



Glowing products for science

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

Biotin Azide

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



Product attributes

Product Description

Biotin azide reacts with alkyne to form 1,2,3-triazole by copper(I) catalyzed 1,3-dipolar Huisgen cycloaddition. It can also undergo Staudinger ligation with phosphine containing compounds.

This datasheet was generated on June 4, 2026 at 08:23:50 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-azide/>