



Glowing products for science

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

## Biotin Picolyl Azide

Biotin picolyl azide reacts with alkyne to form 1,2,3-triazole by 1,3-dipolar Huisgen cycloaddition through the use of a much lower copper (I) concentration without sacrificing reaction efficiency.



### Product attributes

## Product Description

Biotin picolyl azide reacts with alkyne to form 1,2,3-triazole by 1,3-dipolar Huisgen cycloaddition through the use of a much lower copper (I) concentration without sacrificing reaction efficiency.

This datasheet was generated on December 22, 2025 at 10:32:36 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/biotin-picolyl-azide/>