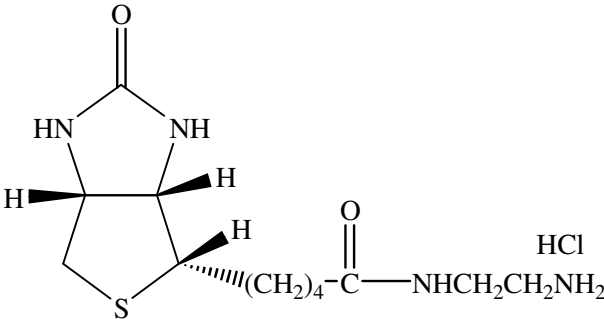


Functionally equivalent to Neurobiotin

## Instruction Manual

Catalog Number	PK-CA707-90075
Description	Biotin-Ethylenediamine, hydrochloride is similar to Biotin-Ethylenediamine, hydrobromide (PK-CA707-90057) and identical to Neurobiotin (a trademark of Vector Laboratories), which is a useful anterograde and transneuronal tracers.
Quantity	25 mg
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	322.5 Da; C <sub>12</sub> H <sub>23</sub> N <sub>4</sub> O <sub>2</sub> S
Purity	>99% (as determined by TLC)
Appearance / Formulation / Solubility	White solid; soluble in DMSO or water.
Storage & Stability	Store at 4°C.
Applications	See Description
References	<ol style="list-style-type: none"> <li>1) J. Neurosci. 14, 3805((1994)</li> <li>2) J. Neurosci. Meth. 41, 31(1992)</li> <li>3) J. Neurosci. Meth. 37, 141(1991)</li> </ol>
Caution	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.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.