

6-TAMRA, SE [6-Carboxytetramethylrhodamine, succinimidyl ester] *CAS#: 150810-69-8*

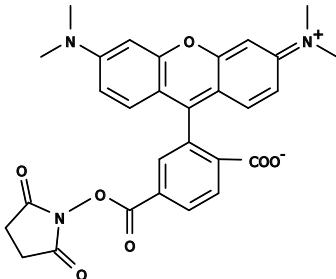
Catalog Number: 376, 377, 378

Unit Size: 5 mg, 100 mg, 1 g

Product Details

| | |
|--------------------|--------------------------------------------|
| Storage Conditions | Freeze (< -15 °C), Minimize light exposure |
| Expiration Date | 12 months upon receiving |

Chemical Properties

| | |
|--------------------|------------------------------------------------------------------------------------|
| Appearance | Solid red |
| Molecular Weight | 527.53 |
| Soluble In | DMSO |
| Chemical Structure |  |

Spectral Properties

| | |
|-----------------------|--------|
| Excitation Wavelength | 552 nm |
| Emission Wavelength | 578 nm |

Applications

The single TAMRA isomers are increasingly preferred for labeling peptides and nucleotides because they give better resolution in HPLC purification that is often required in the conjugation processes. 6-TAMRA, SE is predominantly used for labeling nucleotides and nucleic acids.