



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## MTS-BCN

MTS-BCN is a fast-reacting and highly selective thiol-reactive label for attaching BCN to thiol groups on biomolecules. BCN, an alternative to DIBO and DBCO, reacts with azide to form 1,2,3-triazole by copper free 1,3-dipolar Huisgen cycloaddition.



### Product attributes

Chemical reactivity (reacts with)	Azides/Picolyl azides, Thiols
Functional group	BCN, MTS
Storage Conditions	Store at -10 to -35 °C

## Product Description

BCN (bicyclo[6.1.0]nonyne) with an MTS (methanethiosulfonate) reactive group is a fast-reacting, highly selective thiol-reactive label that can be used to attach BCN moieties to thiol groups on target molecules. BCN, an alternative to DIBO and DBCO, reacts with azide to form 1,2,3-triazole by copper-free 1,3-dipolar Huisgen cycloaddition. This copper-free bioorthogonal reaction allows staining the surface of live cells or in cell extracts when there are concerns about native protein function loss with copper.

- Thiol-reactive BCN derivative.
- BCN, an alternate to DIBO and DBCO, reacts with azide groups via copper-free bioorthogonal reaction.
- Supports copper-free bioorthogonal reaction for labeling the surface of live cells or in cell extracts.

This datasheet was generated on January 7, 2026 at 06:26:04 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/mts-bcn/>