

Z-DEVD-FMKCatalog number: 13303
Unit size: 1 mg**Product Details**

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Soluble In	DMSO

Applications

Z-DEVD-FMK is a specific, irreversible Caspase-3 inhibitor, and also shows potent inhibition on caspase-6, caspase-7, caspase-8, and caspase-10.