

Digitalize Your Bioprocess Control Training with BIOSTAT[®] T CHO



turning science into solutions



Virtual Bioreactor Training Tool

00:00:00 05/25/2017 13:17:55

CS - History

The BIOSTAT[®] T CHO is an interactive software training tool and ideal for the education of students and operators in bioprocess control. It enables users to learn about bioprocess engineering and to practice with menu navigation and control of a bioreactor prior to operating an actual bioreactor.

Integrative Learning

Theoretical principles of bioprocess engineering can immediately be applied and verified by using the BIOSTAT[®] T, ensuring that the principles are really understood and consolidated in the long term.

Close to Reality

The BIOSTAT[®] T CHO is based on the human machine interface of a BIOSTAT STR[®] 200 L single-use bioreactor and contains real cultivation data. As a result, users of the software can be trained realistically ensuring that they are ideally prepared for the actual operation of a bioreactor.

Quality in Hands-On Operation

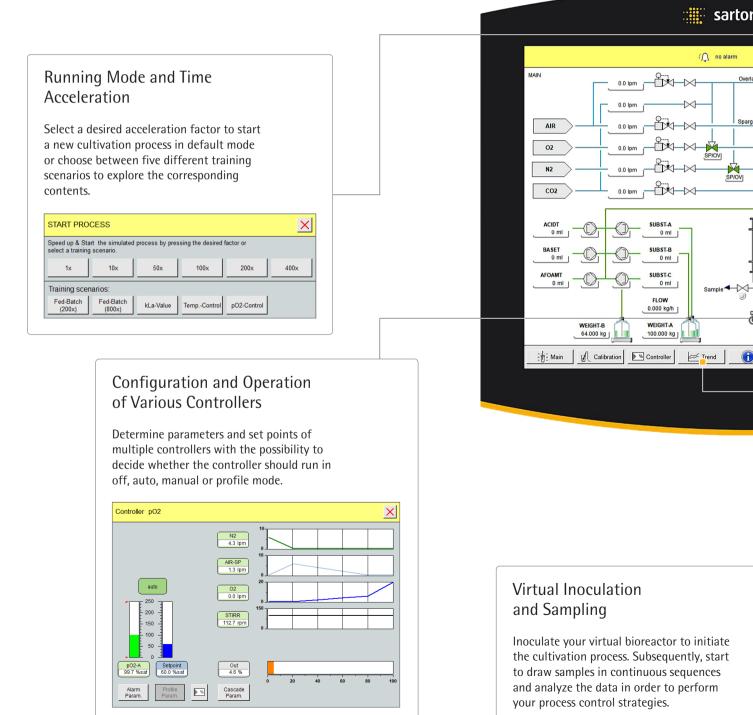
Users of the BIOSTAT[®] T CHO can be introduced to the operation of a bioreactor completely virtually. This enables them to achieve high quality results when using a bioreactor in real life from the very beginning.

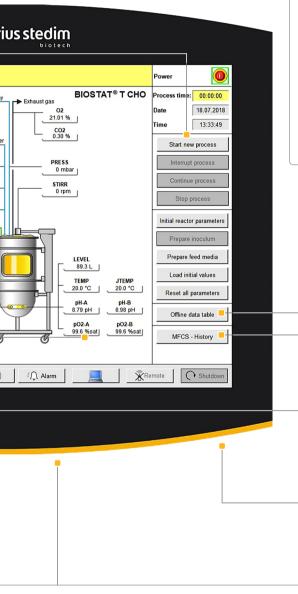
Time and Cost Saving

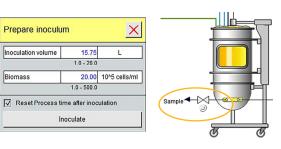
The virtual bioreactor training tool is free-to-use and requires only a computer to be used. Virtual bioprocess control trainings allow the risk mitigation of human error and the prevention of batch losses. Time- and cost-intensive activities can be minimized as using the BIOSTAT[®] T CHO does not require any laboratory clean-ups or procurements of microorganisms, cell cultures or corresponding feed media.

