

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.

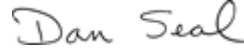
Further clarifications regarding the scope of this certificate and the applicability of ISO 9001:2015 requirements may be obtained by consulting the organization

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

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 US15/842193.02

The management system of

Sartorius Corporation

1173 North Service Road West Unit D-4
Oakville ON L6M 2V9, Canada

has been assessed and certified as meeting the requirements of

ISO 9001:2015

For the following activities:

Sales, support, service, installation, qualification and calibration of laboratory equipment including weighing equipment, pipettes and related peripherals.

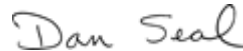
Further clarifications regarding the scope of this certificate and the applicability of ISO 9001:2015 requirements may be obtained by consulting the organization.

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 2: 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

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

Page 1 of 1

