Biotin-PEO4-amine

Biotin-PEO₄-amine is a water soluble biotin derivative that can be readily coupled to activated carboxylic acids.



Call us: 800-304-5357 Product attributes

CAS number

359860-27-8

Product Description

Biotin-PEO₄-amine is a water soluble biotin derivative that can be readily coupled to activated carboxylic acids.

- White solid soluble in water or DMSO
- Store at 4°C
- C₁₈H₃₄N₄O₅S
- MW: 419
- [359860-27-8]

References

1. Biochemistry 21, 978 (1982).

This datasheet was generated on November 28, 2025 at 05:53:10 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-peo4-amine/