



# YOUR AUTOMATED, INTEGRATED BIOPROCESS CULTURE MONITORING SOLUTION



The need for automated bioreactor sampling and sample dilution is being driven by several factors, including increased demand for high-throughput bioprocess development and the progression of high-cell-density cell culture manufacturing processes.

As a result of these factors, industry trends are moving toward the implementation of automated sampling and sample preparation strategies that are designed to bring analytics closer to the operation, eliminate gaps in real-time process monitoring and develop a more rapid, deeper understanding of the bioprocess.

## Vi-CELL BLU Cell Viability Analyzer & Seg-Flow S3 Automated Sampling System

The Vi-CELL BLU cell viability analyzer automates the trypan blue dye exclusion method for cell counting and viability analysis. The Seg-Flow S3 is an automated sampling system from our partners at Flownamics. By integrating them, we've created a precise, rapid cell analysis platform for real-time characterization of bioprocess cell cultures.

This automated integration:

- Removes error-prone manual steps, including sample preparation and sample dilution
- Enables load-and-go sample runs for up to eight bioreactors
- Helps ensure GMP compliance with electronic data management
- Gives you 24/7 control of viable cell density
- Is fully compatible with both automated samples and walkup samples

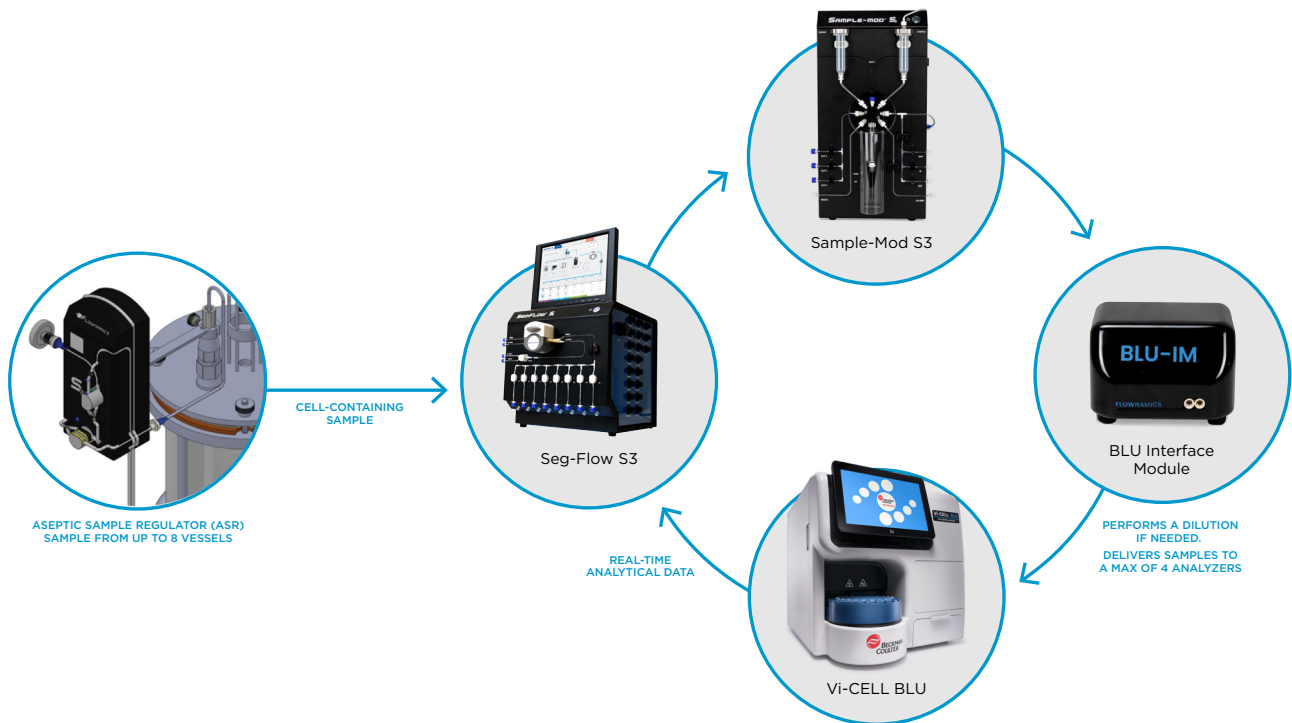
**CULTURE**  
*confidence*

 **BECKMAN  
COULTER**  
*Life Sciences*

# How the Seamless Integration Works

The Seg-Flow S3 pulls continuous and/or scheduled samples from a bioreactor, performs in-line dilution, if needed, and delivers the samples to the Vi-CELL BLU cell viability analyzer. The Vi-CELL BLU provides the analyzed data to the Seg-Flow S3 via the API, which is then sent to the SCADA system. All fluid paths and sample reservoirs are fully cleaned to ensure consistent and accurate measurements for the next sample.

## Sample Flow



To learn more, please visit:

