

FMOC-Asp(TQ2)-OH

Catalog number: 5206 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 761.89

Soluble In DMF

Chemical Structure

Spectral Properties

Excitation Wavelength 516 nm

Emission Wavelength nm

Applications

FMOC-Asp(TQ2)-OH is a building block for in-sequence Lys labeling by Tide Quencher™ 2 (TQ2). TQ2 is the best quencher that is used to develop a FRET probe with fluorescein or other fluorescent dyes that have the similar fluorescence spectra such as Bodipy FL dyes. It can also be used with rhodamine 6G, 6-TET and 6-JOE derivatives.