Revised: February 10, 2022

Product Information

ROX Cadaverine, Trifluoroacetate Salt

Product List

Catalog no.	Product	Unit size
91060	5-ROX Cadaverine, Trifluoroacetate Salt	5 mg
90121	6-ROX Cadaverine, Trifluoroacetate Salt	5 mg
90108	5(6)-ROX Cadaverine, Mixed Isomer	10 mg

Storage and Handling

Store at -20°C desiccated and protected from light. Product is stable for at least 12 months from date of receipt when stored as recommended.

Stock solutions may be prepared in DMSO or DMF. Solutions can be stored protected from light at 4°C, for up to 12 months.

Molecular Information: $C_{42}H_{44}F_6N_4O_8$

Molecular Weight: 847 Color and Form: Dark puple solid Solubility: Soluble in DMSO or DMF Absorption/Emission: 568/595 nm

Figure 1. 5(6)-ROX cadaverine (mixed isomer), trifluoroacetate salt

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.

Related Products

Catalog number	Product
90082	DMSO, Anhydrous, 10 mL
92001	TAMRA Cadaverine (Mixed Isomers)
91055	ROX Maleimide, 5 and 6 Isomers
90098	ROX, SE, 6 Isomer in DMSO
90104	Aminooxy-5(6)-ROX
91037	MTS-ROX, Mixed Isomers
90116	4-Sulforhodamine 101 Cadaverine, Trifluoroacetate Salt
92035-92043, 92065, 92102, 96010	CF® Dye Amine
92050-92059	CF® Dye Aminooxy
59002	EDC (EDAC), 100 mg

Please visit our website at www.biotium.com for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF® Dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.