

# Series S Sensor Chip Protein G

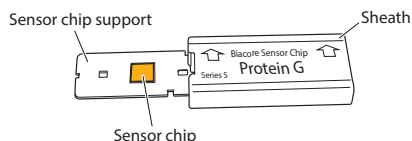
## Instructions for Use

### A Biacore™ Extend product

This product is to be considered a development product and not a standard Biacore consumables product. For more information, see [cytiva.com/biacoreextend](https://www.cytiva.com/biacoreextend).

Order code: 29179315 (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.

Sensor Chip Protein G consists of a carboxymethylated dextran matrix pre-immobilized with a recombinant Protein G - GammaBind G, Type 2.

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

# Application areas

Series S Sensor Chip Protein G is designed to bind antibodies of different species and subclasses for interaction analysis in Biacore systems. Series S Sensor Chip Protein G is a good choice for kinetic characterization and concentration analyses in a wide range of applications.

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

# Surface specificity

The recombinant Protein G binds a broad range of IgG, such as human (including IgG<sub>3</sub>), rat, rabbit, mouse, guinea pig, goat, sheep, and cow.

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

# Analysis temperature

Series S Sensor Chip Protein G 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 30-second injection of 10 mM Glycine-HCl, pH 1.5 (available from Cytiva, product code BR100354). This will remove captured ligand together with any analyte bound to them.

Alternative regeneration procedures are:

- One additional 30 s injection of 0.05-0.5% Sodium dodecyl sulfate (SDS), or
- One 30 s injection of 10 mM Glycine-HCl pH 1.5 with 0.5% Surfactant P20 added (available from Cytiva, product code BR100054).

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)

29186433 AB V:7 07/2020