



sartorius stedim  
biotech

## Environmental impact of single-use technologies



turning science **into solutions**

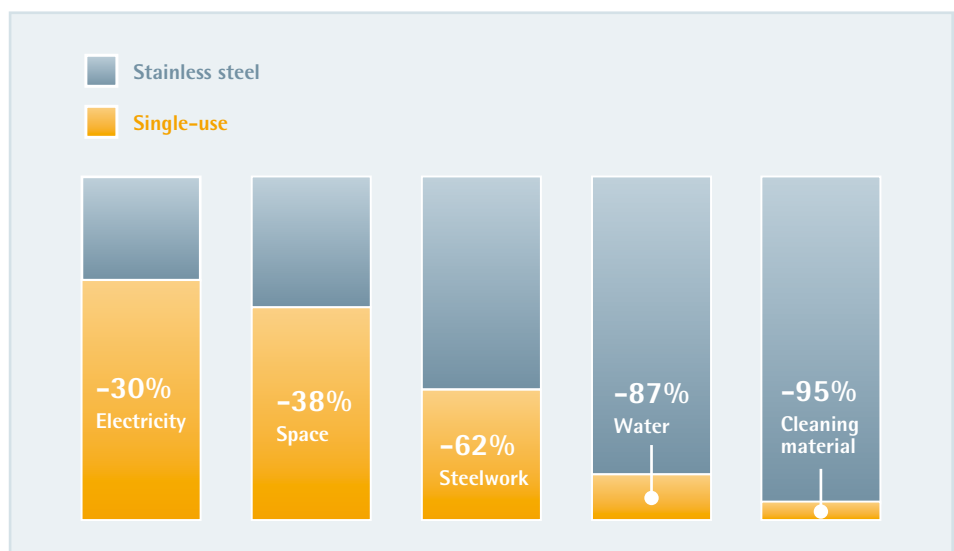
# Implementation of single-use technologies can reduce a facility's environmental footprint.

Utilizing single-use technology can decisively improve the environmental footprint of biomanufacturing processes. Single-use equipment from Sartorius Stedim Biotech help lower energy consumption and reduce climate-changing emissions. In particular, single-use technology can dramatically reduce the usage of water, a resource indispensable for life, by more than 80%.\*



Utilization of single-use products results in significant savings of resources consumed in manufacture. Single-use systems do not require the amount of energy needed for water heating and preparation. It reduces the water and cleaning agent consumption greatly. As a result, single-use technology helps contribute to substantially reducing your company's overall carbon footprint.\*

\* Reference: Sinclair, A.; Leveen, et al.; The Environmental Impact of Disposable Technologies, The Biopharm International Guide, November 2008; Base of the analysis: Typical mAb process at 3 x 2000 L scale





At Sartorius Stedim Biotech, we work together for one primary goal: To provide our customers with the best solutions for their processes. That's why we deliver choices instead of a narrow band of technologies. Our products and services cover the entire process of the biopharmaceutical industry – from development to quality assurance and production.

Single-use technologies and equipment from Sartorius Stedim Biotech enable you to meet the rapidly changing market and regulatory requirements. Our customers rely on Sartorius Stedim Biotech single-use products in all biopharmaceutical process phases to enhance their production capacity and process flexibility, while maintaining the highest quality. Our single-use concepts prevent cross-contamination, eliminate cost-intensive cleaning cycles and validation, and save time when you need to produce different products simultaneously.

# Sales and Service Contacts

For further contacts, visit [www.sartorius-stedim.com](http://www.sartorius-stedim.com)

## Europe

### Germany

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen

Phone +49.551.308.0  
Fax +49.551.308.3289

[www.sartorius-stedim.com](http://www.sartorius-stedim.com)

Sartorius Stedim Systems GmbH  
Schwarzenberger Weg 73-79  
34212 Melsungen

Phone +49.5661.71.3400  
Fax +49.5661.71.3702

[www.sartorius-stedim.com](http://www.sartorius-stedim.com)

### France

Sartorius Stedim Biotech S.A.  
ZI Les Paluds  
Avenue de Jouques – BP 1051  
13781 Aubagne Cedex

Phone +33.442.845600  
Fax +33.442.845619

Sartorius Stedim France SAS  
ZI Les Paluds  
Avenue de Jouques – CS 71058  
13781 Aubagne Cedex

