

Volume 16
April 2012
pp 24–25

3-4

G.I.T.

LABORATORY
JOURNAL

EUROPE

Reprint



GIT VERLAG

A Wiley Brand

Efficient Cell Cultivation

Single-Use Bioreactor for Process Development Enables More Flexibility

Now, biotechnologists are not only able to choose between single-use and reusable bioreactors and fermentors in the production sector, but in process development as well, where single-use systems are making their mark with advantages like lower costs and faster run times.

During the production of biopharmaceutical products like monoclonal antibodies, enzymes or vaccines, the central process step is performed in bioreactors. In a bioreactor, cells are cultivated aseptically under the most optimally controlled conditions (temperature, pH, dissolved oxygen concentration (DO) and culture media components). The most frequently used fermentors are called stirred tank bioreactors. Traditionally, they are made of glass (150-milliliter to 10-liter working volumes) for seed and small-scale cell cultivation and of stainless steel for larger applications (ten to several 10,000 liters).

Since the mid-90s, single-use bioreactors have been increasingly employed for cultivating mammalian cells from laboratory to production scale. The advantages of single-use systems are clear: Easy handling and no sterilization or cleaning times.

The first instrumented single-use bioreactor systems to become established were two-dimensional bioreactor bags (2D) like the Biostat CultiBag RM from Sartorius Stedim Biotech. These systems work with a platform that rocks back and forth; the wavelike, low-shear motion thoroughly mixes the liquid in the bioreactor bag. Due to their design, the working volume is limited to a maximum of several hundred liters. The considerable demand for single-use stirred tank bioreactors has inspired a variety of manufacturers to develop equivalent systems.

These systems employ three-dimensional (3D) stirred tank bioreactor bags. A bag container made of stainless steel not only guarantees stability, but also controls the temperature. Systems are supplied with working volumes ranging from 50 l to a maximum of 2,000 l.

Ideally, these 3D stirred tank bioreactor bags essentially feature the same design criteria as do conventional stainless steel bioreactors.

Beyond this, it is practical to equip these 3D stirred tank bioreactor bags with single-use sensors for pH and DO measurement to fully exploit all advantages of single-use systems.

with 2-liter working volumes have proven to be particularly well-suited and deliver high performance. These models feature reactor geometries comparable to those of the production vessels; their volume capacity is both affordable and sufficiently large for taking analytical samples, and above all, they are easy to handle.



Single-use Stirred Tank Bioreactors

The latest evolution in single-use stirred tank bioreactors are vessels for the process-scale development, providing an alternative to conventional glass bioreactors. This vessel category is used in the biopharmaceutical industry mainly for developing and optimizing cell lines and processes, as well as for scale-up and scale-down experiments. That is why it is of significant importance that these vessels replicate the design criteria of production-scale stirred tank bioreactors for both reusable and single-use applications. Stirred tank bioreactors

The advantages that single-use culture vessels offer on this scale are remarkable. Beyond the criteria mentioned above, autoclaving relatively cumbersome glass vessels is eliminated as are costs for their servicing and maintenance.

Moreover, single-use culture vessels can also be used interchangeably with glass vessels so that during capacity peaks or maintenance of bioreactors, for instance, reusable vessels can be easily exchanged for single-use bioreactors in the interim. This enables downtimes in bioreactor process control to be reduced or even eliminated. For such uninterrupted cultivation, the only other alterna-

tive is to purchase additional, fully equipped glass bioreactor vessels, which entails high capital costs.

The UniVessel SU is a single-use culture vessel that meets the design criteria of conventional glass and production-scale stirred tank bioreactors. The body of the culture vessel is made of polycarbonate and contains graduation marks. The bottom conforms to the toro-spherical base design of stainless steel stirred tank bioreactors. To ensure a stable upright position during operation, the vessel is placed in a holder. Like on glass stirred tank bioreactors, all vessel inlet ports are located on the lid of the device. The lid is firmly assembled onto the body of the vessel and equipped with a number of ports. All inlet and harvesting ports are supplied with thermoweldable tubing and feature MPC or Luer connectors. A needle-free septum port is installed for sampling. The culture vessel is drained completely through the harvesting tube that is channeled to the center of the culture vessel base.

