



Protein-Protein Conjugation Kit

Product Images



Short Description

The Protein-Protein Conjugation Kit is designed to easily and efficiently conjugate any two proteins. This kit is flexible so that researchers with little or no bioconjugation experience can make their own custom protein-protein conjugates to suit their research needs. It includes all of the necessary components, including S-HyNic and S-4FB, for the rapid and specific crosslinking of any two proteins with molecular weights greater than 7 kDa. This kit contains sufficient reagents for the preparation of two protein-protein conjugates, with up to 650 µg of each protein.

Additional Information

Conjugation Targets	Protein
Label/Modifier Type	Protein
Reactivity	Amine
Recommended Storage	2-8°C
Applications	Antibody Labeling, In Situ Proximity Ligation

Products in this set



Protein-Protein Conjugation Kit, 1.00 Kit

