Biotin Ethylenediamine Iodoacetamide

Biotin ethylenediamine iodoacetamide is readily coupled to thiols.



Product attributes

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

Product Description

Biotin ethylenediamine iodoacetamide is readily coupled to thiols.

- Off-white solid soluble in DMF
- Store at -20°C
- C₁₄H₂₃IN₄O₃S
- MW: 454.33

References

- 1. J Biol Chem 274, 4722 (1999).
- 2. Biochem J 251, 935 (1988).

This datasheet was generated on November 8, 2025 at 09:47:43 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-ethylenediamine-iodoacetamide-n-biotinoyl-n-iodoacetylethylenediamine/