

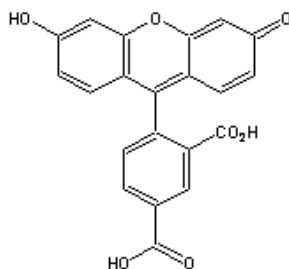
Product Information Sheet

Ordering Information

Product Number:	101
Product Name:	5(6)-FAM [5-(and-6)-Carboxyfluorescein]
Unit Size:	10 g
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	24 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight:	376.32
Appearance:	Orange solid
Chemical Structure:	



Soluble in:	DMSO
Excitation Wavelength:	494 nm
Emission Wavelength:	519 nm

Application Notes

Carboxyfluorescein (commonly called FAM) and its amine-reactive succinimidyl esters are favored over FITC in bioconjugations. FAM reagents give carboxamides that are more resistant to hydrolysis. In addition, FAM reagents require less stringent conjugation conditions and give better conjugation yields, and the resulted conjugates have superior stability. FITC-labeled nucleotides and peptides tend to deteriorate more quickly than the corresponding FAM conjugates. We found that FAM reagents can be used to substitute FITC reagents in most biological applications.