



Monitoring *E. coli* Cultures with the BioLector and Multisizer 4e Instruments

Mariana Gil, Ph.D.

Introduction

Escherichia coli (*E. coli*) is a facultative anaerobic bacterium that lives in the lower intestine of warm-blooded animals, including humans.¹ *E. coli* can be cultured easily and inexpensively in a laboratory setting and has become an important model organism in genetics, microbiology and biotechnology.^{1,2} *E. coli* is the most common organism used for the large-scale production of therapeutic proteins. Indeed, 30% of approved therapeutic proteins are currently being produced using *E. coli*.² This application note will demonstrate how the BioLector microbioreactor and the Multisizer 4e Coulter Counter can be used to optimize *E. coli* culture conditions and characterize cell growth (Figure 1).



FIGURE 1: The BioLector microbioreactor and the Multisizer 4e Coulter Counter can be used to monitor *E. coli* cultures.

Monitoring Optimal Cell Culture Conditions

The BioLector microbioreactor is small-scale, automated system that enables high-throughput screening, cultivation parameter monitoring (e.g., pH, biomass, oxygen saturation, shaking speed, and fluorescence intensity) and feeding strategy optimization. Importantly, all these parameters are monitored online without the necessity to stop shaking, or to take samples. In this example, the BioLector instrument was used to determine the optimal feeding strategy for *E. coli* fed-batch cultures.

Methods

Cultivations of *E. coli* (BL21 wild type) were carried out in BioLector microbioreactor using the microfluidic 32-well FlowerPlate microtiter plate M2P-MTP-MF32-BOH1. This plate is equipped with two standard optodes for the

optical online monitoring of the pH (range 5-7) and dissolved oxygen (DO). A preculture in terrific broth (TB) medium was prepared (initial optical density = 0.1) at 37 °C in a 250 mL shake flask at 250 rpm for 6 hours. In addition, two master mixtures of Wilms MOPS minimal medium with 20 g/L glucose and 10 g/L glycerol were used for the main culture in the microbioreactor (initial optical density = 0.1). The culture volume in the microtiter plate was 800-1200 µL per well. The plate was shaken at 1200 rpm at 37 °C, the humidity level was set at 75 %. The pH-adjusting agent used was 3 M NaOH. The feeding solution was composed of 10 mL glycerol solution (500 g/L glycerol) and 2 mL of nitrogen source (250 g/L $(\text{NH}_4)_2\text{SO}_4$, 25 g/L NH_4Cl). Two different feeding strategies were used: DO-triggered feed and constant feed. In the DO-triggered feed protocol, the MTP was set to deliver a pulse of 4 µL of the feeding solution when the DO dropped below 30 % for the first time for activation, and then whenever the DO was above 70 % during the cultivation process. In the constant feed protocol, the MTP was set to deliver 3 µL/h of the feeding solution during the cultivation process

Results and Discussion

Figure 2 shows the results of the DO-triggered feed experiment. The upper chart shows the biomass gain over 44 hours; the middle chart shows the volume of the pH-adjusting agent (Volume B) and the corresponding pH course; the lower chart shows the DO level and the glycerol feed volume (Volume A). Overall, the results showed that at the beginning of the cultivation, the cell number was low and thus the DO was approximately 100 %. Next, oxygen consumption increased together with cell proliferation, and, consequently, DO levels dropped. The stationary phase started at about 7 hours. When the DO dropped below 30 % for the first time, the system delivered a pulse of 4 µL of the feeding solution for activation. During this phase a slow culture growth rate was observed, which correlated with low oxygen consumption and the consequent rapid increase of the DO. Next, when the DO reached 70 % a new feed pulse was delivered, and, consequently, the growth rate increased again. This cycle was repeated throughout experiment. Therefore, the feed curve showed a step-like increase dependent on the DO fluctuations, whose impact was also observed in the biomass signal. Due to the ongoing addition of glycerol, and the correlated bacterial growth, the culture produced acetate. This led to a growth-dependent drop in the pH value (Figure 2, middle chart). Hence, the volume of the pH-adjusting agent (Volume B in Figure 2) increased according to the culture growth.

