



sartorius stedim
biotech

BIOSTAT STR[®] Selected Bibliography



turning science into solutions



BIOSTAT STR[®]

Selected Bibliography

Journal Articles

(2015) Popova, Daria; Stonier, Adam; Pain, David; Titchener-Hooker, Nigel J.; Farid, Suzanne S.:

Representative mammalian cell culture test materials for assessment of primary recovery technologies: a rapid method with industrial applicability.

🔗 In *Biotechnology journal* 10 (1), pp. 162–170. DOI: 10.1002/biot.201400294.

(2015) Wu, Chia-Ying; Lin, Yi-Wen; Kuo, Chia-Ho; Liu, Wan-Hsin; Tai, Hsiu-Fen; Pan, Chien-Hung et al.:

Inactivated Enterovirus 71 Vaccine Produced by 200-L Scale Serum-Free Microcarrier Bioreactor System Provides Cross-Protective Efficacy in Human SCARB2 Transgenic Mouse.

🔗 In *PLoS one* 10 (8), e0136420. DOI: 10.1371/journal.pone.0136420.

(2014) Blaszczyk, Katharina; Löffelholz, Christian; Eibl, Regine; Eibl, Dieter: Combining single-use stirred bioreactor with standard cross-flow technology in biphasic protein production processes at pilot scale.

🔗 In *Eng. Life Sci.* 14 (3), pp. 327–331. DOI: 10.1002/elsc.201300069.

(2014) Boedeker, Berthold: Disposables for Biomanufacturing. A User's Perspective.

🔗 In *Bioprocess International* 12 (Suppl 5).

(2014) Bou, Jordi Ruano; de Wilde, Davy; c: Comparing Multiuse and Single-Use Bioreactors for Virus Production.

🔗 In *Bioprocess International*.

(2014) de Wilde, Davy; Dreher, Thomas; Zahnow, Christian; Husemann, Ute; Greller, Gerhard; Adams, Thorsten; Fenge, Christel:

Superior Scalability of Single-Use Bioreactors. Innovations in Cell Culture.

🔗 In *Bioprocess International Volume 12* (Suppl. 5), p. 14.

(2014) Dreher, Thomas; Walcarius, Bart; Husemann, Ute; Klingenberg, Franziska; Zahnow, Christian; Adams, Thorsten et al.: Microbial high cell density fermentations in a stirred single-use bioreactor.

🔗 In *Advances in biochemical engineering/biotechnology* 138, pp. 127–147. DOI: 10.1007/10_2013_189.

(2014) Löffelholz, Christian; Kaiser, Stephan C.; Kraume, Matthias; Eibl, Regine; Eibl, Dieter: Dynamic Single-Use Bioreactors Used in Modern Liter- and m(3)- Scale Biotechnological Processes: Engineering Characteristics and Scaling Up.

🔗 In *Advances in biochemical engineering/biotechnology* 138, pp. 1–44. DOI: 10.1007/10_2013_187.

(2014) Reglin, Regina; Ruhl, Sebastian; Weyand, Jörg; Wilde, Davy de; Husemann, Ute; Greller, Gerhard; Fenge, Christel:

Verification of New Flexsafe STR Single-Use Bioreactor Bags. Using a CHO Fed-Batch Monoclonal Antibody Production Process at 1,000-L Scale.

🔗 In *Bioprocess International* 12 (Suppl. 5), p. 53.

(2014) Schirmaier, Carmen; Jossen, Valentin; Kaiser, Stephan C.; Jüngerkes, Frank; Brill, Silke; Safavi-Nab, Agnieszka et al.:

Scale-up of adipose tissue-derived mesenchymal stem cell production in stirred single-use bioreactors under low-serum conditions.

🔗 In *Eng. Life Sci.* 14 (3), pp. 292–303. DOI: 10.1002/elsc.201300134.

(2014) Stering, Magnus; Dahlberg, Martin; Adams, Thorsten; de Wilde, Davy; Fenge, Christel:

Pressure Decay Method for Postinstallation Single-Use Bioreactor Bag Testing.

🔗 In *Bioprocess International* 12 (Suppl. 5).

(2014) Weichert, Henry; Lueders, Julia; Weyand, Joerg; Adams, Thorsten; Becker, Mario: Integrated Optical Single-Use Sensors. Moving Toward a True Single-Use Factory for Biologics and Vaccine Production.

🔗 In *Bioprocess International* 12 (Suppl 5).

(2013) Dreher, Thomas; Husemann, Ute; Ruhl, Sebastian; Greller, Gerhard:

Design space definition for a stirred single-use bioreactor family from 50 to 2000 L scale.