BIOSTAT® T CHO

The BIOSTAT® T CHO mimics the human machine interface and functionalities of a BIOSTAT STR® 200 L single-use bioreactor. The training tool is based on actual cell culture batch data that were collected under normal process conditions. This database enables the software to derive cell growth curves for the mentioned process conditions algorithmically. Users have to control the process conditions to determine the optimal parameters for cell growth and antibody yields utilizing the broad variety of software features.







Three Options to Analyze Your Cultivation Data

Use the software's Trend, Offline data table and MFCS-History feature to evaluate your cultivation progress on the basis of different types of data, from preselected parameters to sample data up to conflated historical data.

Process time	VCD	VIABILITY	Glucose	Glutamine	TAA	pH	Osmolality	Lactate	Ammonium	Antibody	TCD	DCD
Process une	[10'6 Cells/mL]	[%]	[g/L]	[mmoML]	[g/L]	(pH)	[mOsmol/kg]	[g/L]	[mmol/L]	[g/L]	[10^6 Cells/mL]	[10^6 Cells/mL]
00:00:00	0.0	0.0	5.7	7.1	1.0	7.04	320	0.0	0.0	0.0	0.0	0.0
00:01:02	0.3	99.9	4.8	6.0	0.9	7.09	320	0.0	0.0	0.0	0.3	0.0
1d 14:34:22	0.9	99.9	4.6	6.0	0.9	7.19	319	0.1	0.0	0.0	0.9	0.0
3d 00:01:18	2.4	99.7	4.0	6.0	0.8	7.15		0.4	0.0	0.1	2.4	0.0
4d 00:01:34	4.6	99.9	7.2	5.7	1.1	7.09		0.5	0.1	0.1	4.6	0.0
5d 00:01:34	8.3	99.4	9.2	5.5	1.4	7.23	354	0.5	0.1	0.3	8.3	0.0
6d 00:01:34	13.4	99.9		5.2	1.6	7.06	359	0.5		0.6	13.5	0.0
7d 00:01:34	18.9	99.4	9.0	4.9	1.7	7.08	354	0.4	0.4	1.1	18.8	0.0
8d 00:01:34	22.7	99.4	7.3	4.6	1.8	7.19	345	0.4	0.6	1.7	22.9	0.2
9d 00:01:34	24.7	98.2	5.3	4.3	1.8	7.22	334	0.3	0.7	2.3	25.1	0.5
9d 17:13:02	24.6	96.0	5.6	4.1	1.9	7.15		0.3	0.9	2.8	25.6	
10d 00:01:34	24.7	95.1	4.1	4.1	1.8	7.14	326	0.3	0.9	3.0	25.9	
11d 00:01:34	23.4	91.5	4.1	3.8	1.8	7.11	326	0.2	1.1	3.8	25.7	2.2
12d 00:01:34	21.7	86.9	4.2	3.6	1.8	7.12	324	0.2		4.4	25.0	
13d 00:01:34	19.7	82.2	4.2	3.4	1.7	7.24	324	0.2	1.2	5.1	23.9	
14d 00:01:34	17.8	78.8	4.3	3.2	1.7	7.18	323	0.2	1.3	5.6	22.6	4.8
15d 00:01:34	16.0	75.6	4.4	3.0	1.6	7.23	323	0.1	1.3	6.2	21.2	5.2
16d 00:01:34	14.4	72.6	4.4	2.9	1.6	7.27	323	0.1	1.4	6.5	19.8	5.5
17d 00:57:02	13.0	69.9	4.3	2.8	1.6	7.13	321	0.1	1.4	7.1	18.4	5.5

