



Evaluation of BIostat[®] RM for Plant Cell Culture



Application
Note

#06

#07

#08

#09

#10

Introduction

Plants are major sources of natural or secondary products, pharmaceuticals, dyes, or flavours. Over the last decade, plant cell culture was developed as a technique for studying plant metabolism, physiology and development. This new biotechnological technology has reached maturity at industrial scale to offer new perspectives for the production of plant derived active ingredients.

Plant cells were first grown in bioreactors in the 1960s using various systems adapted from animal cell cultures. First cultures were performed in stirred-tank reactor. As shear stress was identified as a critical point for growth, cultures were run afterwards at low mixing. Indeed, plant cells are tolerant or sensitive to mechanical mixing used in classical bioreactors generating high shearing stress that may damage the cells. With the introduction of airlift bioreactor, particularly for biomass production, an alternative to the stirred tank reactor was proposed to overcome the shear stress limitation challenge.

Compared to microbial cells, growth and oxygen demand of plant cells are relatively low, requiring a low $K_L a$ value. Therefore to lead a plant cell culture in bioreactor, mixing has to be efficient with low shear conditions and without excessive aeration.

In general, plant cell cultures in suspensions form aggregates, therefore cells settle rapidly if mixing is stopped. Moreover the sampling could be difficult and non-reproducible. This could be accentuated at the end of growth when the plant cell cultures become viscous forming a kind of meringue. Sampling system has to be adapted to avoid loss of biomass and heterogeneity of samples.

Another characteristic of plant cell cultures, particularly in airlift bioreactors, is foam production. It is a point to consider because it can block air outlet. Antifoam has to be added but it may affect the growth performances of the culture.

Because of the low growth performances of the plant cells and therefore a long period of the culture process there is an increasing contamination risk, hence the control of sterility is an important parameter as well.

All those elements are relevant and in favour to find alternative bioreactor systems for the cultivation of plant cells at industrial scale.

In this application note, BIOSTAT® RM bioreactors were evaluated for the cultivation of plant cells. For this evaluation a suspended cell culture, belonging to the family of the Plantaginaceae was grown in the dark in order to produce cosmetic active ingredients, intended for the Cosmetic industry. We developed a unique range of biochemical actives, with substantiated efficacy for sebum regulation ('adjusting' dry or oily skin).

Our objectives were i) to manage cultivation of plant cells in new devices from small to large scale volume using different BIOSTAT® RM systems and ii) to maintain growth performances and productivity at good level throughout scale-up, as it can be measured in flask cultures.

Materials and methods

Plant cell cultures

Cells were cultivated in Gamborg medium at 25°C. Media were prepared according to internal protocols and added by sterilizing filtration.

Bioreactors

- BIOSTAT® RM20/50
- BIOSTAT® RM50
- BIOSTAT® RM200
- BIOSTAT® RM600

Monitoring

- Growth was characterized by Packed Cell Volume measurement (PCV %)
- Dissolved Oxygen (DO) was monitored by a single use optical sensor installed in the bag
- pH was measured by an external device temperature was set-up and regulated at 25°C

Process scheme

- Feed transfer via connexion bag to the bioreactor bag by using a Biosealer and Biowelder.
- Add fresh MEDIA by filtration (sterilization of filters by autoclaving and sterile connection to the bag)
- Sampling: Connexion to the bag of a sterile vial, filling of the vial realized using a peristaltic pump
- Gas supplied AIR only

This experiment was performed using BIOSTAT® RM20/50, BIOSTAT® RM50, BIOSTAT® RM200 and BIOSTAT® RM600 bioreactors. The two objectives of this study were to demonstrate that we could cultivate the plant cells of interest in these bioreactors and to validate the scale-up. The parameters applied during the different cultures are shown in table 1.

	Erlenmeyerflasks	RM20/50	RM50	RM200	RM600
Culture volume (L)	1	5	25	100	300
Aeration (lpm)	NA	0.6	2	5	10
Shaking (rpm)	100	20–22	22	9–18	13
Angle (°)	NA	8	9	8–9	9

Table 1:
Process parameters for plant cell cultures

The culture cycles were led during 2 to 4 weeks.

The downstream process has been optimized to recover the active ingredient but is not part of this application note.

Results

Plant cell growth profiles according to the type of bioreactor used are shown in figure 1.

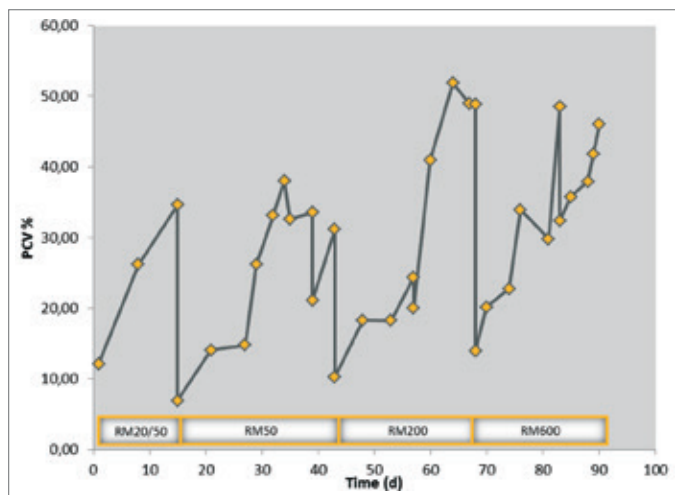


Figure 1:
Growth performances for plant cells production in BIostat® CultiBag bioreactors

A 5 L culture was carried out during 2 weeks in the BIostat® RM20/50, biomass concentration increased from 12% PCV to 35% final PCV value which is in line with flask culture performances (PCV value up to 50% in 2–3 weeks).

Biomass content for 25L culture reached a similar PCV value as in the 5 L culture (38% PCV). For both 5L and 25L cultures, pH values were around 6.3–6.8 and dissolved oxygen (DO) values were between 20 and 30%.

To continue the scale-up, a 100L culture was performed in a BIostat® RM 200. Growth performances reached 50% PCV and pH and DO values were respectively around 6.5 and 10%.

These results were confirmed in 300 L culture which showed satisfactory growth performances of 50% PCV. pH data were collected with an external device and the profile showed values around 6.2–6.4. DO data were collected via the single use optical sensor and remained stable around 10%.

Conclusion

The objective of this study was to demonstrate the ability to cultivate plant cells in BIostat® RM bioreactors and to manage scale-up from Erlenmeyerflasks to industrial production scale.

These objectives were reached and a satisfactory biomass production was achieved of 45–50% PCV during a scale-up of the process from flasks to the BIostat® RM 600 via the BIostat® RM 20, BIostat® RM 50 and BIostat® RM 200. As the evaluation of the active ingredient didn't fall under the scope of this evaluation, further tests are required to also evaluate this important factor of the process.

In terms of practicality, sampling and transfers of the culture were made either under laminar flow hood for the small volume or with Bunsen burner. However using routinely the Biosealer and Biowelder for making the tubing connections can significantly improve handling and eliminate the need for the mentioned alternatives. Even while cells were in aggregates, sampling was rather easy and no losses of biomass was noted.

Thanks to Magella Drouet, Marion Thibault and Frédéric de Baene (Research Manager) for their great work and participation for this study.

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., 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 +65.6778.2494

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



▶ www.sartorius-stedim.com