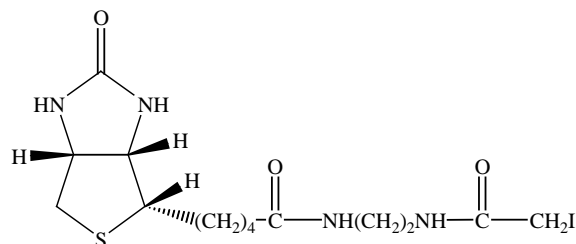


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **Biotin ethylenediamine iodoacetamide** (N-(biotinoyl)-N'-(iodoacetyl)ethylenediamine)

CATALOG #: **90059**

MOLECULAR INFORMATION: $C_{14}H_{23}IN_4O_3S$
MW: 454.33

**PROPERTIES:**

Color & Form Brown solid
Purity ≥ 90% by TLC
Solubility Soluble in DMF or DMSO

STORAGE AND HANDLING:

Store desiccated at -20 °C upon receipt. Protect from light, especially when in solution

APPLICATION:

Biotin ethylenediamine iodoacetamide is readily coupled to thiols.

Ref.: 1) *J Biol Chem* **274**, 4722 (1999); 2) *Biochem. J.* **251**, 935(1988)

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: <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>
