

Biotin Ethylenediamine Iodoacetamide

Biotin ethylenediamine iodoacetamide is readily coupled to thiols.

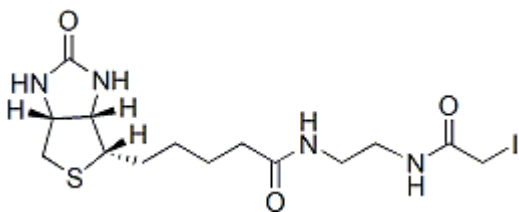


Product attributes

Product Description

Biotin ethylenediamine iodoacetamide is readily coupled to thiols.

- Off-white solid soluble in DMF
- Store at -20 °C
- C₁₄H₂₃N₄O₃S
- MW: 454.33



References

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

This datasheet was generated on June 22, 2026 at 06:22:49 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-ethylenediamine-iodoacetamide-n-biotinoyl-n-iodoacetyleneethylenediamine/>