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

Call us: 800-304-5357

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
- C₂₃H₄₀F₃N₅O₅S
- 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).

I.

- 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 November 28, 2025 at 05:53:10 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-pentylamine-trifluoroacetic-acid-salt/