



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



Application  
Note

#04

#05

#06

#07

#08

By Christel Fenge, Jörg Weyand,  
Gerhard Greller, Thorsten Adams  
Sartorius Stedim Biotech

### Introduction

Perfusion is a well known production approach since many years for a number of biological drugs such as coagulation factors (Factor VIII: Kogenate, Bayer and Refacto, Pfizer; Factor VII: Novoseven, Novo Nordisk), Protein C (Xigris, Eli Lilly) and other enzymes but also monoclonal antibodies (ReoPro and Remicade, both Janssen Biotech; Simulect, Novartis, etc.). Recently, perfusion has been used also to generate high cell density seed cultures to reduce the number of seed steps necessary to reach final production scale [1, 2].

Through the higher cell densities and titers achieved in concentrated fed-batch and perfusion cultures typically smaller culture volumes and bioreactor scales are required to produce the therapeutic protein or antibody. This is a great advantage as the footprint of a production facility can be reduced and scale-up issues are mitigated due to the marginal scale-up factor between clinical and commercial production. Intensified cell culture processes are especially beneficial in the context of single-use facilities as they provide production capacities at 1000L scale that in the past were only achievable with five to ten times larger bioreactors [3, 4]. Furthermore, modern high-end cell culture processes aim to maintain the cells in a defined metabolic state in order to ensure stable product quality through controlling protein folding and glycosylation. In this case, the main aim is not necessarily to reach very high cell densities, but to ensure a steady state of nutrients and metabolites in the bioreactor.

### How to perform concentrated fed-batch or perfusion operation

After inoculation of the bioreactor and an initial 1 – 2 day batch growth phase, the removal of cell free supernatant e.g. with the Refine ATF system is started at a constant harvest flow rate. At the same time, the culture is replenished with fresh medium.

A number of alternative cell retention devices are available such as continuous centrifuges (Centritech Cell from Pneumatic Scale Angelus or kSep from KBI), internal and external spinfilters, settlers, hollow fiber and other membrane based retention systems.

When applying single-use bioreactors such as the BIOSTAT® STR, the addition is controlled via a feed pump that receives a signal from load cells or a platform balance maintaining a defined bioreactor weight. As the cell density grows and the nutrient consumption and metabolite formation increases, the harvest rate is subsequently increased to maintain a certain exchange rate of fresh medium per cell or alternatively a given medium exchange rate per day [2]. On-line biomass measurement, e.g. with the BioPAT® ViaMass probe that will soon be available for single-use BIOSTAT® STR and RM bags, provides an automated option to control the perfusion rate based on cell density. Using at-line glucose and lactate measurement, e.g. with the BioPAT® Trace, an additional concentrated feed can be applied to control the glucose concentration.

Figure 1 provides a schematic depiction of a typical concentrated perfusion or fed-batch set-up based on the BIOSTAT® STR.

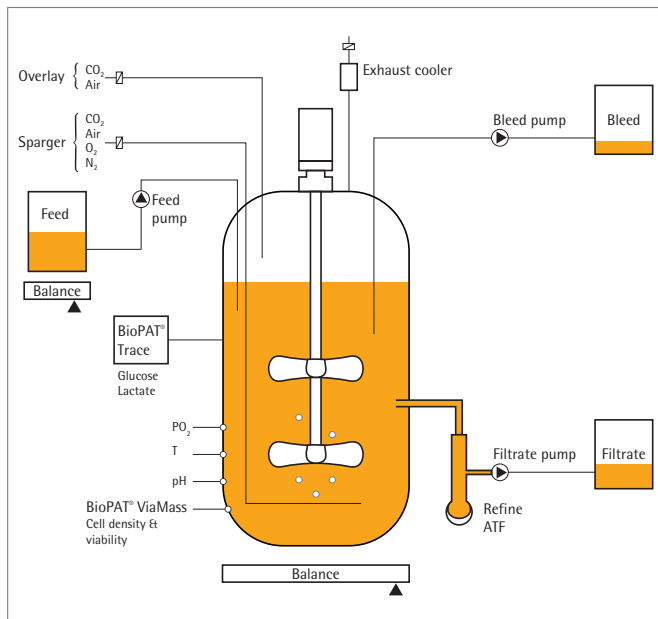


Fig. 1: Set-up of a concentrated fed-batch using the single-use Bioreactor BIOSTAT® STR.

