

# Automated Quantification of Budding *Saccharomyces cerevisiae* using a Novel Image Cytometry Method

Leo L. Chan<sup>1,2</sup>, Daniel J. Laverty<sup>2</sup>, Alexandria L. Kury<sup>2</sup>, Dmitry Kuksin<sup>1,2</sup>, Alnoor Pirani<sup>1</sup>, and Kevin Flanagan<sup>1</sup>

<sup>1</sup>Revvity Health Sciences, Inc., 360 Merrimack St., Suite 200, Lawrence, MA 01843

<sup>2</sup>Center for Biotechnology and Biomedical Sciences, Merrimack College, North Andover, MA 01845

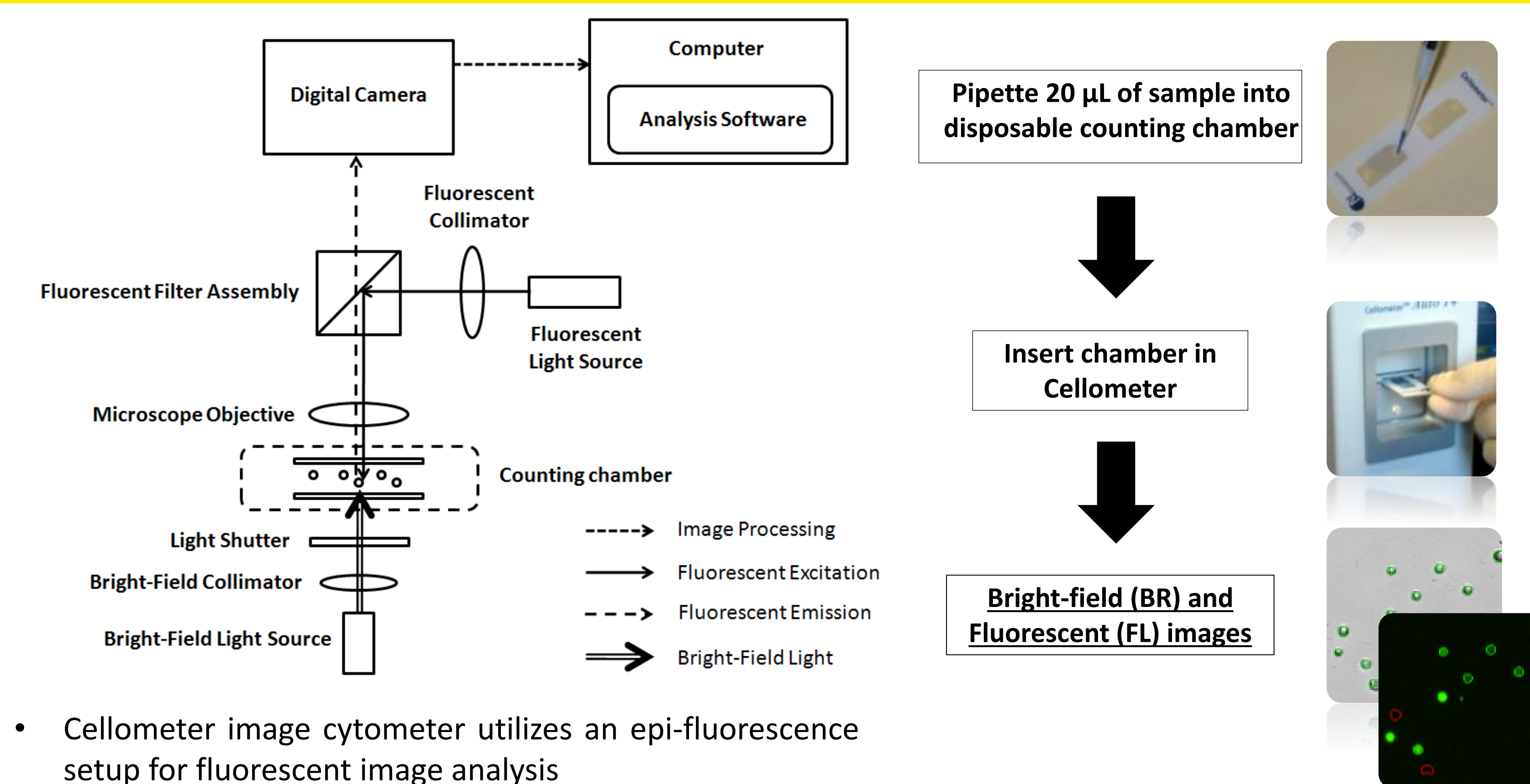
## 1. ABSTRACT

The measurements of concentration, viability, and budding percentages of *Saccharomyces cerevisiae* are performed on a routine basis in the biofuel and brewing industries. Generation of these parameters is of great importance in a manufacturing setting, where they can aid in the estimation of product quality, quantity, and fermentation time of the manufacturing process. Specifically, budding percentages can be used to estimate the reproduction rate of yeast populations, which directly correlates with metabolism of polysaccharides and bioethanol production, and can be monitored to maximize production of bioethanol during fermentation. The traditional method involves manual counting using a hemacytometer, but this is time-consuming and prone to human error. In this study, we developed a novel automated method for the quantification of yeast budding percentages using Cellometer image cytometry. The automated method utilizes a dual-fluorescent nucleic acid dye to specifically stain live cells for imaging analysis of unique morphological characteristics of budding yeast. In addition, cell cycle analysis is performed as an alternative method for budding analysis. We were able to show comparable yeast budding percentages between manual and automated counting, as well as cell cycle analysis. The automated image cytometry method is used to analyze and characterize corn mash samples directly from fermenters during standard fermentation. Since concentration, viability, and budding percentages can be obtained simultaneously, the automated method can be integrated into the fermentation quality assurance protocol, which may improve the quality and efficiency of the bioethanol production process.

