



Glowing products for science

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

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

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 January 9, 2026 at 12:51:33 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-mtsea/>