



Certificate GB18/962158

The management system of

Sartorius Stedim UK Ltd.

Unit 4, Longmead Business Centre, Blenheim Road, Epsom,
Surrey, KT19 9QQ, UK

has been assessed and certified as meeting the requirements of

ISO 9001:2015

For the following activities

The sales, supply, installation, maintenance and after sales support of products and integrated solutions for fermentation, filtration, purification, fluid management and laboratory and process applications.

This certificate is valid from 21 October 2018 until 21 October 2021 and remains valid subject to satisfactory surveillance audits. Recertification audit due a minimum of 60 days before the expiration date. Issue 1. Certified since 03 September 2010



Authorised by



0005

SGS United Kingdom Ltd
Rossmore Business Park Ellesmere Port Cheshire CH65 3EN UK
t +44 (0)151 350-6666 f +44 (0)151 350-6600 www.sgs.com

HC SGS 9001 2015 0818

Page 1 of 1



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.

