

## BioPAT<sup>®</sup> Trace Glucose PID control



#05

Application  
Note

#06

#07

#08

#09

Ajith T George, Anil Kumar Rathod, Dr. Ashok Mundrigi

Sartorius Stedim Biotech, Bangalore, India

Dr. Andreas Prediger, Dr. Stuart Tindal

Sartorius Stedim Biotech GmbH, Goettingen, Germany

## 1. Introduction

Current industry practices for large-scale mammalian cell cultures typically employ a standard platform fed-batch process with fixed volume bolus feeding. Although widely used, these processes are unable to respond to actual nutrient consumption demands from the culture, which can result in accumulation of by-products and depletion of other certain nutrients. This work demonstrates the application of automated glucose control via dynamic feeding with BioPAT® MFCS PID controller. The method is based upon automated glucose measurements obtained from The BioPAT® Trace where cultures were fed to maintain a previously manual achieved target glucose level of 6.0 g/L.

## 2. Material and Methods:

The automated feeding trials were performed in CHO fed batch mode cultivation with following batch parameters:

### Batch Parameters

Cultivation vessel	Biostat® B-DCU II UniVessel® 1L
Temperature set point	36.8°C
pO <sub>2</sub>	Set point 60% (controlled with Air and oxygen)
pH	Set point 7.20 (controlled with CO <sub>2</sub> and 1 M (Na <sub>2</sub> CO <sub>3</sub> , sodium carbonate)
Cultivation media	ActiCHO PM
Feeding quantity	ActiCHO Feed A – 40 mL/L daily on 3 <sup>rd</sup> day onwards ActiCHO Feed B – 4 mL/L on 3 <sup>rd</sup> day onwards
Batch volume	700 mL
Initial cell concentration	0.32 × 10 <sup>6</sup> Cells/mL
Speed	200 rpm – 250 rpm



The BioPAT® Trace system consists of a measurement unit, single-use fluidic system, calibration solutions & transport buffer. The fluidic system assembled with measurement unit was used for calibration and online measurements. The software 'trace\_mon' version 1.3.03 was used to control the device. Calibration solutions with 10 g/L glucose & 5 g/L lactate and 0.5 g/L glucose & 0.25 g/L lactate were used. The BioPAT® Trace system was connected to the bioreactors with a dialysis probe.

The glucose & lactate measurement frequency was set to 20 minutes. Offline reference samples were taken from the bioreactor daily. An internal measurement correction was performed, if there was a deviation from external reference method. The external referencing was performed using spectrophotometric assay methods. For glucose, a test kit from Agappe Diagnostics Ltd. based on glucose oxidase was used. For lactate a test kit from Centronic GmbH based on lactate oxidase was used.

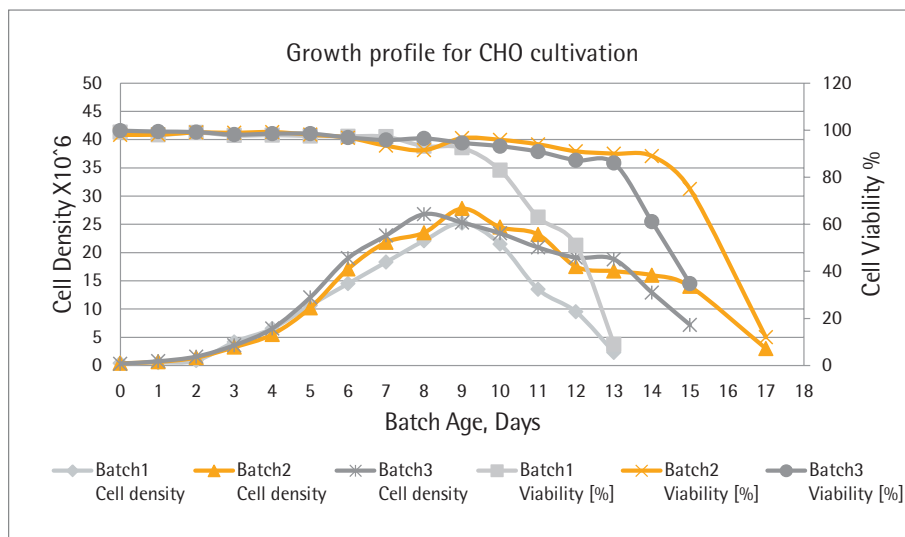
Initial setup trials with water & glucose solution were carried out to optimize the BioPAT® MFCS PID values with different reactor volumes. The optimized PID set up for automatic feeding is shown in Table 1.

