

MoFlo™ Bacterial Proliferation and PHB Production

Introduction

Population dynamics were followed during growth on phenol in response to adaptation processes of *Cupriavidus necator*, bacterial individuals morphologically nearly uniform, yet structurally and functionally diverse. The species belongs to the group of Betaproteobacteria and is known for its capability to degrade numerous toxic substances and therefore is used often as a model organism.¹

Overflow metabolism is induced in cases of imbalanced substrate supplies and temporary changes in nutrient sources, which is thought to be part of a survival strategy. An example is the synthesis of poly- β -hydroxybutyrate (PHB).² Variations were analyzed in bacterial structural cellular parameters like chromosome contents (via 4', 6-diamino-2'-phenylindole -DAPI) and PHB (via Nile red).

Materials and Methods

C. necator JMP 134 was grown aerobically at 30°C and at pH 7.0 in minimal medium in 0.5 l shake flasks for batch experiments and with 0.02 percent phenol as carbon and energy source. The harvested cells were centrifuged at 3,200 x g for six minutes, fixed with 10 percent NaN₃ and stored at 4°C. For flow cytometric measurements³ the preserved cells were centrifuged again, washed in NaCl-phosphate buffer (0.4 M Na₂HPO₄/NaH₂PO₄, 150 mM NaCl, pH 7.2) and resuspended at a concentration of 3 x 10⁸ cells/mL.

DNA was stained with 4', 6-diamino-2'-phenylindole (DAPI) according to Meistrich et al.⁴ The staining is AT-specific and stoichiometric and thus provides quantitative information about the chromosome multiples within single bacterial cells.

To analyse the PHB content of the cells, Nile red was used as follows: 40 μ L of Nile red stock solution (1 mg/mL acetone) was added to 3 x 10⁸ cells/mL DAPI stained cells and incubated for eight minutes.

Flow cytometric measurements were performed using a MoFlo Cell Sorter equipped with two water-cooled argon-ion lasers (Innova 90C and Innova 70C). A wavelength of 488 nm (500 mW) was used to excite Nile red at the first observation point. DAPI was excited by multi-line-UV (333-365 nm, 110 mW) at the second observation point. Forward scatter signals (FSC; cell size related) were analyzed using the 488 nm line and a neutral density filter with an optical density of 2.3. The orthogonal side scatter signal (SSC; cell granularity related) was reflected first by a beam-splitter and then recorded after reflection by a 555 nm long-pass dichroic mirror using a D 488/10. Nile red and DAPI fluorescence was passed through a 580/30 and a 450/65 band-pass filter, respectively. Amplification was linear for all signals and measurement of the events was triggered by the SSC signal. Fluorescent beads (Polybead Microspheres: 0.483 μ m diameter; flow check BB/Green compensation kit, Blue Alignment Grade) were used to align the MoFlo (CV-value about 2 percent). Data were acquired and analyzed using Summit software.

Results

Individual batch grown cells of *C. necator* were characterized with respect to their PHB content, FSC/SSC behavior and proliferation activity. After inoculation, the cells increased in size and an eukaryotic-like cell cycle behavior was observed (Figure 1: 2 hours). PHB production was not yet initiated. The low fluorescence intensity observed is due to unspecific staining of the lipid membranes of the cells. With the decline in nutrient or oxygen concentrations within the cultivation flask, proliferation stopped. The surplus of carbon was stored as PHB in cellular granules (Figure 1: 12 hours). When the carbon and energy sources were depleted, these storage materials were consumed during cell division and maintenance (Figure 1: 16 – 19 hours).