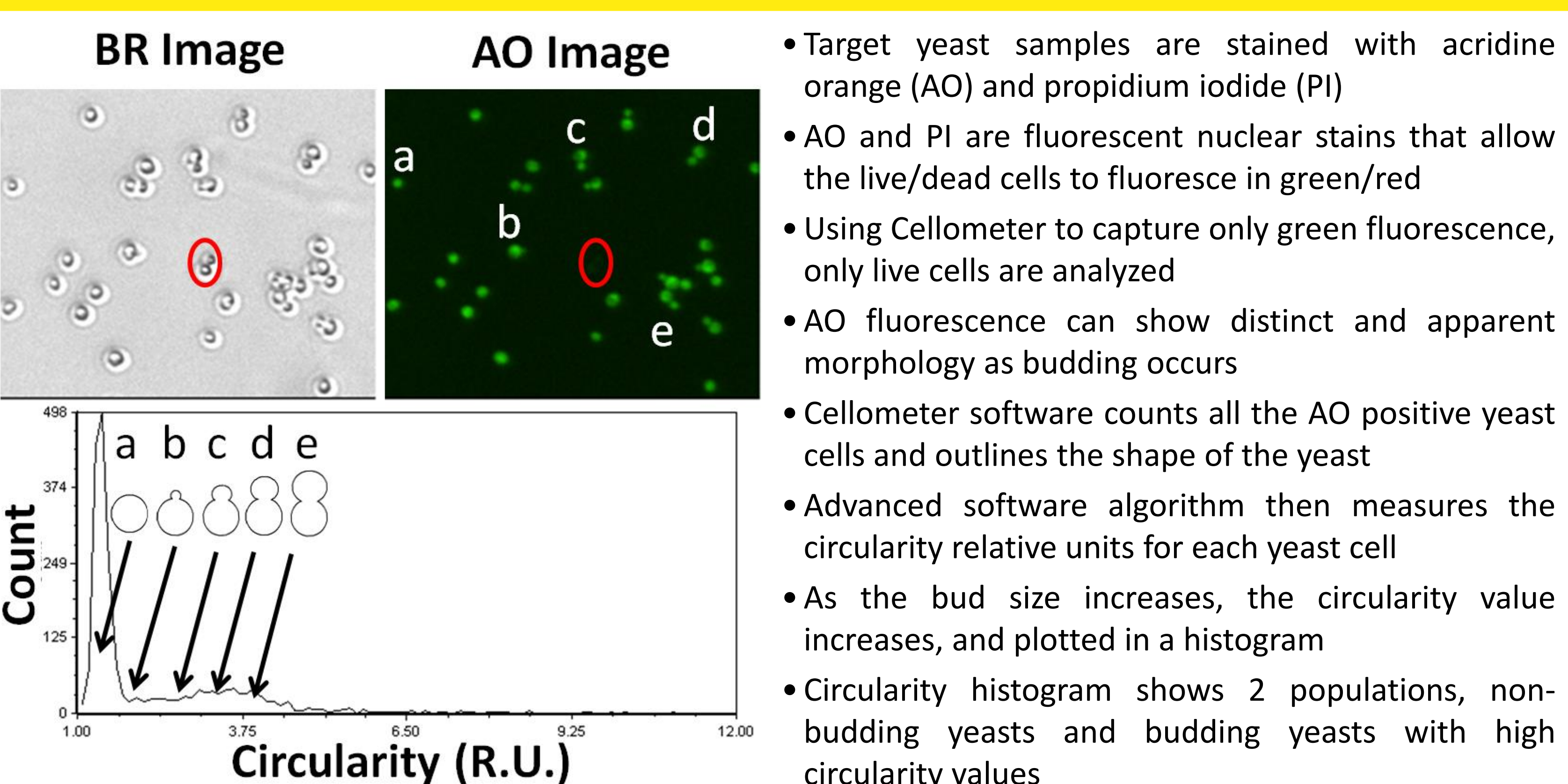
## 2. CURRENT METHODS FOR MEASURING BUDDING YEAST CELLS

Methods	Description	Known Issues
<b>Hemacytometer</b>	Manually counting budding cells	<ul style="list-style-type: none"> <li>Time-consuming and tedious process</li> <li>Requires experienced user for accurate counting</li> </ul>
<b>Fluorescence Microscopy</b>	Visualization of Calcofluor-stained "budding scars"	<ul style="list-style-type: none"> <li>Qualitative observe instead of quantitative analysis</li> <li>Not automated, low throughput</li> </ul>
<b>Flow Cytometry</b>	<ul style="list-style-type: none"> <li>Quantitative cell cycle analysis</li> <li>Automated analysis</li> </ul>	<ul style="list-style-type: none"> <li>Relatively expensive and high maintenance</li> <li>Requires experienced user for proper operation</li> <li>Cannot visually observe budding yeasts</li> </ul>

