



Certificate FI17/5101

The management system of

Sartorius Biohit Liquid Handling Oy

Laippatie 1
FI-00880 Helsinki
Finland



has been assessed and certified as meeting the requirements of

ISO 9001:2015

For the following activities

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

This certificate is valid from 15th of December 2020
until 14th of December 2023
Issue 3. Certified since June 1997

Certified activities performed by additional sites listed on subsequent pages.

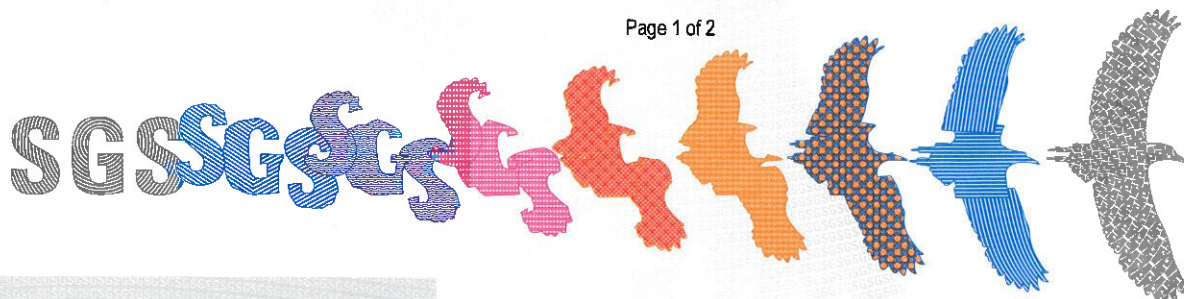
Authorised by

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. 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 9001:2015



Issue 3

Certified activities are performed by the sites on the list.

Sites:

<p>Laippatie 1 00880 Helsinki Finland</p>	<p>Design, development, marketing, training and service of liquid handling products, their disposables for research, healthcare and industrial laboratories</p>
<p>Tulppatie 1 00880 Helsinki Finland</p>	<p>Manufacture of liquid handling and plastic products, design and development of their molds for research, healthcare and industrial laboratories</p>
<p>Tietokatu 4 87400 Kajaani Finland</p>	<p>Design, development and manufacture of liquid handling products, their disposables for research, healthcare and industrial laboratories</p>
<p>Veistäjänkatu 1 87400 Kajaani Finland</p>	<p>Logistic center for liquid handling and their disposables manufactured in Kajaani</p>

