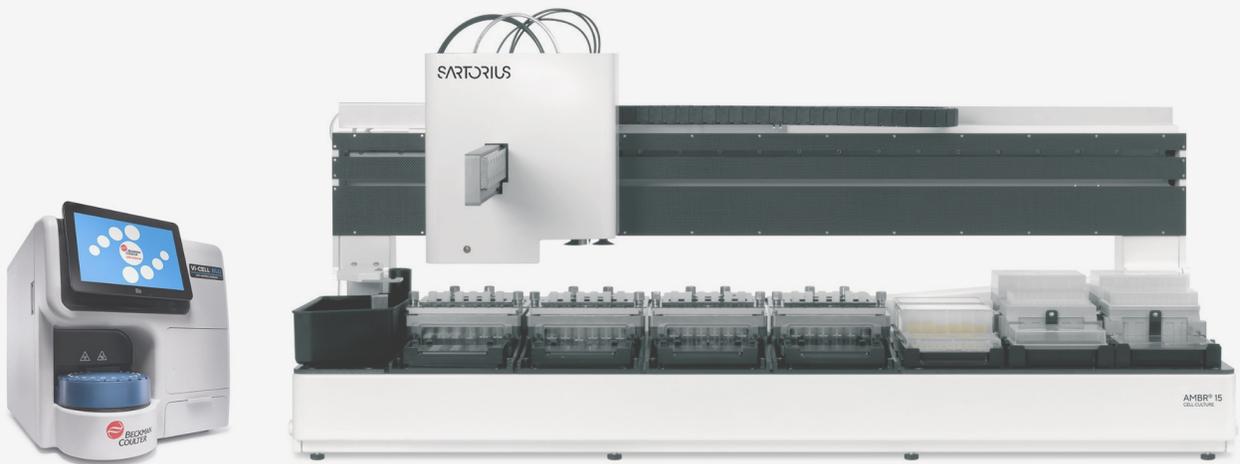


Precision in Every Decision:

Unlock Advanced Cell Culture Analysis with Automation

Integrating the Vi-CELL BLU Cell Viability Analyzer with Sartorius Ambr® 15 Cell Culture Bioreactor System



©Sartorius AG, Image provided by courtesy of Sartorius AG

ACCELERATING
answers

 **BECKMAN
COULTER**
Life Sciences

Introducing a Powerful, Seamless Solution for Cell Culture Monitoring & Process Optimization

With updated software, the Vi-CELL BLU analyzer can now be seamlessly integrated with the Sartorius Ambr[®] Cell Culture Bioreactor System. This integration:

- ✓ Enables walk-away, unattended operation and multiple cell counts 24-hours/day (e.g., overnight and weekends)
- ✓ Improves precision and consistency of results compared to manual sample transfers
- ✓ Controls actions that can be calculated/executed based on cell count (e.g., feed additions)
- ✓ Standardizes sampling protocols between experiments, labs and sites to optimize accuracy and data reproducibility

What's more, the Vi-CELL BLU analyzer can still be used in standalone mode via its 24-position carousel or 96-well plates.

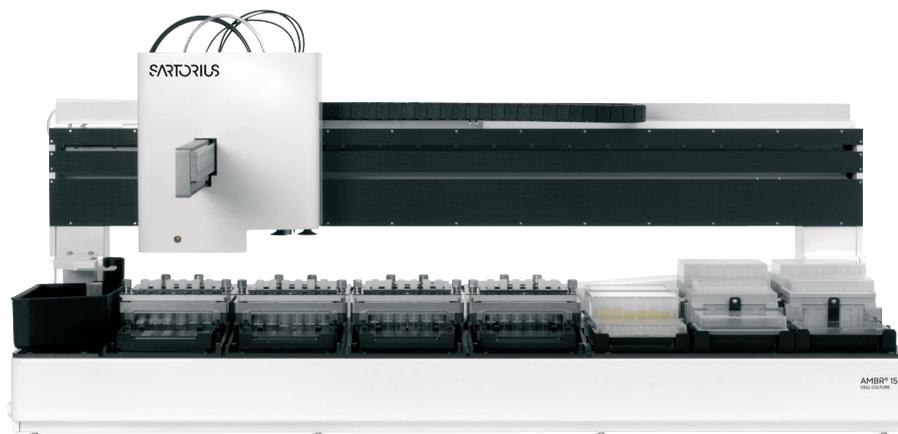
About the Vi-CELL BLU Cell Viability Analyzer

This advanced automated cell viability analyzer is designed to simplify and accelerate cell culture analysis. With its cutting-edge imaging technology and intuitive software, it provides accurate and reliable cell viability and concentration measurements in a matter of minutes.



