

Certificate US19/819943255

SGS

The management system of

Sartorius Stedim Biotech

August-Spindler-Str. 11
37079 Göttingen, Germany



has been assessed and certified as meeting the requirements of

ISO 13485:2016 EN ISO 13485:2016

For the following activities

**Design and development, manufacture and distribution of products
for sterile and particle filtration of liquids and gases.**

This certificate is valid from 02/08/2019 until 16/06/2020
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

Authorised by

Pieter Weterings
Certification Manager
SGS Belgium NV

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

SGS Belgium 13485-2 0308



Accreditation Number

005-CMS
EN ISO/IEC 17021-1:2015

Page 1 of 1



This document is issued by the Company subject to its General Conditions of Certification Services, unless otherwise agreed, 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.