

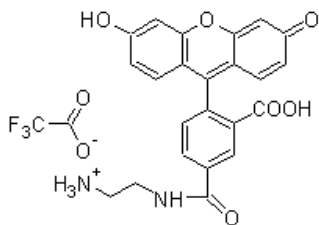
## Product Information Sheet

### Ordering Information

Product Number: 124  
Product Name: 5-FAM ethylenediamine  
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: 532.42  
Appearance: Orange solid  
Chemical Structure:



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

### Application Notes

5-FAM ethylenediamine is an excellent building block for developing fluoresceinated bioconjugates.