



Protein-Oligonucleotide Conjugation Kit

S-9011

[Product Images](#)



Short Description


The Protein-Oligo Conjugation Kit is designed to easily and efficiently prepare two separate protein-oligo conjugates. This kit is flexible so that researchers with little or no conjugation experience can make their own custom protein-oligo 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 protein with any oligo from 20 to 100 nucleotides in length.

The molecular weight range for proteins to be conjugated with this kit is 25,000 to 900,000 Daltons. Protein amounts from 50 – 650 µg at a concentration of 1 – 5 mg/mL in a volume of 50 – 130 µL are suitable for conjugation.

Additional Information

Conjugation Targets	Protein
Label/Modifier Type	Oligonucleotide
Reactivity	Amine
Recommended Storage	2° - 8°C – Do Not Freeze
Applications	Antibody Labeling, In Situ Proximity Ligation

Products in this set



Protein–Oligo Conjugation Kit, 1.00 Kit

