

Biotin-x cadaverine, free base

Biotin-X cadaverine (5-(((N-(Biotinoyl)amino)hexanoyl)amino)pentylamine, free base) can be coupled directly to activated carboxylic acid esters without the use of a base.

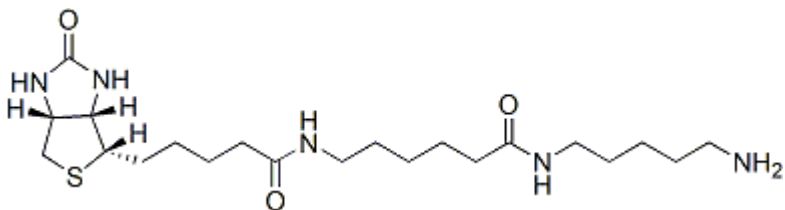


Product attributes

Product Description

Biotin-X cadaverine (5-(((N-(Biotinoyl)amino)hexanoyl)amino)pentylamine, free base) can be coupled directly to activated carboxylic acid esters without the use of a base.

- White solid soluble in DMSO
- Store at 4 °C
- C₂₁H₃₉N₅O₃S
- MW: 441.63



This datasheet was generated on June 3, 2026 at 09:29:55 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-x-cadaverine-free-base-5-n-biotinoylaminohexanoylamino-pentylamine-free-base/>