

PDEA Thiol Coupling Reagent

Instructions for Use

Product description

Product code:	BR100058
Contents:	One bottle 2-(2-pyridinyldithio)ethaneamine hydrochloride, 100 mg
Storage:	2°C to 8°C
Safety:	For use and handling of the product in a safe way, refer to the Safety Data Sheet.

Note: *For in vitro use only.*

Intended use

PDEA Thiol Coupling Reagent is designed to be used for coupling biomolecules on Sensor Chip CM7, Sensor Chip CM5, Sensor Chip CM4, Sensor Chip CM3, and Sensor Chip C1 via thiol/disulphide exchange in Biacore™ systems. The biomolecule or the sensor chip can be modified with PDEA to introduce reactive disulphide groups. Biomolecules are attached via introduced reactive disulphides or native thiol groups.

Formula: $C_7H_{10}N_2S_2 \times HCl$

Chemical structure:



Properties: Off-white to pale yellow crystals.
Molecular weight: 222.7 g/mol.

Handling

For detailed information on immobilization strategies, including preparations of reagent solution and modification of ligand, refer to [cytiva.com/biacore](https://www.cytiva.com/biacore).

Page intentionally left blank

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

© 2020–2022 Cytiva

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

22051359 AE V:6 11/2022