



## Ensuring High Quality and Common Standards

Our centers of excellence in bioprocess engineering ensure highest quality standards through a strong local structure, aligned with the company-wide quality management system:

- We are committed to our customers and to continuous improvement.
- We are regularly audited by certifying agencies and numerous customers.
- Our quality management system is constantly improved and aligned according to internal and external demands.
- Multiple manufacturing plants operating in parallel underline our commitment to flexibility and security for our clients and markets worldwide.



Dr. Klaus J. Walter  
Vice President  
Global Quality

"The harmonization of our quality system guidelines, a robust global change control process and uniform state-of-the-art technical standards guarantee worldwide consistent product quality. Our work is regularly monitored through key performance indicators and assessed by global corporate audits."



# Bioprocess Solutions – Made by Sartorius

## Our Quality Products for Your Success

Since more than 50 years, Sartorius Stedim Biotech serves the biopharma and vaccine industry globally with state of the art cultivation systems.

Our international engineering and manufacturing teams have delivered numerous pilot and production systems worldwide.



Dr. Jens Rupprecht  
Product Manager  
Stainless Steel Fermenter

"Sartorius Stedim Biotech ensures your success thanks to our experience with bioprocess requirements worldwide.

Our long-standing process knowledge combined with the unrivaled quality of our fermenter systems yields to state-of-the-art bioprocess equipment perfectly matching your high demanding applications."



# Engineering & Manufacturing in India

## Securing High Quality of Stainless Steel Products

In India, our center of excellence is located in Bangalore, the home of a newly built, cutting-edge engineering and manufacturing facility opened in 2012.

Since 2008 we forge the stainless steel vessels of our standard instruments (BIOSTAT® Cplus & BIOSTAT® D-DCU) in India. In Germany, where the assembly with the controller takes place, are the vessels integrated and the functionality of the system is being tested according to the customer requirements and our quality directive. Large scale stainless steel projects are typically executed in this facility entirely.

Highest quality across our Bangalore site is ensured by:



Sartorius Stedim India Pvt. Ltd. in Bangalore  
Center of Excellence for Stainless Steel Solutions.

## Excellence in Employee skills

All our engineers and workers are highly qualified. Regular training and qualification sessions ensure highest standards in design and manufacture.

At Sartorius we have exclusively developed a certified welding technique for your high demanding biopharmaceutical applications. Moreover we own welding certifications according to European Standard DIN EN 14 and DIN EN ISO3834-2.





## Certificates & Licenses

Beside global corporate and regular internal audits, certificates from independent testing organisations demonstrate the reliability and high standard of our products. We are DIN EN ISO9001 certified and own manufacturing licenses for pressure vessels according to European, U.S. and Chinese requirements.



## Material Quality & Manufacturing Inspection

At Sartorius we ensure high uniform quality across all our manufacturing sites through e.g.:

- FaroArm coordinate measuring machine to confirm design tolerances in 3D
- Endoscopic inspection of weld seams
- Orbital welding



# 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

Sartorius Stedim Systems GmbH  
Robert-Bosch-Strasse 5 – 7  
34302 Guxhagen

Phone +49.5665.407.0  
Fax +49.5665.407.2200

### France

Sartorius Stedim FMT S.A.S.  
ZI des Paluds  
Avenue de Jouques – CS 91051  
13781 Aubagne Cedex

Phone +33.442.845600  
Fax +33.442.845619

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

Phone +33.442.845600  
Fax +33.442.846545

### Austria

Sartorius Stedim Austria GmbH  
Modcenterstrasse 22  
1030 Vienna

Phone +43.1.7965763.18  
Fax +43.1.796576344

### Belgium

Sartorius Stedim Belgium N.V.  
Rue Colonel Bourg 105  
1030 Bruxelles

Phone +32.2.756.06.80  
Fax +32.2.756.06.81

### Hungary

Sartorius Stedim Hungária Kft.  
Kagyló u. 5  
2092 Budakeszi

Phone +36.23.457.227  
Fax +36.23.457.147

### Italy

Sartorius Stedim Italy S.r.l.  
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.

