

Product Information Sheet

Ordering Information

Product Number: 487

Product Name: 5-TAMRA alkyne

Unit Size: 5 mg

Storage Conditions: Freeze (<-15 °C), Desiccated, Avoid Light

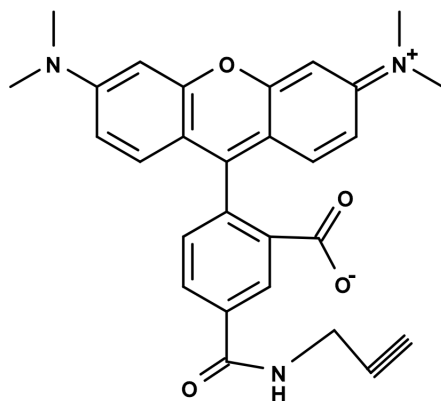
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Appearance: Solid

Molecular Weight: 467.52

Chemical Structure:



Soluble In: DMSO

Excitation Wavelength: 547

Emission Wavelength: 573

Application Notes

5-TAMRA alkyne is an excellent azide-reactive fluorescent reagent for labeling alkyne-containing biological molecules through the well known click chemistry. It has the spectral properties identical to those of TRITC conjugates, thus all the instrument settings of the TRITC conjugates would be applicable to the applications of 5-TAMRA alkyne conjugates.