Glowing products for science

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

Call us: 800-304-5357

Email: btinfo@biotium.com

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 November 3, 2025 at 10:57:44 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-picolyl-azide/