handling products, their disposables for research, health care and industrial laboratories</p> |
| <p>Tietokatu 4<br/>87400 Kajaani<br/>Finland</p>     | <p>Design, development and manufacture of liquid handling products, their disposables for research, health care and industrial laboratories</p>   |
| <p>Veistäjäncatu 1<br/>87400 Kajaani<br/>Finland</p> | <p>Manufacture of liquid handling products, their disposables for research, health care and industrial laboratories</p>   |

