



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Biotin Ethylenediamine, Hydrochloride

Biotin ethylenediamine, hydrochloride is functionally equivalent to Neurobiotin™ (biotin ethylenediamine, hydrobromide, [90057](#)).



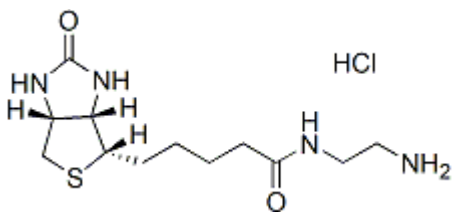
Product attributes

Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Conjugation	Biotin

Product Description

Biotin ethylenediamine, hydrochloride is functionally equivalent to Neurobiotin™ (biotin ethylenediamine, hydrobromide, [90057](#)).

- While solid soluble in DMSO or water
- MW: 322.5



Neurobiotin is a trademark of Vector Laboratories.

This datasheet was generated on June 15, 2026 at 03:33:49 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-ethylenediamine-hydrochloride-functionally-equivalent-to-neurobiotintm/>