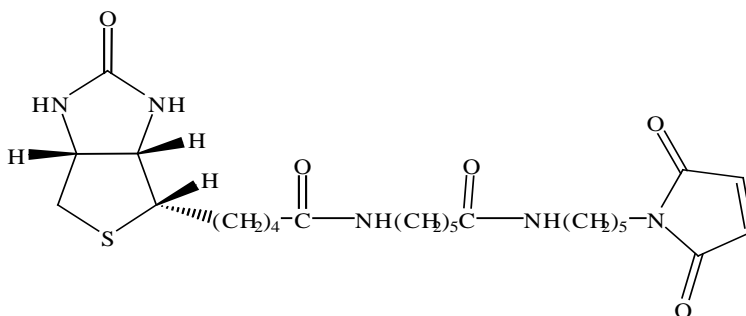


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **Biotin-X-C5-maleimide** (N-(5-(6-((biotinoyl)amino)hexanoyl)amino)pentylmaleimide)

CATALOG #: **90058**

MOLECULAR INFORMATION: C₂₇H₃₉N₅O₇
S
MW: 577

**PROPERTIES:**

Color & Form White solid
Purity ≥ 95% by HPLC
Solubility Soluble in DMSO or DMF

STORAGE AND HANDLING:

Store at -20 °C

APPLICATION:

Biotin-X-C5-maleimide is readily and selectively coupled to thiols under mild conditions. The long linker between biotin and the maleimide group ensures the biotin group to be easily accessible by avidin or streptavidin.

For further information, please see attached “**Protocol for Using Maleimide Products**”

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: <i>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.</i>
--