

FMOC-Lys(5-FITC)-OH

Catalog number: 5070 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 757.81

Soluble In DMF

Chemical Structure

Spectral Properties

Excitation Wavelength 491 nm

Emission Wavelength 516 nm

Applications

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).