



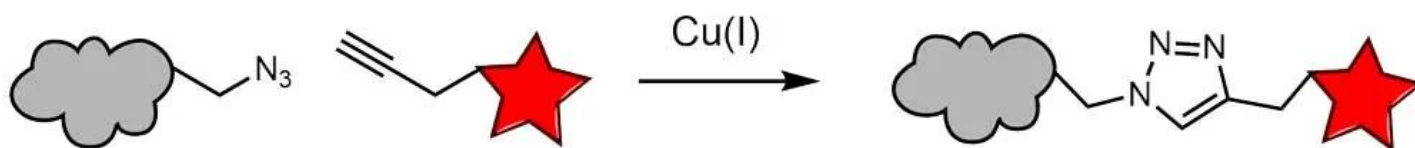
CLICK-&-GO® CELL REACTION BUFFER KIT

SKU: CCT-1263



DESCRIPTION

Click-&-Go™ Cell Reaction Buffer Kit provides researchers everything required to perform the click reaction on cells tagged with an azide or alkyne and with the corresponding click detection reagent for subsequent downstream analysis by flow cytometry, fluorescence microscopy, or high content screening (HCS).



Staining with additional detection reagents can be performed prior or after the click reaction. Each kit includes sufficient material to perform 50 reactions based on 0.5 mL volume.

SPECIFICATIONS

Unit Size	1 Kit
Storage Instructions	4C
Number of Reactions	50
Purchase Note	<p>Add some note</p>
Quantity	50
Shipping Instructions	Ambient temperature

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



SELECTED REFERENCES

1. Costello, A., *et al.* (2021). Directed evolution of rRNA improves translation kinetics and recombinant protein yield. *Nat Commun.*, **12 (1)**, 5638. [[PubMed](#)]

DOCUMENTS

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

GALLERY IMAGES



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