## **Biotin-PEO4-propionate SE**

Biotin-PEO<sub>4</sub>-propionate SE (Biotin-PEO<sub>4</sub>-propionate succinimidyl ester) is a water soluble biotin derivative that can be readily coupled to primary or secondary amines.



## **Product attributes**

Call us: 800-304-5357

## **Product Description**

Biotin-PEO<sub>4</sub>-propionate SE (Biotin-PEO<sub>4</sub>-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<sub>25</sub>H<sub>40</sub>N<sub>4</sub>O<sub>10</sub>S
- MW: 588.7.

This datasheet was generated on November 28, 2025 at 05:58:42 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biotin-peo4-propionate-succinimidyl-ester/">https://biotium.com/product/biotin-peo4-propionate-succinimidyl-ester/</a>