

## Biotin-PEO3-amine

Biotin-PEO<sub>3</sub>-amine is a water soluble biotin derivative that can be readily coupled to activated carboxylic acids.



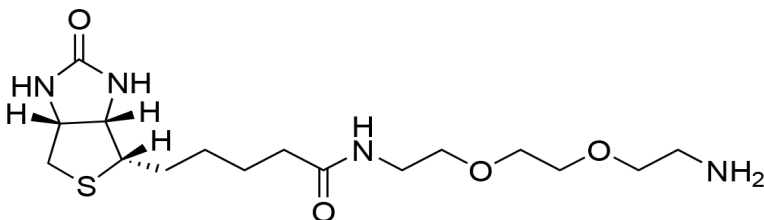
### Product attributes

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

## Product Description

Biotin-PEO<sub>3</sub>-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
- C<sub>16</sub>H<sub>30</sub>N<sub>4</sub>O<sub>4</sub>S
- MW: 375
- [138529-46-1]



## References

1. Biochemistry 21, 978 (1982).

This datasheet was generated on January 7, 2026 at 02:31:21 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/biotin-peo3-amine/>