🔗 In *BMC Proc* 7 (Suppl 6), P55. DOI: 10.1186/1753-6561-7-S6-P55.

(2013) Dreher, Thomas; Husemann, Ute; Zahnow, Christian; de Wilde, Davy; Adams, Thorsten; Greller, Gerhard:

High Cell Density Escherichia coli Cultivation in Different Single-Use Bioreactor Systems.

🔗 In *Chemie Ingenieur Technik* 85 (1–2), pp. 162–171. DOI: 10.1002/cite.201200122.

(2013) Löffelholz, Christian; Husemann, Ute; Greller, Gerhard; Meusel, Wolfram; Kauling, Jörg; Ay, Peter et al.:

Bioengineering Parameters for Single-Use Bioreactors. Overview and Evaluation of Suitable Methods.

🔗 In *Chemie Ingenieur Technik* 85 (1–2), pp. 40–56. DOI: 10.1002/cite.201200125.

(2013) Morrow, K. John:

Scaleup Issues for Single Use and Steel.

🔗 In *Genetic Engineering & Biotechnology News* 33 (18), 32, 34–35. DOI: 10.1089/gen.33.18.13.

(2013) Weber, Anne; Husemann, Ute; Chaussin, Sébastien; Adams, Thorsten; de Wilde, Davy; Gerighausen Susanne et al.: Development and Qualification of a Scalable, Disposable Bioreactor for GMP-Compliant Cell Culture.

🔗 In *Bioprocess International* 11 (Suppl. 2).

(2012) Kusserow, Arne:

Von Klein bis Groß: Fermentationstechnologie entlang der Prozesskette. Interview mit Dr. C. Fenge.

🔗 In *GIT Labor-Fachzeitschrift*, pp. 228–231.

(2010) Dessoy, Sandrine; Battisti, Raphael; Ghislain, Yves; Barbier, Benoit; Gerkens, Pascal; Peeters, Ludovic; Chaubard, Jean-François: Disposable Bioreactors for Viral Vaccine Production: Challenges and Opportunities.

🔗 In *BioPharm International Supplements*.

(2010) de Wilde, Davy; Adams, Thorsten: Improving scalability and comparability by using single-use bioreactors for large-scale manufacturing.

In *European Journal of Parenteral & Pharmaceutical Sciences Volume 15* (Part 2), pp. 41–46.

(2010) Stadler, Ernest L.:

An Exhausting Solution for Fermentors.

In *Pharmaceutical Engineering* 30 (3).

(2010) Ullah, Millie:

Benefits of Disposable Bioreactor Solutions. Novel Configurations Feature Control Capabilities Similar to Traditional Systems.

🔗 In *Genetic Engineering News* 30 (14).

(2010) Ullah, Millie:

Optimizing upstream processes to single use.

Testing a new design for comparability. In *Bioprocess International* 8 (April Suppl).

(2009) de Wilde, Davy; Noack, Ute;

Kahlert, Wolfgang; Barbaroux, Magali: Bridging the gap from reusable to single-use manufacturing with stirred, single-use bioreactors.

🔗 In *Bioprocess International Volume 7* Suppl 4, 46–41.

Conference Presentations

(2016) Bou, Jordi Ruano:

Virus production with single-use bioreactors (SUB).
HIPRA.
Upstream Technology Forum. Göttingen,
Germany, 6/8/2016.

(2016) Toueille, Magali:

Pre-Industrial Manufacturing of Lentiviral Vectors
by Transient Transfection in Single Use Systems.
Genethon.
Cell Therapy Manufacturing & Gene Therapy
Congress. Brussels, Belgium, 2/4/2016.

(2015) Fenge, Christel; Carpio, Melisa:

Continuous culture in the age of single-use.
Integrated Continuous Biomanufacturing II.
Berkeley, USA, 11/2/2015.

(2015) Greller, Gerhard:

Scale up of a CHO fed-batch process from 250 mL
to 2000 L in single-use bioreactors.
ACHEMA 2015. Frankfurt a. M., 6/9/2015.

(2014) Garcia, Lidia:

Veterinary vaccine production using SU
bioreactors: Scale up study from laboratory &
PD up to commercial production.
Zoetis Manufacturing & Research Spain.
Upstream Technology Forum. Göttingen,
Germany, 9/9/2014.

(2014) Greller, Gerhard:

Superior Scalability of Single-Use Bioreactors.
ACHEMA 2014. Frankfurt a. M., 9/9/2014.

