

Product Information Sheet

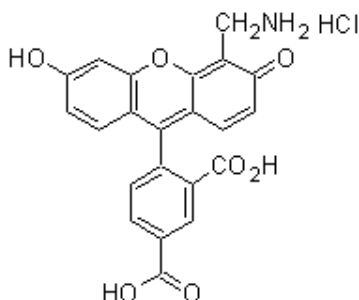
Ordering Information

Product Number: 207
Product Name: 5-Carboxy-4'-aminomethylfluorescein
Unit Size: 10 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 441.82

Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 494 nm
Emission Wavelength: 517 nm

Application Notes

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.