

PRODUCT AND SAFETY DATA SHEET

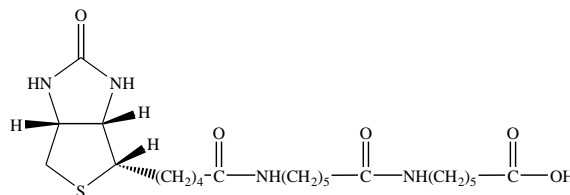
PRODUCT NAME: **Biotin-XX, free acid** (6-((6-((biotinoyl)amino)hexanoyl)amino)hexanoic acid)

CATALOG #: **90054**

MOLECULAR INFORMATION:

C₂₂H₃₈N₄O₅S

MW: 470



PROPERTIES:

Color & Form

White solid

Purity

≥ 95% by HPLC

Solubility

Soluble in DMSO and moderately soluble in DMF

STORAGE AND HANDLING:

Store at 4°C.

APPLICATION:

Biotin-XX free acid can be easily converted to the mixed anhydride form, which is more reactive than the succinimidyl ester and thus can be used to react with aromatic amines or sterically hindered amines.

Ref.: a detailed protocol for making mixed anhydrides is provided upon request with purchase of the product.

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.