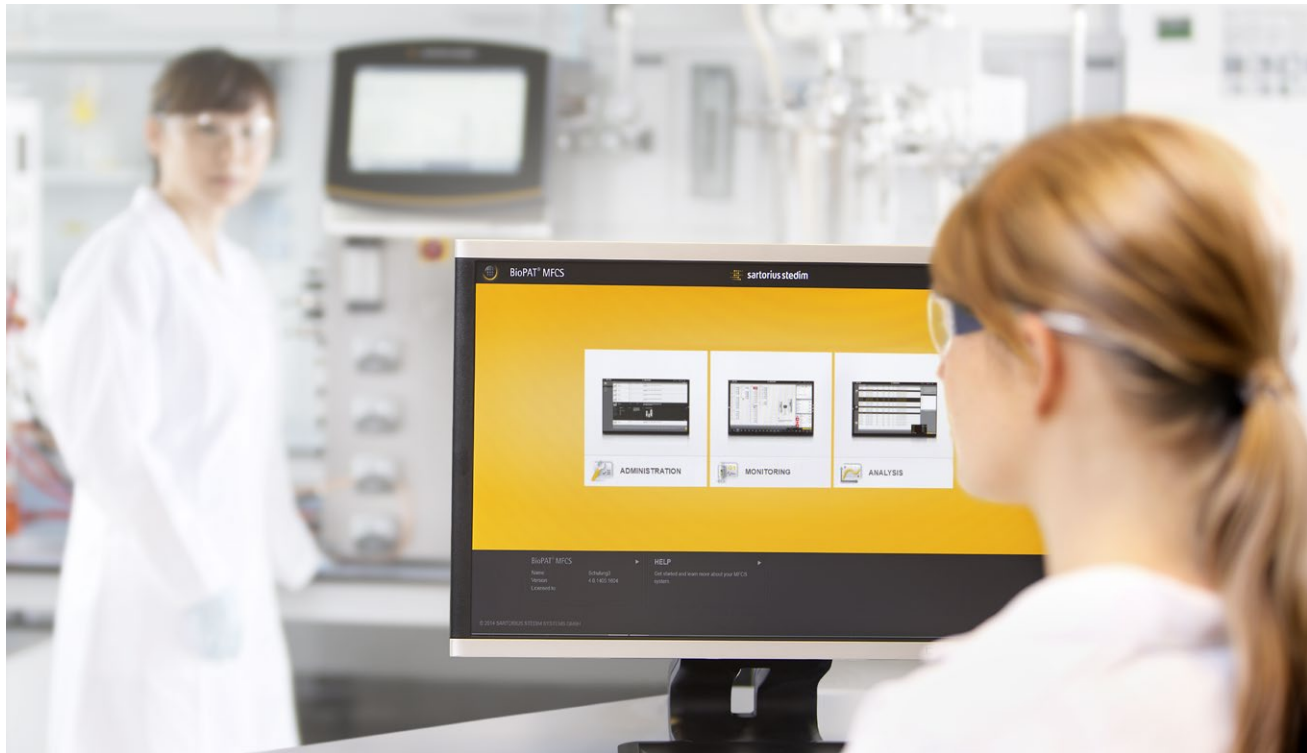




sartorius stedim
biotech

BioPAT[®] MFCS

The New Standard in Bioprocess Data Management



turning science **into solutions**





Developed for the Process – Designed for You!

BioPAT® MFCS enables you to incorporate a new standard in bioprocess data management and automation. Its reliable data acquisition, efficient trend monitoring and advanced recipe control make it an ideal tool for all upstream and downstream processes, no matter if you prefer single-use or reusable systems. The new BioPAT® MFCS is your solution for robust and reproducible processes – all backed by our expertise since 1986.

Clear Composition and Design

Experience the attractive and modern software design of BioPAT® MFCS. The intuitive workflow facilitates your daily tasks and its amazingly simple navigation allows rapid access to all the process-relevant information.

Cutting-edge and Compatible Technology

BioPAT® MFCS is compatible with the latest off-the-shelf hardware and software. Quickly collect your data and store it safely in the searchable SQL-database. You will benefit from high performance, scalability and high availability.

Sartorius Service and Support

The Sartorius Service will help you to get the maximum out of your new BioPAT® MFCS.

Extensive email support and the possibility to download the most recent version with the latest functionality will ensure reliable operation and easy handling.

Intuitive Navigation

Benefit from the easy access to all main operation areas of BioPAT® MFCS.