Figure 3 shows the same variables measured when the culture received a constant glycerol feed (3 µL/h) once the culture reached the stationary phase (at about 7 hours). The constant feeding protocol led to fewer fluctuations in the DO level than the DO-triggered pulse feed protocol. In consequence, both the culture growth and the pH value also showed fewer fluctuations.

It can be concluded that the biomass growth behaved differently depending on the feeding strategy. The final biomass gain was higher when the feeding protocol was constant, even when the culture received a smaller total volume of feeding solution. Moreover, since the constant feed protocol led to almost no fluctuations in the pH level, a smaller total volume of the pH-adjusting agent was required. Thus, a constant feed can be recommended to optimize *E. coli* cultures.

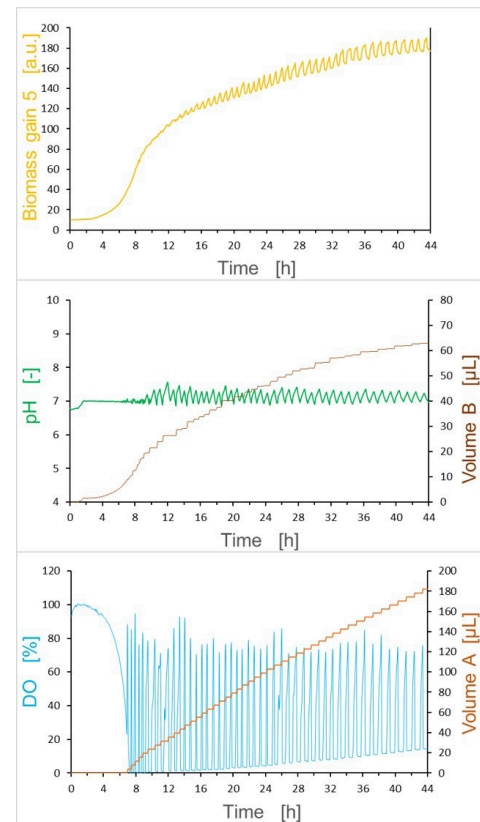


FIGURE 2: Fed-batch cultivation of *E. coli* with a dissolved oxygen (DO)-triggered addition of glycerol (Volume A). The pH was kept constant through addition of NaOH (Volume B).

In conclusion, the BioLector microbioreactor is ideally suited for microbial fed-batch cultivation processes and can be used to optimize bacterial culture conditions. Up to 32 pH-controlled fed-batch cultivations can be performed in one experiment at the same time. Continuous shaking during measurements ensures continuous mass transfer and homogenization of the suspended cells. The BioLector microbioreactor not only allows the online measurement of biomass, pH and dissolved oxygen but also the signal-dependent regulation of the pH value and feeding of nutrient solutions.

Previous studies have shown the efficacy of the BioLector microbioreactor to screen optimal *E. coli* culture conditions prior to scale-up.^{3,4,5} In summary, the system enables the evaluation and discrimination of different cultures improving screening conditions, process development and scale-up procedures.

Monitoring Cell Growth

Cell size is a key parameter used to gain insights into the diversity of cellular mechanisms (e.g., cell cycle, osmotic regulation, cell death, pathogenesis, phagocytosis, species diversity, etc.). The Multisizer 4e instrument uses the Coulter method to detect particles from 200 nm to 1,600 μm regardless of the particle's nature or optical properties. Particles suspended in a 0.9% electrolyte solution are drawn through a small cylindrical aperture. Two submerged electrodes located on each side of the aperture create an electric current. As each particle passes through the aperture, it displaces its own volume of conducting liquid, momentarily increasing the impedance of the aperture. This produces a voltage pulse that is proportional to the volume of the particle. The number of electrical pulses indicates particle count, while the amplitude of the electrical pulse produced depends on the particle's volume.⁶ In the following example, the Multisizer 4e instrument was used to characterize cell sizes in a *E. coli* culture.

