

Product Information

Description	Boc-D-FMK is a Novel inhibitor of Caspase-3
Target:	Apoptosis
Pathway:	Caspase
Chemical Name:	methyl 5-fluoro-3-[(2-methylpropan-2-yl)oxycarbonylamino]-4-oxopentanoate
Molecular Formula:	C ₁₁ H ₁₈ FNO ₅
Supplied As:	Solid
Solubility:	insoluble in H ₂ O; ≥11.65 mg/mL in DMSO; ≥41.65 mg/mL in EtOH. 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:	263.26 Da
Cas No.	187389-53-3, 634911-80-1
Purity:	≥98% by HPLC

Chemical Structure

