

## Biotin Ethylenediamine, Hydrobromide (Neurobiotin™)

Biotin ethylenediamine, hydrobromide (N-(5-(6-((biotinoyl)amino)hexanoyl)amino)pentylmaleimide) is identical to Neurobiotin™, a useful anterograde and transneuronal tracer.



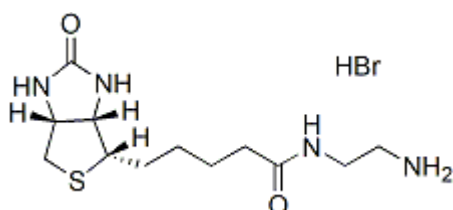
### Product attributes

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

### Product Description

Biotin ethylenediamine, hydrobromide (N-(5-(6-((biotinoyl)amino)hexanoyl)amino)pentylmaleimide) is identical to Neurobiotin™, a useful anterograde and transneuronal tracer.

- White solid soluble in DMSO
- Store at 4°C
- C<sub>12</sub>H<sub>23</sub>BrN<sub>4</sub>O<sub>2</sub>S
- MW: 367



Neurobiotin is a trademark of Vector Laboratories.

### References

1. J Neurosci 14, 3805 (1994).
2. J Neurosci Meth 41, 31 (1992).
3. J Neurosci Meth 37, 141 (1991).