



# Octet® BLI Biosensor Selection Guide

Simplifying Progress

**SARTORIUS**

## Octet® BLI Biosensors: Overview

Biosensor	Description	Intended Use <sup>1</sup>		Application	Specificity	Protein Tag (capture)	Regenerable	Suggested Molecule	
<a href="#">AHC</a>	Anti-Human Fc-Capture			Human IgG or Fc-fusion capture	Human	IgG / Fc Domain			
<a href="#">AHC2</a>	Anti-Human Fc-Capture 2nd Generation			Human IgG or Fc-fusion capture	Human	IgG / Fc Domain			
<a href="#">AHQ</a>	Anti-Human IgG Fc			Human IgG or Fc-fusion capture	Human	IgG / Fc Domain			
<a href="#">AMC</a>	Anti-Mouse Fc-Capture			Mouse IgG or Fc-fusion capture	Mouse	IgG / Fc Domain			
<a href="#">AMC2</a>	Anti-Murine IgG Capture 2nd Generation			Mouse IgG or F(ab')2 capture	Mouse	IgG / F(ab')2			
<a href="#">AMQ</a>	Anti-Murine IgG Fv			Mouse IgG or F(ab')2 capture	Mouse	IgG / F(ab')2			
<a href="#">APS</a>	Aminopropylsilane			Lipids, liposome and hydrophobic proteins capture	Various	N/A			
<a href="#">ARC</a>	Anti-Rabbit Fc-Capture			Rabbit IgG or Fc-fusion capture	Rabbit	IgG / Fc Domain			
<a href="#">AR2G</a>	Amine Reactive 2G			Amine coupling	Various	N/A			
<a href="#">FAB2G</a>	Anti-Human Fab-CH1 2nd Generation			Human IgG or Fab-CH1 capture	Human	CH1			
<a href="#">GST</a>	Anti-GST			GST capture	Various	GST			
<a href="#">HIS1K</a>	Anti-Penta-HIS			His-tagged proteins capture	Various	His tag			
<a href="#">HIS2</a>	Anti-HIS			His-tagged proteins capture	Various	His tag			
<a href="#">NTA</a>	Ni-NTA			His-tagged proteins capture	Various	His tag			
<a href="#">ProA</a>	Protein A			Various species IgG capture	Various	IgG			
<a href="#">ProG</a>	Protein G			Various species IgG capture	Various	IgG			
<a href="#">ProL</a>	Protein L			Various species IgG capture	Various	IgG			
<a href="#">SA</a>	Streptavidin			Immobilization of biotinylated molecules	Various	Biotin, AviTag™			
<a href="#">SAX</a>	High Precision Streptavidin			Immobilization of biotinylated molecules	Various	Biotin, AviTag™			
<a href="#">SAX2</a>	High Precision Streptavidin 2.0			Immobilization of biotinylated molecules	Various	Biotin, AviTag™			
<a href="#">SSA</a>	Super Streptavidin			Immobilization of biotinylated molecules	Various	Biotin, AviTag™			
<a href="#">AAVX</a>	AAV Quantitation			Quantitation of AAV Capsids	Human	N/A			

Kinetics     
 Quantitation     
 Glycan Screening     
 Impurity Testing     
 Yes for Kinetics     
 Yes for Quantitation     
 No for Quantitation     
 Protein and analyte dependent





















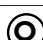

Proteins     
 Antibodies     
 Small Molecules     
 Lipid | liposome     
 DNA

<sup>1</sup> Biosensors are developed, manufactured, and QC is performed for their intended applications; using biosensors outside their intended purpose requires user validation

