

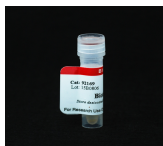


Glowing products for science

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

Biotin BCN

Biotin BCN, an alternative to biotin labeled DIBO and DBCO, reacts with azide to form 1,2,3-triazole by copper free 1,3-dipolar Huisgen cycloaddition.



Product attributes

Chemical reactivity (reacts with)	Azides/Picolyl azides
Functional group	BCN
Storage Conditions	Store at -10 to -35 °C

Product Description

Biotin BCN reacts with azide to form 1,2,3-triazole by copper free 1,3-dipolar Huisgen cycloaddition.

- An alternative to biotin labeled DIBO and DBCO
- Copper free bioorthogonal reaction allows staining the surface of live cells or cell extracts when there are concerns about native protein function loss with copper

Biotin molecules can be detected with any of our [CF@ dye streptavidin conjugates](#). Learn more about [CF@ Dyes](#) or download the [CF@ Dye Brochure](#) for more information.

This datasheet was generated on January 6, 2026 at 05:13:27 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-bcn/>