

Product Information

Description	Z-VEID-FMK is the specific recognition sequence for caspase-6/Mch2.
Target:	Apoptosis
Pathway:	Caspase
Chemical Name:	methyl (4S)-5-[[[(2S)-1-[[[(3S)-5-fluoro-1-methoxy-1,4-dioxopentan-3-yl]amino]-3-methyl-1-oxopentan-2-yl]amino]-4-[[[(2S)-3-methyl-2-(phenylmethoxycarbonylamino)butanoyl]amino]-5-oxopentanoate
Molecular Formula:	C ₃₁ H ₄₅ FN ₄ O ₁₀
Supplied As:	Solid
Solubility:	Insoluble in H ₂ O; ≥113.4 mg/mL in DMSO; ≥3.01 mg/mL in EtOH with gentle warming and ultrasonic For obtaining a higher solubility, please warm the tube at 37°C and shake it in the ultrasonic bath for a while.
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	652.71 Da
Purity:	≥98%

Chemical Structure

