

FMOC-Lys(Cy3)-OH

PRODUCT INFORMATION SHEET

Catalog number: 5054 Unit size: 100 mg

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Red solid
Molecular Weight	1095.21
Soluble In	DMF
Chemical Structure	$ \begin{array}{c} HO \\ O^{*}S^{*}O \\ \downarrow \\ $

Spectral Properties			
Excitation Wavelength	555 nm		
Emission Wavelength	569 nm		

Applications

FMOC-Lys(Cy3)-OH is a building block for in-sequence Lys labeling by Cy3. Cy3 is one of the most common fluorescent dyes for preparing FRET peptides (often paired with Cy5 or TQ3).