



Glowing products for science

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

Biotin TS Ethylenediamine

TS biotin ethylenediamine (Biotin ethylenediamine thiosulfate, sodium salt) is a highly selective thiol labeling probe. Similar to the MTS group, TS (thiosulfate) group is capable of reacting with thiols at nearly neutral pH.



Product attributes

Conjugation
Biotin, Biotin-X, Biotin-XX

Product Description

TS biotin ethylenediamine (Biotin ethylenediamine thiosulfate, sodium salt) is a highly selective thiol labeling probe. Similar to the MTS group, TS (thiosulfate) group is capable of reacting with thiols at nearly neutral pH. However, unlike the MTS group, which is neutral, TS is negatively charged and thus probes bearing a TS group are ideal for specifically labeling thiols exposed at the exterior cell surface because charged probes are less likely to enter cells.

It's available in 3 variations: Biotin, Biotin-X, and Biotin-XX. The X and XX spacer should facilitate the interaction between biotin and avidin or streptavidin.

TS Biotin ethylenediamine

- Biotin ethylenediamine thiosulfate, sodium salt
- $C_{14}H_{23}N_4NaO_6S_3$
- MW: 462.54
- Off-white solid soluble in DMSO or water
- Store desiccated at $-20^{\circ}C$

TS Biotin-X ethylenediamine

- Biotin-X ethylenediamine thiosulfate, sodium salt
- $C_{20}H_{34}N_5NaO_7S_3$
- MW: 575.7
- Off-white solid soluble in DMSO or water
- Store desiccated at $-20^{\circ}C$

TS Biotin-XX ethylenediamine

- Biotin-XX ethylenediamine thiosulfate, sodium salt
- $C_{26}H_{45}N_6NaO_8S_3$
- MW: 688.86
- Off-white solid soluble in DMSO or water
- Store desiccated at $-20^{\circ}C$

This datasheet was generated on January 8, 2026 at 05:10:06 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-ts-ethylenediamine/>