

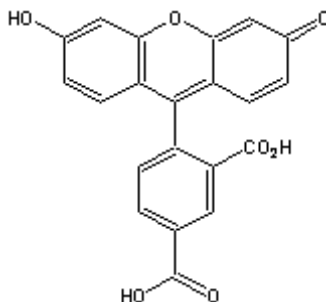
## Product Information Sheet

### Ordering Information

Product Number:	103
Product Name:	5-FAM [5-Carboxyfluorescein] *Single isomer*
Unit Size:	100 mg
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:	492 nm
Emission Wavelength:	518 nm

### Application Notes

5-FAM is the purified single isomer of carboxyfluorescein. It is one of the most popular green fluorescent reagents used for labeling peptides, proteins and nucleotides. It has been predominantly used to develop a variety of green fluorescent peptides that can be excited with the 488 nm line of the Ar laser. It has also been used to prepare various small fluorescent molecules.