

## FMOC-Asp(5-FAM)-OH

Catalog Number: 5004

Unit Size: 100 mg

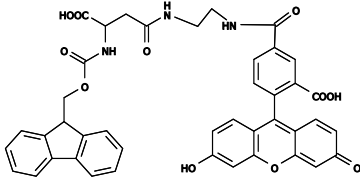
### Product Details

---

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

### Chemical Properties

---

Appearance	Solid
Molecular Weight	755.72
Soluble In	DMF
Chemical Structure	

### Spectral Properties

---

Excitation Wavelength	493 nm
Emission Wavelength	517 nm

### Applications

---

FMOC-Asp(5-FAM)-OH is a building block for in-sequence peptide Asp labeling by FAM. FAM is one of the most common donor dyes for preparing FRET peptides (often paired with TAMRA or TQ2).