

6-HEX acid

Catalog Number: 209

Unit Size: 10 mg

Product Details

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

Expiration Date 12 months upon receiving

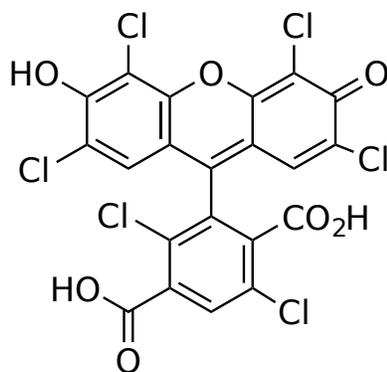
Chemical Properties

Appearance Solid red

Molecular Weight 582.97

Soluble In DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength 533 nm

Emission Wavelength 559 nm

Applications

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