

## Biotin-PEO4-maleimide

Biotin-PEO4-maleimide is a water soluble biotin derivative that can be readily coupled to thiols.

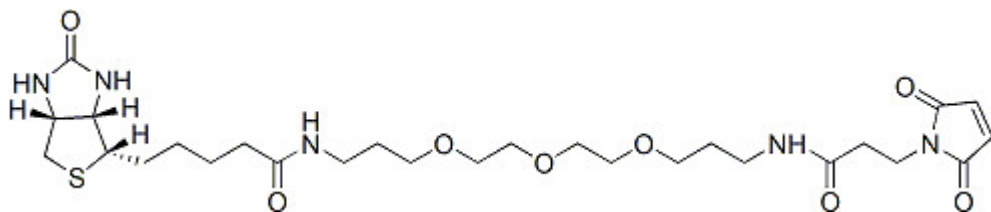


### Product attributes

## Product Description

Biotin-PEO4-maleimide is a water soluble biotin derivative that can be readily coupled to thiols.

- White solid soluble in water or DMSO
- Store at -20 °C
- $C_{27}H_{43}N_5O_8S$
- MW: 597.73



This datasheet was generated on January 26, 2026 at 03:03:18 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/biotin-peo4-maleimide/>