Batch Volume	XP	TI	TD
5 L	200	1000	0
2.5 L	100	1200	0
1 L	100	1200	0

**Table 1:** BioPAT® MFCS PID parameters used for feed control with 20 minute glucose sample frequency

### 3. Experiments and Results:

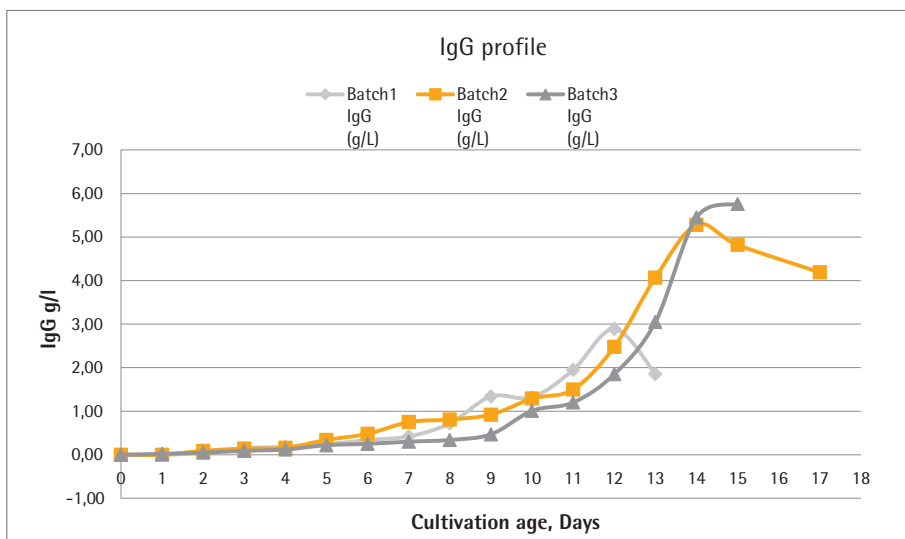
Three replicate CHO cell cultivations in fed batch mode were performed where glucose of 6.0 g/L was maintained with PID & lactate concentrations were monitored with a BioPAT® Trace. During the setup of glucose control loop, feed line was primed and the pump was evaluated for its total flow rate (rpm) capabilities. The starting settings were adjusted to ensure no initial over feeding. An overview of the achieved cell densities and viabilities are presented in figure 2.



**Figure 2:** Overview of the growth patterns of the performed CHO cell cultivations.

During the cultivations, peak cell densities of 25 – 27 million cells/mL were achieved on 8 – 9<sup>th</sup> day of cultivation. During that period the cultivations maintained a viability of over 95%. Afterwards both viability and viable cell density decreased until the end of the cultivation.

Throughout the cultivations, the performance of the batches was monitored by off-line sampling of the IgG media concentration. This data is plotted in figure 3 and shows consistent protein production.



**Figure 3:** Overview of the IgG production of the performed CHO cultivations.

Two different feeding strategies were used for the cultivations and are described below.

Feeding strategy for batch 1 & 2 was

- A constant glucose level of 6.0 g/L was maintained with PID controller from 3<sup>rd</sup> day (end of lag phase) to 9<sup>th</sup> day (peak cell density) by Feed A solution. From 9<sup>th</sup> day onwards, 40% glucose solution replaces feed A in PID controller & 40 mL/L of feed A was added as bolus dosage.

Feeding strategy for batch 3 changed to

- A constant glucose level of 6.0 g/L was maintained with PID controller from 3<sup>rd</sup> day (end of lag phase) to 9<sup>th</sup> day (peak cell density) by Feed A solution. From 9<sup>th</sup> day onwards 40% glucose solution replaces feed A in PID controller & 40 mL/L of feed A was added over a time period of 24 hrs.

In summary, a guided data plot from the BioPAT® Trace and off-line reference measurement points is shown in figure 4.