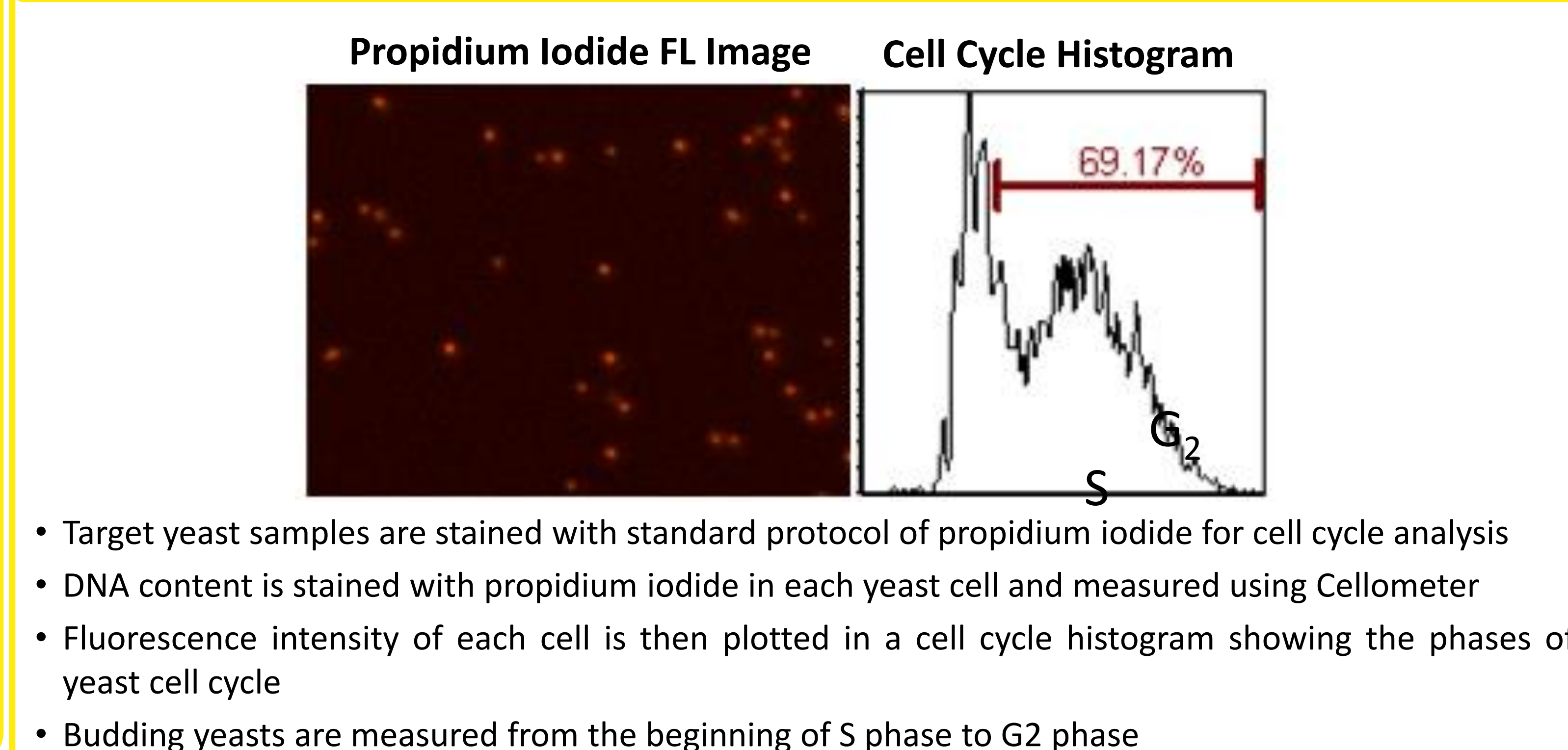
## 3. CELLOMETER IMAGE CYTOMETRY INSTRUMENTATION



