



Certificate GB09/78063

The management system of

Sartorius Stedim Lab Ltd

Sperry Way, Stonehouse Park, Stonehouse,
Gloucestershire, GL10 3UT, UK

has been assessed and certified as meeting the requirements of

ISO 9001:2015



For the following activities

The manufacture of disposable filtration devices.

This certificate is valid from 20 September 2018 until 20 September 2021
and remains valid subject to satisfactory surveillance audits.
Recertification audit due a minimum of 60 days before the expiration date.
Issue 8. Certified since 12 January 2004

Expiry date of last certificate: 15 September 2018
End date of last recertification audit: 23 August 2018



Authorised by

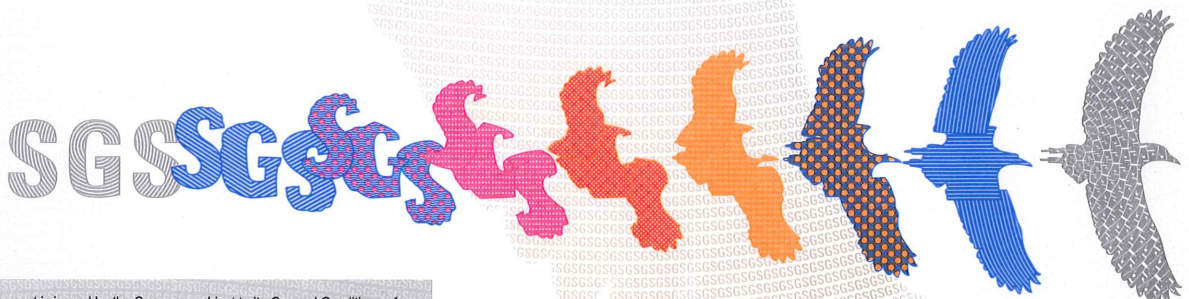
SGS United Kingdom Ltd
Rossmore Business Park Ellesmere Port Cheshire CH65 3EN UK
t +44 (0)151 350-6666 f +44 (0)151 350-6600 www.sgs.com

HC SGS 9001 2015 0818

Page 1 of 1



0005



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.