



Octet® BLI Biosensor Selection Guide

Simplifying Progress

SARTORIUS

| Octet® Consumables | Description | Intended Use ² | Application | Octet® BLI System Quantitation Dynamic Range ¹ | | | Regeneration |
|--|--------------------------------------|---------------------------|--|--|---|---|---|
| | | | | Octet® QK ⁺ , QK384 [‡] , RH96 ≥32 Channel | Octet® RED96e [†] , K2 [‡] , R2, R4, R8, RH16, RH96 8 or 16 Channel | Octet® N1 | |
| 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 | K | 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 | K | 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 assays |
| 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

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

| Octet® Biosensor | Description | Intended Use ² | Application | Octet® BLI System Quantitation Dynamic Range ¹ | | | Regeneration |
|---|---|---------------------------|--|---|---|---|---|
| | | | | Octet® QK [†] , QK384 [†] , RH96 ≥32 Channel | Octet® RED96 [†] , K2 [†] , R2, R4, R8, RH16, RH96 8 or 16 Channel | Octet® N1 | |
| 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 | K | 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 dissociation in buffer since most have fast offrates |
| AAVX (Cat. Nos. 18-5160, 18-5161, 18-5162) | AAV Quantitation | Q | 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 | Yes |
| 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