

# CERTIFICATE OF OCCUPATIONAL HEALTH AND SAFETY MANAGEMENT SYSTEM CERTIFICATION

Certificate No.016ZB21S32535R3M  
Unified social credit code:911101137839553986

This is to Certify that the Occupational Health and Safety Management System of

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

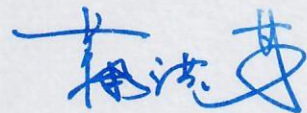
is in conformity with  
GB/T 45001-2020 idt ISO 45001:2018 Standard, applies to

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

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

Date of Initial Issuance:Jun. 12, 2010  
Date of Re-certification Issuance:Dec. 1, 2021  
Date of Expiration:Dec. 16, 2024

BCC Inc.  
President:



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

