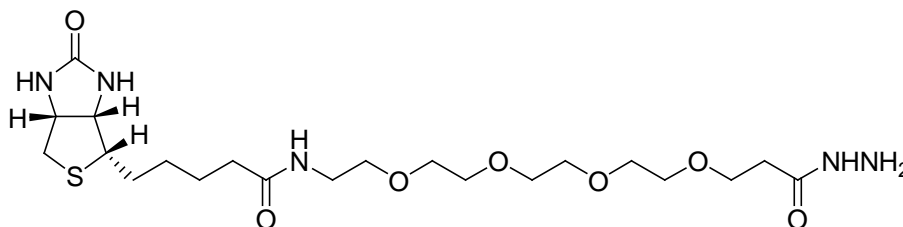


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Biotin-PEO<sub>4</sub>-hydrazide**CATALOG #:** 90111**MOLECULAR INFORMATION:** C<sub>21</sub>H<sub>39</sub>N<sub>5</sub>O<sub>7</sub>S  
Mwt: 505.63**PROPERTIES:****Color & Form**

White solid

**Purity**

≥ 90% by TLC

**Solubility**

Soluble in water or DMSO. If the solution will be stored away for later use, dissolve the solid material in anhydrous DMSO.

**STORAGE AND HANDLING:**

Highly hygroscopic! Store solid desiccated at 4 °C.

**APPLICATION:**Biotin-PEO<sub>4</sub>-hydrazide is a water soluble biotin derivative capable of reacting with aldehydes or ketones to form hydrazones.**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.