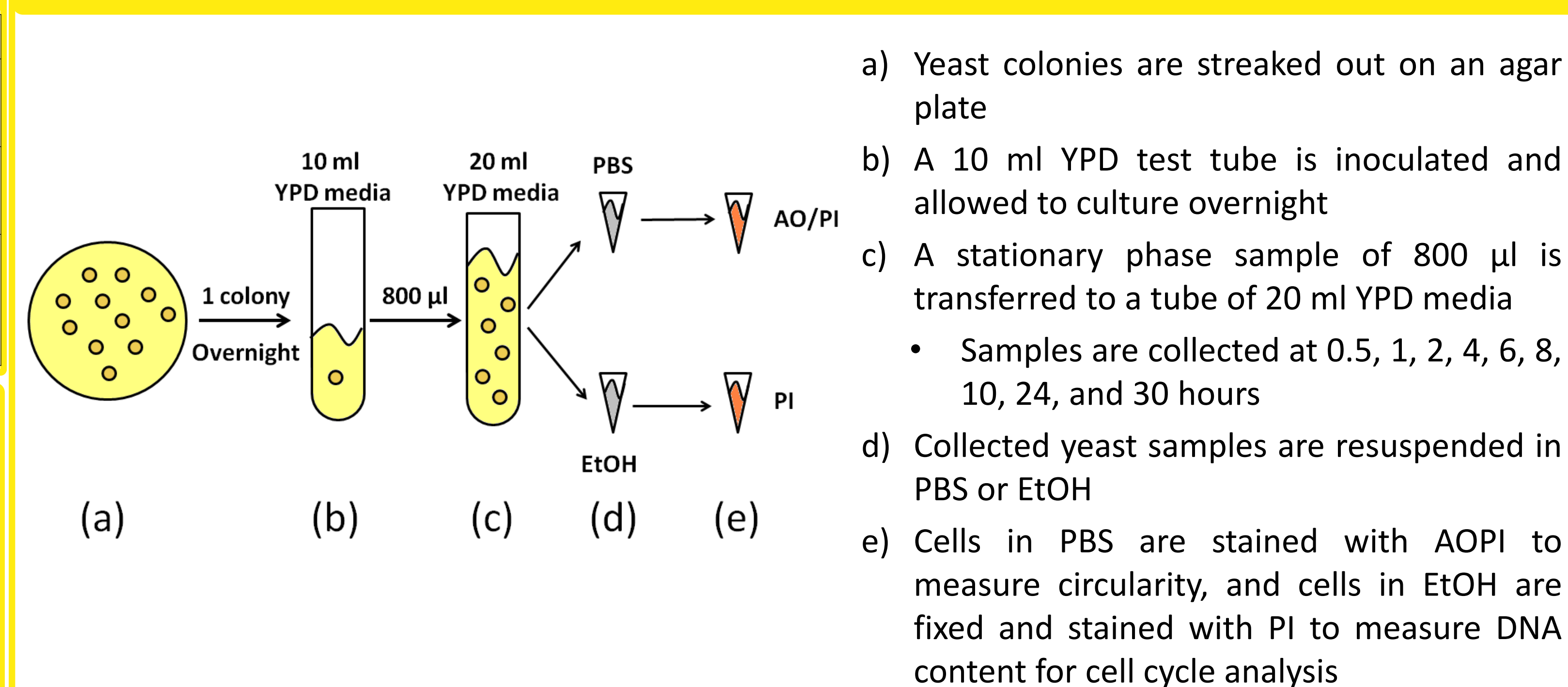
## 4. BUDDING MEASUREMENT VIA MORPHOLOGY USING IMAGE CYTOMETRY



