Glowing products for science

Biotin Azide

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



Product attributes

Call us: 800-304-5357 Email: btinfo@biotium.com

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 November 3, 2025 at 05:13:44 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-azide/