

Product Information Sheet

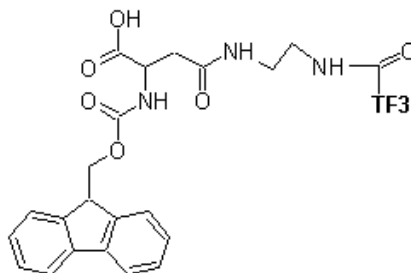
Ordering Information

Product Number: 5007
Product Name: FMOC-Asp(TF3)-OH
Unit Size: 100 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 951.94

Chemical Structure:



Soluble in: DMF
Excitation Wavelength: 554 nm
Emission Wavelength: 583 nm

Application Notes

FMOC-Asp(TF3)-OH is a building block for in-sequence Asp labeling by TF3. TF3 has the spectral properties identical to Cy3®. Compared to Cy3®, TF3 is more stable to acidic and basic deprotection conditions. Our TF3 is also more fluorescent than Cy3® on peptides and oligonucleotides.