

Biotin-PEO4-hydrazide

Biotin-PEO4-hydrazide is a water-soluble biotin derivative capable of reacting with aldehydes or ketones to form hydrazones.

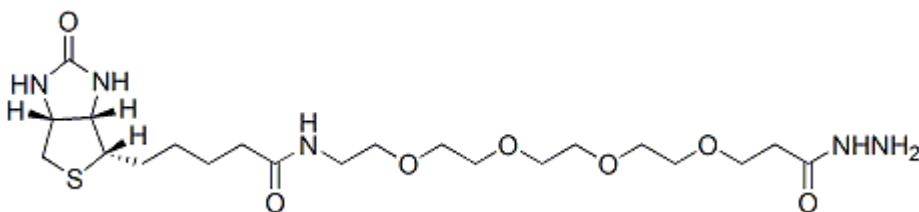


Product attributes

Product Description

Biotin-PEO4-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
- C₂₁H₃₉N₅O₇S
- MW: 505.63



This datasheet was generated on June 22, 2026 at 03:34:46 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-peo4-hydrazide/>