

Biotin-PEO3-amine

Biotin-PEO₃-amine is a water soluble biotin derivative that can be readily coupled to activated carboxylic acids.



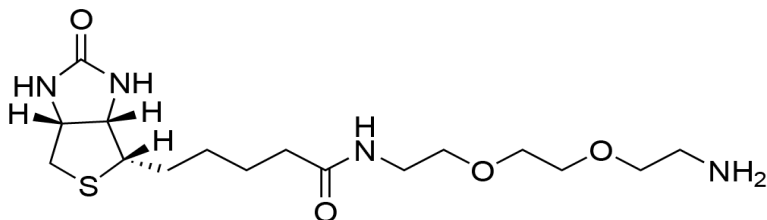
Product attributes

CAS number	138529-46-1
------------	-------------

Product Description

Biotin-PEO₃-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 -20 °C
- C₁₆H₃₀N₄O₄S
- MW: 375
- [138529-46-1]



References

1. Biochemistry 21, 978 (1982).

This datasheet was generated on June 3, 2026 at 08:43:26 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-peo3-amine/>