Glowing products for science

## **Biotin TCO**

Biotin with TCO (trans-cyclooctene) reacts with tetrazine labeled molecule via copper free click chemistry.



## **Product attributes**

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

## **Product Description**

Biotin with TCO (trans-cyclooctene) reacts with tetrazine 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 November 4, 2025 at 10:49:38 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biotin-tco/">https://biotium.com/product/biotin-tco/</a>