Biotin-X, succinimidyl ester

Biotin-X succinimidyl ester (6-((Biotinoyl)amino)hexanoic acid succinimidyl ester) is readily coupled to a primary or secondary amine under mild conditions.



Call us: 800-304-5357

Product attributes

CAS number

72040-63-2

Product Description

Biotin-X succinimidyl ester (6-((Biotinoyl)amino)hexanoic acid succinimidyl ester) is readily coupled to a primary or secondary amine under mild conditions.

- White solid soluble in DMSO and moderately soluble in DMF
- Store at -20°C
- C₂₀H₃₀N₄O₆S
- MW: 454.54
- [72040-63-2]

References

- 1. J Biol Chem 274, 21878 (1999).
- 2. J Biol Chem 273, 793 (1998).
- 3. Meth Mol Biol 80, 173 (1998).

This datasheet was generated on November 28, 2025 at 05:53:10 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-x-succinimidyl-ester-6-biotinoylaminohexanoic-acid-succinimidyl-ester/