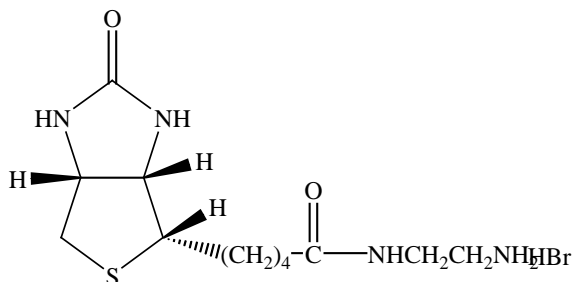


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **Biotin ethylenediamine, hydrobromide** (*N*-(2-aminoethyl)biotinamide, hydrobromide) (equivalent to Neurobiotin)

CATALOG #: **90057**

MOLECULAR INFORMATION: C₁₂H₂₃BrN₄O₂S
MW: 367

**PROPERTIES:**

Color & Form White solid
Purity ≥ 95% by TLC
Solubility Soluble in DMSO or water (solubility in water is at least 100mg/mL)

STORAGE AND HANDLING:

Store at 4°C

APPLICATION:

Biotin ethylenediamine, identical to Neurobiotin (a trademark of Vector Laboratories), 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.

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.