About the Ambr[®] 15 Cell Culture Bioreactor System

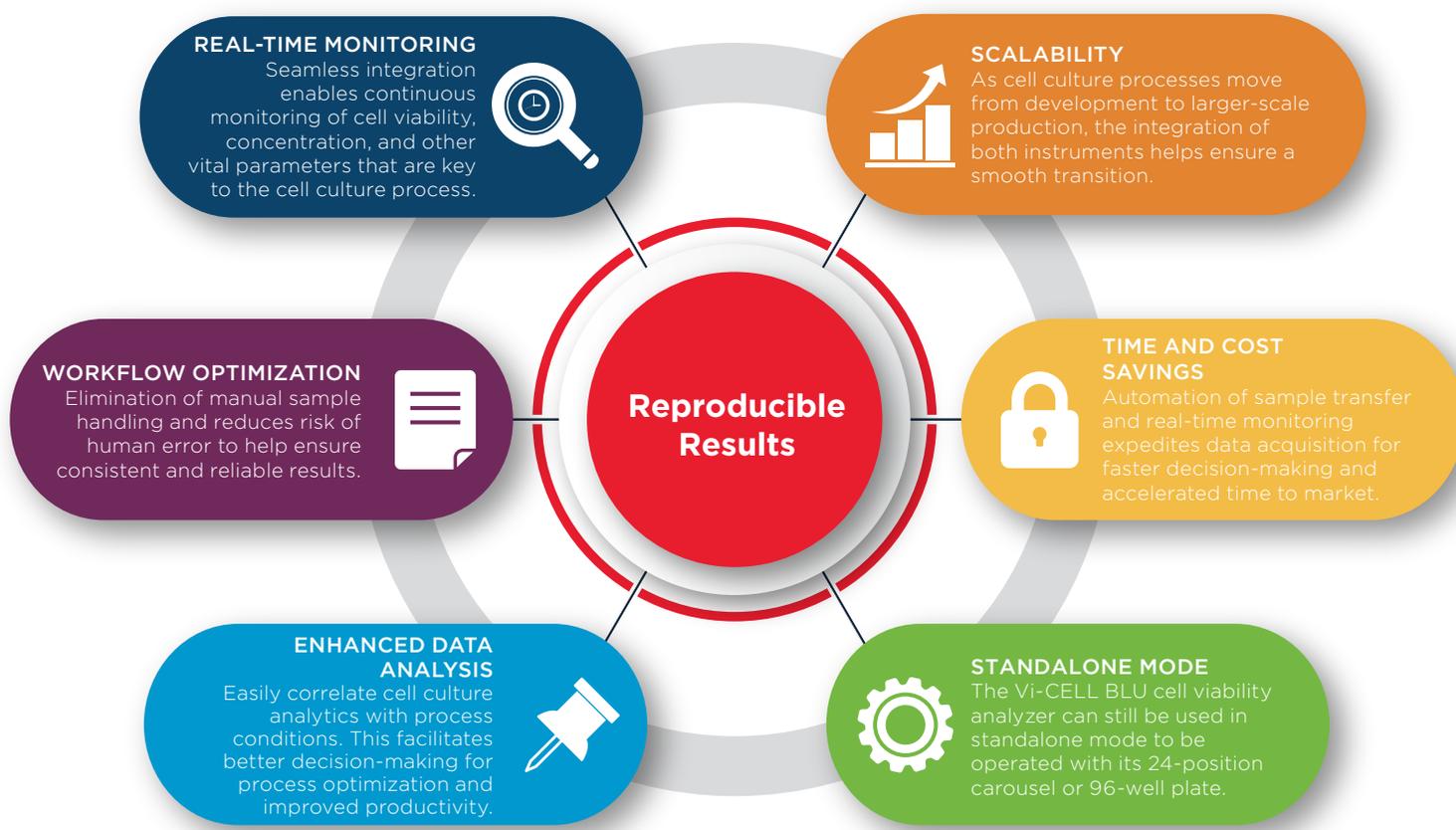
Ambr[®] 15 Cell Culture is a high throughput, automated bioreactor system for 24 or 48 parallel cultivations at the 10 –15 mL microbioreactor scale. This platform enables selection of the best clones, media and feeds under small-scale bioreactor conditions, ensuring a seamless scale-up. Implemented in Cell Line Development (CLD) laboratories worldwide, Ambr[®] 15 Cell Culture makes experimentation more cost effective by saving on labor, facility space, capital, media and consumables.



©Sartorius AG, Image provided by courtesy of Sartorius AG

Unlock Even More Capabilities of your Vi-CELL BLU Analyzer

Integrating Vi-CELL BLU analyzer and Ambr® 15 Cell Culture Bioreactor System leverages the strengths of both instruments to automate cell culture analysis. By combining these two technologies, researchers and bioprocess engineers can unlock benefits such as:



Linking these two robust systems also expands the potential for a wide array of applications, including:

- **BIOPHARMACEUTICAL DEVELOPMENT:** Accelerate development of biopharma processes by optimizing cell culture conditions and ensuring high cell viability and productivity.
- **CELL LINE DEVELOPMENT:** Screen and select high-performing cell lines more efficiently and gain real-time insights into cell growth, viability and productivity to identify the most promising candidates.
- **PROCESS OPTIMIZATION:** Achieve better process control and optimization by leveraging real-time monitoring capabilities of these two instruments. You'll see the impact of process parameters on cell viability, concentration and product quality, which can help improve the effectiveness of your processes.
- **SCALE-UP STRATEGIES:** Seamlessly transition from small-scale experiments with the Ambr® 250 High Throughput system to larger-scale bioreactors by leveraging insights gained with the Vi-CELL BLU analyzer. This enables a smooth scale-up process and the maintenance of process performance at different production scales.

Ordering Information

For current users of the Vi-CELL BLU Analyzer:

Please use the part numbers below to add the Ambr Integration to your existing Vi-CELL BLU instrument.

Part Number	Description
D09687	Sartorius Ambr Automation Integration Kit - Existing Vi-CELL BLU

For NEW users of the Vi-CELL BLU Analyzer:

Part Number	Description
D09686	Sartorius Ambr Automation Integration Kit - New Vi-CELL BLU



For more information

visit beckman.com/cell-counters-and-analyzers/vi-cell-blu/ambrconnector



Product not for use in diagnostic procedures.

© 2024 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 the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at beckman.com
2023-GBL-EN-104632-v1

