

# CERTIFICATE OF REGISTRATION

This is to certify that the management system of:

## Sartorius Stedim North America, Inc.

Main Site: 565 Johnson Avenue, Bohemia, New York,  
11716, United States

Additional Site: 545-1 Johnson Avenue, Bohemia, New York,  
11716, United States

See appendix for additional sites and additional site scopes

has been registered by Intertek as conforming to the requirements of:

## ISO 9001:2015

The management system is applicable to:

Design, Development, Sale and support of products used in bioprocessing.  
Manufacture of membrane filters, single use bags and sampling systems.

*Site: 565 Johnson Avenue, Bohemia, NY 11716, United States*

Management and Sales of bioprocessing components and systems. Management of the design and development of engineered to order bioprocessing systems. Management of field service, repair, and calibration of Sartorius Stedim bioprocessing equipment and installation. Design and development of products used in bioprocessing.

*Site: 545-1 Johnson Avenue, Bohemia, NY 11716, United States*

Management and Sales of bioprocessing components and systems. Management of the design and development of engineered to order bioprocessing systems. Management of field service, repair, and calibration of Sartorius Stedim bioprocessing equipment and installation. Design and development of products used in bioprocessing.

**Certificate Number:**

0096894

**Initial Certification Date:**

September 2002

Certified by SGS before 03/12/2019 under  
Certificate# US09/5619.00

**Date of Certification Decision:**

03 December 2019

**Issuing Date:**

03 December 2019

**Valid Until:**

04 March 2022



TM



**Intertek**

**Calin Moldovean**

President, Business Assurance

Intertek Testing Services NA, Inc. dba Intertek  
900 Chelmsford Street, Lowell, MA, USA



# APPENDIX TO CERTIFICATE OF REGISTRATION

This appendix identifies the locations by the management system of

## Sartorius Stedim North America, Inc.

This appendix is linked to the Main Certificate #0096894 and cannot be shown nor reproduced without it.

Barriada Arturo Lluberas, Road 128, Intersection 376,  
Yauco, 00698, Puerto Rico

**Production of Membrane for filters, production of flat filters, injection molding, product validation laboratory.**

Bo Susua, Susua industrial Park, Road 128 Intersection 368, Km 2.1 Lot 6,  
Yauco, 00698, Puerto Rico

**Production of membranes filters and bag assemblies including the Quality, purchasing, receiving, kitting and shipping for Puerto Rico Operations.**

80 Progress Avenue, New Oxford, Pennsylvania, 17350, United States

**Manufacture of single use sampling systems including the Quality, purchasing, receiving, kitting and shipping operations.**

