

6-TET Acid

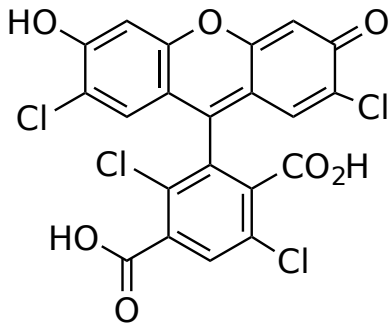
Catalog Number: 210

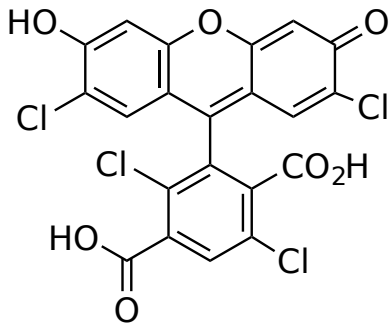
Unit Size: 10 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	514.09
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	521 nm
Emission Wavelength	542 nm

Applications

6-TET is a fluorophore that is widely used in nucleic acid sequencing and related research. 6-TET acid is the most economic form to make a conjugate with amino-containing biomolecules.