## Octet® BLI Consumables: In Depth

Octet® Consumables	Description	Intended Use <sup>2</sup>	Application	Octet® BLI System Quantitation Dynamic Range <sup>1</sup>			Regeneration
				Octet® QK <sup>®</sup> , QK384 <sup>®</sup> , RH96 ≥32 Channel	Octet® RED96 <sup>®</sup> , K2 <sup>®</sup> , R2, R4, R8, RH16, RH96 8 or 16 Channel	Octet® N1	
Biosensors							
<a href="#">AHC</a> (Cat. Nos. 18-5060, 18-5063, 18-5064)	Anti-Human Fc-Capture		Capturing human IgG's or human Fc-fusion proteins for kinetic analysis with various analytes	N/A	N/A	N/A	
<a href="#">AHC2</a> (Cat. Nos. 18-5142, 18-5143, 18-5144)	Anti-Human Fc-Capture 2nd Generation		Capturing human IgG's or human Fc-fusion proteins for both kinetic and quantitation analysis	0.5-2000 µg/mL	0.1-2000 µg/mL	0.5-4000 µg/mL	
<a href="#">AHQ</a> (Cat. Nos. 18-5001, 18-5004, 18-5005)	Anti-Human IgG Fc		Quantitation measurements of human IgG's or human Fc-fusion proteins	0.025-200 µg/mL	0.01-200 µg/mL	0.25-500 µg/mL	
<a href="#">AMC</a> (Cat. Nos. 18-5088, 18-5089, 18-5090)	Anti-Mouse Fc-Capture		Capturing mouse IgG's or mouse Fc-fusion proteins for kinetic analysis with various analytes	N/A	N/A	N/A	
<a href="#">AMC2</a> (Cat. Nos. 18-5163, 18-5164, 18-5165)	Anti-Murine IgG Capture 2nd Generation		Capture of murine IgG's or F(ab'2) for both kinetic and quantitation analysis	0.025-8000 µg/mL	0.025-8000 µg/mL	0.025-8000 µg/mL	
<a href="#">AMQ</a> (Cat. Nos. 18-5022, 18-5023, 18-5024)	Anti-Murine IgG Fv		Quantitation measurements of mouse IgG's or mouse F (ab')2	0.05-200 µg/mL	0.025-200 µg/mL	0.5-500 µg/mL	
<a href="#">APS</a> (Cat. Nos. 18-5045, 18-5046, 18-5047)	Aminopropylsilane		Binding measurement of lipids, liposomes, hydrophobic proteins that don't have other methods of surface attachment	N/A	N/A	N/A	
<a href="#">ARC</a> (Cat. Nos. 18-5168, 18-5169, 18-5170)	Anti-Rabbit Fc-Capture		Capturing rabbit IgG's or rabbit Fc-fusion proteins for both kinetic and quantitation analysis	0.05-4000 µg/mL	0.05-4000 µg/mL	0.05-4000 µg/mL	
<a href="#">AR2G</a> (Cat. Nos. 18-5092, 18-5093, 18-5094)	Amine Reactive 2G		Covalently immobilizing any molecule with a terminal amine group for all kinetic analyses	N/A	N/A	N/A	
<a href="#">FAB2G</a> (Cat. Nos. 18-5125, 18-5126, 18-5127)	Anti-Human Fab-CH1 2nd Generation		Kinetic analysis of human Fab fragments and IgG with target antigen, Fc receptors, or other analytes. Quantitation of Fab and IgG.	Analyte dependent, typically 0.5-1000 µg/mL	Analyte dependent, typically 0.5-1000 µg/mL	Analyte dependent, typically 0.5-1000 µg/mL	
<a href="#">GST</a> (Cat. Nos. 18-5096, 18-5097, 18-5098)	Anti-GST		Quantitation of GST-tagged proteins, direct capturing of GST-tagged proteins for kinetic analyses with analytes	Protein dependent, typically 0.1-2000 µg/mL	Protein dependent, typically 0.1-2000 µg/mL	Protein dependent, typically 0.5-1000 µg/mL**	
<a href="#">HIS1K</a> (Cat. Nos. 18-5120, 18-5121, 18-5122)	Anti-Penta-HIS		Capture of His-tagged proteins for kinetic analysis with target analytes. Quantitation of His-tagged proteins in buffer, media or diluted lysate. Biosensor is pre-coated with Penta-His antibody from Qiagen.	Protein dependent, typically 0.25-200 µg/mL*	Protein dependent, typically 0.25-200 µg/mL*	Protein dependent, typically 10-200 µg/mL*	
<a href="#">HIS2</a> (Cat. Nos. 18-5114, 18-5115, 18-5116)	Anti-HIS		Quantitation of HIS-tagged proteins in crude matrices or buffer or column eluent (pre-coated with anti-His Ab from MBS)	Protein and protocol (time and rpm) dependent, 0.1-200 µg/mL**	Protein and protocol (time and rpm) dependent, 0.1-200 µg/mL**	Protein dependent, typically 0.1-200 µg/mL**	

