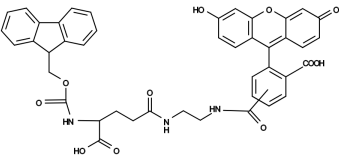


FMOC-Glu(5/6-FAM)-OHCatalog number: 5012
Unit size: 100 mg**Product Details**

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	769.75
Soluble In	DMF
Chemical Structure	

Spectral Properties

Excitation Wavelength	493 nm
Emission Wavelength	517 nm

Applications

FMOC-Glu(5/6-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).