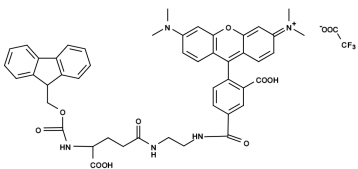


Fmoc-Glu(5-TAMRA)-OHCatalog number: 5015
Unit size: 100 mg**Product Details**

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

Chemical Properties

Appearance	Red solid
Molecular Weight	937.91
Soluble In	DMF
Chemical Structure	

Spectral Properties

Excitation Wavelength	552 nm
Emission Wavelength	578 nm

Applications

Fmoc-Glu(5-TAMRA)-OH is a building block for in-sequence peptide Glu labeling by TAMRA. TAMRA is one of the most common donor dyes for preparing FRET peptides (often paired with FAM or TQ3).