

ROX Cadaverine, Trifluoroacetate Salt

ROX cadaverine can be coupled to activated carboxy groups or sulfonyl chlorides.



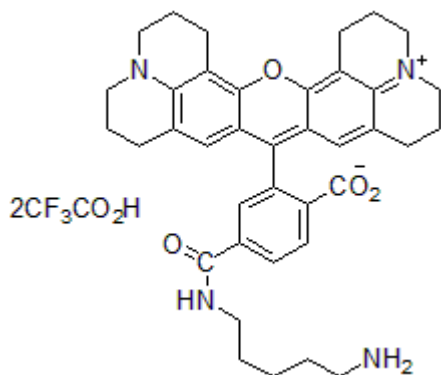
Product attributes

Excitation/Emission	568/595 nm
Isomer	5-ROX Cadaverine, 5(6)-ROX Cadaverine, 6-ROX Cadaverine

Product Description

5-ROX Cadaverine (carboxy-x-rhodamine, cadaverine), 6-ROX Cadaverine, and 5(6)-ROX Cadaverine mixed isomers can be coupled to activated carboxy groups or sulfonyl chlorides.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 568/595 nm
- Dark purple solid soluble in DMSO or DMF
- Store at -20°C and protect from light
- $C_{42}H_{44}F_6N_4O_8$
- MW: 847



Product	Catalog Number	Unit Size
5-ROX Cadaverine, Trifluoroacetate Salt	91060	5 mg
6-ROX Cadaverine, Trifluoroacetate Salt	90121	5 mg
5(6)-ROX Cadaverine, Trifluoroacetate Salt	90108	10 mg

This datasheet was generated on June 3, 2026 at 10:26:54 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/rox-cadaverine-trifluoroacetate-salt/>