

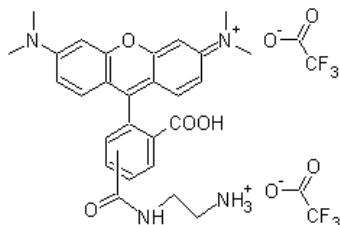
Product Information Sheet

Ordering Information

Product Number: 354
Product Name: 5(6)-TAMRA ethylenediamine
Unit Size: 25 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight: 700.58
Appearance: Purple solid
Chemical Structure:



Soluble in: DMSO or DMF
Excitation Wavelength: 544 nm
Emission Wavelength: 575 nm

Application Notes

5(6)-TAMRA ethylenediamine is a building block for developing TAMRA bioconjugates. TAMRA is one of the most popular bright orange fluorophores used in various bioconjugations. 5(6)-TAMRA is the mixture of two carboxy tetramethylrhodamine (TMR) isomers.