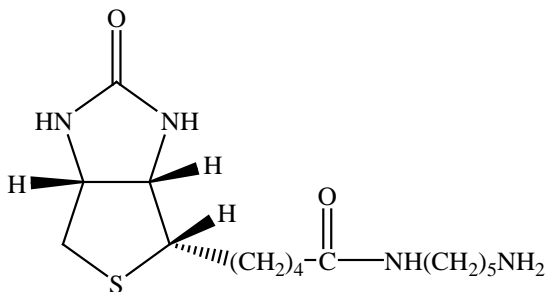


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Biotin cadaverine, free base (5-((N-(biotinoyl)amino) pentylamine)**CATALOG #:** 90079**MOLECULAR INFORMATION:** C<sub>15</sub>H<sub>28</sub>N<sub>4</sub>O<sub>2</sub>S  
MW: 328.47**PROPERTIES:****Color & Form** Tan solid  
**Purity** ≥ 99% by TLC  
**Solubility** Soluble in DMSO**STORAGE AND HANDLING:** Store at 4 °C**APPLICATION:** Biotin cadaverine, free base can be coupled directly to activated carboxylic acid groups such as succinimidyl ester without adding an additional base.**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.