### Key considerations

Typical perfusion rates are in the range of 1–2 bioreactor volumes per day. Applying a small cell bleed stream enables the establishment of a defined cell growth rate and by that a high viability can be maintained which in turn mitigates clogging of the cell retention membrane, either the product is recovered in the cell free harvest (concentrated perfusion) or in the bioreactor content (concentrated fed-batch). As most antibodies are rather stable, concentrated fed-batch with accumulation of the product in the bioreactor is a simple and straight forward approach to increase space time yields of a given facility. Concentrated perfusion is the method of choice for recombinant proteins that in many cases are prone to degradation or might show feedback inhibition and should therefore be removed from the cell culture into a chilled harvest tank and subsequently purified.

### Media logistics

In large-scale continuous processing, media and harvest logistics needs specific attention. Dependent on the bioreactor scale and chosen perfusion rate, 500 L to 2000 L and more fresh medium have to be provided daily. The FlexAct® MP system is an easy to use single-use system for automated media preparation (Fig. 2) especially designed to reduce effort and increase convenience associated with large-scale continuous operation. Its design encompasses dust free transfer of powder media into a mixing bag, a single use sensor for pH measurement and adjustment and a controller that manages automatic sterile filtration and transfer into a storage bag. In case of perfusion, equal volumes have to be collected and processed in subsequent purification steps. Our range of chilled Palletank provide a proven single-use option for storage of large harvest volumes in disposable bags. In summary, the FlexAct® MP system together with Palletank and Flexel® bags manage the challenging fluid handling of large scale, single-use continuous processing.



Fig. 2: FlexAct® MP – Disposable solution for media preparation.

### Scaling up continuous processing

Concentrated fed-batch and perfusion processes can be developed and successfully implemented at production scale using Sartorius stirred tank single-use bioreactors in combination with different sizes of the Refine ATF module. At the 2L bench scale, e.g. our UniVessel® in combination with our BIOSTAT® B or B-DCU controller provides a fully scalable development system. Subsequent scale-up from 50L and 200L can be achieved to the 1000L scale in the BIOSTAT® STR. At large-scale, the ATF modules might be connected via side ports of the single-use bioreactor bag using up to two 1" sterile connectors and operated in an external loop of the bioreactor (Fig. 7). It is critical that this external loop is as short as possible to avoid that the cell culture is exposed to uncontrolled conditions, e.g. different temperature and potential oxygen limitations.

### Single-use bioreactor configurations suitable for intensified cell cultures

Key to successful concentrated fed-batch and perfusion operation is an efficient aeration system that provides  $k_L a$  values above  $10 - 15 \text{ h}^{-1}$  to supply the culture with sufficient oxygen (Fig. 3). At the same time, excessive  $\text{CO}_2$  is formed in the intensified culture which needs to be removed to avoid any inhibitory effect on productivity or even product quality.

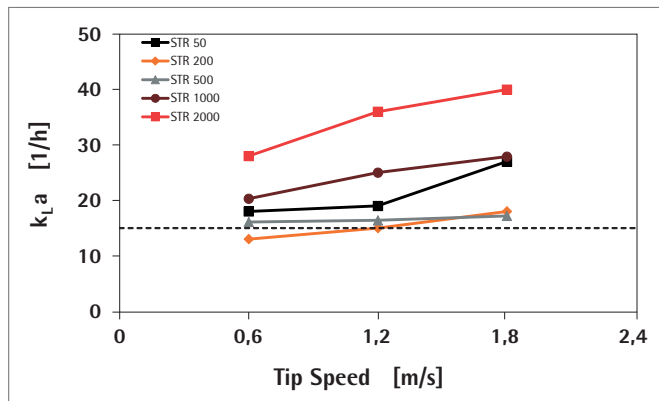


Fig. 3:  $k_L a$  values determined in different single-use bioreactor bag volumes of the BIOSTAT® STR, equipped with  $2 \times 3$  blade segment impellers, using the gassing out method in phosphate buffered saline, aeration rate  $0.1 \text{ vvm}$ ,  $150 \mu\text{m}$  holes of combisparger, temperature  $25^\circ\text{C}$ .

