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

Call us: 800-304-5357

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
- C₂₄H₃₇N₅O₇S
- MW: 539.64

This datasheet was generated on November 4, 2025 at 10:35:57 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-12-se-4-6-biotinoylaminohexanoylaminobutyric-acid-succinimidyl-ester/