

# Series S Sensor Chip Protein L

## Instructions for Use

### A Biacore Extend product

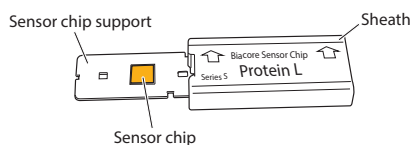
This product is part of the Biacore™ Extend product line and is not a regular Biacore consumables product. For more information,

see [cytiva.com/biacoreextend](https://www.cytiva.com/biacoreextend).

### Product description

Order code: 29205138 (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.

Series S Sensor Chip Protein L consists of a carboxymethylated dextran matrix pre-immobilized with a recombinant Protein L.

**Note:** For *in vitro* use only.

# Application areas

Series S Sensor Chip Protein L is suitable for concentration analysis and yes/no screening. Series S Sensor Chip Protein L is designed to bind antibodies and antibody fragments of various species containing a kappa light chain for interaction analysis in Biacore systems.

Refer to [cytiva.com/biacore](http://cytiva.com/biacore) for updates on applications and scientific publications, and to *Handbook* (29019400).

# Surface specificity

Series S Sensor Chip Protein L is suitable for the capture of a wide range of antibody fragments such as Fab, single-chain variable fragments (scFv), and domain antibodies (Dab).

The recombinant Protein L binds to representatives of all human antibody classes, including IgG, IgM, IgA, IgE and IgD, and to antibodies from some other mammalian species, including mouse. It binds to the variable region of the kappa light chain (subtypes 1, 3 and 4 in humans, and subtype 1 in mice). No part of the heavy chain is involved in the binding interaction. Species such as bovine, goat, sheep, and horse whose antibodies contain almost exclusively lambda chains will not bind well, if at all, to Protein L.

# Preparations for use

Step	Action
1	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.

**Note:** *Biacore 4000 requires a modified procedure for hydrodynamic addressing. Please refer to the instructions called "Hydrodynamic addressing in Biacore 4000 when using pre-immobilized sensor chips" available at [cytiva.com/proteinlsensorchip](http://cytiva.com/proteinlsensorchip).*

**Note:** *Storage stability is affected by exposure to air. Keep sensor chip in unopened pouch until use.*

## Analysis temperature

Series S Sensor Chip Protein L is designed for use at 25°C.

## Start-up cycles

For best assay performance, run at least one start-up cycle using sample or buffer as analyte and identical settings as for the analysis cycles.

## Regeneration

Regenerate the surface with one 120 s injection of 10 mM Glycine-HCl, pH 1.7. This will remove captured ligand together with any analyte bound to them. Some ligands may require one additional 120 s injection of 10 mM Glycine-HCl, pH 1.7.

10 mM Glycine-HCl pH 1.7 can be prepared from 10 mM Glycine-HCl pH 1.5 and pH 2.0 (available from Cytiva, product code BR100354 and BR100355). Example: Mix 4.6 mL 10 mM Glycine-HCl pH 1.5 with 5.4 mL 10 mM Glycine-HCl pH 2.0 to obtain pH 1.7.

Avoid using basic regeneration solutions with a pH >10



## [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)

29212800 AB V:5 07/2020