## Octet® BLI Consumables: In Depth (con't)

Octet® Consumables	Description	Intended Use <sup>2</sup>	Application	Octet® BLI System Quantitation Dynamic Range <sup>1</sup>			Regeneration	
				Octet® QK <sup>®</sup> , QK384 <sup>†</sup> , RH96 ≥32 Channel	Octet® RED96e <sup>†</sup> , K2 <sup>†</sup> , R2, R4, R8, RH16, RH96 8 or 16 Channel	Octet® N1		
Biosensors								
<a href="#">NTA</a> (Cat. Nos. 18-5101, 18-5102, 18-5103)	Ni-NTA	 	Quantitation of HIS-tagged proteins in buffer or diluted matrix, capturing of HIS-tagged proteins for kinetic analyses with various analytes	Protein dependent, typically 0.5–1000 µg/mL	Protein dependent, typically 0.5–1000 µg/mL	Protein dependent, typically 0.5–1000 µg/mL	 	
<a href="#">ProA</a> (Cat. Nos. 18-5010, 18-5012, 18-5013)	Protein A		Quantitation of IgG’s of various species including human	0.1–700 µg/mL	0.025–2000 µg/mL	0.5–4000 µg/mL		
<a href="#">ProG</a> (Cat. Nos. 18-5082, 18-5083, 18-5084)	Protein G		Quantitation of IgG’s of various species including human	0.1–700 µg/mL	0.025–2000 µg/mL	0.5–4000 µg/mL		
<a href="#">ProL</a> (Cat. Nos. 18-5085, 18-5086, 18-5087)	Protein L		Quantitation of IgG’s of various species via the kappa light chain	0.1–700 µg/mL	0.05–2000 µg/mL	0.5–2000 µg/mL		
<a href="#">SA</a> (Cat. Nos. 18-5019, 18-5020, 18-5021)	Streptavidin		Immobilizing biotinylated molecules for all kinetic analyses	N/A	N/A	N/A		
<a href="#">SAX</a> (Cat. Nos. 18-5117, 18-5118, 18-5119)	High Precision Streptavidin	 	Immobilizing biotinylated molecules for high precision quantitation and kinetic measurements	Protein dependent	Protein dependent	Protein dependent		
<a href="#">SAX2</a> (Cat. Nos. 18-5136, 18-5137, 18-5138)	High Precision Streptavidin 2.0	 	Immobilizing biotinylated molecules for high precision and reproducible kinetic characterization and custom quantitation	Protein dependent	Protein dependent	Protein dependent		
<a href="#">SSA</a> (Cat. Nos. 18-5057, 18-5065, 18-5070)	Super Streptavidin		Small molecule and fragment analyses only, should not be used for large molecule measurements	N/A	N/A	N/A		
<a href="#">AAVX</a> (Cat. Nos. 18-5160, 18-5161, 18-5162)	AAV Quantitation		Quantitation of AAV Capsids for various AAV serotypes, including AAV1 -AAV9 and AAVrh10	AAV serotype and sample dependent, typically 8.5E8–1.0E13 vp/mL	AAV serotype and sample dependent, typically 8.5E8–1.0E13 vp/mL	AAV serotype and sample dependent, typically 8.5E8–1.0E13 vp/mL		

 Kinetics  
  Quantitation  
  Glycan Screening  
  Impurity Testing  
  Yes for Kinetics  
  Yes for Quantitation  
  No for Quantitation  
  Protein and analyte dependent

<sup>1</sup> Dynamic range might vary for different background conditions, numbers listed are guidelines only and are based on testing of intended analyte molecules, users should validate range for their own samples

<sup>2</sup> Biosensors are developed, manufactured, and QC is performed for their intended applications; using biosensors outside their intended purpose requires user validation

