

质量管理体系认证证书

注册号: 016ZB19Q30195R0S
统一社会信用代码: 91110113677401025F

兹证明

赛多利斯斯泰迪生物设备（北京）有限公司

质量管理体系符合
GB/T19001-2016 idt ISO9001:2015 标准，适用于

生产应用于生物制药的一次性容器（塑料类）

注册地址: 北京市顺义区临空经济核心区裕安路33号二楼西侧
经营地址: 北京市顺义区临空经济核心区裕安路33号二楼西侧

初次发证日期: 2019年01月30日
证书有效期至: 2022年01月29日

新世纪检验认证股份有限公司
总裁:

尚志强



BCC地址: 北京市西城区国英园1号楼11层1101室
本证书在国家规定的各行政许可、资质许可有效期内使用有效
获证组织必须定期接受监督审核并经审核合格, 此证书方继续有效。
证书有效性可通过网站: www.bcc.com.cn 查询, 也可二维码查询
本证书信息可在国家认监委网站www.cnca.gov.cn 查询



中国认可
国际互认
管理体系
MANAGEMENT SYSTEM
CNAS C016-M



CERTIFICATE OF QUALITY MANAGEMENT SYSTEM CERTIFICATION

Certificate No.016ZB19Q30195R0S
Unified social credit code:91110113677401025F

This is to Certify that the Quality Management System of

SARTORIUS STEDIM BIOTECH (BEIJING)CO., LIMITED

is in conformity with
GB/T19001-2016 idt ISO9001:2015 Standard, applies to

MANUFACTURE OF SINGLE USE CONTAINERS FOR BIOPHARMACENEICAL APPLICATIONS
(PLASTIC)

REGISTERED ADDRESS:WEST SIDE OF THE 2ND/F, NO.33 YU'AN RD., AIRPORT ECONOMIC
ZONE, SHUNYI DISTRICT, BEIJING, P.R.C.
OPERATION ADDRESS:WEST SIDE OF THE 2ND/F, NO.33 YU'AN RD., AIRPORT ECONOMIC
ZONE, SHUNYI DISTRICT, BEIJING, P.R.C.

Date of Initial Issuance: Jan. 30, 2019
Date of Expiration: Jan. 29, 2022

BCC Inc.
President:

尚志强



BCC Address: Room 1101, Floor 11, Building 1, Guoyingyuan, Xicheng District, Beijing, P.R.C.
This certificate is valid within state - specified validities of administrative and qualification licensing.
The effectiveness of this certificate shall be maintained by regular surveillance audit,
the validity of the certificate can be inquired through www.bcc.com.cn or by QR code.
The information of this certificate available for inquiry on CNCA's website: www.cnca.gov.cn.



中国认可
国际互认
管理体系
MANAGEMENT SYSTEM
CNAS C016-M

