



# High Precision Streptavidin (SAX) Biosensors

*For high precision quantitation and kinetic characterization*

## Key features

- Rapid and stable capture of biotinylated molecules
- High precision quantitation and kinetic analyses
- QC-tested to meet precision-controlled CV specification of 4%

## Overview

Streptavidin-coated surfaces are widely used as a simple and straightforward method of molecular immobilization. Utilized with Bio-Layer Interferometry (BLI), Streptavidin biosensors enable quick and easy modification and customization of the biosensor with any biotin-tagged molecule for quantitative and kinetic measurements. The Dip and Read™ High Precision Streptavidin Biosensor (SAX) was specifically developed and qualified for applications in downstream drug discovery and regulated environments that have more stringent assay precision requirements. SAX biosensors are QC-tested at ForteBio to meet our precision-controlled coefficient of variation (CV) specification of <4%.

## High precision assays

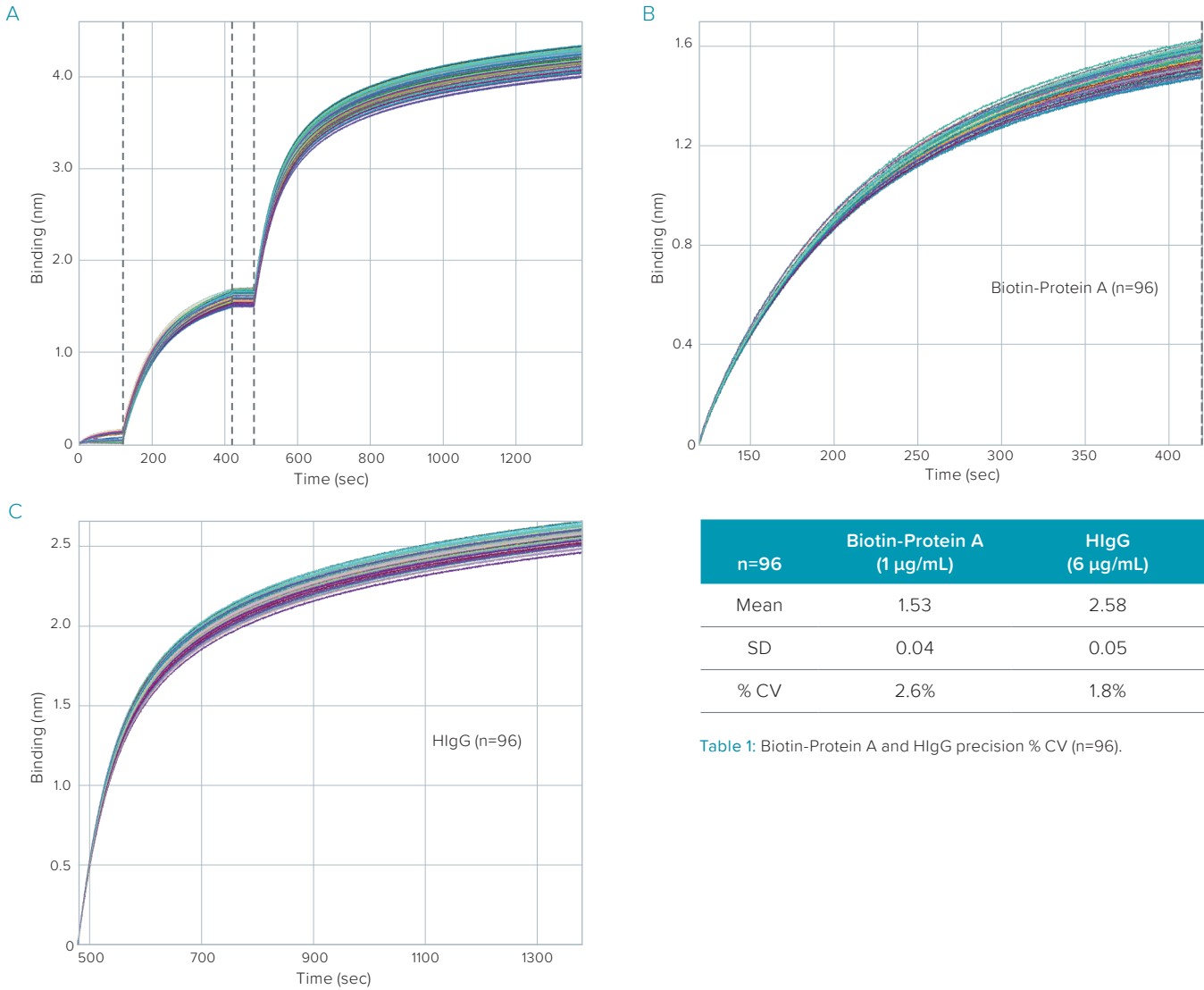
Characterization of protein interactions provides valuable information in every stage of a drug candidate's development and processing. The High Precision Streptavidin Biosensor (SAX) enables researchers to perform direct immobilization of biotin-labeled proteins for both kinetics characterization and quantitation measurements on Octet® systems. Intra-assay precision is significantly improved. Researchers can now develop and perform custom quantitation or stringent kinetic assays with high confidence in the reliability and consistency of results.