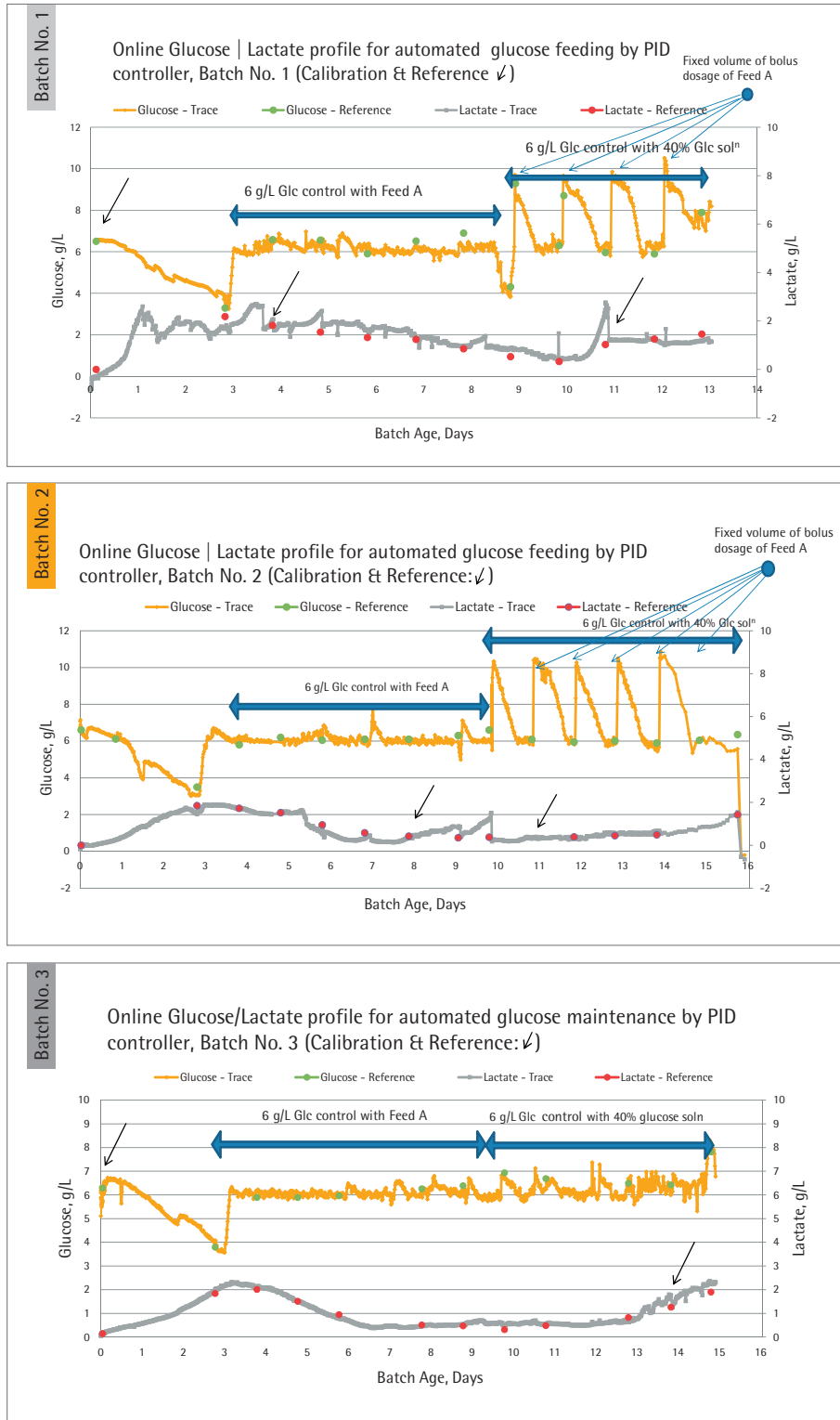
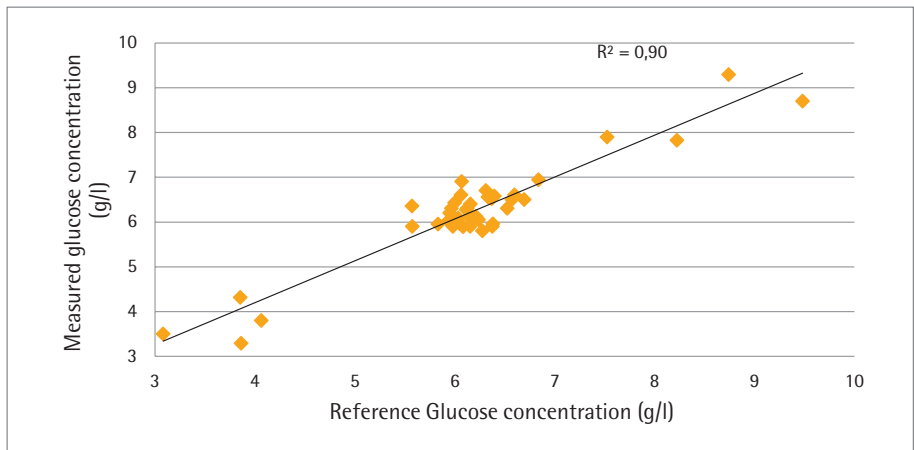


