

# Top Ten Reasons to Choose Sartorius Octet® BLI Technology to Accelerate Your Research, Development and Manufacturing

1

## Sartorius—A Trusted Partner for the Biopharmaceutical Industry and Research Laboratories



With a track record of over 150 years of success, we enable pharmaceutical researchers and lab personnel to simplify workflows, minimize human errors, and speed up their work to help bring medications safely and quickly to the patient.

2

## An Eniable Range of Applications



Octet® BLI provides the highest flexibility for biophysical characterization in biopharma, industrial, and academic research. The applications of this industry standard for label-free, real-time analysis reach from discovery, through development, to manufacturing, to final QC.

3

## Measure Protein-Protein and Protein-Drug Interactions Label-Free Under Native Conditions



Octet® BLI uses optical biosensors to measure molecular interactions in parallel, without the need for expensive detection reagents, and with temperature control, the interactions can also be measured at physiological temperature.

4

## Experimental Simplicity to De-Risk Your Journey and Simplify Progress



A microfluidic-free design allows for unprecedented time and cost savings for binding kinetics, affinity and analyte concentration measurement with high precision, accuracy, and a broad dynamic range.

5

## ISO 9000 Certified Manufacturing



Rigorous manufacturing standards combined with robust microplate-based design and disposable biosensors reduces the risk of instrument downtime.

**6**

## Boost Your Productivity by Speeding up Assay Performance



Intuitive workflows, walkaway automation, no system maintenance between experiments, and no sample loading time provides the fastest assay run times and minimizes hands-on time to really generate data.

**7**

## Perfect for Multi-User Environments



Octet® BLI systems are known for their reliable, user-friendly, trouble-free operation, and have quickly become the go-to systems in both academic and industrial labs because of their ease of use, robustness, and application flexibility.

**8**

## Analyze Your Toughest Samples Faster



Octet® BLI allows you to analyze complex and unpurified samples in just 2 minutes. Parallel processing coupled with minimal sample preparation provides unprecedented throughput. Process up to 96 crude lysates simultaneously for analyte quantitation in just 2 minutes.

**9**

## Comprehensive Insight into Expression Levels, Stability, Potency, and Glycosylation from One Sample



Combine orthogonal data, such as titer data with glycosylation data, to make more informed decisions and facilitate the selection of clones that produce proteins with favorable expression profiles during cell line development.

**10**

## Walk-In Ready



Assays performed on Octet® BLI systems do not require prior preparation such as cleaning, priming, or equilibration. Simply design the experiment, load the samples and run protein stability, purity, and functionality assays on one system. Enabling you to be confident in handing over high-quality biologics faster.

### Germany

Sartorius Lab Instruments  
Phone +49 551 308 0

### USA

Sartorius Corporation  
Phone +1 631 254 4249  
Toll-free +1 800 635 2906