

Biotin-PEO4-amine

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



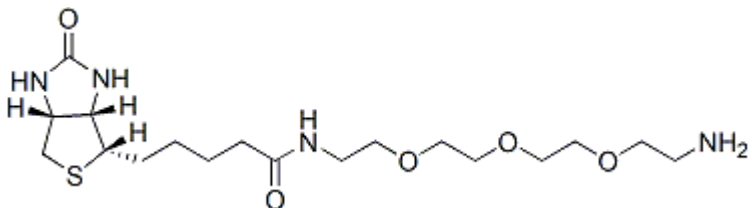
Product attributes

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 June 3, 2026 at 08:43:59 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-peo4-amine/>