# SVISCISVS

## Product Datasheet

## Octet<sup>®</sup> High Precision Streptavidin 2.0 (SAX2) Biosensors

For High Precision and Reproducible Kinetic Characterization and Custom Quantitation

### Key Features

- QC-tested to limit biotinylated ligand loading variance, providing lot-to-lot biosensor consistency
- Minimal lot-to-lot biosensor variance enables confident detection of variance from samples
- Rapid and stable capture of biotinylated molecules



## 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 2.0 (SAX2) is specifically developed and qualified for applications in downstream drug discovery and regulated environments that have stringent assay precision requirements.

## Quick Facts

- Within lot CV (binding of biotin-tagged molecule): < 4%
- Lot-to-lot CV (binding of biotin-tagged molecule): < 20%</p>

SAX2 Biosensors are QC-tested at Sartorius to meet our precision-controlled lot-to-lot coefficient of variation (CV) specification of < 20% for binding of biotin-tagged molecule onto the biosensor. Within a lot, SAX2 Biosensors are QC-tested to meet a CV of < 4%.

## Example Data

#### А Loading Association Dissociation 2.0 1.8 1.6 1.4 -1.2 -Binding (nm) 0.8 0.6 0.4 0.2 400 800 1600 2000 2400 2800 1200 Time (sec) В Global fitting for kinetics 1.0 0.8 Binding (nm) 0. 0.4 0.2 0 200 400 600 1000 1200 1600 1800 800 1400 Time (sec)

#### Kinetics Assay Using 3 Different SAX2 Lots

Table 1: CV calculations for loading of biotin-Fcy receptor onto SAX2 Biosensor shown in Figure 1.

	Lot 1	Lot 2	Lot 3	SAX2 Specification	
Loading CV within lot	1.5%	1.3%	1.3%	< 4%	
Loading CV across lots		1.5%		< 20%	
$K_{\rm D}$ CV within lot	3.4%	2.7%	4.1%	Dependent on the	
$K_{\rm D}$ range across lots		4%	4% sample pair and		

#### **Custom Quantitation Assay**



Figure 1: (A) Kinetic analysis of the interaction between a ligand biotin-Fcy Receptor (~55 kDa) and analyte hIgG (150 kDa) with SAX2 Biosensors, overlaying three biosensor lots and raw data aligned at the ligand loading step. (B) Data was processed and curve fitted using a 1:1 binding model. Yellow lines represent fitted curves; other colors represent raw curves.

Figure 2: Detection of PSA standards and unknowns after Biotin-anti-PSA loading (30 µg/mL, 400 seconds) using SAX2 on the Octet® RED384 system with assay parameters (1000 rpm, 120 seconds). (A) PSA dose response and unknown response. (B) Representative resulting calibration curves and unknowns prediction from A.

PSA	Expected concentration (µg/mL)	Expected concentration (µg/mL) (n=3)	% CV (n=3)	% Recovery (n=3)
Standards	30	30.2	4.5%	100.6%
	20	19.9	6.5%	99.6%
	10	10.0	0.8%	100.4%
	5	5.0	2.6%	99.7%
	2.5	2.5	4.5%	100.6%
	1.25	1.2	4.3%	99.6%
	0.625	0.6	6.3%	100.0%
Unknowns	15	16.1	3.1%	107.5%
	3	3.1	0.9%	104.3%

Table 2: Average calculated concentration, % CV and Recovery for PSA binding curves in Figure 2.

### What SA Biosensor is Right for Me?



## Range of Applications

SAX2 Biosensors are designed and optimized for applications that require minimal variation from consumables to allow for accurate measurements of variance from samples. They are best suited for:

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

## Ordering Information

Part No.	UOM	Description
18-5136	Tray	One tray of 96 Octet® High Precision Streptavidin 2.0 (SAX2) Biosensors
18-5137	Pack	Five trays of 96 Octet® High Precision Streptavidin 2.0 (SAX2) Biosensors
18-5138	Case	Twenty trays of 96 Octet® High Precision Streptavidin 2.0 (SAX2) Biosensors

#### Germany

USA

Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Strasse 20 37079 Goettingen Phone +49 551 308 0 Sartorius Corporation 565 Johnson Avenue Bohemia, NY 11716 Phone +1 888 OCTET 75 Or +1 650 322 1360

For further contacts, visit www.sartorius.com/octet-support