



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Biotin cadaverine, free base

Biotin cadaverine free base (*N*-(5-Aminopentyl)biotinamide) can be directly coupled to activated carboxylic acid groups without using an additional base.

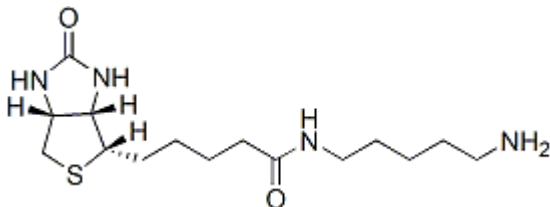


Product attributes

Product Description

Biotin cadaverine free base (*N*-(5-Aminopentyl)biotinamide) can be directly coupled to activated carboxylic acid groups without using an additional base. Also see biotin-X cadaverine free base ([90080](#)).

- Tan solid soluble in DMSO
- MW: 328.47



This datasheet was generated on June 22, 2026 at 06:23:57 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-cadaverine-free-base-n-5-aminopentylbiotinamide/>