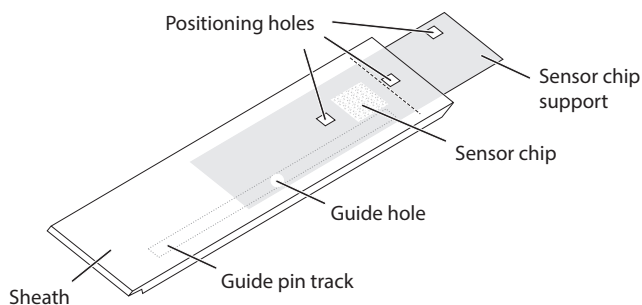


Sensor Chip CM7

Instructions for Use

Product description

Order code:	28957332 (Package of one sensor chip) 29147017 (Package of three sensor chips)
Content:	Sensor Chip CM7
Storage:	The use-before date applies to chips stored at 4°C to 8°C in unopened pouches.



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

Note: For *in vitro* use only.

Application areas

Sensor Chip CM7 is intended for interaction analysis in Biacore systems. The surface has a carboxymethylated dextran matrix covalently attached to a gold film, with similar chemical properties to Sensor Chip CM5 but prepared to provide a higher immobilization capacity.

Sensor Chip CM7 is designed primarily for work with low molecular weight analytes. With macromolecular analytes, the analyte binding capacity (R_{max}) is lower than expected on the basis of the amount of ligand immobilized, possibly because of multi-site attachment of the ligand and matrix crosslinking on the higher capacity surface.

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. If the chip is not inserted into the instrument immediately, it should be stored in a plastic bag.
4	Dock the sensor chip in the instrument as described in the Instrument Handbook.

Immobilizing the ligand

The ligand 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 $-\text{NH}_2$, $-\text{SH}$, $-\text{CHO}$, $-\text{OH}$ and $-\text{COOH}$.

For more detailed information on immobilization strategies and procedures, refer to the *Biacore Sensor Surface Handbook*.

Interaction analysis

Interaction analysis is performed by injection of samples over the sensor chip surface. Analyte molecules in the injected sample bind directly to the covalently immobilized ligand.

Refer to Biacore handbooks and [cytiva.com/biacore](https://www.cytiva.com/biacore) for details on experimental protocols and methodology.

Regeneration

Regeneration of the immobilized ligand is performed by 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 Sensor Chip CM7 is resistant to a wide range of agents for this purpose (see the following chapter). The choice of regeneration procedure may be limited by the stability of the ligand. Do not use harsher conditions than required to remove the analyte.

For more detailed information on regeneration strategies, refer to the *Biacore Sensor Surface Handbook*.

Chemical resistance

The surface of Sensor Chip CM7 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	40%
Ethanolamine	1 M
Ethylene glycol	100%
Formamide	40%
Formic acid	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
HCl	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

Page intentionally left blank



[cytiva.com/biacore](https://www.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](https://www.cytiva.com/contact)

28957862 AC V:4 07/2020