

Octet® BLI Biosensor Selection Guide Simplifying Progress

SARTURIUS

Octet [®] Consumables			Application	Octet® S			
	Description	Intended Use²		Octet® QK ^{e‡} , QK384 [‡] , RH96 ≥32 Channel	Octet® RED96e [‡] , K2 [‡] , R2, R4, R8, RH16, RH96 8 or 16 Channel	Octet [®] N1	Regeneration
Biosensors							_
AHC (Cat. Nos. 18-5060, 18-5063, 18-5064)	Anti-Human Fc-Capture	K	Capturing human IgG's or human Fc-fusion proteins for kinetic analysis with various analytes	N/A	N/A	N/A	Yes for K
AHQ (Cat. Nos. 18-5001, 18-5004, 18-5005)	Anti-Human IgG Fc	Q	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	No for Q
AHC2 (Cat. Nos. 18-5142, 18-5143, 18-5144)	Anti-Human Fc-Capture 2nd Generation	Q and K	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	Yes for Q and K
AMC (Cat. Nos. 18-5088, 18-5089, 18-5090)	Anti-Mouse Fc-Capture	К	Capturing mouse IgG's or mouse Fc-fusion proteins for kinetic analysis with various analytes	N/A	N/A	N/A	Yes for K
AMQ (Cat. Nos. 18-5022, 18-5023, 18-5024)	Anti-Murine IgG Fv	Q	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	No for Q
APS (Cat. Nos. 18-5045, 18-5046, 18-5047)	Aminopropylsilane	К	Binding measurement of lipids, liposomes, hydrophobic proteins that don't have other methods of surface attachment	N/A	N/A	N/A	Protein and analyte dependent, users should validate their own assay
AR2G (Cat. Nos. 18-5092, 18-5093, 18-5094)	Amine Reactive 2G	K	Covalently immobilizing any molecule with a terminal amine group for all kinetic analyses	N/A	N/A	N/A	Protein dependent
FAB2G (Cat. Nos. 18-5125, 18-5126, 18-5127)	Anti-Human Fab-CH1 2nd Generation	Q and K	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	Yes for Q and K
GST (Cat. Nos. 18-5096, 18-5097, 18-5098)	Anti-GST	Q and K	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**	Yes for K, no for Q
HIS1K (Cat. Nos. 18-5120, 18-5121, 18-5122)	Anti-Penta-HIS	Q and K	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*	Yes for K
HIS2 (Cat. Nos. 18-5114, 18-5115, 18-5116)	Anti-HIS	Q	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**	Protein dependent

¹Dynamic 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 ² Biosensors are developed, manufactured, and QC is performed for their intended applications; using biosensors outside their intended purpose requires user validation

Octet® Biosensor	Description	Intended Use²	Application	Octet® System Quantitation Dynamic Range¹			
				Octet® QK° [‡] , QK384 [‡] , RH96 ≥32 Channel	Octet® RED96e [‡] , K2 [‡] , R2, R4, R8, RH16, RH96 8 or 16 Channel	Octet [®] N1	Regeneration
NTA (Cat. Nos. 18-5101, 18-5102, 18-5103)	Ni-NTA	Q and K	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	Yes for K, no for Q
ProA (Cat. Nos. 18-5010, 18-5012, 18-5013)	Protein A	Q	Quantitation of IgG's of various species including human	0.1-700 μg/mL	0.025-2000 μg/mL	0.5-4000 μg/mL	Yes
ProG (Cat. Nos. 18-5082, 18-5083, 18-5084)	Protein G	Q	Quantitation of IgG's of various species including human	0.1-700 μg/mL	0.025-2000 μg/mL	0.5-4000 μg/mL	Yes
ProL (Cat. Nos. 18-5085, 18-5086, 18-5087)	Protein L	Q	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	Yes
SA (Cat. Nos. 18-5019, 18-5020, 18-5021)	Streptavidin	K	Immobilizing biotinylated molecules for all kinetic analyses	N/A	N/A	N/A	Protein dependent
SAX (Cat. Nos. 18-5117, 18-5118, 18-5119)	High Precision Streptavidin	Q and K	Immobilizing biotinylated molecules for high precision quantitation and kinetic measurements	Protein dependent	Protein dependent	Protein dependent	Protein dependent
SAX2 (Cat. Nos. 18-5136, 18-5137, 18-5138)	High Precision Streptavidin 2.0	Q and K	Immobilizing biotinylated molecules for high precision and reproducible kinetic characterization and custom quantitation	Protein dependent	Protein dependent	Protein dependent	Protein dependent
SSA (Cat. Nos. 18-5057, 18-5065, 18-5070)	Super Streptavidin	К	Small molecule and fragment analyses only, should not be used for large molecule measurements	N/A	N/A	N/A	Analyte compounds can be washed off during dis- sociation in buffer since most have fast offrates
Kits							
AR2G (Cat. No. 18-5095)	Amine Coupling 2nd Generation Reagent Kit	K	Reagent kit for immobilizing any molecule with a terminal amine group onto Octet® AR2G biosensors	N/A	N/A	N/A	Protein dependent
GlyM (Cat. No. 18-5139)	Mannose Screening Kit	Q	Relative screening of Mannose glycans in crude or purified cell culture samples	Sample dependent	Sample dependent	Sample dependent	No for Q
GlyS (Cat. No. 18-5135)	Sialic Acid Screening Kit	Q	Relative screening of sialic acid in crude or purified cell culture samples	Sample dependent	Sample dependent	Sample dependent	No for Q
HCP (Cat. Nos. 18-5141, 18-5158)	Anti-CHO HCP Detection Kit	Q	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	Sample dependent, typically 0.5-200 ng/mL	No for Q
RPA (Cat. No. 18-5128)	Residual Protein A Detection Kit	Q	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	Sample dependent, typically 0.1-25 ng/mL	No for Q

 $^{^*}$ Assay conditions and dynamic range should be validated ** Users should validate their assay ‡ Discontinued model

Germany

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

USA

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