

Desorb Kit

Instructions for Use

Product description

Product code: BR100823

Content: Desorb Kit contains the following products.

Content	Description
One bottle of BIAdesorb solution 1, 500 mL	0.5% (w/v) sodium dodecyl sulphate (SDS)
One bottle of BIAdesorb solution 2, 500 mL	50 mM glycine-NaOH, pH 9.5

Additional materials required:

- Series S Maintenance Chip (BR100562).
- HBS-EP+ 10× buffer (BR100826 or BR100669). Dilute 1+9 (v/v) with deionized, 0.2 µm filtered water to make the running buffer.
- Sodium hypochlorite solution (10% to 15%).

Storage on arrival:

Store BIAdesorb solution 1 at room temperature (SDS precipitates at low temperature).

Store BIAdesorb solution 2 at 2°C to 8°C.

Kit capacity:

The solutions in the kit are sufficient for 2 to 6 months of normal use depending on the Biacore™ system used.

Safety:

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

Intended use

Desorb Kit is intended to be used with the maintenance application in any Biacore system. For more information, refer to the Biacore system specific *Operating Instructions*.

Software procedures are identified by **bold italic** text.

The table below shows the procedures to use with Desorb Kit and their recommended frequency.

Procedure	Purpose	Frequency
<i>Desorb</i>	Clean the flow system	Weekly
<i>Desorb and sanitize</i>	Disinfect the flow system	Monthly

Note: *For research use only.*

Clean the flow system

Make sure that the running buffer is compatible with BIAdesorb solution 1 before cleaning the flow system. Do not use running buffer with additives containing metal ions as SDS in BIAdesorb solution 1 will precipitate with some common metal ions including potassium and magnesium.

Follow the steps below to clean the flow system.

Step	Action
1	Set the analysis and sample compartment temperatures to 25°C. Note: <i>SDS in BIAdesorb solution 1 can precipitate at low temperature.</i>
2	Make sure that the surface of the Series S Maintenance Chip is clean before use. If needed, rinse with water. Wipe the glass and support dry with a lint-free tissue. Note: <i>Store Series S Maintenance Chip in the enclosed plastic bag when not inserted in the system.</i>

Step	Action
3	Remove the used sensor chip from the docking station and insert Series S Maintenance Chip. Note: <i>The ligands and capture molecules on the sensor chip can be damaged by the Desorb procedure. If required, store used sensor chip in suitable storage buffer at 2°C to 8°C. For reuse, refer to the sensor chip Instructions for Use.</i>
4	Open Desorb and follow the instructions on the screen to place the tubing for running buffer, water, and reagent in the relevant solutions.
5	Prepare BIAdsorb solutions 1 and 2 in vials, bottles, or plates according to the Desorb instructions, and position these within the Biacore instrument.
6	Follow the instructions on the screen to run Desorb .

Disinfect the flow system

Sodium hypochlorite solution

Sodium hypochlorite, 10% to 15% is used to prepare the disinfectant solution used in **Desorb and sanitize**. The following products are recommended.

Product name	Contents	Manufacturer	Product Code
Sodium hypochlorite, 10% to 15%	500 mL	Acros Organics, part of Thermo Fisher Scientific	219255000
Sodium hypochlorite, 10% to 15%	500 mL	Alfa Aesar, part of Thermo Fisher Scientific	33369 ¹

¹ Contact the manufacturer and ask for the date of quality control of a certain lot.

Note: *The active chlorine concentration decreases with time. Do not use hypochlorite product longer than 12 months after the date of quality control to make sure that the concentration is sufficient for disinfection. The quality control date is stated in the Certificate of Analysis for the product.*

Disinfectant solution

The disinfectant solution used in **Desorb and sanitize** is sodium hypochlorite solution, 0.6% to 1.0%. This is prepared from sodium hypochlorite, 10% to 15% described above.

Prepare the disinfectant solution according to the following table. Mix and use within 12 hours.

Solution	Volume (mL)
Sodium hypochlorite, 10% to 15%	27
Distilled and filtered water	360
Total volume	387

The total consumption of sodium hypochlorite solution can be found in the Biacore system specific *Operating Instructions* and is stated in a software dialog box when starting the procedure.

Disinfect the flow system

Make sure that the running buffer is compatible with BIAdesorb solution 1 before disinfecting the flow system. Do not use running buffer with additives containing metal ions as SDS in BIAdesorb solution 1 will precipitate with some common metal ions including potassium and magnesium.

Do not use plain water as running buffer. Water rinses sodium hypochlorite insufficiently compared to running buffer. Residues of sodium hypochlorite can damage sensor chips.

Follow the steps below to disinfect the flow system.

Step	Action
1	Set the analysis and sample compartment temperatures to 25°C. Note: <i>SDS in BIAdesorb solution 1 can precipitate at low temperature.</i>
2	Make sure that the surface of the Series S Maintenance Chip is clean before use. If needed, rinse with water. Wipe the glass and support dry with a lint-free tissue. Note: <i>Store Series S Maintenance Chip in the enclosed plastic bag when not inserted in the system.</i>

Step	Action
3	<p>Remove used sensor chip from the docking station and insert Series S Maintenance Chip.</p> <p>Note: <i>The ligands and capture molecules on the sensor chip can be damaged by Desorb and sanitize procedure. If required, store used sensor chip in suitable storage buffer at 2°C to 8°C. For reuse, refer to the sensor chip Instructions for Use.</i></p>
4	<p>Open Desorb and sanitize and prepare recommended volumes of disinfectant solution, and BIA desorb solutions 1 and 2 in vials, bottles, or plates according to the software instruction.</p>
5	<p>Follow the instructions on the screen to place the specified tubing for buffer, water, and reagent in the solutions during each step of Desorb and sanitize.</p> <p>Note: <i>Be careful when removing flow tubing from the disinfectant solution to avoid splashing.</i></p>
6	<p>After completion of the procedure, run running buffer in standby mode for at least 3 to 4 hours or overnight before docking a new sensor chip.</p>

**Give feedback on this document**

Visit cytiva.com/techdocfeedback or scan the QR code.



cytiva.com/biacore

Cytiva and the Drop logo are trademarks of Life Sciences IP Holdings Corp. or an affiliate doing business as Cytiva.

Biacore is a trademark of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Any third-party trademarks are the property of their respective owners.

© 2020–2022 Cytiva

For local office contact information, visit cytiva.com/contact

22064879 AD V:8 09/2022