

## FMOC-Asp(DABCYL)-OH

**PRODUCT INFORMATION SHEET** 

Catalog number: 5204 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	648.71
Soluble In	DMF
Chemical Structure	
Creatural Dranautica	

Spectral	Properties
----------	------------

Excitation Wavelength	454 nm
Emission Wavelength	nm

## Applications

A building block for in-sequence Asp labeling by DABCYL.