Phone +33.442.845600  
Fax +33.442.846545

### Austria

Sartorius Stedim Austria GmbH  
Franzosengraben 12  
A-1030 Vienna

Phone +43.1.7965763.18  
Fax +43.1.796576344

### Belgium

Sartorius Stedim Belgium N.V.  
Leuvensesteenweg, 248/B  
1800 Vilvoorde

Phone +32.2.756.06.80  
Fax +32.2.756.06.81

### Denmark

Sartorius Stedim Nordic A/S  
Hoerskaetten 6D, 1.  
DK-2630 Taastrup

Phone +45.7023.4400  
Fax +45.4630.4030

### Italy

Sartorius Stedim Italy S.p.A.  
Via dell'Antella, 76/A  
50012 Antella-Bagno a Ripoli (FI)

Phone +39.055.63.40.41  
Fax +39.055.63.40.526

### Netherlands

Sartorius Stedim Netherlands B.V.  
Edisonbaan 24  
3439 MN Nieuwegein

Phone +31.30.6025080  
Fax +31.30.6025099

### Spain

Sartorius Stedim Spain SA  
C/Isabel Colbrand 10,  
Oficina 70  
Polígono Industrial de Fuencarral  
28050 Madrid

Phone +34.90.2110935  
Fax +34.91.3589623

### Switzerland

Sartorius Stedim Switzerland GmbH  
Lerzenstrasse 21  
8953 Dietikon

Phone +41.44.741.05.00  
Fax +41.44.741.05.09

### U.K.

Sartorius Stedim UK Limited  
Longmead Business Park  
Blenheim Road, Epsom  
Surrey KT19 9 QQ

Phone +44.1372.737159  
Fax +44.1372.726171

## America

### USA

Sartorius Stedim North America Inc.  
5 Orville Drive  
Bohemia, NY 11716

Toll-Free +1.800.368.7178  
Fax +1.631.254.4253

Sartorius Stedim SUS Inc.  
1910 Mark Court  
Concord, CA 94520

Phone +1.925.689.6650  
Toll Free +1.800.914.6644  
Fax +1.925.689.6988

### Sartorius Stedim Systems Inc.

201 South Ingram Mill Road  
Springfield, MO 65802

Phone +1.417.873.9636  
Fax +1.417.873.9275

### Argentina

Sartorius Argentina S.A.  
Int. A. Avalos 4251  
B1605ECS Munro  
Buenos Aires

Phone +54.11.4721.0505  
Fax +54.11.4762.2333

### Brazil

Sartorius do Brasil Ltda  
Av. Dom Pedro I, 241  
Bairro Vila Pires  
Santo André  
São Paulo  
Cep 09110-001

Phone +55.11.4451.6226  
Fax +55.11.4451.4369

### Mexico

Sartorius de México S.A. de C.V.  
Circuito Circunvalación Poniente No. 149  
Ciudad Satélite  
53100 Naucalpan, Estado de México

Phone +52.5555.62.1102  
Fax +52.5555.62.2942

## Asia | Pacific

### China

Sartorius Stedim Beijing  
Representative Office  
No. 33, Yu'an Road,  
Airport Industrial Zone B, Shunyi District  
Beijing 101300

Phone +86.10.80426516  
Fax +86.10.80426580

### Sartorius Stedim Shanghai

Representative Office  
Room 618, Tower 1, German Centre,  
Shanghai, PRC., 201203

Phone +86.21.28986393  
Fax +86.21.28986392.11

### Sartorius Stedim Guangzhou Office

Room 704, Broadway Plaza,  
No. 233-234 Dong Feng West Road  
Guangzhou 510180

Phone +86.20.8351.7921  
Fax +86.20.8351.7931

### India

Sartorius Stedim India Pvt. Ltd.  
10, 6th Main, 3rd Phase Peenya  
KIADB Industrial Area  
Bangalore – 560 058

Phone +91.80.2839.1963|0461  
Fax +91.80.2839.8262

### Japan

Sartorius Stedim Japan K.K.  
KY Building, 8-11  
Kita Shinagawa 1-chome  
Shinagawa-ku  
Tokyo 140-0001

Phone +81.3.3740.5407  
Fax +81.3.3740.5406

### Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.  
Lot L3-E-3B, Enterprise 4  
Technology Park Malaysia  
Bukit Jalil  
57000 Kuala Lumpur

Phone +60.3.8996.0622  
Fax +60.3.8996.0755

### Singapore

Sartorius Stedim Singapore Pte. Ltd.  
10, Science Park Road, The Alpha  
#02-25, Singapore Science Park 2  
Singapore 117684

Phone +65.6872.3966  
Fax +65.6778.2494

### Australia

Sartorius Stedim Australia Pty. Ltd.  
Unit 5, 7-11 Rodeo Drive  
Dandenong South Vic 3175

Phone +61.3.8762.1800  
Fax +61.3.8762.1828