

## Biotin Cadaverine, Trifluoroacetate Salt

Biotin cadaverine, biotin-X cadaverine ([90061](#)), and biotin ethylenediamine ([90057](#)) can be coupled to carboxylic acids, DNA and other biomolecules.

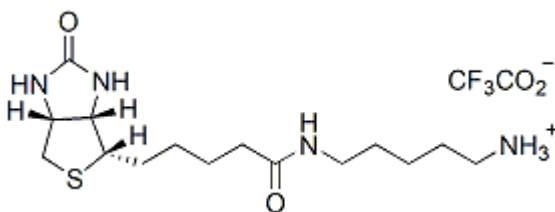


### Product attributes

### Product Description

Biotin cadaverine, biotin-X cadaverine ([90061](#)), and biotin ethylenediamine ([90057](#)) can be coupled to carboxylic acids, DNA and other biomolecules.

- White solid soluble in DMSO
- Store at -20°C or 4°C
- C<sub>17</sub>H<sub>29</sub>F<sub>3</sub>N<sub>4</sub>O<sub>4</sub>S
- MW: 443



### References

1. J Histochem Cytochem 38, 377 (1990).
2. Ann NY Acad Sci 463, 214 (1984).
3. Biochem J 251,935 (1988).

This datasheet was generated on June 3, 2026 at 08:49:14 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-cadaverine-trifluoroacetate-salt-n-5-aminopentylbiotinamide-trifluoroacetic-acid-salt/>