

5(6)-TAMRA, SE [5-(and-6)-Carboxytetramethylrhodamine, succinimidyl ester] *CAS 246256-50-8*

Catalog number: 370, 371, 372 Unit size: 25 mg, 100 mg, 1 g

Product Details

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Red purple solid

Molecular Weight 527.53

Soluble In DMSO

Chemical Structure

Spectral Properties

Excitation Wavelength 552 nm

Emission Wavelength 578 nm

Applications

The succinimidyl esters of 5-TAMRA, 6-TAMRA or the mixed isomers are the primary labeling reagents for the preparation of orange fluorescent bioconjugates, including peptide, protein, nucleotide and nucleic acid conjugates, especially fluorescent antibodies and avidin derivatives used in immunochemistry. TMR dyes have also been widely used as acceptors for FAM fluorophores in a variety of FRET studies.