

## Product Information Sheet

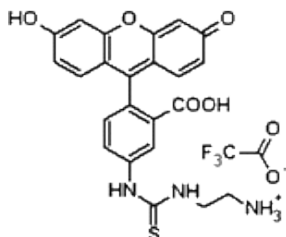
### Ordering Information

Product Number:	126
Product Name:	5-FITC ethylenediamine [Fluorescein thiocarbamylethylenediamine]
Unit Size:	100 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

### Chemical and Spectral Properties

Molecular Weight: 563.5

Chemical Structure:



Soluble in:	DMSO
Excitation Wavelength:	494 nm
Emission Wavelength:	521 nm

### Application Notes

5-FITC ethylenediamine is an excellent building block for developing fluoresceinated bioconjugates. It has been used for labeling peptides and small biomolecules that contain a carboxy group. It can also be used for labeling carbohydrates or glycoproteins via reductive amination.