Calculation Capabilities

Calculate important bioprocess parameters and non-direct measurable variables with the calculation module.

Connectivity

Integrate PAT sensors and analyzers via OPC for enhanced process knowledge.



Quick Overview

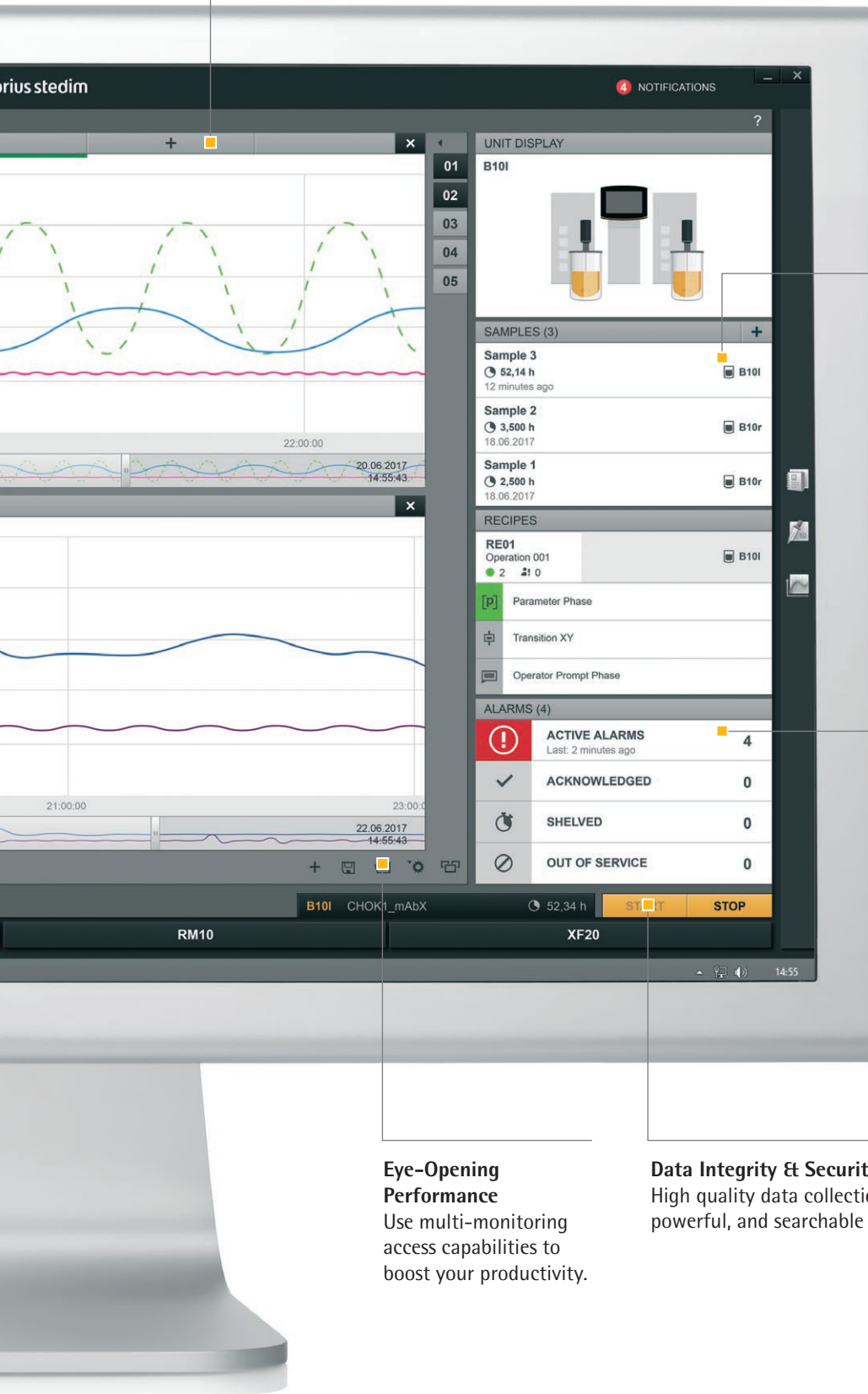
Quickly switch between up to 24 Units and get all relevant information at a glance.

Focus on Essentials

Get the right focus for a fast and clear data inspection with the slidable time bar.

Easy Comparison

Compare up to 12 real-time process values in an online configurable trend.



Centralized Storage

Enter off-line process data with the sample data module.

Process Safety

Let the alarm management system take care of process limits.

