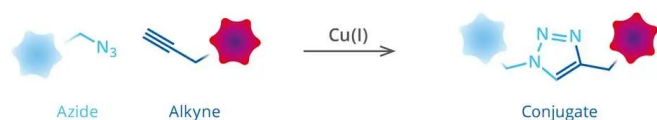




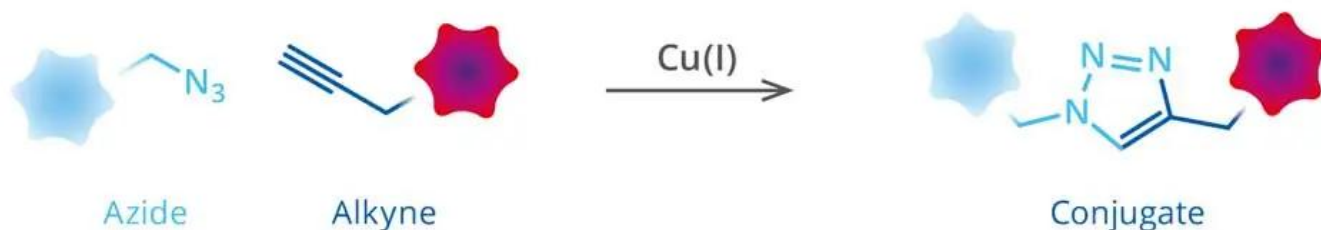
CLICK-&-GO® PROTEIN REACTION BUFFER KIT

SKU: CCT-1262



DESCRIPTION

Click-&-Go Click Chemistry Reaction Buffer Kit provides researchers who have biomolecules labeled with an azide or alkyne and the corresponding click detection reagent with all the necessary reagents to perform an up to 25 copper-catalyzed click reactions for subsequent analysis by gel electrophoresis, western blot or mass spectrometry.



The kit utilizes a copper chelating reagent to stabilize a catalytically active Cu(I) to accelerate the rate of CuAAC. The use of copper-chelating ligand also minimizes amount of free copper in solution, thus limiting the generation of reactive oxygen species (ROS) and protecting against undesired oxidative damage side reactions to proteins (e.g., GFP, RPE), nucleic acids (e.g., RNA, oligos), and other biomolecules.

SPECIFICATIONS

Unit Size	1 Kit
Storage Instructions	4C
Number of Reactions	25

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



Quantity

25

Shipping Instructions

Ambient temperature

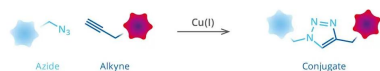
SELECTED REFERENCES

1. Bruna, R. E., *et al.* (2021). Limitation of phosphate assimilation maintains cytoplasmic magnesium homeostasis. *Proc Natl Acad Sci U S A.*, **118 (11)**, e2021370118. [[PubMed](#)]
2. Kwong, J. Q., *et al.* (2021). Mitochondrial functional resilience after TFAM ablation in the adult heart. *Am J Physiol Cell Physiol.*, **320 (6)**, C929-C942. [[PubMed](#)]
3. Coates, H. W. *et al.* (2021). The mammalian cholesterol synthesis enzyme squalene monooxygenase is proteasomally truncated to a constitutively active form. *J Biol Chem.*, Online ahead of print. [[PubMed](#)]

DOCUMENTS

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

GALLERY IMAGES



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