

## TAMRA cadaverine (mixed isomers)

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



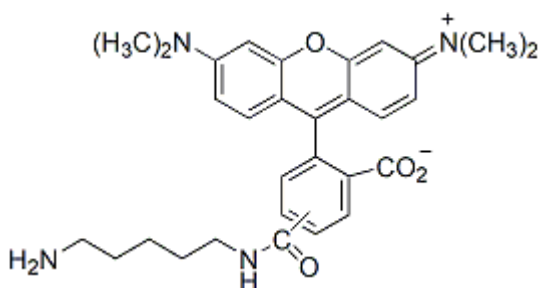
### Product attributes

Excitation/Emission	544/571 nm
---------------------	------------

## Product Description

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

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 544/571 nm
- Dark purple solid soluble in DMSO or DMF
- Store at -20°C and protect from light
- $C_{30}H_{34}N_4O_4$
- MW: 514.62



This datasheet was generated on June 22, 2026 at 11:17:24 AM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/tetramethylrhodamine-cadaverine-mixed-isomers-5-and-6-n-5-aminopentylaminocarbonyltetramethylrhodamine/>