



Propionic acid production of *Propionibacterium acidipropionici* in a BIOSTAT® A



Application
Note

#09

#10

#11

#12

#13

Carlos H Luna-Flores, Lars Nielsen
and Esteban Marcellin*

Australian Institute for Bioengineering and
Nanotechnology (AIBN), The University of
Queensland, Brisbane, Qld 4072, Australia

* Corresponding author
Dr. Esteban Marcellin, Australian Institute
for Bioengineering and Nanotechnology,
The University of Queensland,
Brisbane, QLD 4072, Australia.
Phone: +61 7 334 64298,
Fax: +61 7 3346 3973,
Email: e.marcellin@uq.edu.au

Introduction

Traditionally derived from fossil fuels, propionic acid (PA) is widely used in the food and pharmaceutical industries. PA is used as a preservative and increasingly, in the synthesis of monomers. Recently, the market for PA has grown to over 410,000 tonnes per year with a steady 3% increase per annum. Mounting environmental concerns have encouraged end-users to look for sustainable PA alternatives, opening a niche market for biological production of PA (Liu et al., 2012). Bio-production of PA addresses many environmental concerns and offers a sustainable alternative for the production of C3 chemicals including propylene, propanol and vinyl propionate. *Propionibacterium acidipropionici* (*P. acidipropionici*) is a gram-positive, anaerobic, rod bacteria that naturally produce PA as the main fermentation product through the Wood-Werckman cycle (Liu et al., 2012; Parizzi et al., 2012). Natively, PA is produced along other organic acids such as lactate, succinate and acetate.

The anaerobic fermentation of *P. acidipropionici* is sensitive to environmental, physicochemical, and hydrodynamic conditions, including temperature, pH, dissolved oxygen, and shear stress. These conditions need to be tightly controlled during the fermentation. *P. acidipropionici* uses a complex mixture of nutrients to produce optimally PA. This generates a dynamic metabolism through the different growth phases. In this study, we demonstrated the robustness of the BIOSTAT® A to ferment *P. acidipropionici* cells. The equipment was able to control the pH successfully despite the large production of PA. The pH control parameters had to be adjusted to have an adequate pH control. Nitrogen was constantly fed into the reactor to ensure anaerobic conditions. The temperature was controlled with the thermo-jacket surrounding the vessel and the recirculation chiller. The chiller also helped to control evaporation by cooling the exhaust gases. The hydrodynamic conditions were maintained using two Rushton impellers.

1. Material and Methods

1.1 Bacteria and media

P. acidipropionici ATCC 55737 selected from a collection of 17 strains (Stowers, Cox, & Rodriguez, 2014). The strain was kept at -80°C using glycerol (20%) as cryoprotector. The culture media (PAM) was made from (g/L): yeast extract (10), trypticase soy (5), K₂HPO₄ (0.05), MnSO₄ (0.05), and glucose (75). Media components and the carbon source were sterilized separately for 20 min at 121°C.

1.2 Cultivation

Glycerol stock cultures were revived in 1.5 mL Eppendorf tubes containing 1 mL of PAM media inoculated with 0.8% (v/v) of the glycerol stock. This culture grew for 24 hours at 32°C. The culture was transferred to a 15 mL Falcon tube containing 14 mL of PAM media and allowed to grow for an additional 24 hours. 5% (v/v) of this culture was used to inoculate a 250 mL serum bottles containing 100 mL of PAM media which grew for an additional 24 hours. Cells from the serum bottles in mid-exponential phase were used to inoculate the fermenter at an initial OD_{600nm} of 0.3. The fermentation was performed using a 1 L BIOSTAT® A fermenter. The fermenter was equipped with standard probes and standard controllers, controlling pH, dissolved oxygen, temperature, and agitation. The agitation rate was kept constant at 300 rpm. The pH was controlled at 6.4 using 10 M NaOH. The temperature of the culture was maintained at 32°C using an electric jacket and a recirculation chiller. The exhaust gases condenser was set to 20% to avoid evaporation media. Prior to inoculation, the fermenter was sparged with N₂ for at least 15 minutes. A constant N₂ flow was kept for the entire fermentation at a flow rate of 300 ccm using a correction factor for the mass flow controller of 0.992.

1.3. Analytical methods

Optical density of the culture was measured at 600 nm using a Biochrom Libra S12 UV/Vis Spectrophotometer. Organic acids and carbohydrates were quantified by ion-exclusion chromatography using an Agilent 1200 HPLC system and an Agilent Hiplax H column (300 × 7.7 mm, PL1170-6830) with guard column (SecurityGuard Carbo-H, Phenomenex PN: AJO-4490). Sugars were monitored using a refractive index detector (Agilent RID, G1362A) set on positive polarity and optical unit temperature of 40°C, while organic acids were monitored at 210 nm (Agilent MWD, G1365B). 30 µL of sample was injected onto the column using an auto-sampler (Agilent HiP-ALS, G1367B) and column temperature kept at 65°C [70°C] using a thermostatted column compartment (Agilent TCC, G1316A). Analytes were eluted isocratically with 4 mM H₂SO₄ at 0.6 mL/min for 26 min. Chromatograms were integrated using ChemStation (Rev B.03.02[341]).

2. Results

P. acidipropionici produces organic acids which generates a strong pH dynamic which needs to be controlled. Figure 1 shows the control parameters in the bioreactor using the BIOSTAT[®] A. As illustrated, the pH was maintained constant via the addition of base using the integrated peristaltic pump which supplied the required 10 M NaOH. The PID control parameters were adjusted to P: 300%, I: 0%, and D: 0%. N₂ was continuously supplied to maintained the anaerobic environment in the fermenter. The temperature and agitation were also controlled and monitored.

