



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Biotin-PEO4-propionate SE

Biotin-PEO4-propionate SE (Biotin-PEO4-propionate succinimidyl ester) is a water soluble biotin derivative that can be readily coupled to primary or secondary amines.

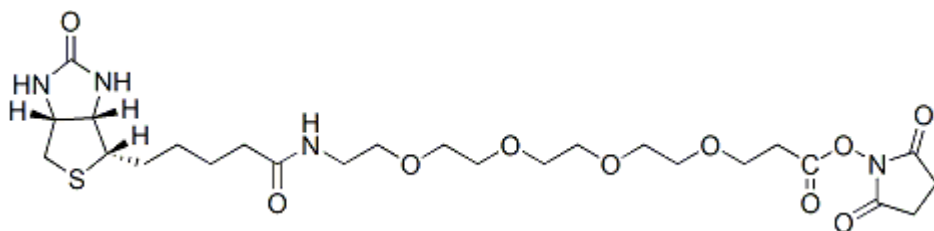


Product attributes

Product Description

Biotin-PEO4-propionate SE (Biotin-PEO4-propionate succinimidyl ester) is a water soluble biotin derivative that can be readily coupled to primary or secondary amines.

- White solid soluble in water or DMSO
- Store at -20 °C
- $C_{25}H_{40}N_4O_{10}S$
- MW: 588.7.



This datasheet was generated on January 21, 2026 at 08:27:50 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-peo4-propionate-succinimidyl-ester/>