\* Assay conditions and dynamic range should be validated

\*\* Users should validate their assay

† Discontinued model

## Octet® BLI Consumables: In Depth (con't)

Octet® Consumables	Description	Intended Use	Application	Octet® BLI System Quantitation Dynamic Range¹			Regeneration
				Octet® QK™, QK384™, RH96 ≥32 Channel	Octet® RED96e™, K2™, R2, R4, R8, RH16, RH96 8 or 16 Channel	Octet® N1	
Kits and Reagents							
<a href="#">AR2G</a> (Cat. No. 18-5095)	Amine Coupling 2nd Generation Reagent Kit	Ⓚ	Reagent kit for immobilizing any molecule with a terminal amine group onto Octet® AR2G biosensors	N/A	N/A	N/A	◀
<a href="#">GlyM</a> (Cat. No. 18-5139)	Mannose Screening Kit	ⓐ	Relative screening of Mannose glycans in crude or purified cell culture samples	Sample dependent	Sample dependent	N/A	Ⓚ
<a href="#">GlyS</a> (Cat. No. 18-5135)	Sialic Acid Screening Kit	ⓐ	Relative screening of sialic acid in crude or purified cell culture samples	Sample dependent	Sample dependent	N/A	Ⓚ
<a href="#">HCP</a> (Cat. Nos. 18-5141, 18-5158)	Anti-CHO HCP Detection Kit	Ⓛ	High sensitivity assay kit for generic analyses of CHO HCP	Sample dependent, typically 0.5–200 ng/mL	Sample dependent, typically 0.5–200 ng/mL	N/A	Ⓚ
<a href="#">RPA</a> (Cat. No. 18-5128)	Residual Protein A Detection Kit	Ⓛ	High sensitivity assay kit for analyses of residual Protein A	Sample dependent, typically 0.1–25 ng/mL	Sample dependent, typically 0.1–25 ng/mL	N/A	Ⓚ
<a href="#">Regeneration Buffer Kit</a> (Cat. No. 18-5171)	Regeneration Buffer Kit	Ⓚ ⓐ	Set of four ready-to-use buffers for screening regeneration conditions for various Octet® Biosensors.	N/A	N/A	N/A	◀
<a href="#">Kinetics Buffer 10X</a> (Cat. No. 18-1105)	Optimized buffer matrix to be used in kinetics assays	N/A	Sartorius’ Octet® Kinetics Buffer 10X (10x KB) is essential for kinetics applications performed on the Octet® platform with Octet® biosensors.	N/A	N/A	N/A	N/A
<a href="#">ProA Calibrator Set</a> (Cat. No. 18-1118)	Calibration of the Octet® ProA Biosensors	N/A	Octet® ProA Calibrator Set is intended for the calibration of the Octet® ProA Biosensors and generation of a standard curve for IgG titer measurement. The set includes 8 calibrators with IgG concentrations ranging from 1 – 700 µg/mL.	N/A	N/A	N/A	N/A
<a href="#">Octet® Sample Diluent</a> (Cat. No. 18-1104)	Octet® sample dilution buffer for quantitation assays	N/A	Octet® sample dilution buffer for quantitation assays, 50mL. Contains Kathon.	N/A	N/A	N/A	N/A
Accessories							
<a href="#">Octet® AT</a> (Cat. No. 18-5159)	Biosensor Transfer Tool	N/A	The Octet®AT is a tool for Octet® BLI biosensor transfer. Its ergonomic design delivers exceptional comfort and makes biosensor pickup and release quick and easy.				
<a href="#">Octet® AS</a> (Cat. No. OCTET-AS)	Offline Biosensor Immobilization Station	N/A	Simultaneous and Uniform reagent loading capable of simultaneously and uniformly loading reagents onto all 96 biosensors in a biosensor tray				
<a href="#">Octet® AC</a> (Cat. No. 18-5133)	Biosensor mount cleaning tray	N/A	Octet® AC is a biosensor mount cleaning tray for regular automated cleaning of metal biosensor mounts on Octet® RH96 and RH16 instruments.				

 Kinetics  
  Quantitation  
  Glycan Screening  
  Impurity Testing  
  Yes for Kinetics  
  Yes for Quantitation  
  No for Quantitation  
  Protein and analyte dependent

## Germany

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49 551 308 0

## USA

Sartorius Corporation  
3874 Research Park Dr.  
Ann Arbor, MI 48108  
Phone +1 734 769 16006



For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)