6-TAMRA Cadaverine, Trifluoroacetate Salt

stictum

Product attributes

Excitation/Emission 544/571 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

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

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
- \bullet C₃₀H₃₄N₄O₄
- MW: 743

This datasheet was generated on November 2, 2025 at 06:18:54 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/6-tamra-cadaverine-trifluoroacetate-salt/