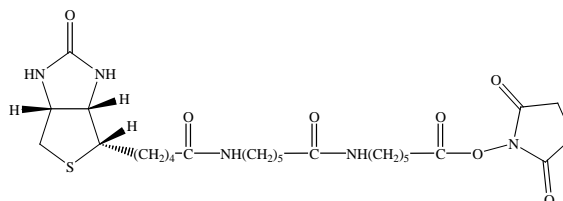


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **Biotin-XX, succinimidyl ester** (6-((6-((biotinoyl)amino)hexanoyl)amino)hexanoic acid succinimidyl ester)

CATALOG #: **90052**

MOLECULAR INFORMATION: C₂₆H₄₁N₅O₇S
MW: 567.7

**PROPERTIES:**

Color & Form White solid
Purity ≥ 95% by TLC
Solubility Soluble in anhydrous DMSO

STORAGE AND HANDLING:

Store desiccated at -20 °C. Keep away from moisture.

APPLICATION:

Biotin-XX succinimidyl ester is readily coupled to a primary or secondary amine under mild conditions. The longer XX arm facilitates the interaction between biotin and avidin or streptavidin.

Ref.: 1) *Bioconjugate Chem* **10**, 32 (1999); 2) *J Cell Biol* **144**, 99 (1999); 3) *Anal. Chem.* **70**, 1242 (1998); 4) *Meth Mol Biol* **80**, 173 (1998).

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.