

The management system of

Sartorius Stedim Biotech

August-Spindler-Str. 11
37079 Göttingen, Germany

has been assessed and certified as meeting the requirements of

Directive 93/42/EEC

on medical devices, Annex II (excluding Section 4)

For the following products

Sterile and non-sterile syringe filters as injection filters and vent filters for removal of particles from sterile liquids or from sterile dry air and gases or for ultracleaning of air and gases or liquids for eventual infusion, administration or introduction into the human body.

This certificate is valid from 02/08/2019 until 16/06/2022 and remains valid subject to satisfactory surveillance audits.

Issue 1. Certified since 16/06/2008 and first certified by SGS Belgium on 02/08/2019
Re certification audit due before 05/05/2020

Certification is based on reports numbered WW/PCI 614440

Authorised by

Pieter Weterings
Certification Manager

SGS Belgium NV, Notified Body 1639

SGS House Noorderlaan 87 2030 Antwerp Belgium
t +32 (0)3 545-48-48 f +32 (0)3 545-48-49 www.sgs.com

LPMD5007 - Certificate CE1639 Annex II-4_EN rev. 02

Page 1 of 1