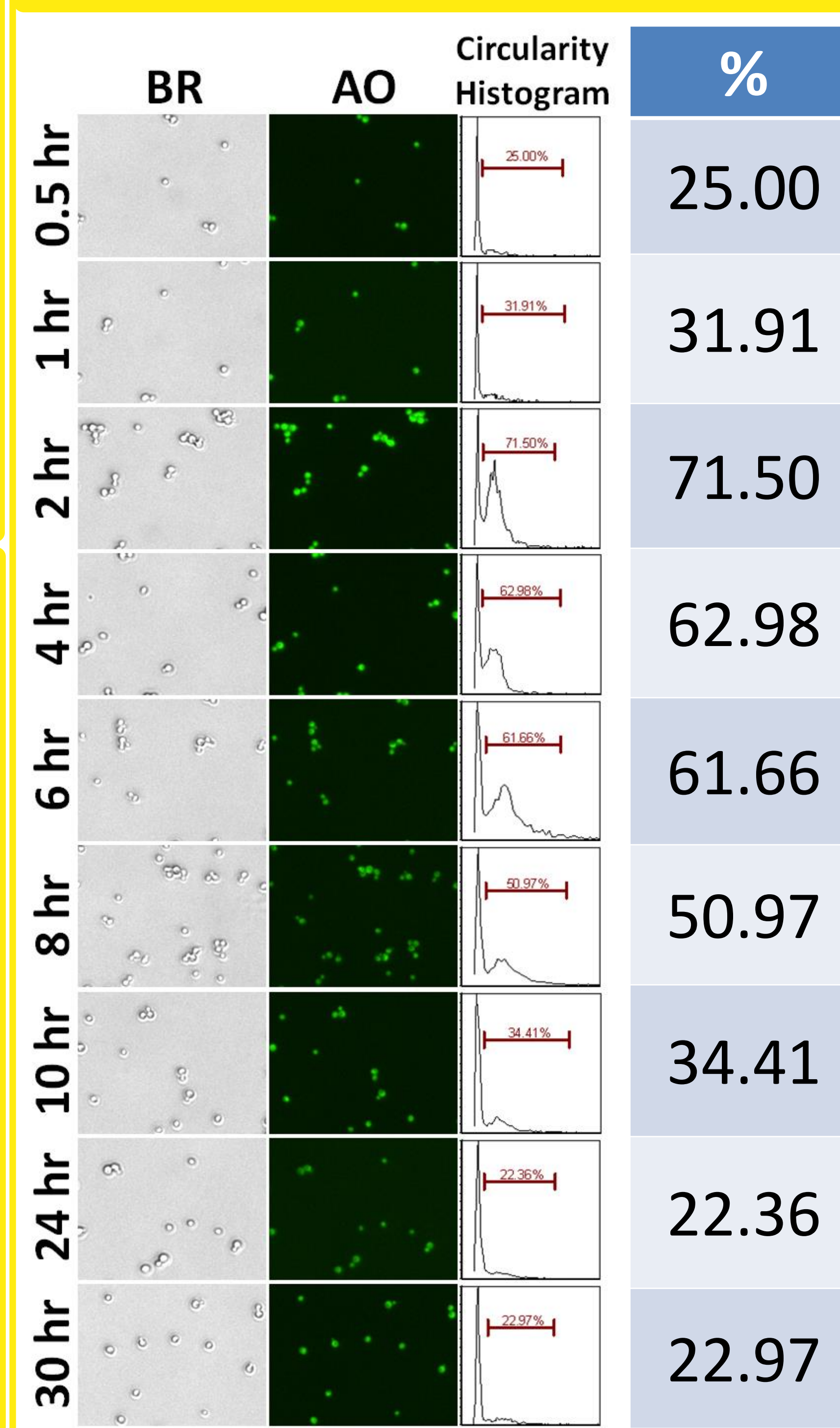
## 5. BUDDING MEASUREMENT VIA CELL CYCLE USING IMAGE