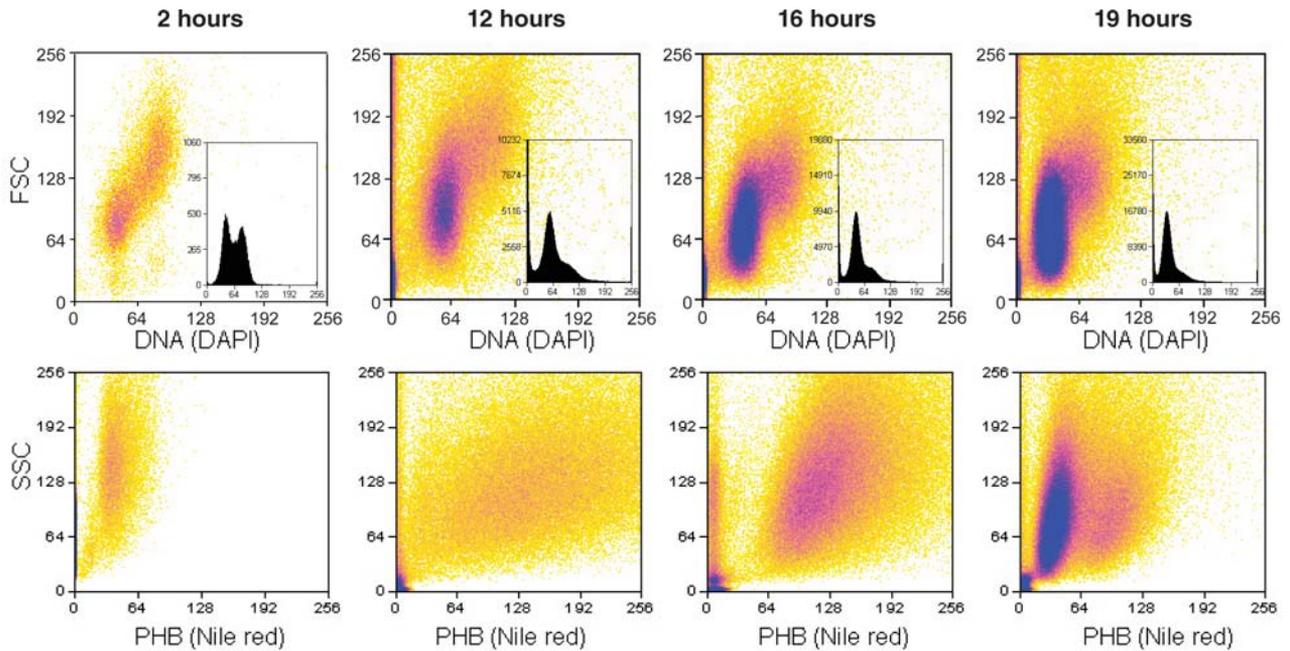


Figure 1a

DNA content (vs. FSC, upper row) and PHB content (vs. SSC, lower row) of batch cultivated *C. necator* JMP134 cells grown on 0.02% phenol. Sampling times are indicated. The insets represent the chromosome contents.

Conclusion

The cell cycle distribution profiles of *C. necator*, represent an eukaryotic-like cell cycle behavior when grown on phenol under batch conditions. Since the species is usually able to induce uncoupled DNA synthesis on other substrates, these patterns provide information on phenol being a good but suboptimal growth substrate. Also the strain is able to use PHB to survive disturbed growth conditions and to use it for maintaining turnover processes and cell-homeostasis.

References

- Clément P, Matus V, Cárdenas L, González B. Degradation of trichlorophenols by *Alcaligenes eutrophus* JMP134. *FEMS Microbiol. Lett.* 1995; 127: 51-55.
- Srienc F, Arnold B, and Bailey JE. Characterization of intracellular accumulation of poly-β-hydroxybutyrate (PHB) in individual cells of *Alcaligenes eutrophus* H16 by flow cytometry. *Biotechnol. Bioeng.* 1984; 26: 982-987.
- Müller S, Bley Th, Babel W. Adaptive responses of *Ralstonia eutropha* to master feast and famine conditions analysed by flow cytometry. *J. Biotechnol.* 1999; 75: 81-97.
- Meistrich ML, Göhde W, White RA, Schuhmann J. Resolution of X and Y spermatids by pulse cytophotometry, *Nature* 1978; 274: 821-823.
- Maskow T, Müller S, Lösche A, Kemp R. Appropriateness of different heat measurement principles and flow cytometry used to control the continuous synthesis of polyhydroxyalcanoates. *Biotechnol. Bioeng.* 2006; 93 (3): 541-552.

For Research Use Only. Not for use in Diagnostic Procedures.

Australia, Gladesville (61) 2 9844-6000 Canada, Mississauga (1) 905 819 1234 China, Beijing (86) 10 6515 6028
 Czech Republic, Prague (420) 272 01 73 32 Eastern Europe, Middle East, North Africa, South West Asia: Switzerland, Nyon (41) 22 365 3707
 France, Villepinte (33) 1 49 90 90 00 Germany, Krefeld (49) 2151 33 35 Hong Kong (852) 2814 7431
 India, Mumbai (91) 22 3080 5101 Italy, Cassina de' Pecchi, Milan (39) 02 953921 Japan, Tokyo (81) 3 5530 8500 Korea, Seoul (82) 2 404 2146
 Latin America (1) (305) 380 4709 Mexico, Mexico City (001) 52 55 9183 2800 Netherlands, Mijdrecht (31) 297 230630 Puerto Rico (787) 747 3335
 Singapore (65) 6339 3633 South Africa/Sub-Saharan Africa, Johannesburg (27) 11 564 3203 Spain, Madrid (34) 91 3836080
 Sweden, Bromma (46) 8 564 85 900 Switzerland, Nyon (41) 0800 850 810 Taiwan, Taipei (886) 2 2378 3456 Turkey, Istanbul (90) 216 309 1900
 UK, High Wycombe (44) 01494 441181 USA, Fullerton, CA (1) 800 742 2345

B2008-8956

www.beckmancoulter.com

© 2008 Beckman Coulter, Inc.

BMR-PRINTED IN U.S.A.



Simplify • Automate • Innovate