Figure 2 shows the growth kinetic profiles of *P. acidipropionici*. PA was the main fermentation product along with minor quantities of acetic acid and succinic acid. As can be observed in the figure, the growth stopped after 40 h. However, PA the production continued until the end of the fermentation, resulting in an increase in the addition of NaOH until the 96th h.

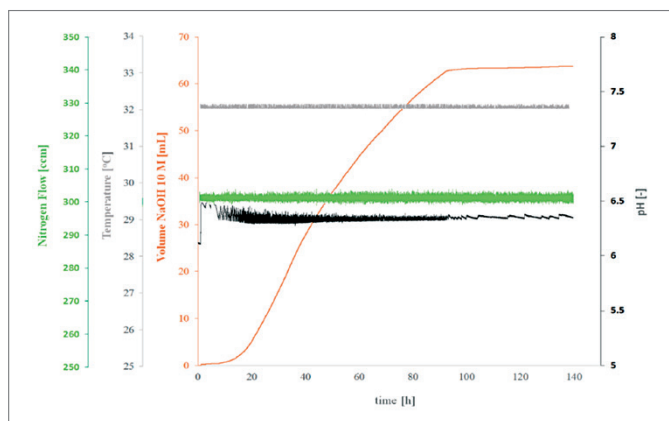


Figure 1: BIOSTAT[®] A control parameters during the fermentation of *P. acidipropionici*. Nitrogen flow, temperature, pH, and volume of NaOH 10 M are shown.

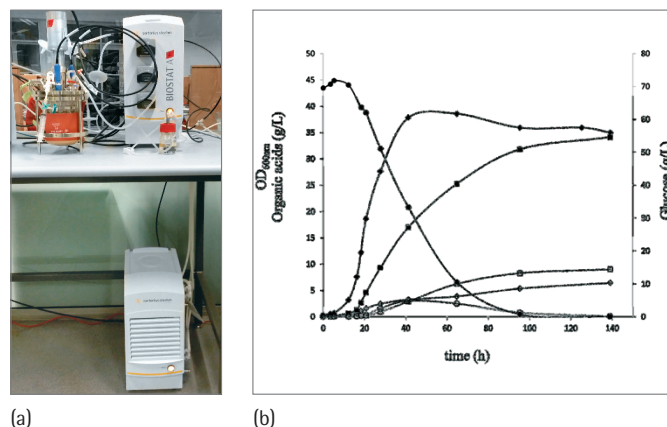


Figure 2: Growth kinetic of *P. acidipropionici* fermentation in a BIOSTAT[®] A.

- (a) BIOSTAT[®] A equipment.
 (b) Growth and production kinetic.
 Optical Density: Black line and ◆;
 Propionic Acid: Black line and ■;
 Glucose: Black line and ●;
 Succinic Acid: Grey line and □;
 Acetic Acid: Grey line and ◇;
 Pyruvate: Grey line and ○.

3. Conclusion

P. acidipropionici fermentation for the production of PA can be performed in a BIOSTAT[®] A. The most important parameter to control was the pH, which was controlled by changing the PID. In conclusion, BIOSTAT[®] A can be used to ferment anaerobic microorganisms for the production of more than 30 g/L of propionic acid.

References

- Liu, L., Zhu, Y., Li, J., Wang, M., Lee, P., Du, G., & Chen, J. (2012). Microbial production of propionic acid from propionibacteria: current state, challenges and perspectives. *Critical Reviews in Biotechnology*, 32(4), 374–81. doi:10.3109/07388551.2011.651428
- Parizzi, L. P., Grassi, M. C. B., Llerena, L. a, Carazzolle, M. F., Queiroz, V. L., Lunardi, I., ... Pereira, G. a G. (2012). The genome sequence of *Propionibacterium acidipropionici* provides insights into its biotechnological and industrial potential. *BMC Genomics*, 13, 562. doi:10.1186/1471-2164-13-562
- Stowers, C. C., Cox, B. M., & Rodriguez, B. a. (2014). Development of an industrializable fermentation process for propionic acid production. *Journal of Industrial Microbiology & Biotechnology*, 41(5), 837–52. doi:10.1007/s10295-014-1423-6

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 FMT S.A.S.
ZI des Paluds
Avenue de Jouques – CS 91051
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.845619

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques – CS 71058
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.846545

Austria

Sartorius Stedim Austria GmbH
Modcenterstrasse 22
1030 Vienna

Phone +43.1.7965763.18
Fax +43.1.796576344

Belgium

Sartorius Stedim Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles

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

Netherlands

Sartorius Stedim Netherlands B.V.

Phone +31.30.60.25.080
Fax +31.30.60.25.099

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
Fax +48.61.879.25.04

Russian Federation

LLC "Sartorius Stedim RUS"
Uralskaya str. 4, Lit. B
199155 St. Petersburg

Phone +7.812.327.53.27
Fax +7.812.327.53.23

Spain

Sartorius Stedim Spain, S.A.U.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)

Phone +34.902.110.935
Fax +34.91.358.96.23

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

Ukraine

LLC "Biohit"
Post Box 440 "B"
01001 Kiev, Ukraine

Phone +380.44.411.4918
Fax +380.50.623.3162

Americas

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
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP - Brasil

Phone +55.11.4362.8900
Fax +55.11.4362.8901

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

Peru

Sartorius Peru S.A.C.
Av. Emilio Cavenecia 264 San Isidro
15073 Lima, Perú

Phone +51.1.441 0158
Fax +51.1.422 6100

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 (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
Fax +86.21.6878.2882

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