Methods

A sample of the *E. coli* culture was pre-diluted 1:1000 with Isoton 2 in a 10 mL Accuvette ST. From this dilution, 50 μL was used for each measurement. The aperture used was 20 μm and each measurement lasted 36 seconds. The complete list of settings is shown in Table 1. The generated data were processed using Digital Pulse Processing (DPP) technology. DPP enables the acquisition, storage and display of each individual pulse. This means that individual areas of the pulse spectrum can be evaluated separately at a later date.

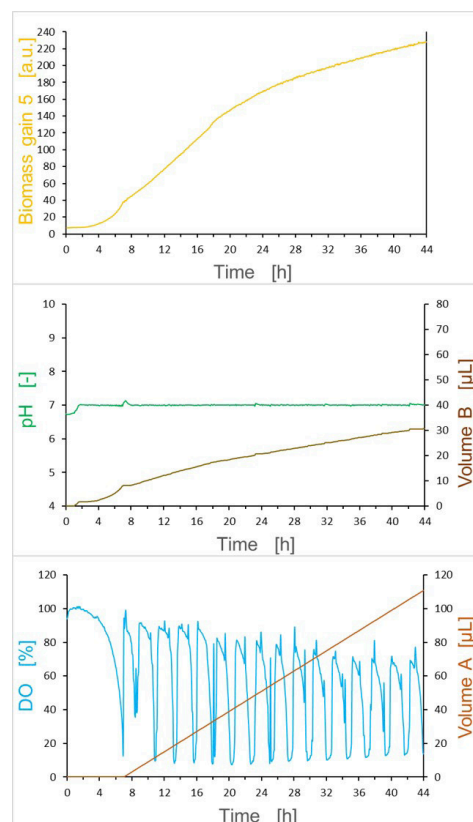


FIGURE 3: Fed-batch cultivation of *E. coli* with a constant glycerol feed (Volume A). The pH was kept constant through addition of NaOH (Volume B).

Table 1: Settings used in the experiment

System	Multisizer 4e
Aperture	20 μm
Current/Gain	200 $\mu\text{A}/16$
SOM	Volumetric 50 μL
Vessel	Accuvette ST 10 mL
Threshold	0,4 μm
Measuring Time	36 sec
Size bins	400
Dilution	1:100
Counts	140474

Results and Discussion

The DPP data of the experiment indicates that the size of the *E. coli* size in this culture ranged between 0.4 to 2.2 μm (Figure 4).

