

BIOSERO 

SARTORIUS

 **GoSimple**  
**Workcell for**  
**Sartorius Octet<sup>®</sup>**



# GoSimple Workcell for Sartorius Octet®



The GoSimple Workcell for the Sartorius Octet® is a compact, integrated lab automation solution for increasing the throughput and walkaway time of your Octet® BLI platform. The GoSimple Workcell comes with predefined hardware, a benchtop table setup, and established methods for quick setup and simple operator use.

## WHAT IT CAN DO FOR YOU



### Deploy Quickly with Pre-Configured Methods and Coordinated Support

Installs on an existing lab bench with minimal time and effort. Pre-programmed methods require no development. Get 12 months of GoCare for technical support and access to our online training academy.



### Extend Walkaway Time to Maximize Instrument Use

Load up to 44 MTP plates to extend the walkaway time for Octet® BLI systems — enabling longer, unattended runs and reducing costly instrument idle time.



### Improve Data Security and Traceability

Ensure high quality results with 21CFR11 Manager compliance module — supporting secure data handling and traceability.



### Increase Operator Efficiency

Offload tedious, repetitive tasks, and leave more time for knowledge work and creative thinking.

**NOTE:** The GoSimple Workcell is available for the Octet® RH96 and RH16.

## BIOSERO'S PROVEN PROCESS SUPPORTS INDUSTRY LEADERS

1,000+ Laboratories

Serving 20 of the 20 Top Pharma

Providing GBG-Scheduled Workcells Since 2012

Supporting Scientists with Automation Since 2003

Providing Orchestrated Solutions Since 2018

3,000+ GBG Installs

# What's Included in the GoSimple Workcell for Sartorius Octet® BLI

**SUPPORTS 1X OCTET® RH96 OR RH16, PURCHASED SEPARATELY FROM SARTORIUS**

## 1. Hardware

- 1 2x Plate Hotels
- 2 1x Robotic Arm
- 3 1x E-Stop
- 4 1x Shaker
- 5 1x Regrip Nest
- 6 1x Biosensor Tray Hotel
- 7 1x Baseplate with System Controller

Optional: Computer

## 2. Integration Software

- Green Button Go Scheduler Basic
- Preconfigured methods for quick setup and simple operator use
- 21CFR11 Manager Extension

## 3. Services

- Installation and setup are provided at lab site
- Biosero GoCare Plus Support Subscription for 12 months from date of shipping, which includes access to Biosero Academy



# Do More with Automation

## GETTING STARTED IS A BREEZE WITH GUIDED USER INPUT SCREENS



Wizard built with existing Green Button Go tools



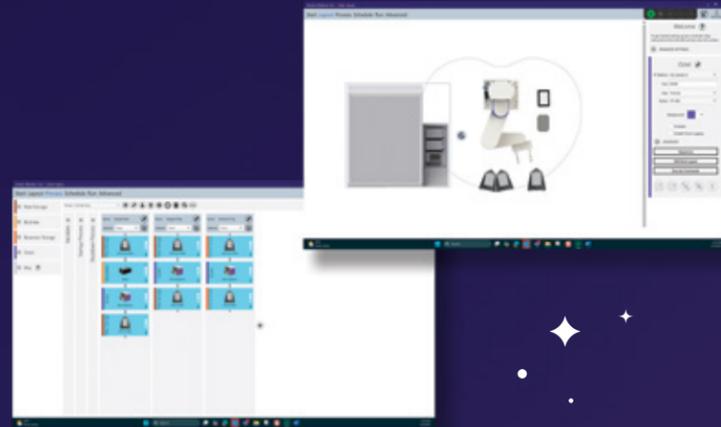
No prior or specialized knowledge required to use



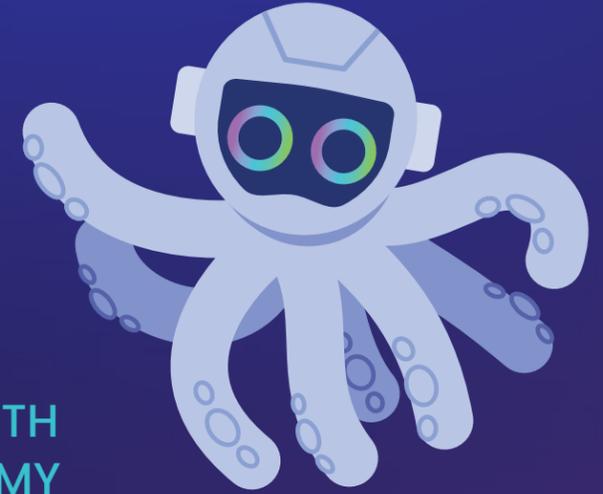
Accepts operator input with parametrized fields



Simple entry point into integrations



## BUILD YOUR TEAM'S CAPABILITIES WITH BIOSERO'S ONLINE TRAINING ACADEMY



Learn how to modify your pre-configured Green Button Go Methods

Learn about advanced uses of Green Button Go to plan future automation expansions

Learn about how to teach the robotic arm for placing additional instruments

Get access to wiki articles on Frequently Asked Questions (FAQs)



# What GoSimple Can Do For You

## GOSIMPLE WORKCELL FOR THE SARTORIUS OCTET® BLI SUPPORTS A WIDE RANGE OF APPLICATION AREAS

The Biosero GoSimple Workcell for Octet® BLI systems supports increased throughput across many application areas.



