

FMOC-Asp(5/6-FAM)-OH

Catalog Number: 5003

Unit Size: 100 mg

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

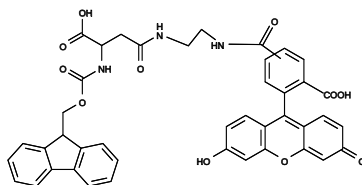
Chemical Properties

Appearance Solid orange

Molecular Weight 755.72

Soluble In DMF

Chemical Structure



Spectral Properties

Excitation Wavelength 493 nm

Emission Wavelength 517 nm

Applications

FMOC-Asp(5/6-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).