Glowing products for science

## **Biotin MTSEA**

MTSEA biotin selectively labels thiols under mild conditions. It's available in 3 variations: Biotin, Biotin-X, and Biotin-XX, all of which allow selective labeling of cell surface thiols.



## **Product attributes**

Conjugation

Biotin, Biotin-X, Biotin-XX

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

## **Product Description**

MTSEA biotin selectively labels thiols under mild conditions. It's available in 3 variations with different linker lengths, all of which allow selective labeling of cell surface thiols. Longer linkers should facilitate interaction between biotin and streptavidin.

## References

- 1. J Membr Biol 180 (1), 83 (2001).

- 2. J Neurosci 21 (1), 67 (2001). 3. J Neurosci 19(12), 4847 (1999) 4. J Biol Chem 273(20), 12675 (1998).

This datasheet was generated on November 8, 2025 at 04:52:04 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-mtsea/