Quantification



Analyte Screening



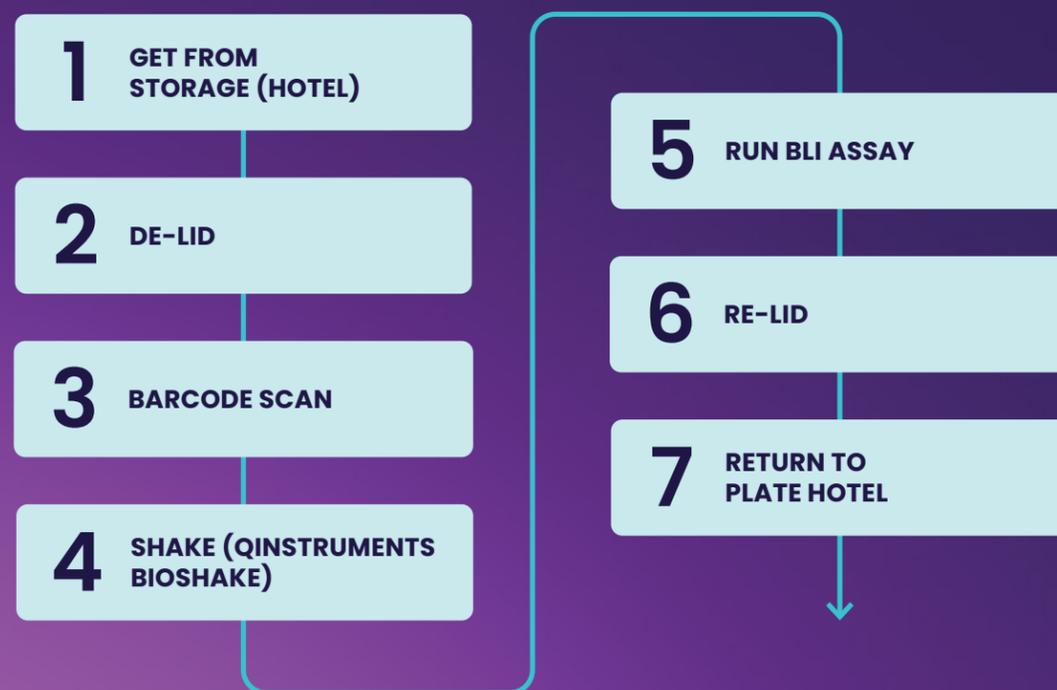
Clone Screening



Cell Line Development

## GoSimple Workcell for the Sartorius Octet® BLI Automation Workflow

The GoSimple Workcell for Octet® BLI supports the following workflow steps. These steps can be scheduled to run in continuous cycles.



## GOSIMPLE WORKCELL FOR SARTORIUS OCTET® COMPATIBILITY



### OCTET® RH96

#### 96-Channel Ultra High Throughput

Fluidic-free biolayer interferometry (BLI) platform that monitors 96 biosensors simultaneously



### OCTET® RH16

#### High Throughput Platform for Biologics Development

Fluidic-free biolayer interferometry (BLI) platform ideal for the rapid characterization and development of biologics drug molecules

## SUPPORT FOR YOUR AUTOMATION SOLUTION



### Biosero GoCare Subscription

Biosero ensures that you have visibility and control at your fingertips to maximize your investment.

#### Training at Your Own Pace

Self-serve training courses and certification to learn at your own pace

#### Virtual Troubleshooting

For help via email or phone with challenges in your automation

#### 24/7 Access to Biosero Portal

View and manage your users, organizations, licenses, and subscriptions any time of day

#### Easy to Upgrade and Patch

Easy access to software upgrades, patches, and driver updates

## TESTIMONIAL

“The [Octet] is already automated by itself to a certain extent, but what we’re seeing is a lot of labs and companies that need some more. They want to analyze thousands of samples instead of hundreds.

So this automation will allow that, bringing efficiency into these labs.

And then Biosero is a great partner to do that, because a lot of our customers have mentioned they like the Biosero products as well.

– **NILSHAD SALIM,**  
Product Manager,  
Sartorius

## Contact Biosero Today to Start Your Automation Journey

Biosero and Green Button Go are trademarks owned and/or registered by Biosero in the U.S. and/or other countries. Sartorius®, the Sartorius Logo, and Octet® are registered trademarks of Sartorius AG or any of its affiliates.