## Range of applications

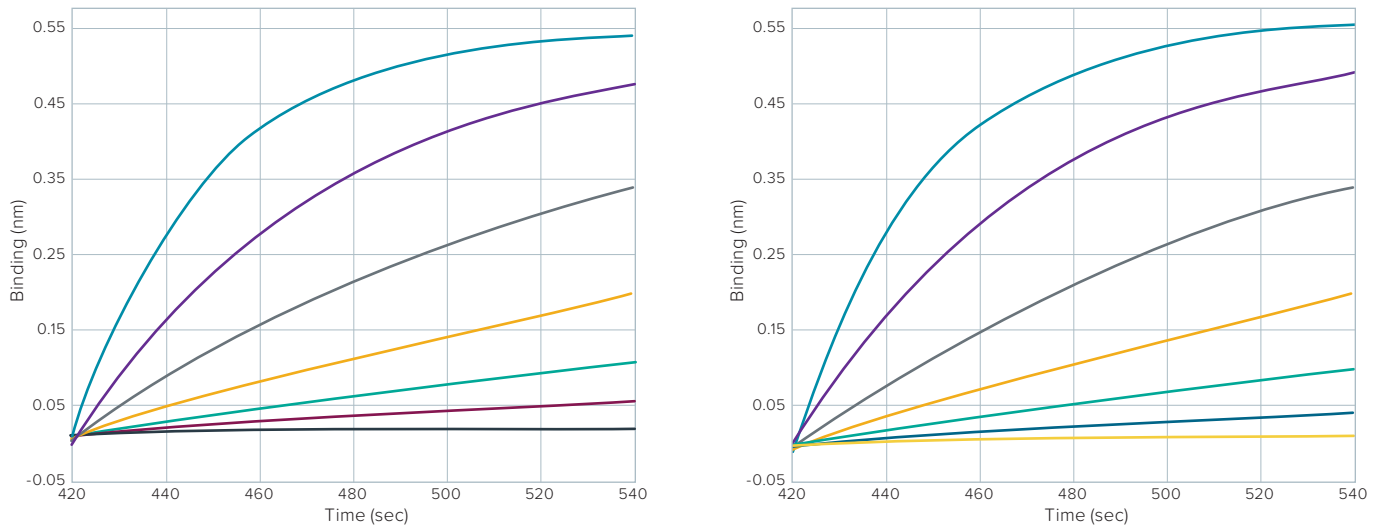
The High Precision Streptavidin Biosensor offers researchers unparalleled ease of use for applications such as:

- Highly precise titer determination
- Activity assays in manufacturing
- Product release QC





**Figure 1:** Data showing an experiment performed with High Precision Streptavidin biosensors (SAX), n=96. (A) Raw data showing the consecutive steps in the experiment where the 96 SAX biosensors are dipped into buffer for a baseline, followed by loading of the biotinylated ligand (biotinylated Protein A), buffer again for a second baseline step, and binding of the analyte (human IgG) to the surface-immobilized ligand (biotinylated Protein A). (B) Data showing very consistent loading of biotinylated Protein A onto all 96 SAX biosensors. (C) Data showing very consistent binding of human IgG onto all 96 biotinylated Protein A-coated SAX biosensors.



**Figure 2:** Binding experiment of prostate-specific antigen (PSA) showing high precision data with High Precision (SAX) Streptavidin biosensors. Biotinylated anti-PSA was loaded onto four different sets of High Precision Streptavidin biosensors, and the subsequent binding of PSA was measured on each set of anti-PSA coated biosensors, showing data consistency and precision.

## Ordering information

Part No.	UOM	Description
18-5117	Tray	One tray of 96 High Precision Streptavidin (SAX) Biosensors
18-5118	Pack	Five trays of 96 High Precision Streptavidin (SAX) Biosensors
18-5119	Case	Twenty trays of 96 High Precision Streptavidin (SAX) Biosensors

For more information about the Octet and BLItz platforms for label-free, real-time detection of biomolecular interactions, applications, and services, visit [www.fortebio.com](http://www.fortebio.com) or contact us directly.



**FortéBio**  
47661 Fremont Boulevard  
Fremont, CA 94538  
888.OCTET-75 or 650.322.1360  
[fortebio.info@moldev.com](mailto:fortebio.info@moldev.com)

**FortéBio Analytics (Shanghai) Co., Ltd.**  
No. 88 Shang Ke Road  
Zhangjiang Hi-tech Park  
Shanghai, China 201210  
[salesops.china@moldev.com](mailto:salesops.china@moldev.com)

**Molecular Devices (UK) Ltd.**  
660-665 Eskdale  
Winnersh Triangle  
Wokingham, Berkshire  
RG41 5TS, United Kingdom  
+44 118 944 8000  
[uk@moldev.com](mailto:uk@moldev.com)

**Molecular Devices (Germany) GmbH**  
Bismarckring 39  
88400 Biberach an der Riss  
Germany  
+ 00800 665 32860