## **Biotin-PEO4-maleimide**

Biotin-PEO<sub>4</sub>-maleimide is a water soluble biotin derivative that can be readily coupled to thiols.



## **Product attributes**

Call us: 800-304-5357

## **Product Description**

Biotin-PEO<sub>4</sub>-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<sub>27</sub>H<sub>43</sub>N<sub>5</sub>O<sub>8</sub>S
- MW: 597.73

This datasheet was generated on November 27, 2025 at 12:56:48 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biotin-peo4-maleimide/">https://biotium.com/product/biotin-peo4-maleimide/</a>