

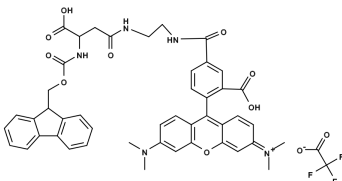
**FMOC-Asp(5-TAMRA)-OH**Catalog number: 5006  
Unit size: 100 mg**Product Details**

---

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	923.89
Soluble In	DMF
Chemical Structure	

**Spectral Properties**

---

Excitation Wavelength	552 nm
Emission Wavelength	578 nm

**Applications**

---

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