(2013) Chaubard, Jean-François:

Single-Use Bioreactors: how to integrate them into
your Process Development platform. GSK Vaccines.
Upstream Technology Forum. Göttingen,
Germany, 9/9/2013.

(2012) Peeters, Ludovic:

Implementation of Single-Use Bioreactors for
Scale-Up of Vaccine Processes. GSK Vaccines.
Upstream Technology Forum. Basel, Switzerland,
5/10/2012.



Poster Presentations

(2016) Ruhl, Sebastian; Husemann, Ute; Zahnow, Christian; Zehe, Christoph; Greller, Gerhard:

Establishing a new CHO Fed-Batch Process in a Bioreactor Family from Single-Use Process Development to Production Scale. BioTech, Single-Use Technology in Biopharmaceutical Manufacturing, Wädenswil, Switzerland.

(2016) Zahnow, Christian; Bargh, Neil; Dreher, Thomas; Stacey, Adrian J.; Ruhl, Sebastian; Husemann, Ute; Greller, Gerhard:

Superior Scalability of a Single-Use Bioreactor Family from 0.25 to 2000 L. BioTech, Single-Use Technology in Biopharmaceutical Manufacturing, Wädenswil, Switzerland.

(2015) Ruhl, Sebastian; Husemann, Ute; Dreher, Thomas; Zahnow, Christian; Greller, Gerhard:
Challenge of Process Transfer from Single-Use Production to a Process Development Bioreactor. European Society for Animal Cell Technology (ESACT) Meeting, Barcelona, Spain.

(2013) Garcia, Lidia; Mouriño, Mercedes; Urniza, Alicia:

Comparative study of bluetongue virus serotype 8 production on BHK-21 cells in a 50L BIOSTAT® STR single-use bioreactors vs roller bottles. Zoetis Manufacturing & Research Spain, S.L Pfizer Olot S.L.U.

(2013) Husemann, Ute; Dreher, Thomas; Greller, Gerhard:

Design space definition for a stirred single-use bioreactor family from 50 to 2000 L scale. Sartorius Stedim Biotech. European Society for Animal Cell Technology (ESACT) Meeting, Lille, France.

(2013) Husemann, Ute; Ruhl, Sebastian; Jurkiewicz, Elke; Dreher, Thomas; Greller, Gerhard:

The successful transfer of a modern CHO fed-batch process to different single-use bioreactors. Sartorius Stedim Biotech. European Society for Animal Cell Technology (ESACT) Meeting, Lille, France.

(2013) Jossen, Valentin; Kaiser, Stephan C.; Schirmaier, Carmen; van den Bos, Christian; Eibl, Dieter; Eibl, Regine:

Investigations of fluid flow and microcarrier suspension for hMSC expansion in single-use stirred bioreactors. BioTech, Single-Use Technology in Biopharmaceutical Manufacturing, Wädenswil, Switzerland.

(2013) Reglin, Regina; Husemann, Ute; Dreher, Thomas; Greller, Gerhard:

Scale-up of a modern CHO process to a 1000 L stirred single-use bioreactor. Sartorius Stedim Biotech GmbH. Upstream Technology Forum, Göttingen, Germany.

(2013) Schirmaier, Carmen; Kaiser, Stephan C.; Jossen, Valentin; Brill, Silke; Jüngerkes, Frank; van den Bos, Christian et al.:

Expansion of mesenchymal adipose-tissue derived stem cells in a stirred single-use bioreactor under low-serum conditions. 23rd European Society for Animal Cell Technology (ESACT) Meeting: Better Cells for Better Health.

(2013) Steiger, Nina; Imseng, Nicole; Schirmaier, Carmen; Jurkiewicz, Elke; Husemann, Ute; Eibl, Regine; Greller, Gerhard:
Towards a standardized biological test for the identification of cell growth influencing leachables. Zurich University of Applied Sciences, Sartorius Stedim Biotech GmbH. European Society for Animal Cell Technology (ESACT) Meeting, Lille, France.

(2012) Dreher, Thomas; Zahnow, Christian; Husemann, Ute; Greller, Gerhard:
Towards a complete single-use upstream process for Escherichia coli high cell density fermentations. Sartorius Stedim Biotech GmbH. BioProScale, Berlin, Germany.

(2011) Hundt, Boris; Molle, Noreen; Stefaniak, Sabine; Durrwald, Ralf; Weyand, Joerg:
Large pilot scale cultivation process study of adherent MDBK cells for porcine Influenza A virus propagation using a novel disposable stirred-tank bioreactor. IDT Biologika GmbH, Sartorius Stedim Biotech GmbH.