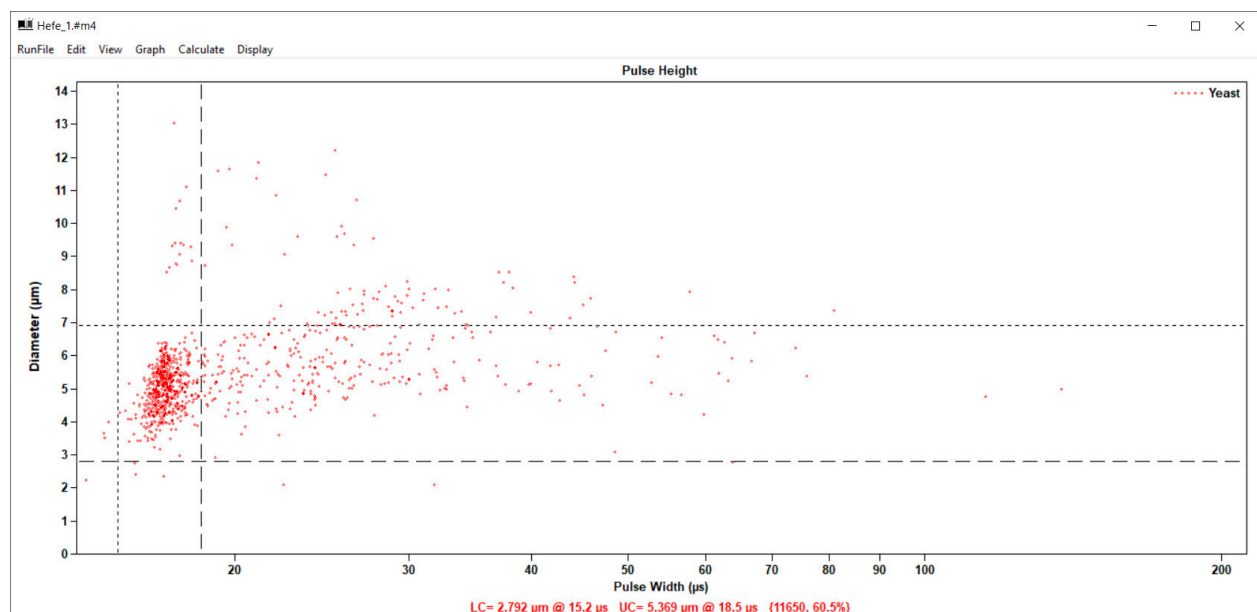


FIGURE 4: Digital Pulse data. Impulse width vs. impulse height. The dotted lines indicate the data points selected for further analysis.

From this raw data it is possible to select a range of interest (dotted lines in figure 4) and create different frequency histograms (i.e. the number of cells for each volume). For example, figure 5 shows the resulting histogram if a range between 0.8 and 1.5 μm is selected.

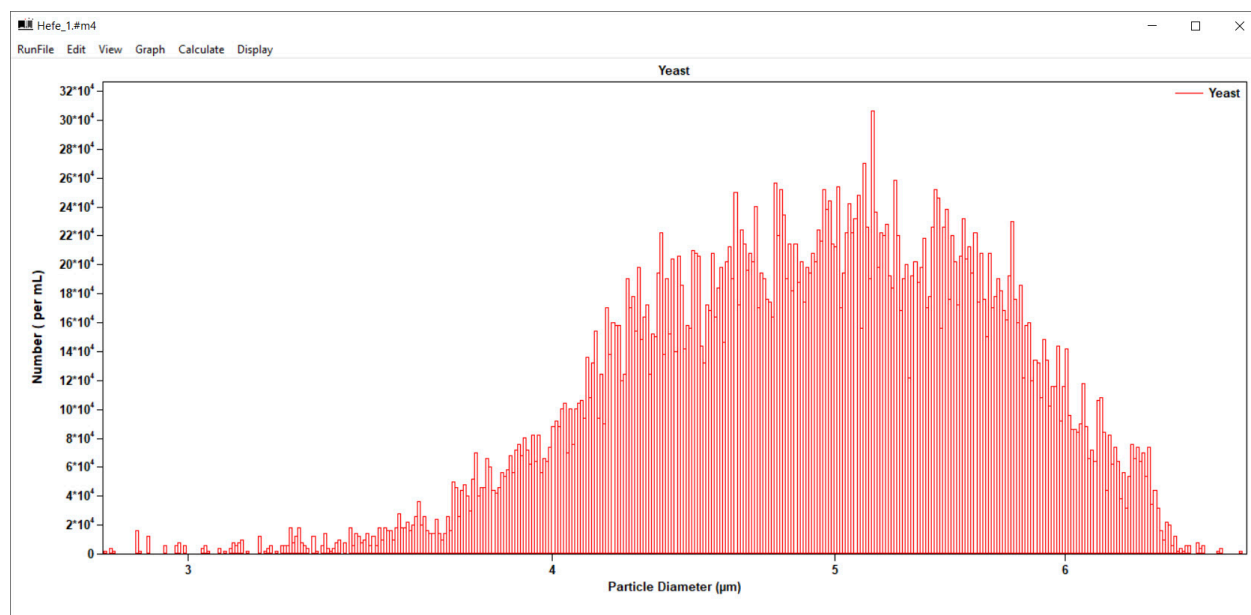


FIGURE 5: Distribution of cell sizes across the sample (diameter, in μm).

