

Revised: September 10, 2020

Product Information

TAMRA Cadaverine

Product List

Catalog no.	Product	Unit size	MW	
90122	6-TAMRA Cadaverine, trifluoroacetate	5 mg	743	
92001	5(6)-TAMRA Cadaverine, mixed isomer	10 mg	514	

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 aliquoted and stored with desiccant and protected from light at 4° C, for up to 12 months.

Molecular Information: $C_{30}H_{34}N_4O_4$ Color and Form: Dark pupile solid Solubility: Soluble in DMSO or DMF Absorption/Emission: 544/571 nm

$$\begin{array}{c|c} (H_3C)_2N & O & \stackrel{+}{N}(CH_3)_2 \\ \hline \\ H_2N & O & \\ \end{array}$$

Figure 1. 5(6)-tetramethylrhodamine (TAMRA) cadaverine (mixed isomer)

Product Description

Tetramethylrhodamine (TAMRA) cadaverine (5-(and-6-)-((N-(5-Aminopentyl)amino) carbonyl)tetramethylrhodamine) can be coupled to activated carboxy groups or sulfonyl chlorides.

Related Products

Catalog number	Product
90082	DMSO, Anhydrous, 10 mL
90108	5(6)- ROX Cadaverine
90018	TRITC, 5 isomer
90019	TRITC, 6 isomer
90020	FITC (Isomer I)
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.