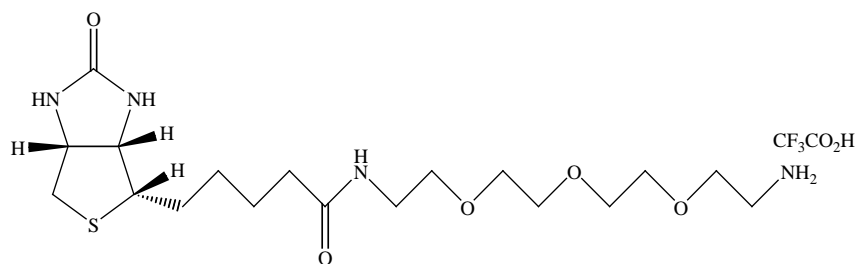


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Biotin-PEO₄-amine**CATALOG #:** 90068**MOLECULAR INFORMATION:** C₂₀H₃₅F₃N₄O₇S
Mwt: 532.58**PROPERTIES:**

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

STORAGE AND HANDLING: Highly hygroscopic! Store desiccated at 4 °C.

APPLICATION: Biotin-PEO₄-amine is a water soluble biotin derivative that can be readily coupled to activated carboxylic acids

Ref.: *Biochemistry* 21, 978(1982).

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.