

Product Information Sheet

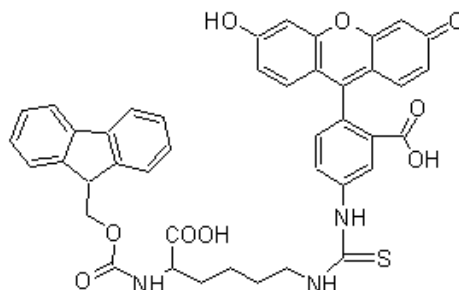
Ordering Information

Product Number: 5070
Product Name: FMOC-Lys(5-FITC)-OH
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: 757.81

Chemical Structure:



Soluble in: DMF
Excitation Wavelength: 492 nm
Emission Wavelength: 515 nm

Application Notes

FMOC-Lys(5-FITC)-OH is a building block for in-sequence Lys labeling by FITC. FITC is one of the most common donor dyes for preparing FRET peptides (often paired with TAMRA or TQ2).