## **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

Call us: 800-304-5357 Email: btinfo@biotium.com

## **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 November 5, 2025 at 02:59:13 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biotin-peo3-amine/">https://biotium.com/product/biotin-peo3-amine/</a>