This can be achieved with the Combisparger that microsparges compressed air or pure oxygen through defined  $150 \mu\text{m}$  holes and provides a stripping gas flow through  $0.8 \text{ mm}$  holes at the same time (Fig. 4). This single-use sparger design emulates a successful aeration strategy applied since many years in conventional stainless steel bioreactors.



Fig. 4: Combisparger with  $150 \mu\text{m}$  defined micro-holes and  $0.8 \text{ mm}$  holes.

A problem that should not be underestimated is excessive aerosol formation in the exhaust gas due to high gas flow rates and high protein content in the concentrated cell cultures. A specifically developed single-use exhaust cooler design based on the well-known principle of plate heat exchangers (Fig. 5) mitigates the risk of blocked filters and increases process reliability dramatically [6]. Additional safety locks in the bioreactor control software prevent bioreactor overflow in case of clogging of the cell retention device. As a worst case safety lock, all feed pumps and gas flows are interrupted if the pressure in the bioreactor exceeds the maximum defined operating pressure.



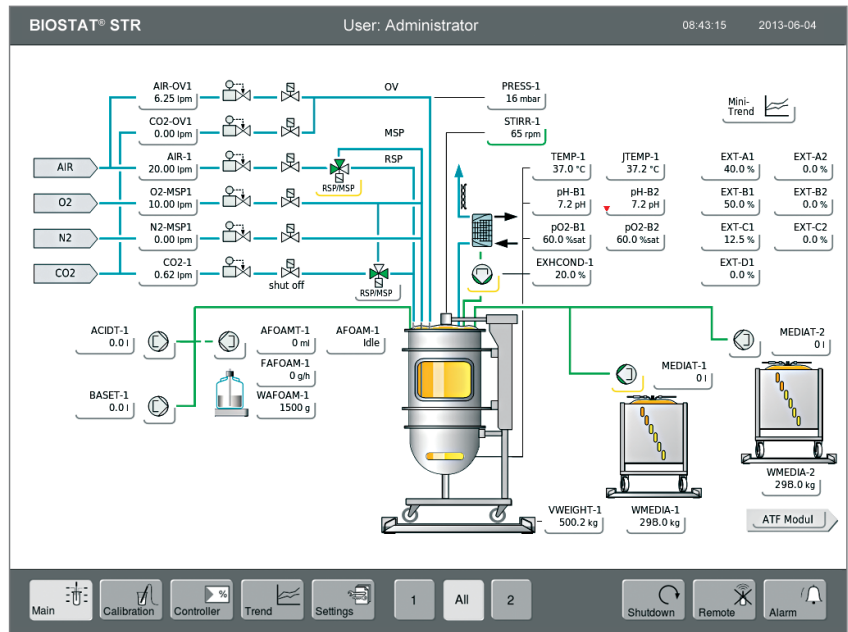
Fig. 5: Exhaust Cooler

**Conclusion**

Modern single-use bioreactor designs such as of the BIOSTAT® STR allow advanced, intensified cultivation strategies whilst providing tools to mitigate operational risks associated to a complex bioprocessing strategy and thus enabling robust single-use production for clinical trials and commercial drug manufacturing.

