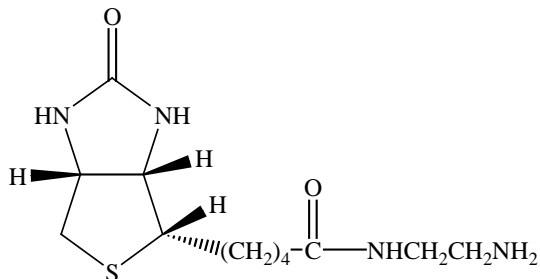


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Biotin ethylenediamine (free base form)**CATALOG #:** 90071**MOLECULAR INFORMATION:** C<sub>12</sub>H<sub>22</sub>N<sub>4</sub>O<sub>2</sub>S  
Mwt: 286**PROPERTIES:**

<b>Color &amp; Form</b>	White solid
<b>Purity</b>	≥ 95% by TLC
<b>Solubility</b>	Soluble in DMSO

**STORAGE AND HANDLING:** Store desiccated at ≤4°C.**APPLICATION:** Similar to Biotin ethylenediamine hydrobromide (**90057**), which is identical to Neurobiotin ( a trademark of Vector Laboratories), and is a useful anterograde and transneuronal tracers.Ref.: 1) *J. Neurosci.* **14**, 3805(1994); 2) *J. Neurosci. Meth.* **41**, 31(1992); 3) *J. Neurosci. Meth.* **37**, 141(1991).**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>
--