

6-TAMRA Cadaverine, Trifluoroacetate Salt

6-Tetramethylrhodamine (TAMRA) cadaverine can be coupled to activated carboxy groups or sulfonyl chlorides.



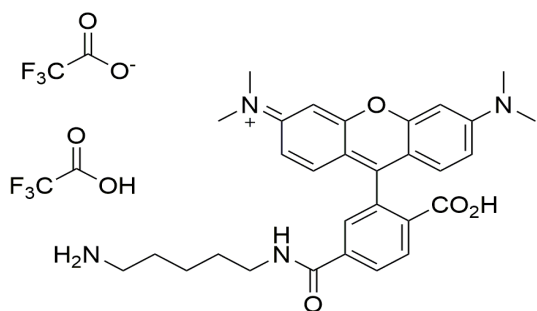
Product attributes

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

Product Description

6-Tetramethylrhodamine (TAMRA) cadaverine (full name: 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: 743



This datasheet was generated on June 3, 2026 at 10:32:16 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/6-tamra-cadaverine-trifluoroacetate-salt/>