

FMOC-Glu(5-FAM)-OH

Catalog number: 5013 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 769.75

Soluble In DMF

Chemical Structure

Spectral Properties

Excitation Wavelength 493 nm

Emission Wavelength 517 nm

Applications

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