



CLICK-&-GO® PLUS 546 IMAGING KIT

SKU: CCT-1316



DESCRIPTION

Click-&-Go® Plus 546 Imaging Kit is an all-inclusive kit optimized for imaging alkyne-tagged biopolymers with orange-fluorescent AZDye 546 Azide Plus. Azide Plus reagents contain a complete copper-chelating system in their structure, allowing for the formation of strong, active copper complexes that act simultaneously as both reactant and catalyst in the CuAAC reaction. This azide-copper complex reacts almost instantaneously with alkynes under diluted conditions. This unprecedented reactivity in the CuAAC reaction is of special value for the detection of low abundance targets, improving biocompatibility, and is also valuable for any other application where greatly improved S/N ratio is highly desired. Sufficient materials is supplied to perform at least 30 cell or tissue assays based on a total reaction volume of 500 µL.

SPECIFICATIONS

Dye Type	AZDye 546
Unit Size	1 Kit
Storage Instructions	4 C
Laser Line	532 nm

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



Excitation/Emission Maximum	543/563
Number of Reactions	30
Purchase Note	Add some note
Shipping Instructions	Ambient temperature

SELECTED REFERENCES

1. Jiang, H., *et al.* (2014). Monitoring Dynamic Glycosylation in Vivo Using Supersensitive Click Chemistry. *Bioconjugate Chem.*, **25**, 698-706. [[PubMed](#)]
2. Uttamapinant, C., *et al.* (2012). Fast, Cell-Compatible Click Chemistry with Copper-Chelating Azides for Biomolecular Labeling. *Angew. Chem. Int. Ed.*, **51**, 5852-56. [[PubMed](#)]
3. Gaebler, A., *et al.* (2016). A highly sensitive protocol for microscopy of alkyne lipids and fluorescently tagged or immunostained proteins. *J. Lipid. Res.*, **57**, 1934-47. [[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.