

## Biotin-X cadaverine, trifluoroacetate salt

Biotin-X cadaverine trifluoroacetic acid salt (5-(((N-(biotinoyl)amino)hexanoyl)amino) pentylamine, trifluoroacetic acid salt) can be coupled to activated carboxylic acid esters or sulfonyl chlorides in the presence of a base or in a basic buffer. Also see biotin-X cadaverine free base ([90080](#)).

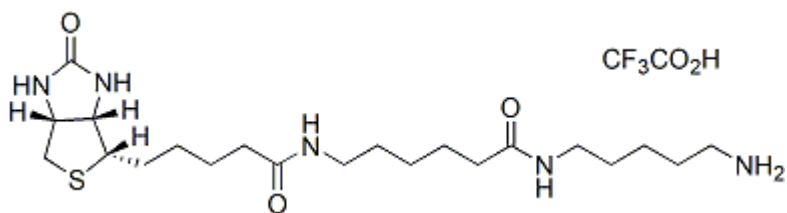


### Product attributes

## Product Description

Biotin-X cadaverine trifluoroacetic acid salt (5-(((N-(biotinoyl)amino)hexanoyl)amino) pentylamine, trifluoroacetic acid salt) can be coupled to activated carboxylic acid esters or sulfonyl chlorides in the presence of a base or in a basic buffer. Also see biotin-X cadaverine free base ([90080](#)).

- Tan solid soluble in DMSO
- Store at -20 °C or 4 °C
- C<sub>23</sub>H<sub>40</sub>F<sub>3</sub>N<sub>5</sub>O<sub>5</sub>S
- MW: 555.65



## References

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

|

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 January 18, 2026 at 07:43:26 PM. Visit product page to check for updated information before use.

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