



Glowing products for science

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

Biotin Methyltetrazine

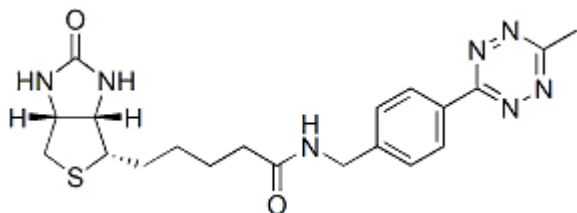
Biotin with methyltetrazine reacts with cyclooctene (TCO, methylcyclopropene) labeled molecule via copper free click chemistry.



Product attributes

Product Description

Biotin with methyltetrazine reacts with cyclooctene (TCO, methylcyclopropene) containing molecules via copper free click chemistry. This copper free bioorthogonal reaction is useful for labeling live cells, or when there are concerns about native protein function loss with copper in cell extracts.



This datasheet was generated on January 8, 2026 at 10:04:09 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-methyltetrazine/>