(2011) Zijlstra, Gerben; Noack, Ute; Weisshaar, Stefan; de Wilde, Davy; Greller, Gerhard; Fenge, Christel; Douwenga, Rolf:
High cell density XD® cultivation of CHO cells in the BIOSTAT® CultiBag STR 50 L single use bioreactor with novel micro-sparger and single-use exhaust cooler. DSM Biologics; Sartorius Stedim Biotech GmbH; Sartorius Stedim Belgium NV. European Society for Animal Cell Technology (ESACT) Meeting, Vienna, Austria.

(2010) Noack, Ute; Brandt, Sophie; de Wilde, Davy; Greller, Gerhard:
Evaluation of the BIOSTAT® CultiBag STR family in terms of scalability. Sartorius Stedim Biotech. BioTech, Single-Use Technology in Biopharmaceutical Manufacturing, Wädenswil, Switzerland.



Book Chapters

(2015) Kaiser, Stephan C.; Kraume, Matthias; Eibl, Dieter; Eibl, Regine:

Single-Use Bioreactors for Animal and Human Cells.

🔗 In Mohamed Al-Rubeai (Ed.): *Animal Cell Culture*. Cham: Springer International Publishing, pp. 445–500.

(2014) Jossen, Valentin; Pörtner, Ralf; Kaiser, Stephan C.; Kraume, Matthias; Eibl, Dieter; Eibl, Regine:

Mass Production of Mesenchymal Stem Cells – Impact of Bioreactor Design and Flow Conditions on Proliferation and Differentiation.

🔗 In Daniel Eberli (Ed.): *Cells and biomaterials in regenerative medicine*. [Elektronische Ressource]. Rijeka, Croatia: InTech Europe.

(2012) Greller, Gerhard; Noack, Ute:

Evaluation of Disposable Bioreactor Design Compared to Reusable Stainless Steel Reactors.

🔗 In Nigel Jenkins, Niall Barron, Paula Alves (Eds.): *Proceedings of the 21st Annual Meeting of the European Society for Animal Cell Technology (ESACT)*, Dublin, Ireland, June 7–10, 2009. Dordrecht: Springer Netherlands, pp. 359–362.

(2011) Löffelholz, Christian; Kaiser, Stephan C.; Werner, Sören; Eibl, Dieter:

CFD as a Tool to Characterize Single-Use Bioreactors.

🔗 In Regine Eibl, Dieter Eibl (Eds.): *Single-use technology in biopharmaceutical manufacture*. [Online-Ausg.]. Chichester: Wiley-Blackwell, pp. 263–279.

(2011) Noack, Ute; Wilde, Davy de;

Verhoeve, Francis; Balbirnie, Edward;

Kahlert, Wolfgang; Adams, Thorsten et al.:

Single-Use Stirred Tank Reactor BIOSTAT® CultiBag STR: Characterization and Applications.

🔗 In Regine Eibl, Dieter Eibl (Eds.): *Single-use technology in biopharmaceutical manufacture*. [Online-Ausg.]. Chichester: Wiley-Blackwell, pp. 225–240.

(2009) Eibl, Regine; Catapano, Gerardo;

Czermak, Peter; Eibl, Dieter; Pörtner, Ralf (Eds.): *Cell and Tissue Reaction Engineering*.

🔗 With a Contribution by Martin Fussenegger and Wilfried Weber. Berlin, Heidelberg: Springer (Principles and Practice).

Application Notes

(2015) Mueller, Miriam; Ruhl, Sebastian;

Noack, Ute; Greller, Gerhard; de Wilde, Davy;

Adams, Thorsten:

Large scale cultivation of CHO cells in the stirred single-use bioreactor BIOSTAT® CultiBag STR.

🔗 Sartorius Stedim Biotech.

(2013) Fenge, Christel; Weyand, Joerg;

Greller, Gerhard; Adams, Thorsten:

Large-Scale Perfusion and Concentrated Fed-Batch Operation of BIOSTAT® STR Single-Use Bioreactor.

🔗 Sartorius Stedim Biotech.

(2011) Dreher, Thomas; Husemann, Ute;

Greller, Gerhard; Weyand, Joerg:

High cell density Escherichia coli cultivation in the stirred single use bioreactor BIOSTAT® CultiBag STR 50 plus MO.

🔗 Sartorius Stedim Biotech.





Sales and Service Contacts

For further contacts, visit 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
Modectcenterstrasse 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

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
5th 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. Avalos 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

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