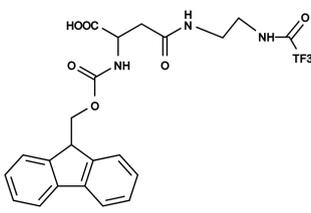
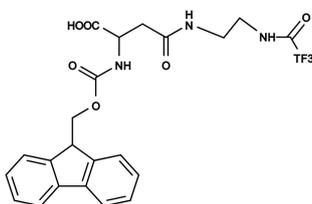


FMOC-Asp(TF3)-OHCatalog number: 5007
Unit size: 100 mg**Product Details**

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

Chemical Properties

Appearance	Solid
Molecular Weight	951.94
Soluble In	DMF
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	546 nm
Emission Wavelength	571 nm

Applications

FMOC-Asp(TF3)-OH is a building block for in-sequence Asp labeling by TF3. TF3 has the spectral properties identical to Cy3. Compared to Cy3, TF3 is more stable to acidic and basic deprotection conditions. Our TF3 is also more fluorescent than Cy3 on peptides and oligonucleotides.