

## Biotin-12 SE

Biotin-12 SE is an amine-reactive biotin derivative with a long spacer group. It is useful for making biotin labeled nucleoside triphosphates among other applications.

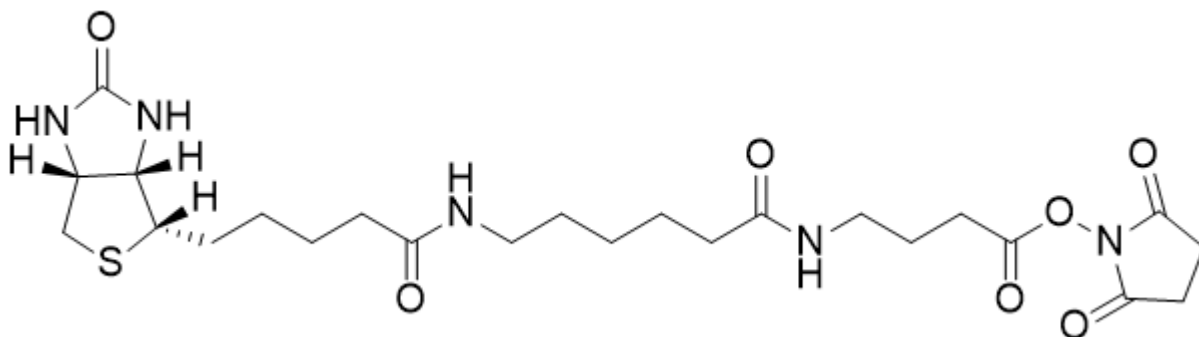


### Product attributes

## Product Description

Biotin-12 SE is an amine-reactive biotin derivative with a long spacer group. It is useful for making biotin labeled nucleoside triphosphates among other applications.

- White solid soluble in DMSO or DMF
- Store at -20 °C
- $C_{24}H_{37}N_5O_7S$
- MW: 539.64



This datasheet was generated on August 10, 2026 at 02:29:40 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-12-se-4-6-biotinoylamino-hexanoylamino-butyl-ester/>