

6-JOE acid

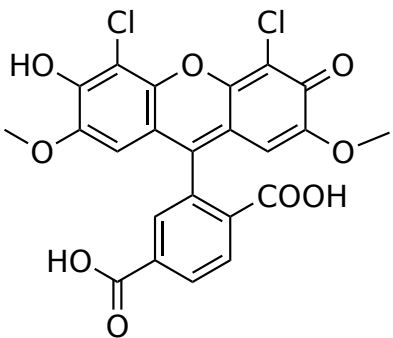
Catalog Number: 204

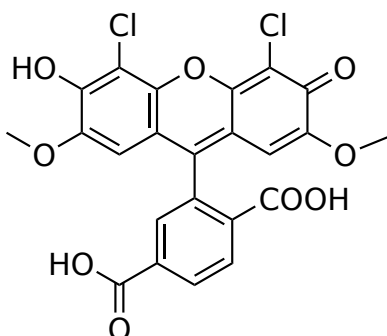
Unit Size: 10 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	505.26
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	520 nm
Emission Wavelength	545 nm

Applications

6-JOE has similar absorption and emission spectra to those of rhodamine 6G, and its SE is predominantly used in automated DNA sequencing. It is also used to label nucleotides. 6-JOE acid is the most economic form to make a conjugate with amino-containing biomolecules.