The DPP technology provides ultra-high resolution and accuracy (i.e. detection of 1 particle in 1 mL of a sample), which is unattainable using other current technologies. Thus, it can be used to accurately determine the volume, number, and cell concentration and detect real-time size changes during the cultivation process. For example, different studies used the instrument to characterize *E.coli* growth dynamics² and growth fitness.⁸ The Multisizer instrument also has been used to investigate the growth rate dependence of cell size of *E. coli* cultures under different nutrient conditions. Under normal conditions, *E. coli* growth rate positively correlated with cell size (i.e., large cells grow faster). However, this study found that when nutrients are scarce, protein overexpression inverts this relationship and leads to large cells with a low growing rate. This finding reveals an important role of protein synthesis in cell division control in *E. coli* cultures.⁹

Conclusion

The BioLector microbioreactor and the Multisizer 4e instruments are bench-top devices with an intuitive user interface that can be integrated to any laboratory to optimize and monitor *E. coli* cultures for a broad range of applications.

References

1. Tenaillon O, Skurnik D, Picard B, *et al.* The population genetics of commensal *Escherichia coli*. *Nat. Rev. Microbiol.* 2010;8:207-217. doi: [10.1038/nrmicro2298](https://doi.org/10.1038/nrmicro2298).
2. Baeshen MN, Hejin AMA, Bora RS, *et al.* Production of biopharmaceuticals in *E. coli*: current scenario and future perspectives. *J. Microbiol. Biotechnol.* 2015;25:953-962. doi: [10.4014/jmb.1412.12079](https://doi.org/10.4014/jmb.1412.12079).
3. Toeroek C, Cserjan-Puschmann M, Bayer K, Striedner G. Fed-batch like cultivation in a micro-bioreactor: screening conditions relevant for *Escherichia coli* based production processes. *Springerplus.* 2015;4:490. doi: [10.1186/s40064-015-1313-z](https://doi.org/10.1186/s40064-015-1313-z).
4. Ladner T, Mühlmann M, Schulte A, *et al.* Prediction of *Escherichia coli* expression performance in microtiter plates by analyzing only the temporal development of scattered light during culture. *J. Biol. Eng.* 2017;11:20. doi: [10.1186/s13036-017-0064-5](https://doi.org/10.1186/s13036-017-0064-5).
5. Fink M, Cserjan-Puschmann M, Reinisch D, *et al.* High-throughput microbioreactor provides a capable tool for early stage bioprocess development. *Sci. Rep.* 2021;11:2056. doi: [10.1038/s41598-021-81633-6](https://doi.org/10.1038/s41598-021-81633-6).
6. Rhyner M, Prestigiacomo G, Kumar K, Lee L. Cellular analysis using the Coulter principle. measurement with great accuracy and speed. Available from: <https://www.mybeckman.uk/resources/reading-material/application-notes/cellular->

[analysis-using-the-coulter-principle](#).

7. Tsuchiya K, Cao YY, Kurokawa M, *et al.* A decay effect of the growth rate associated with genome reduction in *Escherichia coli*. *BMC Microbiol.* 2018;18:101. doi: [10.1186/s12866-018-1242-4](#).
8. Ying BW, Yama K, Kitahara K, Yomo T. The *Escherichia coli* transcriptome linked to growth fitness. *Genom. Data.* 2015;7:1-3. doi: [10.1016/j.gdata.2015.11.011](#).
9. Basan M, Zhu M, Dai X, *et al.* Inflating bacterial cells by increased protein synthesis. *Mol. Syst. Biol.* 2015;11(10):836. doi: [10.15252/msb.20156178](#).

*The data shown in this Application Note was generated on BioLector II and BioLector Pro microbioreactors. The BioLector model shown in this Application Note is the latest model of the series, a BioLector XT microbioreactor.



©2022 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit [beckman.com](#)

22.04.4672.PCC