

Sartorius Biohit Liquid Handling Oy

Laippatie 1
FI-00880 Helsinki
Finland



has been assessed and certified as meeting the requirements of

ISO 14001:2015

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 15th of December 2017
until 14th of December 2020
Issue 2. Certified since February 2003

This is a multi-site certification
Additional site details are listed on the subsequent page

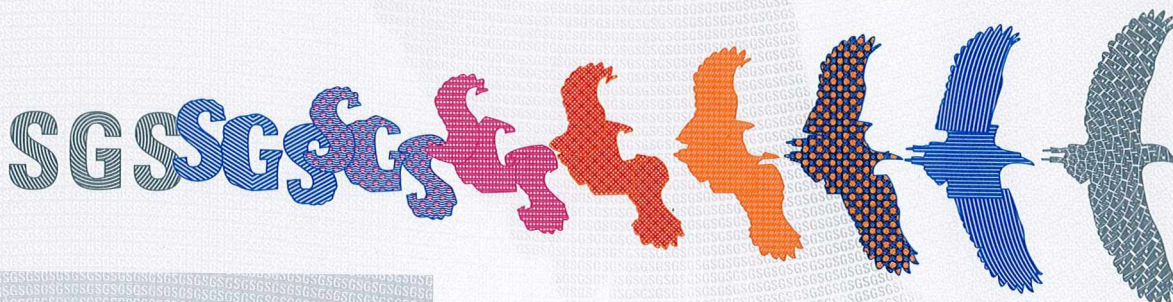
Authorised by



Harri Kaartinen
Manager



SGS Fimko Ltd Certification and Business Enhancement
Särkiniementie 3, 00210 Helsinki, Finland
t +358 (0)9 6963 61 - f +358 (0)9 692 5474 www.sgs.fi



Sartorius Biohit Liquid Handling Oy

ISO 14001:2015



Issue 2.

Detailed scope

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

Additional facilities

Laippatie1
FI-00880 Helsinki
Finland

Tietokatu 4
FI-87400 Kajaani
Finland

