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

Call us: 800-304-5357 Email: btinfo@biotium.com

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 November 6, 2025 at 09:18:14 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biotin-x-cadaverine-free-base-5-n-biotinoylamino-pentylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotinoylamino-pentylamine-free-base-5-n-biotino-free-ba