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

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 November 3, 2025 at 11:11:22 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-cadaverine-free-base-n-5-aminopentylbiotinamide/