

# Ambr® 250 High Throughput

Fully Automated for Accelerated Process Development Simplifying Progress

### **SVISCISVS**

# Ambr® 250 High Throughput

A Fully Automated Bioreactor System for Parallel Fermentation or Cell Culture

Ambr<sup>®</sup> 250 High Throughput is a enhanced parallel microbial or cell culture system using 100 - 250 mL single-use bioreactors with a fully automated liquid handling platform. The system provides a high precision environment that supports the demands of microbial fermentation and mammalian cultures. Ambr<sup>®</sup> 250 High Throughput provides an efficient and rapid scaledown system to explore a wide range of conditions and strains with scalability to both bench top and larger stirred tank bioreactors.

#### Productivity

Ambr<sup>®</sup> 250 High Throughput dramatically improves productivity and enables full DoE experiments to be performed at a fraction of the cost and with higher throughput than can be done in traditional bench top reactors.

#### Wide Range of Cell | Strain Conditions and Performance

Ambr<sup>®</sup> 250 High Throughput provides a highly parallel system to rapidly develop processes for clones or strains in bioreactors that have full individual control of culture conditions.

#### Scalability

Because the bioreactor vessels are geometrically similar to larger bioreactors, all processes on the system correlate empirically to those of larger bioreactors making for optimum scalability.

#### Convenience

With a fully automated liquid handler together with single-use bioreactors that are fully integrated to sensors, users no longer need to spend long hours in the lab, manipulating, cleaning or setting up.



# System Features

Ambr<sup>®</sup> 250 High Throughput System Combines 12 or 24 "Easy Connect" Single-Use Bioreactors, Fully Automated Platform, Bioreactor Controller and Flexible System Control Software



SVIECTEVS

#### Ambr<sup>®</sup> 250 High Throughput platform

Integrates bioreactor stations, sensor readers, pumps and a liquid handler resulting in an efficient, high throughput scale-down model for process development and optimization.

### 01

#### Automated liquid handler

Fully automated liquid handler, for all liquid samples and additions, significantly reducing manual operator interactions.

#### 02

### Bioreactor station array

12 or 24 bioreactor stations in parallel with full individual process control, offers convenient set-up and simplifies running of full DOE experiments.

#### 03 Integrated biological safety cabinet

A class II biological safety cabinet designed to maintain an aseptic environment even during robot manipulation of bioreactors, significantly reducing any risk of vessel contamination



## Single-Use Bioreactor Vessels

Each bioreactor is equipped with impellers for stirring, a pH and DO sensor and ports to connect to the system's liquid and gas lines. The single-use technology means there is no need to clean bioreactors and sensors between runs.

#### Mammalian or microbial vessels

- 100 250 mL working volume with baffles
- Dual 20 mm pitched-blade or Rushton impeller
- Spot based DO sensor
- Disposable pH electrode
- Integrated gas and liquid inlet filters
- Robotic compatible cap for sampling
- Sparge and overlay gassing options
- Integrated condenser for output to exhaust gas analysis (OUR | CER)



#### 01 Septum cap Allows for rapid liquid

additions with a syringe.

#### 02 Gas tube

Gases can either be delivered into the headspace or sparged into the media. These delivery systems are independent and can function in parallel.

#### 03

### Integrated pH and DO sensors

vessel design incorporates disposable sensors to simplify the process of system preparation.

### 04

Double impeller – Rushton or pitch – blade For microbial or mammalian vessels respectively.



# Functions

Ambr® 250 High Throughput automatically controls, feeds and samples 12 or 24 bioreactor | fermenter vessels in parallel

### **Bioreactor controller**

- Three gasses per bioreactor with mass flow sensor:
- Mammalian
  - O<sub>2</sub>
  - CO<sub>2</sub>
  - N<sub>2</sub> | air
- Microbial
  - O<sub>2</sub>
  - Air
- N<sub>2</sub>
- Four positive displacement liquid pumps per bioreactor for high precision at low flow rates
- Individual bioreactor temperature control with heating or cooling
- Individual impeller speed control per bioreactor
- Optional exhaust gas analysis for OUR | CER
- Integrated CIP | SIP for pumps and liquid lines.

### Control software

- Fully flexible set up and control interface with simple, dialogue based interaction for quick and easy process creation
- Fine tuning options including programmable PID control loops and other advanced features
- Ambr<sup>®</sup> 250 High Throughput software is fully integrated with Umetrics MODDE Design of Experiment (DoE) application, enabling streamlined operation in execution of large DoE studies.





## **Applications**

Ambr<sup>®</sup> 250 High Throughput is configurable for microbial fermentation or mammalian cell culture and able to model a wide range of requirements in a variety of applications across biopharm as well as industrial biotech.

- Process optimization
- Process characterization
- Process robustness experimentation in support of QbD studies.
- Process scale-down model.

# Scalability

### Single-Use from Cell Line and Process Development to Production Scale

- Geometrical similarity of vessel design
- Consistent mixing and gassing strategies
- High performance gas transfer and mixing
- Reliable single-use platforms

#### **BioPAT® Process Insights**

Included with the Ambr<sup>®</sup> 250 High Throughput system is a 1 year license of BioPAT<sup>®</sup> Process Insights software application.

BioPAT® Process Insights enables predictive scale conversion and low risk process transfer between supported Sartorius bioreactors: Ambr® 15 and 250, Univessel® SU and Glass and Biostat STR® 50-2000 L.

BioPAT® Process Insights is a standalone, on-premise software application. Each license permits up to 10 concurrent seats. A renewal license (not included) is required after 1 year to continue usage of the software. Please check the technical specification for minimum computer hardware and software requirements.



Ambr® 250 High Throughput



Biostat® B Univessel® SU 2L



Biostat STR® 50



Biostat STR® 200





Ambr<sup>®</sup> 250 High Throughput



Biostat® B-DCU with Univessel® Glass 1 - 10 L



Biostat® D-DCU 10 - 200 L

Also scalable to multi-use technologies



#### Germany

Sartorius Stedim Biotech GmbH August-Spindler-Strasse 11 37079 Goettingen Phone +49 551 308 0

For further contacts, visit www.sartorius.com

**USA** Sartorius Stedim North America Inc. 565 Johnson Avenue

Bohemia, NY 11716 Toll-Free +1 800 368 7178