Figure 4: Graph of online values for glucose & lactate for 3 batches & external reference values

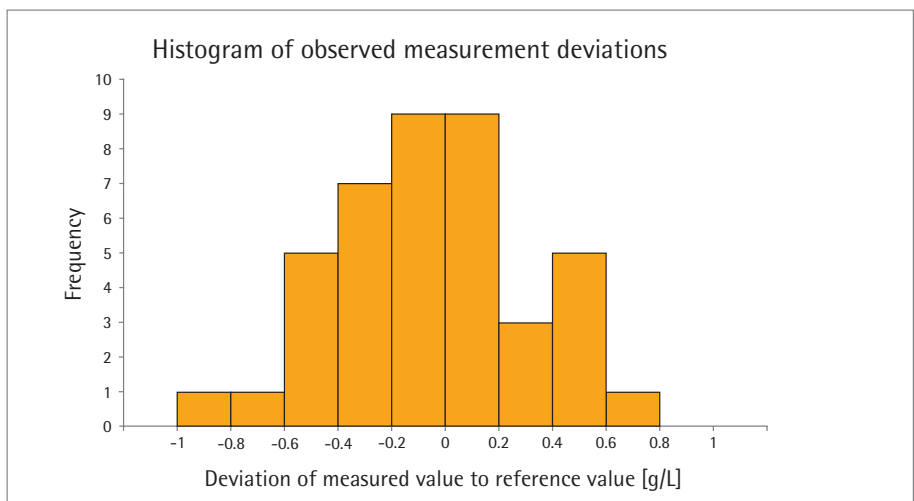
With the optimized PID values, the constant glucose level was maintained with feed A up to 9th day (Peak cell density day). The average consumption of the feed A per day was 27 – 30 mL for 0.75 L of working volume to maintain this constant glucose level. Afterwards 40% glucose solution replaced feed A in PID to maintain the glucose level. Approximately 4 – 5 mL of 40% glucose solution was consumed daily to maintain the set-point.

A calibration and referencing procedure of the BioPAT® Trace was performed at the start of the cultivations and during cultivation when the estimation of the device deviated from the reference measurements (indicated with black arrows in figure 4). Any referencing is needed if any basic process parameters are changed like temperature, feeding profiles or measurement ranges of the system. In figure 5 the measured concentration values from the BioPAT® Trace are compared to the reference values and the accuracy displayed.



**Figure 5:** Measured values from BioPAT® Trace vs values from reference method.

Shown are only values inside the calibration range between 1 – 10 g/L. It shows that the values generated by the BioPAT® Trace correlate well with the off-line glucose concentration measurement. A histogram of the observed deviations of the BioPAT® Trace measurements from the reference values is presented in figure 6.



**Figure 6:** Histogram of the observed deviations of the BioPAT® Trace measurements from the reference values.

It can be seen that the observed deviations are normally distributed. The standard deviation of the observed deviations is 0.36 g/L and all deviations are within a range of  $\pm 1$  g/L.

#### **4. Conclusion**

The BioPAT® Trace ensures a high degree of accurate results for online glucose measurements and dynamic feed control. The standard deviation of the observed error of the BioPAT® Trace values in comparison to the reference values is 0.36 g/L.

BioPAT® MFCS PID controls the glucose level of 6.0 g/L throughout the cultivation very efficiently. The dynamic feeding has the ability to automatically adjust feed rates according to cultures dynamic behavior.

High levels of glucose can result in high levels of lactate through glycolysis. Lactate accumulation can reduce the pH and the low pH can be detrimental to cell viability and productivity.

Traditionally, the control of glucose is being done by taking samples and measuring the sample with an external system such as spectrophotometry. This is a time consuming and operator dependent task that has inherent manual inconsistency.

The BioPAT® Trace is automated handle free equipment after installing the consumables, which provides actual monitoring of glucose level & 100% direct control on the entire cell cultivation process.

# 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  
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.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.  
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"  
Uralskaya str. 4, Lit. B  
199155 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 +34.902.110.935  
Fax +34.91.358.96.23

**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 "Biohit"  
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.  
Circuito Circunvalación Poniente  
No. 149  
Ciudad Satélite  
53100, Estado de México  
México  
Phone +52.5555.62.1102  
Fax +52.5555.62.2942

## 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.  
1 Science Park Road,  
The Capricorn, #05-08A,  
Singapore Science Park II  
Singapore 117528  
Phone +65.6872.3966  
Fax +65.6872.2494

**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

