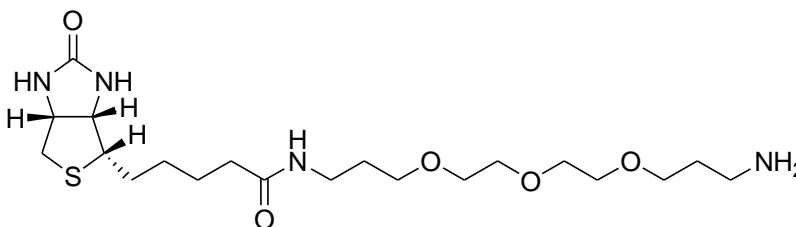


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Biotin-PEO<sub>2</sub>-PPO<sub>2</sub>-amine, trifluoroacetic acid salt

**CATALOG #:** 90078

**MOLECULAR INFORMATION:** C<sub>22</sub>H<sub>39</sub>F<sub>3</sub>N<sub>4</sub>O<sub>7</sub>S  
Mwt: 560.63



**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-PEO<sub>2</sub>-PPO<sub>2</sub>-amine is a water soluble biotin derivative with an amine group.

**TOXICITY:** Unknown.

<b>FIRST AID:</b>	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.
-------------------	---

<b>Disclaimer:</b> <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>
--