

# Product Information

## 5-TAMRA-PEO3-Amine, TFA salt

**Catalog Number:** 90107

**Unit Size:** 5 mg

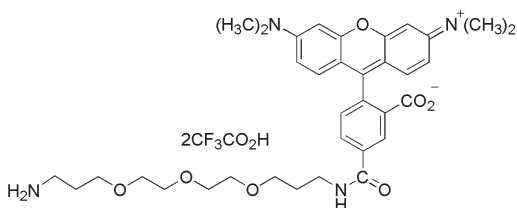
### Storage and Handling

Store at 4°C, desiccated and protected from light. Product is stable for at least 12 months from date of receipt when stored as recommended.

### Molecular Information

$C_{39}H_{46}F_6N_4O_{11}$

MW: 860.74



### Properties

**Color & Form:** Dark purple solid

**Solubility:** Soluble in DMSO, DMF, or water

**Abs/Em maxima:** 544/571 nm (see Fig. 1)

### Product Description

5-TAMRA-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.

Please visit our website at [www.biotium.com](http://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.

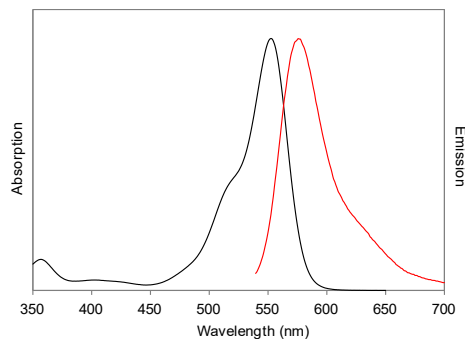


Figure 1. Absorbance and emission spectra of TAMRA.