

## Product Information Sheet

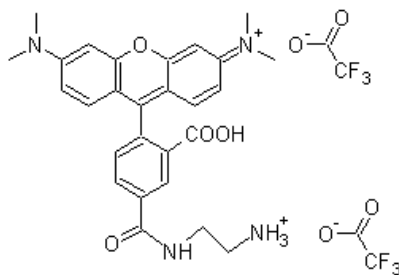
### Ordering Information

Product Number: 358  
Product Name: 5-TAMRA ethylenediamine  
Unit Size: 5 mg  
Storage Conditions: <-15 °C and kept from light and moisture  
Expiration Date: 12 months upon receiving

### Chemical and Spectral Properties

Molecular Weight: 700.58

Chemical Structure:



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

### Application Notes

5-TAMRA ethylenediamine is a building block for developing TAMRA bioconjugates. TAMRA is one of the most popular bright orange fluorophores used in various bioconjugations.