

# 质量管理体系认证证书

注册号: 016ZB18Q35060R3M

统一社会信用代码: 911101137839553986

兹证明

赛多利斯科学仪器(北京)有限公司

质量管理体系符合

GB/T19001-2016/ISO9001:2015标准, 适用于

电子天平、酸度计、热重水分测定仪和称重模块的组装、销售和售后服务

注册地址: 北京市顺义区天竺空港工业区B区裕安路33号厂房办公楼

经营地址: 北京市顺义区天竺空港工业区B区裕安路33号厂房办公楼

初次发证日期: 2010年06月12日

再认证日期: 2018年12月19日

证书有效期至: 2021年12月16日

(本次再认证审核日期: 2018年11月26日到2018年11月28日, 上一认证周期截止时间: 2018年12月16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](http://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](http://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.016ZB18Q35060R3M  
Unified social credit code:911101137839553986

This is to Certify that the Quality Management System of

**SARTORIUS SCIENTIFIC INSTRUMENTS (BEIJING) CO.,  
LIMITED**

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

ASSEMBLY, SALES AND SERVICE OF ELECTRONIC BALANCE, ACIDITY METER,  
THERMOGRAVIMETRIC MOISTURE METER AND WEIGHING MODULE

REGISTERED ADDRESS:WORKSHOP BULIDING, NO.33, YU'AN RD., SHUNYI DISTRICT, ZONE B,  
TIANZHU AIRPORT INDUSTRIAL PARK, BEIJING, P.R.C.  
OPERATION ADDRESS:WORKSHOP BULIDING, NO.33, YU'AN RD., SHUNYI DISTRICT, ZONE B,  
TIANZHU AIRPORT INDUSTRIAL PARK, BEIJING, P.R.C.

Date of Initial Issuance: Jun. 12, 2010

Date of Re-Certification: Dec. 19, 2018

Date of Expiration: Dec. 16, 2021

(Date of Recertification Audit:Nov. 26, 2018 To Nov. 28, 2018,expiry date of the last certification cycle:Dec. 16, 2018)

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](http://www.bcc.com.cn) or by QR code.  
The information of this certificate available for inquiry on CNCA's website: [www.cnca.gov.cn](http://www.cnca.gov.cn).



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

