



Certificate GB12/86848

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 13485:2016 EN ISO 13485:2016



For the following activities

The manufacture of non-sterile single use filtration devices for the concentration of macromolecules and biological samples for analysis in IVD applications.

This certificate is valid from 20 October 2018 until 20 October 2021 and remains valid subject to satisfactory surveillance audits.

Re certification audit due before 05 October 2021

Issue 4. Certified since 17 October 2012

Authorised by

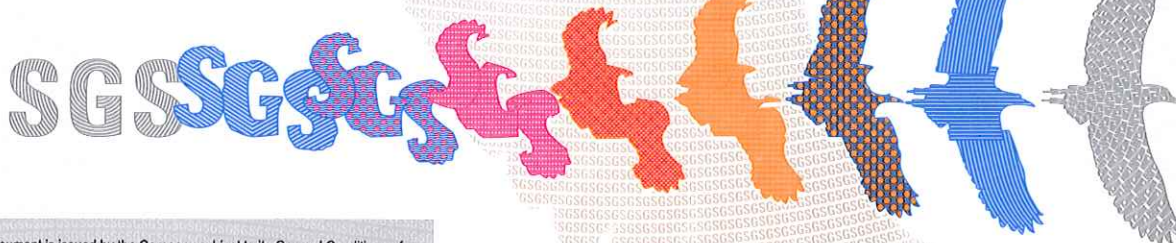
SGS United Kingdom Ltd
Rossmore Business Park Ellesmere Port Cheshire CH65 3EN UK
t +44 (0)151 350-6666 f +44 (0)151 350-6600 www.sgs.com

HC SGS 13485 2016 0118

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.