



Product Information Sheet

Ordering Information

Product Number: 5310

Product Name: BOC-Asp(OMe)-FMK [Boc-D(OMe)-FMK]

Unit Size: 1 mg

Storage Conditions: Freeze (<-15 °C), Desiccated, Avoid Light

Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Appearance: Solid

Molecular Weight: 263.27

Soluble In: DMF

Excitation Wavelength: N/A

Emission Wavelength: N/A

Application Notes

The FMK peptides are potent, cell permeable, irreversible capase inhibitors. It is useful for detecting and monitoring apoptosis in vitro and in vivo applications. BOC-ASP(OMe)-FMK is the key intermediate for developing and preparing FMK-based capase inhibitors and probes. The Asp(OMe)-FMK-based caspase inhibitors are cell permeable and irreversibly binds to activated caspases to block apoptosis. Caspases, which act as molecular chainsaw are the central components of apoptosis response. Caspases are cysteine proteases that cleave after aspartate residue in their substrates. At least 7 of 14 known mammalian caspases have important role in apoptosis.