

Certificate US15/842193.00

The management system of

# Sartorius Corporation

565 Johnson Avenue  
Bohemia, NY 11716, United States

has been assessed and certified as meeting the requirements of

## ISO 9001:2015

For the following activities:

**Sales, light manufacturing, support, distribution, in-house service and calibration of laboratory equipment including weighing equipment, pipettes and related peripherals, and moisture lab applications including testing. Management of field service, repair, installation, qualification and calibration of divisional laboratory products.**

Further clarifications regarding the scope of this certificate and the applicability of ISO 9001:2015 requirements may be obtained by consulting the organization

This certificate is valid from 9 June 2019 until 9 June 2021 and remains valid subject to satisfactory surveillance audits. Recertification audit due a minimum of 60 days before the expiration date. Issue 6: 16 September 2019. Certified since June 2015.

The audit leading to this certificate commenced on 26/07/2019. Previous issue certificate validity date was until 09/06/2021.

This is a multi-site certification.  
Additional site details are listed on subsequent pages.

Authorized by:



Dan Seal

Technical Accreditation Manager, Certification &  
Business Enhancement North America  
SGS North America, Inc.

201 Route 17 North, Rutherford, NJ 07070, USA  
t (201) 508-3000 f (201) 935-4555 [www.us.sgs.com](http://www.us.sgs.com)

This certificate remains the property of SGS and shall be returned upon request

Page 1 of 2



# Sartorius Corporation

## ISO 9001:2015

Issue 6:16 September 2019



Additional facilities:

**565 Johnson Avenue, Bohemia, NY 11716, United States**  
**Scope: Senior management, sales, order processing, support.**

**1173 North Service Road West, Unit D-4, Oakville, ON L6M 2V9, Canada**  
**Scope: Sales, support, service, installation, qualification and calibration of laboratory equipment including weighing equipment, pipettes and related peripherals.**

**545-1 Johnson Avenue, Bohemia, NY 11716, United States**  
**Scope: Sales, support, Technical service, in-house service and calibration of laboratory equipment including weighing equipment, pipettes and related peripherals, and moisture lab applications including testing. Management of field service, repair, installation, qualification and calibration of divisional laboratory products.**

**6542 Fig Street, Arvada, CO 80004, United States**  
**Scope: Product distribution of regulated and non-regulated products, Light manufacturing, final assembly, configuration management and in-house servicing of non-regulated products.**



Certificate US15/842193.01

The management system of

# Sartorius Corporation

6542 Fig St.  
Arvada, CO 80004 United States

has been assessed and certified as meeting the requirements of

## ISO 9001:2015

For the following activities:

**Product distribution of regulated and non-regulated products, Light manufacturing, final assembly, configuration management and in-house servicing of non-regulated products.**

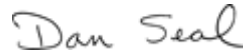
Further clarifications regarding the scope of this certificate and the applicability of ISO 9001:2015 requirements may be obtained by consulting the organization.

This certificate is valid from 9 June 2018 until 9 June 2021 and remains valid subject to satisfactory surveillance audits. Recertification audit due a minimum of 60 days before the expiration date. Issue 3: 16 September 2019. Certified since June 2015.

Multiple certificates have been issued for this scope. The main certificate is numbered US15/842193.00

The audit leading to this certificate commenced on 07/05/2019. Previous issue certificate validity date was until 09/06/2021.

Authorized by:



Dan Seal

Technical Accreditation Manager, Certification & Business Enhancement North America  
SGS North America, Inc.

201 Route 17 North, Rutherford, NJ 07070, USA  
t (201) 508-3000 f (201) 935-4555 [www.us.sgs.com](http://www.us.sgs.com)

This certificate remains the property of SGS and shall be returned upon request

Page 1 of 1

