## Biotin-PEO4-hydrazide

Biotin-PEO<sub>4</sub>-hydrazide is a water-soluble biotin derivative capable of reacting with aldehydes or ketones to form hydrazones.



## **Product attributes**

Call us: 800-304-5357

## **Product Description**

Biotin-PEO<sub>4</sub>-hydrazide is a water-soluble biotin derivative capable of reacting with aldehydes or ketones to form hydrazones.

- White solid soluble in water
- Store at 4°C
- C<sub>21</sub>H<sub>39</sub>N<sub>5</sub>O<sub>7</sub>S
- MW: 505.63

This datasheet was generated on November 28, 2025 at 05:55:37 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biotin-peo4-hydrazide/">https://biotium.com/product/biotin-peo4-hydrazide/</a>