

Series S Sensor Chip CM5 Instructions for Use

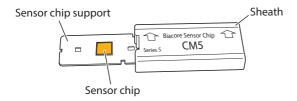
Product description

Order code: 29149603 (package of ten sensor chips)

BR100530 (package of three sensor chips) 29104988 (package of one sensor chip)

Storage: The use-before date applies to chips stored at 2°C to 8°C in unopened

pouches.



The sensor chip is fixed to a polystyrene support sheath. Each cassette, consisting of a sensor chip and sheath assembly, is individually packed under a nitrogen atmosphere in a sealed pouch.

Note: For in vitro use only.

cytiva.com 22064823 AF

Application areas

Series S Sensor Chip CM5 is designed for interaction analysis in Biacore™ systems. The surface has a carboxymethylated dextran matrix covalently coupled to a surface coating on a gold film. Ligands ranging from small molecules to proteins, nucleic acids and carbohydrates may be attached via covalent coupling using well-established chemistries. Alternatively, a capturing molecule may be coupled for subsequent capture. Series S Sensor Chip CM5 is suitable for most general applications.

Refer to cytiva.com/biacore for updates on applications and scientific publications.

Preparations for use

Step Action 1 If you are working in a humid environment, allow the sealed sensor chip pouch to equilibrate at room temperature for 15 to 30 minutes in order to prevent condensation on the chip surface. 2 Prepare the Biacore instrument with running buffer. The buffer should be filtered (0.22 µm), and degassed for systems that do not have an integrated buffer degasser. 3 Open the sensor chip pouch. Make sure that the sensor chip support remains fully inserted into the sheath at all times to protect the chip from dust particles. Dock the sensor chip in the instrument as described in the instrument 4 handbook. Note: Sensor chips that are not docked in the instrument should be stored in closed containers.

Immobilizing the ligand

The ligand or capturing molecule is covalently bound to the sensor chip surface via carboxyl groups on the dextran. Functional groups on the molecule that can be used for coupling include -NH $_2$, -SH, -CHO, -OH, and -COOH.

Refer to $\it Biacore Sensor Surface Handbook\ for\ more\ information\ on\ immobilization\ strategies\ and\ procedures.$

Interaction analysis

Interaction analysis is performed by injection of samples over the sensor chip surface. The covalently immobilized molecule either participates directly in the interaction under study or is used for affinity capture of one of the interacting molecules.

Refer to Biacore handbooks and *cytiva.com/biacore* for details on experimental protocols and methodology.

Regeneration

Regeneration of the immobilized ligand may be performed by selective dissociation of the bound analyte. Conditions should be chosen to achieve complete dissociation of the analyte without affecting the binding characteristics of the ligand. The surface of Series S Sensor Chip CM5 is resistant to a wide range of agents for this purpose (for more information see *Chapter Chemical resistance*, on page 7). The choice of regeneration procedure may be limited by the stability of the ligand.

Refer to $Biacore\ Sensor\ Surface\ Handbook\ for\ more\ detailed\ information\ on\ regeneration\ strategies.$

Chemical resistance

The surface of Series S Sensor Chip CM5 is resistant to 1-minute pulses of many commonly used agents.

Agent	Concentration
Acetonitrile	30%
DMSO	10%
DTE	0.1 M
EDTA	0.35 M
Ethanol	70%
Ethanolamine	1 M
Ethylene glycol	100%
Formamide	40%
Formicacid	20%
Glycine-HCl pH 1.5 to 3.0	100 mM
Glycine-NaOH pH 9.5 (BIAdesorb Solution 2)	50 mM
Guanidine hydrochloride	6 M
HCI	100 mM
Imidazole	300 mM
MgCl ₂	4 M
NaOH	100 mM
NaCl	5 M
SDS (BIAdesorb Solution 1)	0.5%
Surfactant P20	5%
Urea	8 M



cytiva.com/biacore

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

Biacore is a trademark of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

All other third-party trademarks are the property of their respective owners.

© 2020 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit cytiva.com/contact

22064823 AF V:4 06/2020