



Glowing products for science

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

## Biotin Alkyne

Biotin alkyne reacts with azide to form 1,2,3-triazole by copper(I) catalyzed 1,3-dipolar Huisgen cycloaddition.



### Product attributes

## Product Description

Biotin alkyne reacts with azide to form 1,2,3-triazole by copper(I) catalyzed 1,3-dipolar Huisgen cycloaddition. It can also be used as a monomer to copolymerize with other monomers to form biotin labeled polymers.

This datasheet was generated on June 4, 2026 at 08:27:41 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/biotin-alkyne/>