Biotin Ethylenediamine (free base)

Biotin ethylenediamine in free base form. White solid soluble in $\ensuremath{\mathsf{DMSO}}$.



Product attributes

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

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

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

- Store at 4°C
- C₁₂H₂₂N₄O₂S
- MW: 286

This datasheet was generated on November 7, 2025 at 07:38:43 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-ethylenediamine-free-base-form/