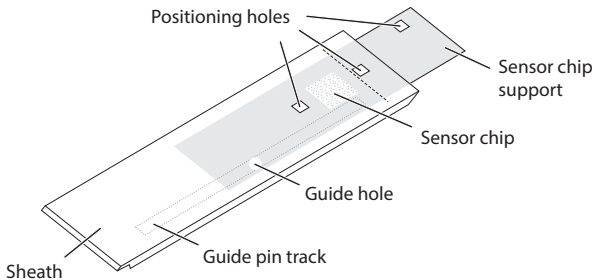


# Sensor Chip CM3

## Instructions for Use

### Product description

Order code:	BR-1005-41 (package of three sensor chips)
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.

## Application areas

Sensor Chip CM3 is designed for interaction analysis in Biacore™ systems. The surface has a carboxymethylated dextran matrix covalently attached to 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 can be coupled for subsequent capture.

Sensor Chip CM3 has a shorter dextran matrix than Sensor Chip CM5. This may be an advantage for example in work with large molecules, such as virus particles or cells. Non-specific binding may also be reduced for some samples.

Due to the shorter matrix, immobilization levels will typically be around 30% of those obtained on Sensor Chip CM5 under comparable conditions.

Refer to [cytiva.com/biacore](http://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.
4	Dock the sensor chip in the instrument as described in the instrument handbook.
	<b>Note:</b> <i>Sensor chips that are not docked in the instrument should be stored in closed containers.</i>

## Immobilizing the ligand

The ligand or capturing molecule is covalently bound to the sensor chip surface via carboxyl moieties on the dextran. Functional groups on the molecule that can be used for coupling include -NH<sub>2</sub>, -SH, -CHO, -OH, and -COOH.

Refer to *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 ligand 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](http://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 Sensor Chip CM3 is resistant to a wide range of agents for this purpose (for more information see the following chapter). The choice of regeneration procedure may be limited by the stability of the ligand.

Refer to *Biacore Sensor Surface Handbook* for more information on regeneration strategies.

# Chemical resistance

The surface of Sensor Chip CM3 is resistant to one-minute pulses of many commonly used agents. See table below for information of common agents compatible with Sensor Chip CM3.

Agent	Concentration
Acetonitrile	30%
DMSO	10%
DTE	0.1 M
EDTA	0.35 M
Ethanol	70%
Ethanolamine	1 M
Ethylene glycol	100%
Formamide	40%
Formic acid	20%
Glycine pH 1.5 to 3.0	100 mM

Agent	Concentration
HCl	100 mM
Imidazole	300 mM
MgCl <sub>2</sub>	4 M
NaOH	100 mM
NaCl	5 M
SDS	0.5%
Surfactant P20	5%
Urea	8 M

Page intentionally left blank



[www.cytiva.com/sampleprep](http://www.cytiva.com/sampleprep)

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](http://cytiva.com/contact)

22060728 AD V:7 07/2020