

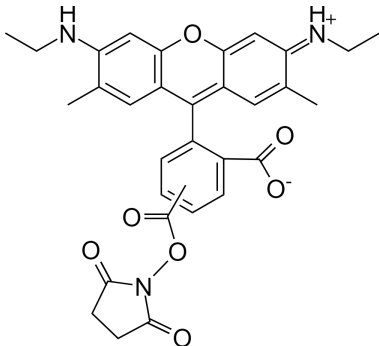
5(6)-CR6G, SE [5-(and 6)-Carboxyrhodamine 6G, succinimidyl ester] *Mixed isomers*

Catalog number: 340
Unit size: 10 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	555.59
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	522 nm
Emission Wavelength	546 nm

Applications

The absorption maxima of CR6G conjugates match well with the 514 nm spectral line of the argon-ion laser that is used in most of fluorescence instruments. In addition, the excitation and emission peaks of rhodamine 6G conjugates falling between those of fluorescein and tetramethylrhodamine provide another color choice for the multicolor imaging applications. 5(6)-CR6G, SE is the amine-reactive form of 5(6)-CR6G. Although 5(6)-CR6G, SE has excellent fluorescence properties, it has very poor solubility in buffers, DMSO or DMF, limiting its utility. iFluor™ 514 has the essentially same spectral properties to 5(6)-CR6G, making iFluor™ 514 an excellent replacement for 5(6)-CR6G.