CYTOMETRY



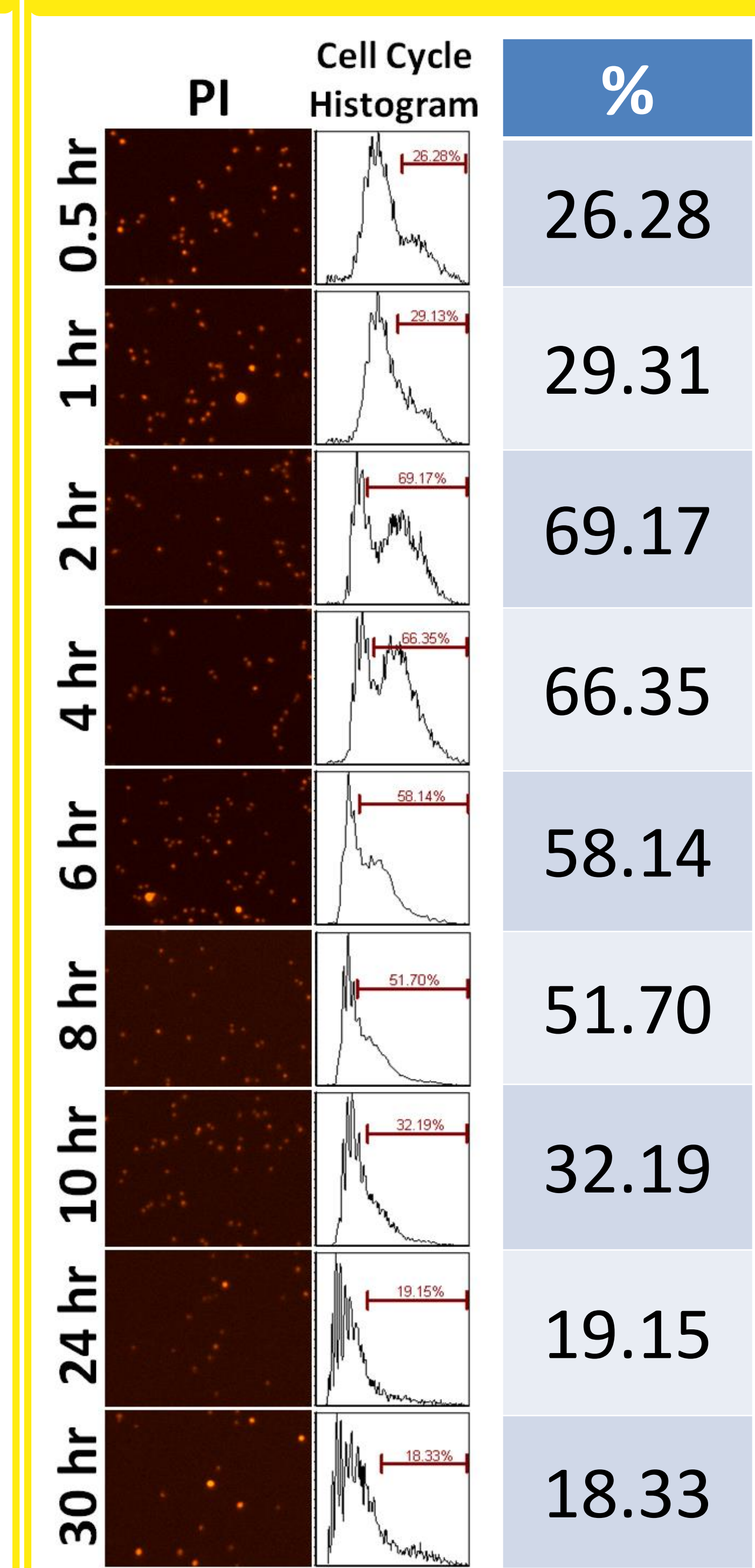
## 6. EXPERIMENTAL PROTOCOL FOR BUDDING MEASUREMENT



