

Product Information

Description	Caspase-3 inhibitor. Inhibits MG 132-induced small cell lung cancer cell death in vitro.
Target:	Apoptosis
Pathway:	Caspase
Chemical Name:	methyl (3S)-3-[[[(2S)-2-[[[(2S)-5-amino-2-[[[(2S)-4-methoxy-4-oxo-2-(phenylmethoxycarbonylamino)butanoyl]amino]-5-oxopentanoyl]amino]-4-methylsulfanylbutanoyl]amino]-5-fluoro-4-oxopentanoate
Molecular Formula:	C ₂₉ H ₄₀ FN ₅ O ₁₁ S
Supplied As:	Solid
Solubility:	≥29.2 mg/mL in DMSO; insoluble in EtOH; insoluble in H ₂ O. 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:	685.72 Da
Cas No.	767287-99-0
Purity:	≥98%

Chemical Structure

