

Biotin-XX, free acid

Biotin-XX free acid

(6-((6-((Biotinoyl)amino)hexanoyl)amino)hexanoic acid) can be easily converted to the mixed anhydride form, which is more reactive than the succinimidyl ester.

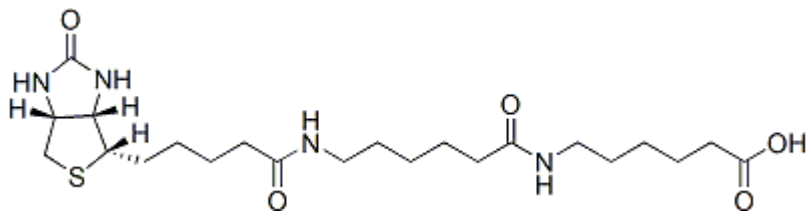


Product attributes

Product Description

Biotin-XX free acid (6-((6-((Biotinoyl)amino)hexanoyl)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. A detailed protocol for making mixed anhydrides is provided upon request with purchase of the product.

- White solid soluble in DMSO (moderately soluble in DMF)
- MW: 470



This datasheet was generated on January 28, 2026 at 06:37:47 PM. Visit product page to check for updated information before use.

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