

The management system of

Sartorius Stedim India Pvt. Ltd.

No.-69/2 & 69/3, Jakkasandra, Kunigal Road, Nelamangala Taluk,
Bangalore, 562123, India

has been assessed and certified as meeting the requirements of

ISO 9001:2015

For the following activities

- (1) Design, Manufacture, Installation and Service of Products and Systems for Bioprocess Applications
- (2) Product Development, Validation Projects and Services for Pharma and Bio-Pharmaceuticals industry

This certificate is valid from 26 June 2019 until 18 April 2022
and remains valid subject to satisfactory surveillance audits.
Recertification audit due before 17 February 2022
Issue 5. Certified since 01 July 2010

The audit leading to this certificate commenced on 18 March 2019
Previous issue certificate validity date was until 17 April 2019

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

