

## 5-Carboxy-4'-aminomethylfluorescein

Catalog number: 207

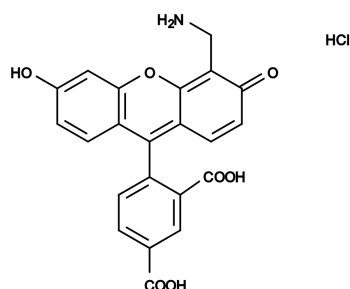
Unit size: 10 mg

### Product Details

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

### Chemical Properties

Appearance	Orange solid
Molecular Weight	441.82
Soluble In	DMSO
Chemical Structure	



### Spectral Properties

Excitation Wavelength	498 nm
Emission Wavelength	525 nm

### Applications

5-Carboxy-4'-aminomethylfluorescein is a key building block for developing Big Dye DNA sequencing terminators. It can also be incorporated into a peptide sequence to give fluoresceinated peptides.