Aeration can take place either in a submerged configuration and/or through the headspace; both air inlet and exhaust channels are equipped with sterilizing-grade filters. Submerged aeration takes place through an L-sparger with tiny holes. Moreover, the device features three threaded ports (steel conduit thread PG 13.5) for attaching standard 12 mm sensors. The stirrer shaft has two ball bearings that ensure stability and a lip seal that maintains sterility. Thorough, low-shear mixing of the cell culture medium is accomplished by two 3-blade segment impellers (down-flow).

Additionally, the device includes non-invasive, optical, single-use pH and DO sensors. These sensor patches contain special dyes that are excited to fluorescence when exposed to light emitted at a characteristic wavelength through the culture vessel wall.

The properties of the longer wavelength fluorescent light emitted by the sensor patches are re-

	2L- reusable glass bioreactor *	UniVessel SU 2L	500 L reusable stainless steel bioreactor	CultiBag STR 200L
Stirred-tank bioreactor design / process parameters				
Working volume [L]	2	2	500	200
Stirrer speed [rpm]	200	212	65	100
Aeration rate [vvm]	0,1	0,1	0,16	0,1
Height-to-diameter ratio	1.5 :1	1:4 :1	1:1 1	1:3 : 1
Impeller design	3-blade segment impeller (standard stirrer)	3-blade segment impeller (standard stirrer)	Lightnin A320 (3-blade segment impeller)	3-blade segment impeller (standard stirrer)
Number of impellers	1	2	1	2
Sparger design	L-sparger with holes	L-sparger with holes	2-line aeration system/open pipe sparger with microsparger	Ring sparger
Process-related parameters				
Mixing time [s]	20	< 3	48	16
Specific power input [W/m ³]	2.8	1,4	13,6	17.5
kLa [1/h]	10	6,9	8	6

* Data from A-mab: A Case Study in Bioprocess Development; CMC Biotech Working Group

Tab.1: Comparison of reusable vs. single-use bioreactors

spectively influenced by the parameters of pH or dissolved oxygen saturation of the culture medium in contact with the sensor patches. The Sensolux optoelectronic components built into the culture vessel holder are used both to excite fluorescence and to detect the fluorescent light emitted (see Fig. 2).

Scale-up and Comparability

The scale-up or scale-down of fermentation processes is not an exact science; rather it combines experience with the application of the equivalent process parameters and design criteria. The key scaling criteria include bioreactor design along with process-related parameters,

such as mixing behavior and gas transfer. Design criteria essentially comprise the bioreactor's height-to-diameter ratio, impeller design, ratio of the impeller to bioreactor diameter as well as the construction of the sparger design.

With respect to mixing behavior, the mixing times and the specific power input are the parameters considered, while the volumetric oxygen transfer coefficient is important for gas transfer (see Table 1).

The single-use culture vessel can be easily integrated into both new as well as existing bioreactor control units. All that is needed for this purpose is a heating jacket for regulating the temperature. A reusable motor adapter enables

connection of the device to the existing impeller motor; a variety of adapters are supplied for the various stirred tank bioreactors. The presterilized pH and DO sensors are attached to the vessel lid under a sterile workbench. The temperature sensor is simply inserted into the open port in the vessel lid. For enhanced time efficiency, a barcode scanner can be used to read in the calibration parameters of the precalibrated sensor patches. As a result, the labor-intensive steps involving autoclaving, calibrating and attaching the pH and DO sensors are eliminated, which means users can fully benefit from all the advantages of single-use vessels.

The UniVessel SU provides the flexibility required for connection to

existing bioreactor control units. Further culture vessel components, such as an exhaust cooler, various impeller geometries, etc., are in the development pipeline and, moving forward, will allow microbial high cell-density cultivation with instrumented single-use culture vessels. Beyond all this, more vessel sizes will be added to the product family.

