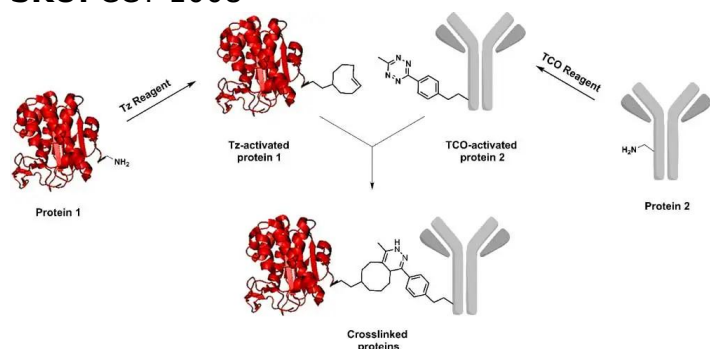




CLICK-&-GO® LYS-TO-LYS PROTEIN-PROTEIN CONJUGATION KIT

SKU: CCT-1008



DESCRIPTION

The Lys-to-Lys Protein-Protein Conjugation Kit (Cat. # 1008) provides a sufficient amount of reagents to perform two protein-protein conjugation reactions. Any two lysine containing proteins (100-500 µg) in a volume of 100 µL (1-5 mg/mL) can be efficiently conjugated in less than 3 hours, start-to-finish.

The Lys-to-Lys Protein-Protein Conjugation kit provides all of the necessary reagents to perform two protein-protein conjugation reactions. Conjugates are formed by the very efficient, catalyst-free bio-orthogonal ligation reaction between trans-cyclooctene (TCO) and tetrazine (Tz) functional groups. In the protocol provided, two proteins (or any other amine-containing biopolymers) are labeled separately with TCO and Tetrazine reagents. The labeled molecules are then mixed for 60 minutes in any suitable buffer to form a conjugate.

The extremely fast kinetics of the TCO/Tz ligation reaction enables rapid conjugation (30-60 min) of lysine containing proteins to each other at low concentrations (e.g. 5-10 µM), with >99% conversion of the limiting protein in mild buffered media (e.g. PBS pH 7.5). Another benefit is the long-term stability of TCO and Tz functional groups on modified proteins stored in aqueous buffered media (e.g. maintaining > 90% reactivity after 1 month at 4°C, pH 7.5). This stability allows for worry-free conjugation results without the sensitive timing of reactions associated with classical thiol/maleimide chemistry.

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.

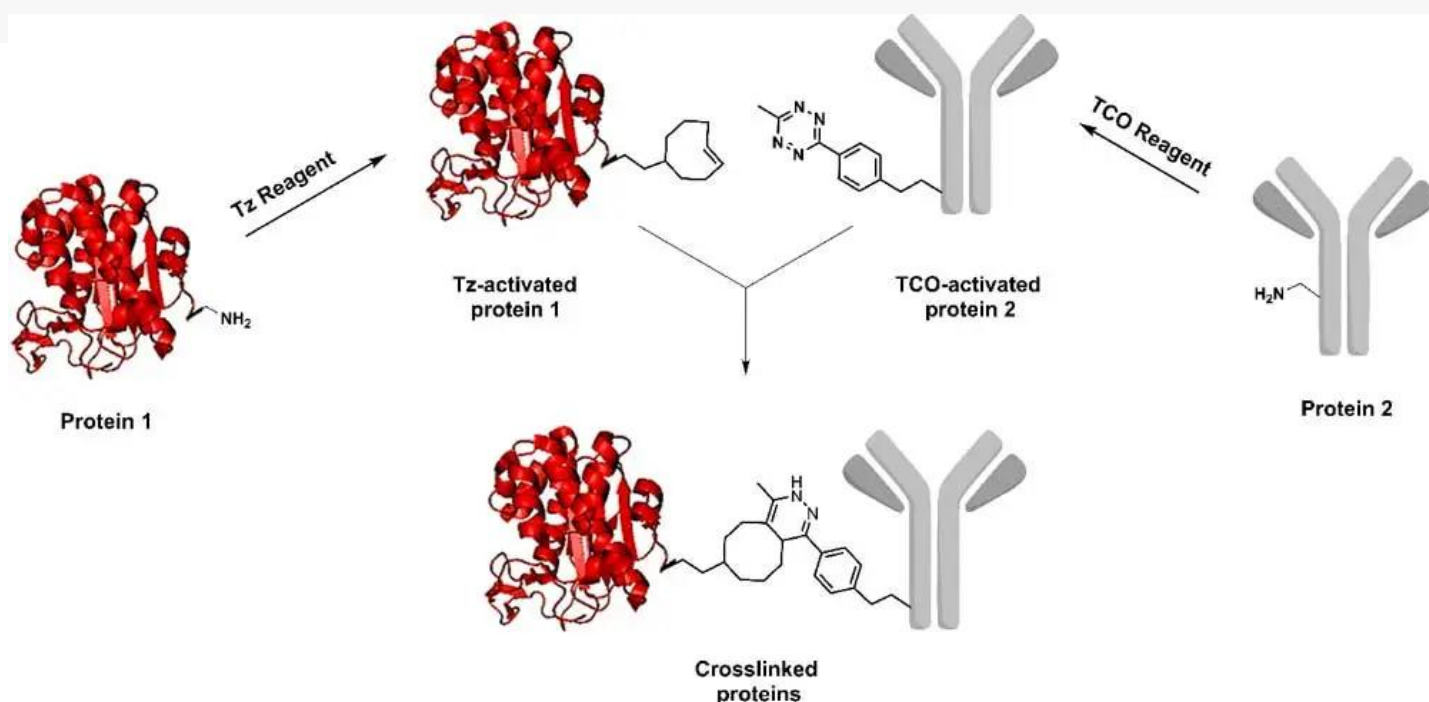


Figure 1. Schematic representation of protein-protein conjugation.

SPECIFICATIONS

Unit Size	1 Kit
Storage Instructions	4C
Number of Reactions	2
Quantity	2
Shipping Instructions	Ambient temperature

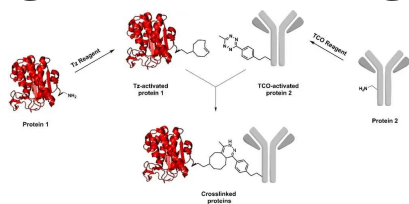
DOCUMENTS

- [Safety Data Sheet](#)
- [Download CoA](#)
- [User Guide](#)
- [Datasheet](#)

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



GALLERY IMAGES



For research use only. Not intended for therapeutic or diagnostic use in animals or humans.