## 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**

Call us: 800-304-5357

## **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 November 28, 2025 at 05:53:06 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biotin-xx-free-acid-6-6-biotinoylaminohexanoy