

5-(and-6)-Carboxy SNARF-5F AM Ester

Catalog Number: 21248

Unit Size: 10x50 ug

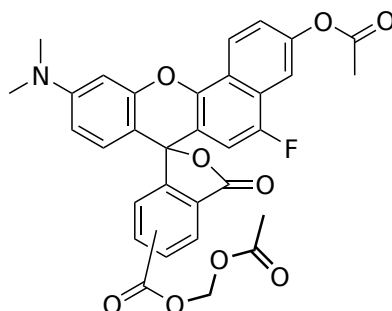
Product Details

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

Chemical Properties

| | |
|------------------|-----------------|
| Appearance | Solid off white |
| Molecular Weight | 585.54 |
| Soluble In | DMSO |

Chemical Structure



Spectral Properties

| | |
|-----------------------|-----|
| Excitation Wavelength | N/A |
| Emission Wavelength | N/A |

Applications

5-(and-6)-Carboxy SNARF-5F AM Ester is an excellent cell-permeable pH probes for ratiometrically monitoring pH changes in live cells. SNARF-5F has pKa of ~7.2. 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.