**References**

1. R. Heidemann et al., Cytotechnol. 38 (1-3), 99-108 (2002)
2. T. Adams et al., BioPharm Int. Supplement May 2011, S4-S11
3. J. Lim, et al., BioPharm Int. 24 (2), 54-60 (2011)
4. J. Pollock et al., Biotechnol. and Bioeng. 110 (1), 206-219 (2013)
5. C. Fenge et al., in "Animal cell technology: Development, processes and products", Butterworth-Heinemann, 365-370 (1992)
6. Zijlstra, G. et al., poster presented on 22. ESACT meeting, Vienna (2011)



**Fig. 6:** Screenshot of BIOSTAT® STR controller configured for perfusion operation using ATF module.



**Fig. 7:** BIOSTAT® STR 200L with connected Refine ATF system

# 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

Sartorius Stedim Systems GmbH  
Robert-Bosch-Strasse 5-7  
34302 Guxhagen  
Phone +49.5665.407.0

**France**  
Sartorius Stedim FMT S.A.S.  
ZI des Paluds  
Avenue de Jouques - CS 91051  
13781 Aubagne Cedex  
Phone +33.442.845600

Sartorius Stedim France SAS  
ZI des Paluds  
Avenue de Jouques - CS 71058  
13781 Aubagne Cedex  
Phone +33.442.845600

**Austria**  
Sartorius Stedim Austria GmbH  
Modecenterstrasse 22  
1030 Vienna  
Phone +43.1.7965763.18

**Belgium**  
Sartorius Stedim Belgium N.V.  
Rue Colonel Bourg 105  
1030 Bruxelles  
Phone +32.2.756.06.80

**Hungary**  
Sartorius Stedim Hungária Kft.  
Kagyló u. 5  
2092 Budakeszi  
Phone +36.23.457.227

**Italy**  
Sartorius Stedim Italy S.r.l.  
Via dell'Antella, 76/A  
50012 Antella-Bagno a Ripoli (FI)  
Phone +39.055.63.40.41

**Netherlands**  
Sartorius Stedim Netherlands B.V.  
Phone +31.30.60.25.080  
[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

**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

**Spain**  
Sartorius Stedim Spain, S.A.U.  
Avda. de la Industria, 32  
Edificio PAYMA  
28108 Alcobendas (Madrid)  
Phone +34.913.586.098

**Switzerland**  
Sartorius Stedim Switzerland AG  
Ringstrasse 24 a  
8317 Tagelswangen  
Phone +41.52.354.36.36

**U.K.**  
Sartorius Stedim UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9 QQ  
Phone +44.1372.737159

**Ukraine**  
LLC "Sartorius Stedim RUS"  
Post Box 440 "B"  
01001 Kiev, Ukraine  
Phone +380.44.411.4918

## Americas

**USA**  
Sartorius Stedim North America Inc.  
5 Orville Drive, Suite 200  
Bohemia, NY 11716  
Toll-Free +1.800.368.7178

**Argentina**  
Sartorius Argentina S.A.  
Int. A. Ávalos 4251  
B1605ECS Munro  
Buenos Aires  
Phone +54.11.4721.0505

**Brazil**  
Sartorius do Brasil Ltda  
Avenida Senador Vergueiro 2962  
São Bernardo do Campo  
CEP 09600-000 - SP- Brasil  
Phone +55.11.4362.8900

**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  
[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

## Asia | Pacific

**Australia**  
Sartorius Stedim Australia Pty. Ltd.  
Unit 5, 7-11 Rodeo Drive  
Dandenong South Vic 3175  
Phone +61.3.8762.1800

**China**  
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

Sartorius Stedim (Shanghai)  
Trading Co., Ltd.  
Beijing Branch Office  
No. 33 Yu'an Road  
Airport Industrial Park Zone B  
Shunyi District, Beijing 101300  
Phone +86.10.8042.6501

Sartorius Stedim (Shanghai)  
Trading Co., Ltd.  
Guangzhou Branch Office  
Room 1105  
Xing Guang Ying Jing Building  
No. 119, Shui Yin Road  
Yue Xiu District, Guangzhou 510075  
Phone +86.20.3836.4193

**India**  
Sartorius Stedim India Pvt. Ltd.  
#69/2-69/3, NH 48, Jakkasandra  
Nelamangala Tq  
562 123 Bangalore, India  
Phone +91.80.4350.5250

**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

**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

**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

**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



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