Eye-Opening Performance

Use multi-monitoring access capabilities to boost your productivity.

Data Integrity & Security

High quality data collection in a fast, powerful, and searchable SQL-database.

Turn Advanced Process Strategies into Reality

Easy Configuration
Get started immediately and implement your process strategies.

Quick Access
Pre-defined functions are right where you need them.

Features

- Intuitive graphically guided configuration
- Drag & drop operations and phases
- Sequential, parallel and repeated execution of phases
- State and time-dependent transitions
- Pre-defined phase types for setpoints, feeding profiles or timers

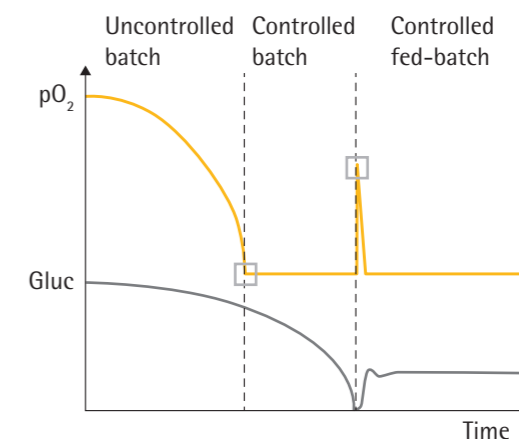
Benefits

- Improved batch-to-batch consistency
- Decreased risk of errors
- Minimized number of rejected lots
- Automated processing to free up operator time
- ANSI-88 compliant standardized automation

Application Example

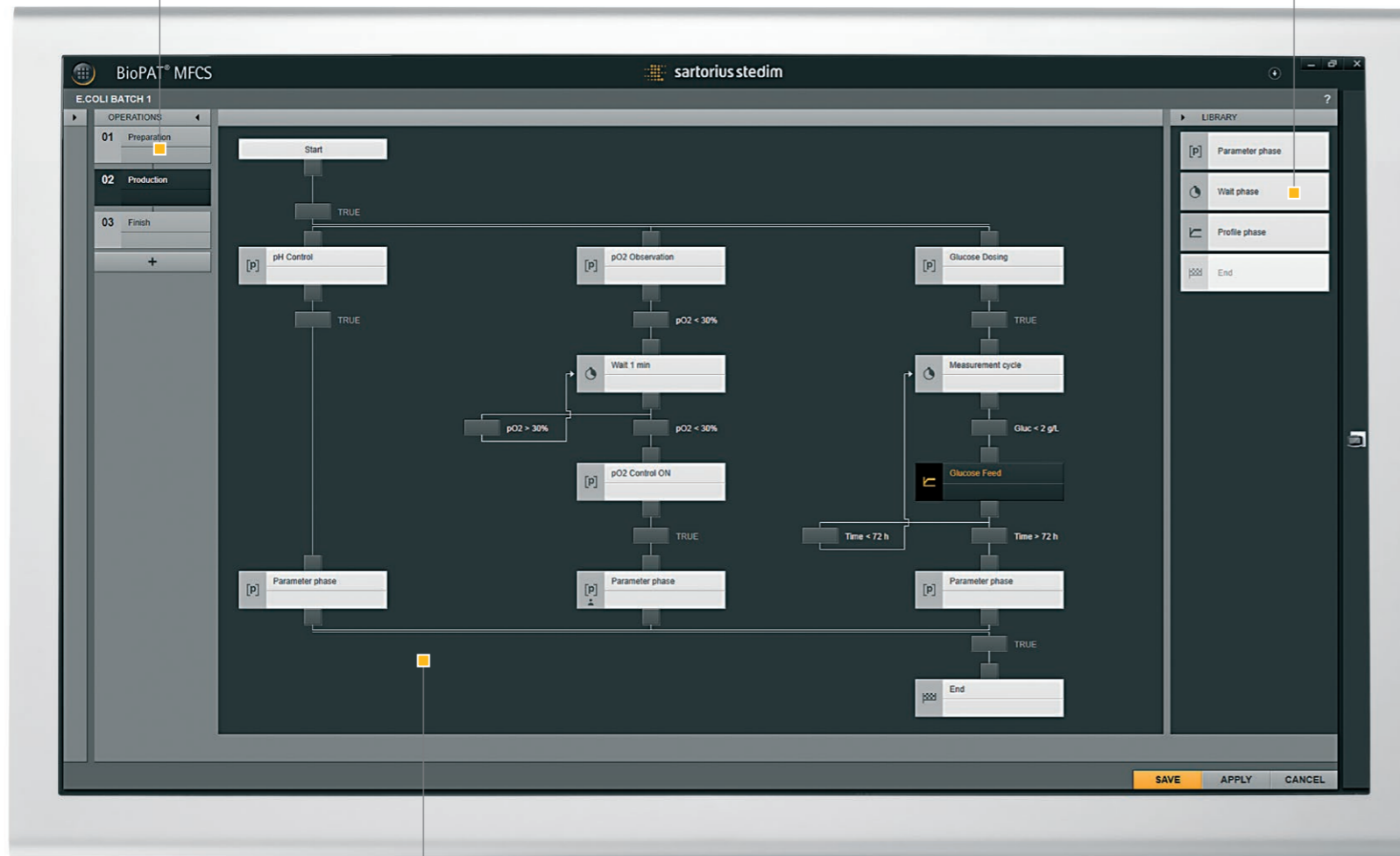
- Event-based activation of control loops
- Automatic batch end detection, e.g. based on pO_2 | offgas measurements
- Implement advanced feeding strategies for fed-batch processes

