

Biotin-XX, succinimidyl ester

Biotin-XX succinimidyl ester (6-((6-((Biotinoyl)amino)hexanoyl)amino)hexanoic acid succinimidyl ester) is readily coupled to a primary or secondary amine under mild conditions. The longer XX arm facilitates the interaction between biotin and avidin or streptavidin.

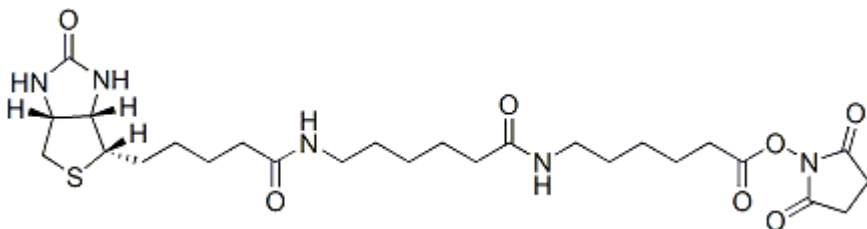


Product attributes

Product Description

Biotin-XX succinimidyl ester (6-((6-((Biotinoyl)amino)hexanoyl)amino)hexanoic acid succinimidyl ester) is readily coupled to a primary or secondary amine under mild conditions. The longer XX arm facilitates the interaction between biotin and avidin or streptavidin.

- White solid soluble in DMSO and moderately soluble in DMF
- Store at -20 °C
- C₂₆H₄₁N₅O₇S
- MW: 567.7
- [89889-52-1]



References

1. Bioconjugate Chem 10, 32 (1999).
2. J Cell Biol 144, 99 (1999).
3. Anal Chem 70, 1242 (1998).
4. Meth Mol Biol 80, 173 (1998).

This datasheet was generated on June 3, 2026 at 07:14:30 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-xx-succinimidyl-ester-6-6-biotinoylaminohexanoylaminohexanoic-acid-succinimidyl-ester/>