

Product Information Sheet

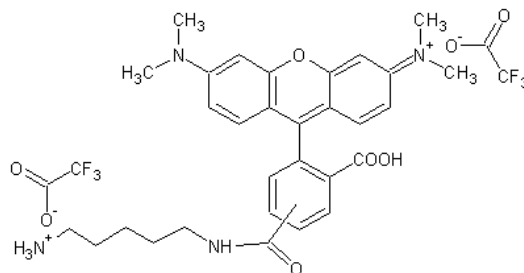
Ordering Information

Product Number: 355
Product Name: 5(6)-TAMRA cadaverine
Unit Size: 25 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 742.66

Chemical Structure:



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

Application Notes

5(6)-TAMRA cadaverine is a building block for developing fluoresceinated bioconjugates. It is also a fluorescent transglutaminase substrate to label proteins by transamination. TAMRA is one of the most popular fluorophores used in various bioconjugations. TMR is a bright orange fluorophore. 5(6)-TAMRA is the mixture of two carboxy tetramethylrhodamine (TMR) isomers.