Automatically Switch between Process Phases



pO₂ control

Glucose control



BioPAT® MFCS Recipe Control Module

Intuitive Use
Clear workflows for maximum productivity.

Extend the Possibilities of BioPAT[®] MFCS



Recipe Control Module

ANSI-88 compliant batch control for robust process automation



OPC DA Client Module

Connect additional equipment, e.g. BioPAT[®] Trace, BioPAT[®] Xgas



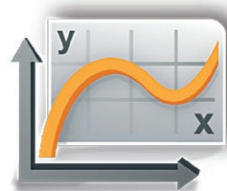
Sample Data Module

Integrate and manage off-line sample data



Calculation Module

Create soft sensors to access non-directly measurable quantities, e.g. OUR, RQ



Charting Module

Advanced visualization capabilities to compare live and historical batch processes



Multi User Option

Monitor and control your process from up to 30 BioPAT[®] MFCS clients simultaneously (RDP supported)

Minimum System Requirements

Server & Client standalone Installation

4-core 2.5 GHz processor
8 GB memory (RAM)
DirectX 10 graphics device (1 GB)
Network card (1 GB)
50 GB hard disk space
15.6" monitor (1366 × 768)

Multi User Installation

Client

4-core 2.5 GHz processor
4 GB memory (RAM)
DirectX 10 graphics device (1 GB)
Network card (1 GB)
50 GB hard disk space
15.6" monitor (1366 × 768)

Server

8-core 2.5 GHz processor
16 GB memory (RAM) +
2GB per Terminal Server session
Network card (1 GB)
250 GB hard disk space RAID 1 or 5

Supported Operating Systems (64-bit)

Windows® 8.1
Windows® 10

Supported Virtual Environments

VMware® Workstation
Hyper-V®

Ordering Information

Software

BioPAT® MFCS 4	Standard Package	1 Unit
----------------	------------------	--------

Additional Software Modules and Options

BioPAT® MFCS 4	Recipe Control Module
BioPAT® MFCS 4	OPC DA Client Module
BioPAT® MFCS 4	Sample Data Module
BioPAT® MFCS 4	Calculation Module
BioPAT® MFCS 4	Charting Module
BioPAT® MFCS 4	Multi User Option
BioPAT® MFCS 4	Microsoft® SQL Server 2014 Standard Runtime (Server, 64-bit)
BioPAT® MFCS 4	Microsoft® SQL Server 2014 Standard Runtime (Server, 64-bit) University

BioPAT® MFCS is available for up to 24 Units.
Please contact your local sales representative regarding a license for more than 1 Unit.

Upgradable with Upcoming Modules

Additional modules (e.g. User Management and OPC UA Server) for the adaptation of BioPAT® MFCS to your individual process requirements will be available soon.



Learn more about BioPAT® MFCS and download the demo version. Stay up-to-date about the latest product and security updates and visit BioPAT® MFCS website for more information.
◀ www.sartorius-stedim.com/biopatmfcs

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

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.
Kagyó 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

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
5th 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

Africa

Republic of South Africa

Sartorius South Africa (Pty) Ltd
Unit 4, Alphen Square South
853 16th Road
Midrand 1685, RSA

Phone +27.11.315.5444

Americas

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
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

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

Microsoft® and Windows® are registered trademarks of Microsoft Corporation, USA.