## 7. CIRCULARITY ANALYSIS RESULTS



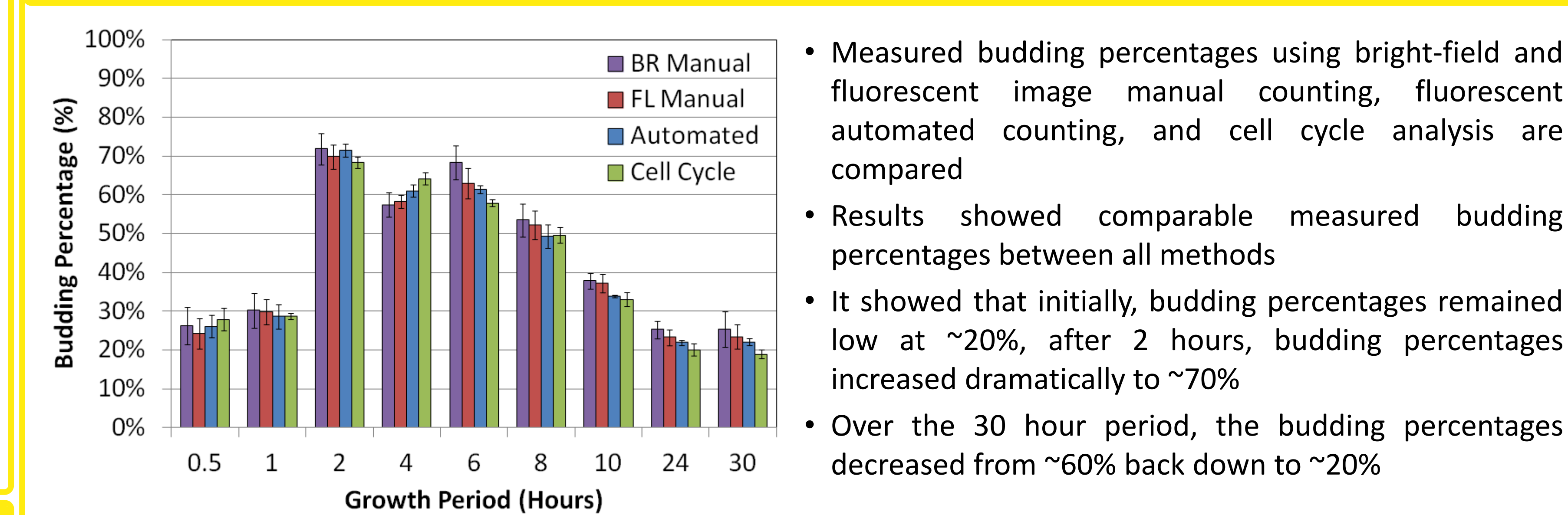
## 8. CELL CYCLE ANALYSIS RESULTS



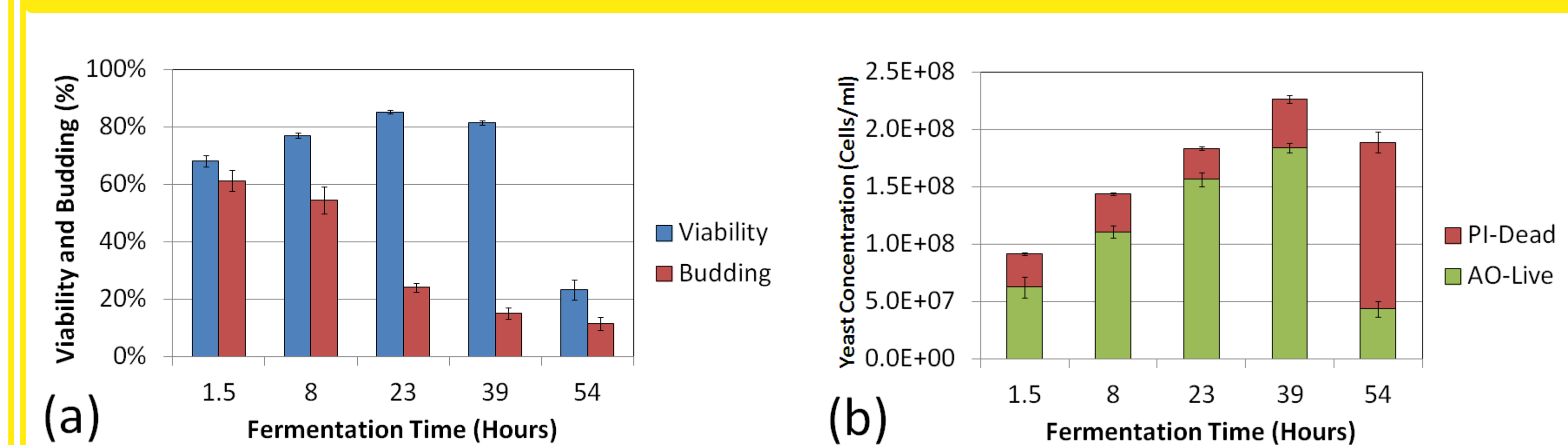
## 9. COMPARISON OF BUDDING MEASUREMENT METHODS

- Bright-field image manual counting**
  - Bright-field images are captured by Cellometer image cytometer and the number of budding yeasts and total yeasts are manually counted
  - Percent budding yeasts are then calculated
- Fluorescence image manual counting**
  - Fluorescent AO images are captured by Cellometer image cytometer and the number of budding and total yeasts are manually counted
  - Percent budding yeast are then calculated
- Fluorescence automated counting via morphology analysis**
  - Fluorescent AO images are automatically counted and the circularity data is exported to FCS Express (De Novo Software)
  - Yeast population with large circularity values are gated to measure the budding percentages
- Cell cycle analysis**
  - Fluorescent PI images are automatically counted and the fluorescence intensity data is exported to FCS Express (De Novo Software)
  - Yeast population with fluorescence intensity values that indicates S phase to G2 phase are gated to measure the budding percentages

## 10. COMPARISON OF TIME-COURSE BUDDING MEASUREMENT RESULTS



## 11. BUDDING MEASUREMENT OF BIOETHANOL FERMENTATION SAMPLES



## 12. CONCLUSION

The development of a fast, accurate, and simple yeast analysis method can improve the current industry standard method, which relies mainly on manual counting using a hemacytometer. With the combination of these three parameters (concentration, viability, budding percentage), the fluorescence-based image cytometry method can be used to easily monitor yeast population characteristics during fermentation, which can allow researchers in the biofuel or brewing industry to improve their fermentation process, as well as improve the efficiency of quality assurance protocols. Future work may also involve supplementing the detection process with a yeast vitality parameter to complete the characterization of yeast during fermentation. We have demonstrated the capability of the image cytometry method for quantifying yeast budding via morphology and DNA content. This automated method can reduce the time required to obtain yeast characteristics in an industry setting, which is of great importance for the optimization of the fermentation process.