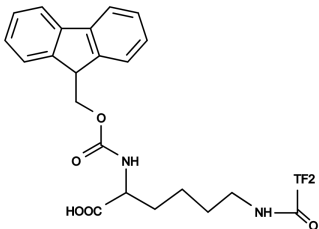
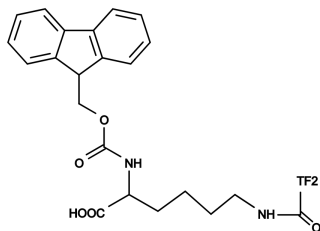


FMOC-Lys(TF2-Boc)-OHCatalog number: 5050
Unit size: 100 mg**Product Details**

Storage Conditions	Freeze (<math>< -15\text{ }^{\circ}\text{C}</math>), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	924.99
Soluble In	DMF
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	503 nm
Emission Wavelength	525 nm

Applications

FMOC-Lys(TF2-Boc)-OH is a building block for in-sequence Lys labeling by TF2, an excellent and more affordable replacement for the least affordable Alexa Fluor 488. TF2 is one of the most common donor dyes for preparing FRET peptides (often paired with TAMRA or TQ2) since TF2 and Alexa Fluor 488 share the same R110 fluorophore.