Phone +31.30.60.25.080  
Fax +31.30.60.25.099

[filtratie.nederland@sartorius-stedim.com](mailto:filtratie.nederland@sartorius-stedim.com)

### Poland

Sartorius Stedim Poland Sp. z o.o.  
ul. Wrzesinska 70  
62-025 Kostrzyn

Phone +48.61.647.38.40  
Fax +48.61.879.25.04

### Russian Federation

LLC "Sartorius Stedim RUS"  
Vasilyevsky Island  
5<sup>th</sup> line 70, Lit. A  
199178 St. Petersburg

Phone +7.812.327.53.27  
Fax +7.812.327.53.23

### Spain

Sartorius Stedim Spain, S.A.U.  
Avda. de la Industria, 32  
Edificio PAYMA  
28108 Alcobendas (Madrid)

Phone +34.913.586.098  
Fax +34.913.589.623

### Switzerland

Sartorius Stedim Switzerland AG  
Ringstrasse 24 a  
8317 Tagelswangen

Phone +41.52.354.36.36  
Fax +41.52.354.36.46

### U.K.

Sartorius Stedim UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9 QQ

Phone +44.1372.737159  
Fax +44.1372.726171

### Ukraine

LLC "Sartorius Stedim RUS"  
Post Box 440 "B"  
01001 Kiev, Ukraine

Phone +380.44.411.4918  
Fax +380.50.623.3162

## Americas

### USA

Sartorius Stedim North America Inc.  
5 Orville Drive, Suite 200  
Bohemia, NY 11716

Toll-Free +1.800.368.7178  
Fax +1.631.254.4253

### Argentina

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

Phone +54.11.4721.0505  
Fax +54.11.4762.2333

### Brazil

Sartorius do Brasil Ltda  
Avenida Senador Vergueiro 2962  
São Bernardo do Campo  
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900  
Fax +55.11.4362.8901

### Mexico

Sartorius de México, S.A. de C.V.  
Libramiento Norte de Tepotzotlan s/n,  
Colonia Barrio Tlacateco,  
Municipio de Tepotzotlan,  
Estado de México,  
C.P. 54605

Phone +52.55.5562.1102  
Fax +52.55.5562.2942

[leadsmex@sartorius.com](mailto:leadsmex@sartorius.com)

### Peru

Sartorius Peru S.A.C.  
Avenue Alberto del Campo 411  
Floor 12 – The Office  
15076 – San Isidro, Lima

Phone +51.1.441 0158  
Fax +51.1.422 6100

## Asia | Pacific

### 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

### China

Sartorius Stedim Biotech (Beijing) Co. Ltd.  
No. 33 Yu'an Road  
Airport Industrial Park Zone B  
Shunyi District, Beijing 101300

Phone +86.10.80426516  
Fax +86.10.80426580

Sartorius Stedim (Shanghai)  
Trading Co., Ltd.

3rd Floor, North Wing, Tower 1  
No. 4560 Jinke Road  
Zhangjiang Hi-Tech Park  
Pudong District  
Shanghai 201210, P.R. China

Phone +86.21.6878.2300  
Fax +86.21.6878.2882

Sartorius Stedim Biotech (Beijing) Co. Ltd.  
Guangzhou Representative Office

Unit K, Building 23  
Huihua Commerce & Trade Building  
No. 80 Xianlie Middle Road  
Guangzhou 510070

Phone +86.20.37618687 | 37618651  
Fax +86.20.37619051

### India

Sartorius Stedim India Pvt. Ltd.  
#69/2-69/3, NH 48, Jakkasandra  
Nelamangala Tq  
562 123 Bangalore, India

Phone +91.80.4350.5250  
Fax +91.80.4350.5253

### Japan

Sartorius Stedim Japan K.K.  
4th Fl., Daiwa Shinagawa North Bldg.  
8-11, Kita-Shinagawa 1-chome  
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.4331.4300  
Fax +81.3.4331.4301

### Malaysia

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

57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622  
Fax +60.3.8996.0755

### Singapore

Sartorius Stedim Singapore Pte. Ltd.  
10 Science Park Rd  
The Alpha #02-13/14  
Singapore Science Park II  
Singapore 117684

Phone +65.6872.3966  
Fax +65.6778.2494

### South Korea

Sartorius Korea Biotech Co., Ltd.  
8th Floor, Solid Space B/D,  
PanGyoYeok-Ro 220, BunDang-Gu  
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700  
Fax +82.31.622.5799



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