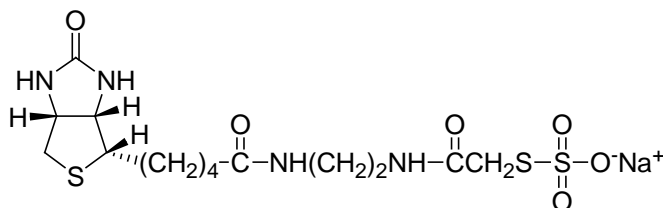


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: TS Biotin ethylenediamine (biotin ethylenediamine thiosulfate, sodium salt)**CATALOG #:** 92010**MOLECULAR INFORMATION:** C₁₄H₂₃N₄NaO₆S₃
MW: 462.54**PROPERTIES:**

Color & Form Off-white solid
Purity ≥ 95% by TLC
Solubility Soluble in DMSO or water

STORAGE AND HANDLING: Store desiccated at -20°C

APPLICATION: TS biotin ethylenediamine 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 neutral MTS group, 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.

RELATED PRODUCTS *TS Bioti-X ethylenediamine (cat# 92011); TS Biotin-XX ethylenediamine (cat# 92012)*

TOXICITY: Unknown.

FIRST AID: Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.