

Fmoc-Glu(TQ3)-OH

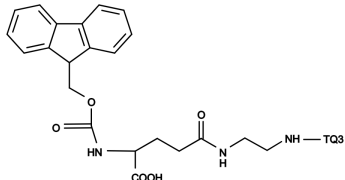
Catalog number: 5238

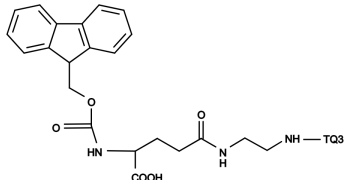
Unit size: 100 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	846.95
Soluble In	DMF
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	573 nm
Emission Wavelength	nm

Applications

Fmoc-Glu(TQ3)-OH is a building block for in-sequence Lys labeling by Tide Quencher™ 3 (TQ3). TQ3 is the best quencher that is used to develop a FRET probe with TAMRA and Cy3 or other fluorescent dyes that have the similar fluorescence spectra (to TAMRA and Cy3) such as Bodipy TMR dyes. It can also be used with 6-HEX derivatives.