Calibration SUBST-B

Totalize

30.00 ml/mir

Mode

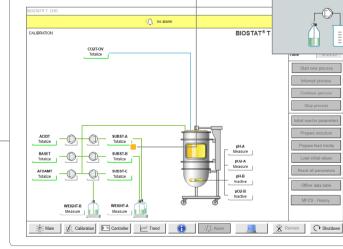
SUBST-B

Flow

×

All Calibration Functions of a Real Bioreactor

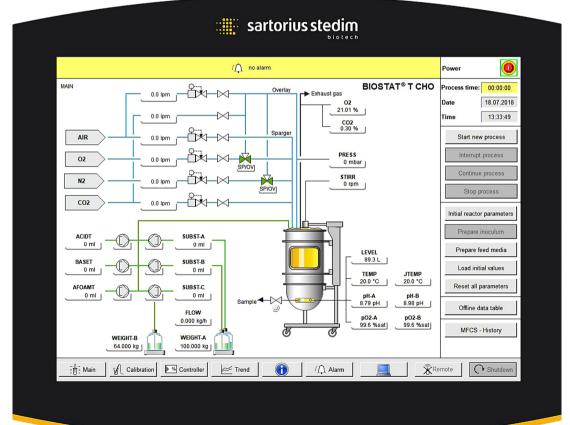
Calibrate the probes, pumps and totalizers of your virtual bioreactor to ensure high accuracy during your cultivation processes – like in reality.



Ready To Take The Next Step

After successfully accomplishing your virtual training with BIOSTAT[®] T you have acquired the knowledge and skills to start your cultivation in a real BIOSTAT[®] bioreactor.

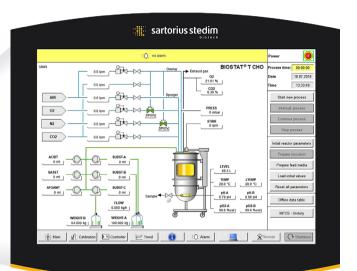




BIOSTAT STR®

Virtual training with the BIOSTAT® T CHO facilitates you to easily start operating a BIOSTAT STR® single-use bioreactor for commercial manufacturing. The BIOSTAT® T CHO mimics BIOSTAT STR® control software with it's user interface and process control tools.





Experiencing behavior of controllers and processes by using the BIOSTAT[®] T CHO reduces training time for new BIOSTAT STR[®] users and prevents human error. There is no faster knowledge transfer to the real world than with the BIOSTAT[®] T CHO.

Technical Data BIOSTAT® T CHO

	Minimum	Recommended		
	System Requirements	System Requirements		
Operating system	Windows [®] XP or newer	Windows [®] 7 or newer		
Processor	1 GHz Dual-Core or faster	2.5 GHz Dual-Core or faster		
RAM	2 GB	4 GB		
Available hard disk space	1000 MB	1000 MB		

For additional information and software access, please contact your local Sartorius representative or FRT.Support@Sartorius.com.

Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Italy

Sartorius Stedim Italy S.r.l.

50012 Antella-Bagno a Ripoli (FI)

Sartorius Stedim Netherlands B.V.

Sartorius Stedim Poland Sp. z o.o.

filtratie.nederland@sartorius-stedim.com

Via dell'Antella, 76/A

Netherlands

Poland

ul Wrzesinska 70

62-025 Kostrzyn

Phone +39.055.63.40.41

Phone +31.30.60.25.080

Phone +48.61.647.38.40

LLC "Sartorius Stedim RUS"

Russian Federation

199178 St. Petersburg

Phone +7.812.327.53.27

Avda. de la Industria, 32

Phone +34.913.586.098

28108 Alcobendas (Madrid)

Sartorius Stedim Spain, S.A.U.

Vasilvevsky Island

5th line 70, Lit. A

Edificio PAYMA

Spain

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

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

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

Specifications subject to change without notice. Copyright Sartorius Stedim Biotech GmbH. Printed in the EU on paper bleached without chlorine. Publication No.: SB11573-e180901 - Order No.: 88037-561-75 - Ver. 09 [2018



Switzerland Sartorius Stedim Switzerland AG Ringstrasse 24 a 8317 Tagelswangen Phone +41.52.354.36.36 U.K. Sartorius Stedim UK Ltd.

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