



Glowing products for science

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

Biotin Ethylenediamine (free base)

Biotin ethylenediamine in free base form. White solid soluble in DMSO.



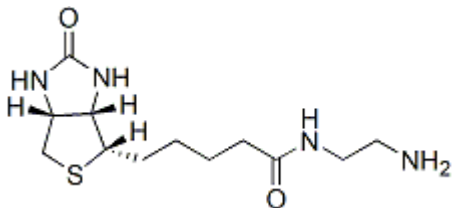
Product attributes

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

Product Description

Biotin ethylenediamine in free base form. White solid soluble in DMSO.

- Store at 4 °C
- $C_{12}H_{22}N_4O_2S$
- MW: 286



This datasheet was generated on June 22, 2026 at 03:02:40 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-ethylenediamine-free-base-form/>