

5-(and-6)-Carboxy SNARF-4F Maleimide

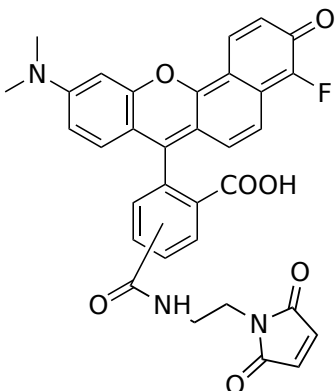
Catalog Number: 21172

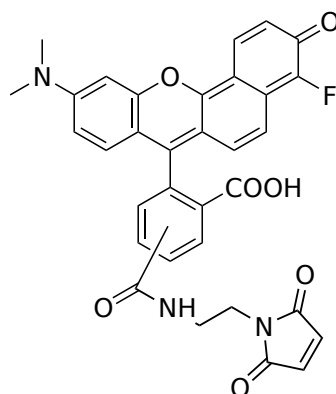
Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	593.57
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

5-(and-6)-Carboxy SNARF-4F Maleimide is an excellent thiol-reactive building block that can be readily used for developing SNARF-4F pH probes. It can be readily conjugated to peptides, antibodies and other thiol-containing biomolecules such as thiol-modified oligos. SNARF-4F has pKa of ~6.4. It exhibits a significant pH-dependent emission shift from yellow orange to deep red fluorescence under acidic and basic conditions, respectively. The ratioable pH dependent fluorescence intensity change allows the ratio of the fluorescence intensities from the dye at two emission wavelengths - typically 580 nm and 640 nm - to be used for quantitative determinations of pH in cells and other biological media.