

5-Carboxyrhodamine 6G maleimide

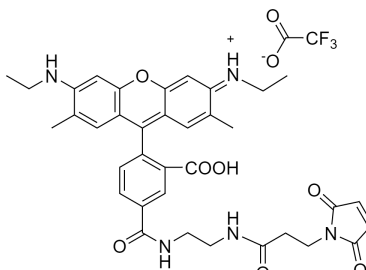
Catalog number: 353

Unit size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	765.74
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	525 nm
Emission Wavelength	548 nm

Applications

5-CR6G maleimide is a thiol-reactive derivative of rhodamine 6G. It is preferred choice (than the mixed isomers) for some complicated biological applications where reproducibility is more critical than material cost since the minor positional difference between 5-CR6G and 6-CR6G might significantly affect some biological properties of the underlying conjugates. Rhodamine 6G is one of the brightest rhodamine dyes, but most of the rhodamine 6G dyes have much lower solubility than other rhodamine dyes. For similar or equivalent spectral properties, iFluor dyes have much better water solubilities.