

www.aatbio.com; E-Mail: info@aatbio.com

Product Information Sheet

Ordering Information

Product Number: 5000

Product Name: FMOC-Asp(EDANS)-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: 603.64

Chemical Structure:

Soluble in: DMF
Excitation Wavelength: 335 nm
Emission Wavelength: 493 nm

Application Notes

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