

# Amine Coupling Kit, type 2

## Instructions for Use

### Product description

Order code:	BR-1006-33
Contents:	<ul style="list-style-type: none"><li>• N-Hydroxysuccinimide (NHS), 115 mg</li><li>• 1-Ethyl-3-(3-dimethylaminopropyl) carbodiimide hydrochloride (EDC), 750 mg</li><li>• 1 M Ethanolamine hydrochloride-NaOH pH 8.5, 2 x 10.5 ml</li></ul>
Storage:	+2 to 8°C
For use with:	Biacore™ 4000, Biacore A100 and Biacore S51
Capacity:	The kit contains reagents sufficient for approximately 60 immobilizations in Biacore S51, and 80 immobilizations in Biacore 4000 and Biacore A100.
Safety:	For use and handling of the product in a safe way, please refer to the Safety Data Sheet.

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

**Note:** *For other Biacore systems, use Amine Coupling Kit (BR-1000-50).*

### Description

Amine Coupling Kit, type 2 is designed to be used with carboxylated sensor surfaces from Cytiva. Ligands are coupled via primary amine groups to the sensor surface for subsequent analysis.

# Reagent preparation

Step	Action
1	Dissolve the EDC and NHS by adding 10.0 ml of filtered, deionized water to each vial.
2	Cap vials tightly and agitate until the solids are completely dissolved.
3	Dispense EDC and NHS solutions separately in aliquots for storage.
4	Store the aliquots of EDC and NHS frozen at -18°C or below. Ethanolamine may be stored at +2 to 8°C.

**Note:**

*Use aliquots within two months.*

**Note:** *Suitable vials for preparing aliquots for storage and use directly in Biacore systems can be ordered from Cytiva, for example Plastic Vials, 7 mm (BR-1002-12) with snap caps (BR-1002-13).*

# Handling

Step	Action
1	Thaw frozen aliquots and agitate them gently to make sure that the solutions are homogeneous.
2	Place the aliquots separately in the autosampler as instructed in the software.

Page intentionally left blank



## [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)

22060777 AD V:4 06/2020