

Biotin-X, free acid

Biotin-X free acid (6-((Biotinoyl)amino)hexanoic acid) can be easily converted to the mixed anhydride form, which is more reactive than the succinimidyl ester and thus can be used to react with aromatic amines or sterically hindered amines.

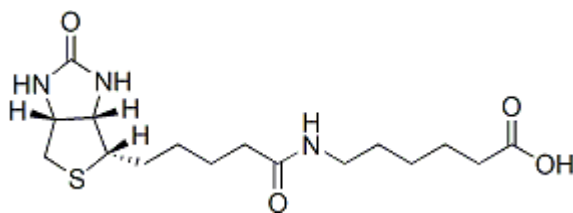


Product attributes

Product Description

Biotin-X free acid (6-((Biotinoyl)amino)hexanoic acid) can be easily converted to the mixed anhydride form, which is more reactive than the succinimidyl ester and thus can be used to react with aromatic amines or sterically hindered amines.

- White solid soluble in DMSO and moderately soluble in DMF
- Store at 4 °C
- $C_{16}H_{27}N_3O_4S$
- MW: 357.47
- [72040-64-3]



References

1. Anal Chem 67, 1014 (1995).

This datasheet was generated on January 28, 2026 at 06:35:57 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-x-free-acid-6-biotinoylaminohexanoic-acid/>