Authors:

A. Grebe, J. Weyand:
Sartorius Stedim Biotech

Contact:

Sartorius Stedim Biotech
Goettingen, Germany
info@sartorius-stedim.com
www.sartorius-stedim.com

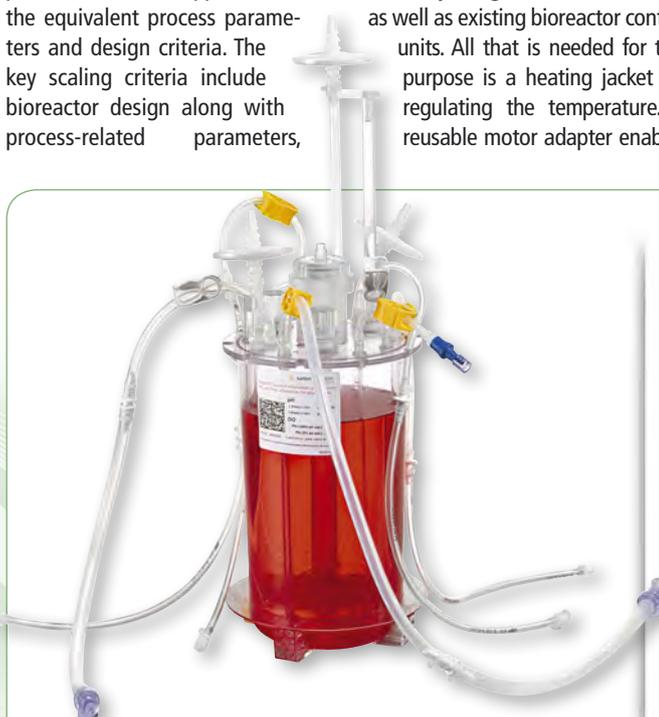


Fig. 1: The UniVessel SU is a novel, single-use culture vessel specially designed for process development.

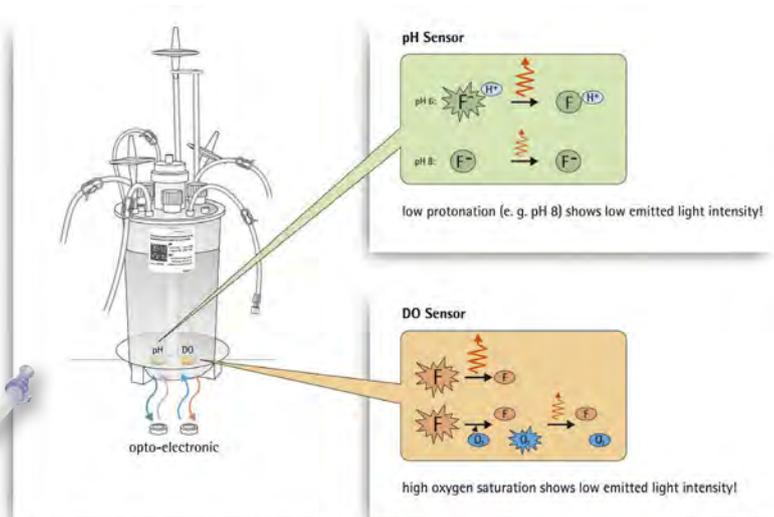


Fig. 2: Functional Diagram of the UniVessel's Optical pH and DO Sensors

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
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 Biotech S.A.
ZI Les Paluds
Avenue de Jouques - CS 91051
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.845619

Sartorius Stedim France SAS
ZI Les Paluds
Avenue de Jouques - CS 71058
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.846545

Austria

Sartorius Stedim Austria GmbH
Franzosengraben 12
1030 Vienna

Phone +43.1.7965763.18
Fax +43.1.796576344

Belgium

Sartorius Stedim Belgium N.V.
Leuvensesteenweg, 248/B
1800 Vilvoorde

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

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 ICR"
Uralskaya str. 4, Lit. B
199155, Saint-Petersburg

Phone +7.812.327.5.327
Fax +7.812.327.5.323

Spain

Sartorius Stedim Spain SA
C/Isabel Colbrand 10,
Oficina 70
Polígono Industrial de Fuencarral
28050 Madrid

Phone +34.90.2110935
Fax +34.91.3589623

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

America

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
Av. Dom Pedro I, 241
Bairro Vila Pires
Santo André
São Paulo
Cep 09110-001

Phone +55.11.4451.6226
Fax +55.11.4451.4369

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 Biotech (Beijing) Co. Ltd.
Shanghai Branch Office
3rd Floor, North Wing, Tower 1
No. 4560 Jin Ke Road
Pudong District, Shanghai 201210

Phone +86.21.68782300
Fax +86.21.68782332 | 68782882

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., Diawa 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 +65.6778.2494

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

