

TAMRA-PEO3-amine, TFA salt, 5 isomer

5-TAMRA-PEO3-amine (Tetramethylrhodamine-PEO3-amine) has a long water-soluble PEG spacer separating the dye and the amine functional group. The dye can be coupled to activated carboxy groups or sulfonyl chlorides.



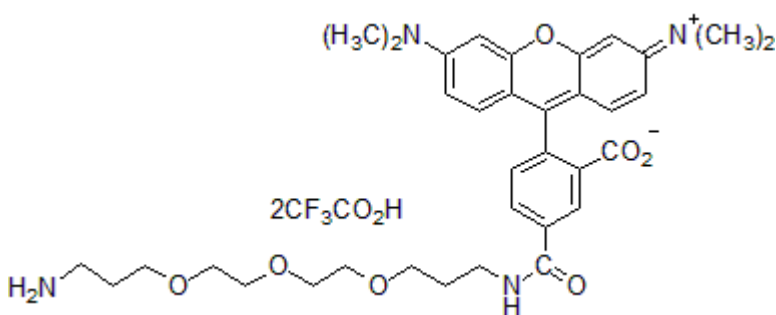
Product attributes

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

Product Description

5-TAMRA-PEO3-amine (Tetramethylrhodamine-PEO3-amine) has a long water-soluble PEG spacer separating the dye and the amine functional group. The dye can be coupled to activated carboxy groups or sulfonyl chlorides.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 544/571 nm
- Dark purple solid soluble in DMSO, DMF, or H₂O
- Store desiccated at 4 °C and protect from light
- C₃₉H₄₆F₆N₄O₁₁
- MW: 860.74



This datasheet was generated on June 22, 2026 at 03:31:18 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/5-tamra-peo3-amine-tfa-salt-tetramethylrhodamine-peo3-amine/>