



Octet® SPR Sensor Chip Selection Guide

Simplifying Progress

SARTORIUS

Octet® SPR Sensor Chip Selection Guide

Sensor Chip	Part No.	Surface Chemistry	Capacity	Immobilization Method	Recommended Use	Advantages
COOH1	19-0053	Planar carboxylated oligoethylene oxide	Low	Amine coupling	Protein-protein or other large molecule kinetic assays when lectin binding may occur, or dextran alternative is desired.	<ul style="list-style-type: none"> Immobilization of targets without derivatization or tags Produces a highly stable covalent bond Can be used to immobilize affinity ligands to create additional capture chemistries (i.e. Protein A, antibodies, etc.) Effective over a wide pH range
CDL	19-0127	Thin, low density carboxymethyl dextran layer	Medium	Amine coupling	Measuring accurate kinetics of protein-protein or other intermediate (>1 kDa) to large (>25 kDa) molecule interactions.	<ul style="list-style-type: none"> Immobilization of targets without derivatization or tags Biocompatible with a range of molecules Produces a highly stable covalent bond Can be used to immobilize affinity ligands to create additional capture chemistries (i.e. Protein A, antibodies, etc.) Effective over a wide pH range
CDH	19-0128	Carboxymethyl dextran three-dimensional hydrogel surface with carboxylic acids	High	Amine coupling	Small molecule-protein interactions, fragment screening, or other target interactions with low binding activity. Assays where high capacity is needed.	<ul style="list-style-type: none"> Immobilization of targets without derivatization or tags High-capacity carboxymethyl dextran surface, ideal for immobilizing a large amount of target for small molecule and fragment analysis Biocompatible with a wide range of molecules Produces a highly stable covalent bond Effective over a wide pH range
PCH	19-0129	Non-dextran polycarboxylate hydrogel surface	High+	Amine coupling	Small molecule (<1 kDa)-protein interactions, fragment screening, or other target interactions with low binding activity. Assays where high capacity is needed and/or dextran alternative is desired.	<ul style="list-style-type: none"> Immobilization of targets without derivatization or tags Highest capacity surface, ideal for immobilizing targets for small molecule and fragment analysis Provides an alternative to dextran surface Produces a highly stable covalent bond Effective over a wide pH range
HisCap	19-0058	Non-dextran polysaccharide three-dimensional surface with carboxylic acids pre-immobilized nitrilotriacetic acid (NTA), regenerable with imidazole, EDTA	High	Capture via His tag	Small molecule or peptide kinetics with His-tagged protein ligands. Large molecule kinetic assays with His-tagged proteins.	<ul style="list-style-type: none"> High ligand capture capacity for low molecular weight analytes such as fragments Stable baseline, important for accurate kinetic analysis Provides a convenient means of directed capture of His-tagged proteins Can be regenerated using a variety of conditions, such as imidazole, SDS, or EDTA Suitable alternative for proteins that are not amenable to amine coupling
SADH	19-0130	Streptavidin immobilized in three-dimensional carboxymethyl dextran hydrogel	Medium-high	Capture via biotin	Intermediate (>1 kDa) to large (>25 kDa) molecule kinetics with biotinylated ligands.	<ul style="list-style-type: none"> Highly efficient capture over a wide pH range Requires low quantities of ligand (nanomolar concentrations) Single step capture Surface has lower electrostatic charge compared to amine coupling sensors



SARTORIUS

Maintenance Chip for Octet® SPR systems

Part Number: 19-0XXX, Lot: 03142021

Store at Room Temperature. For Research Use Only

Safety information for this product is provided on the outside package

sartorius.com/Octet-support, Emergency 800-424-9300

WARNING

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