

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.



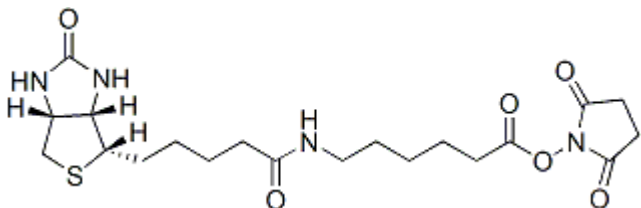
Product attributes

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
- $C_{20}H_{30}N_4O_6S$
- 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 January 28, 2026 at 08:20:55 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-x-succinimidyl-ester-6-biotinoylamino-hexanoic-acid-succinimidyl-ester/>