Glowing products for science

Biotin Alkyne

Biotin alkyne reacts with azide 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 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 November 3, 2025 at 11:20:04 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-alkyne/