

Amine Coupling Kit

Instructions for Use

Product description

Order code:	BR-1000-50
Contents:	<ul style="list-style-type: none">• N-Hydroxysuccinimide (NHS), 115 mg• 1-Ethyl-3-(3-dimethylaminopropyl) carbodiimide hydrochloride (EDC), 750 mg• 1 M Ethanolamine hydrochloride-NaOH pH 8.5, 10.5 mL
Storage:	2°C to 8°C
For use with:	All Biacore™ systems except Biacore 4000, Biacore A100 and Biacore S51.
Capacity:	The kit contains reagents sufficient for approximately 80 immobilizations.
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 Biacore 4000, Biacore A100 and Biacore S51, use Amine Coupling Kit, type 2 (BR-1006-33).*

Description

Amine Coupling Kit 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 the 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°C to 8°C.
	Note: <i>Use aliquots within two months.</i>

Note: Suitable vials for preparing aliquots for storage can be ordered from Cytiva, for example, Plastic Vials 7 mm (BR-1002-12) and Caps (BR-1002-13), or Plastic Vials and Caps 11 mm (BR-1002-14).

Handling

Step	Action
1	Thaw frozen aliquots and agitate them gently to make sure that the solutions are homogeneous.
2	For Biacore systems with autosampler: Place the aliquots separately in the autosampler as instructed in the software.
	For Biacore systems without autosampler: Mix equal volumes of EDC and NHS immediately before use.

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)

22051062 AI V:6 07/2020