

# Biacore Maintenance Kit, type 2

## Instructions for Use

### Product description

Order code:	29394519
Contents:	<ul style="list-style-type: none"><li>• One bottle of BIAtest solution, 65 mL (15% (w/w) sucrose in HBS-N buffer with 3.4 mM EDTA)</li><li>• One bottle of BIAnormalizing solution, 90 mL (70% (w/w) glycerol)</li><li>• Two bottles of BIAdesorb solution 1, 95 mL (0.5% (w/v) sodium dodecyl sulphate (SDS))</li><li>• Two bottles of BIAdesorb solution 2, 95 mL (50 mM glycine-NaOH pH 9.5)</li><li>• One bottle of HBS-N buffer, 10x solution, 50 mL (0.1M HEPES, 1.5M NaCl)</li><li>• One Sensor Chip Maintenance</li></ul>
Storage:	Store at 2°C to 8°C. BIAdesorb solution 1, should be stored at room temperature (SDS precipitates at low temperature).
For use with:	Biacore™ Maintenance Kit, type 2 is designed to be used with the <b>Maintenance tools</b> and <b>Test tools</b> in Biacore 4000, Biacore T200, and Biacore S200 SPR systems. For more information, see the respective instrument handbook.

Kit capacity: With the exception of the HBS-N 10x buffer, the solutions in the kit are sufficient for 3 to 4 months (Biacore T200 and Biacore S200 SPR systems) or 1 to 2 months (Biacore 4000 SPR system) of normal use. Additional buffer packs can be purchased from Cytiva (order code BR-1006-70).

Safety: For use and handling of the product in a safe way, please refer to the Safety Data Sheet.

**Note:** *For research use only.*

**Note:** *Not included in the kit: Sodium hypochlorite with 8% to 15% active chlorine*

# Reagent preparation

## Sodium hypochlorite solution

Depending on the consumption multiples of 21.5 mL are prepared according to the pipetting scheme given in the following table. Mix and use within 12 hours. The total consumption of Sodium hypochlorite solution can be found in the respective instrument handbook and is stated in a software dialog box when starting the procedure.

**Note:** *Prepare solutions just prior to use.*

Sodium hypochlorite solution (conc.) (mL)	Distilled and filtered water (mL)	Total volume (mL)
1.5	20	21.5
3.0	40	43
6.0	80	86
15	200	215

**Note:** *Take care when removing flow tubings from the disinfectant solution to avoid splashing.*

## HBS-N buffer

Dilute the 10x solution 1+9 (v/v) with deionized, 0.2 µm filtered water, for example by mixing 10 mL 10x solution and 90 mL water.

# Handling

Make sure that the running buffer is compatible with BIAdesorb solution 1 before running the **Desorb** and **Desorb and sanitize** procedures (SDS will precipitate with some common metal ions including potassium and magnesium).

Make sure to prime the system with running buffer, not water, before docking pre-immobilized sensor chip for example sensor chip SA or CAP as these surfaces are sensitive to sodium hypochlorite residues.

**Note:** See the respective instrument handbook for details.

## Maintenance intervals

Recommended maintenance intervals are weekly for **Desorb** and monthly for **Desorb and sanitize**. Depending on the instrument usage, maintenance intervals can be shorter.

## General recommendations

Use Sensor Chip Maintenance for **Desorb** and **Desorb and sanitize** procedures.

Store Sensor Chip Maintenance in the enclosed plastic bag when not inserted in the system. Inspect the surface before use. If the surface is not clean, it should be rinsed with water. Wipe the glass and support dry with a lint-free tissue before docking.

## Flow system cleaning

**Desorb** removes adsorbed proteins from the autosampler and IFC. Use BIAdesorb solutions 1 and 2 for the **Desorb** procedure. Make sure that the analysis and sample compartment temperatures are set to 25°C, since the SDS in BIAdesorb solution 1 may precipitate at low temperature.

## Flow system disinfection

**Desorb and sanitize** cleans the flow system of microorganisms. Following completion of the procedure, an instrument stabilization period of 3 to 4 hours is recommended to achieve a stable baseline and particularly, if maximum sensitivity in measurements is required.

## Instrument normalizing

**Normalize** normalizes the resonance signal in all flow cells to compensate for small differences between individual sensor chips.

Perform **Normalize** when docking a new sensor chip or when changing analysis temperature and if maximum sensitivity in measurements is required.

The BIAnormalizing solution should be at room temperature for this procedure.

## Instrument performance testing - system check

**System check** is used to perform a comprehensive check of system performance using a standard sucrose solution. Use a non-immobilized sensor chip surface for **System check**.

**Note:** *Do not use Sensor Chip Maintenance for System check.*

Use BIAtest solution for instrument performance testing with **Test tools**. Use HBS-N buffer as running buffer. For **System check**, a newly opened HBS-N buffer package or a fresh dilution of concentrated HBS-N buffer must be used as running buffer to achieve correct results.

**Note:** *An accurate concentration of the HBS-N buffer is imperative for correct test results.*



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

29396777 AB V:4 10/2020