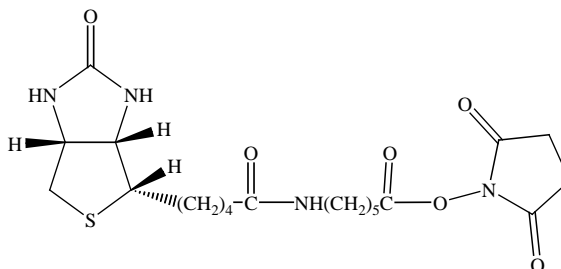


PRODUCT AND SAFETY DATA SHEET

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

CATALOG #: **90051**

MOLECULAR INFORMATION: $C_{20}H_{30}N_4O_6S$
MW: 454.54

**PROPERTIES:**

Color & Form

Off-white to pinkish solid

Purity

≥ 95% by TLC

Solubility

Soluble in DMSO and moderately soluble in DMF

STORAGE AND HANDLING:

Store desiccated at -20 °C.

APPLICATION:

Biotin-X succinimidyl ester is readily coupled to a primary or secondary amine under mild conditions.

Ref.: 1) *J Biol Chem* **274**, 21878 (1999); 2) *J Biol Chem* **273**, 793 (1998); 3) *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.