

## 5(6)-FAM ethylenediamine

Catalog number: 123 Unit size: 100 mg

**Product Details** 

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

**Chemical Properties** 

Appearance Orange solid

Molecular Weight 532.42

Soluble In DMSO

Chemical Structure

**Spectral Properties** 

Excitation Wavelength 493 nm

Emission Wavelength 517 nm

## **Applications**

5(6)-FAM 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.