

## Product Information

<b>Description:</b>	Second generation, irreversible, broad-spectrum caspase inhibitor that is non-toxic, cell permeable, and penetrates the blood-brain barrier.
<b>Target:</b>	Apoptosis
<b>Chemical Name:</b>	(S)-methyl 5-(2,6-difluorophenoxy)-3-((S)-3-methyl-2-(quinoline-2-carboxamido)butanamido)-4-oxopentanoate
<b>Molecular Formula:</b>	C <sub>26</sub> H <sub>25</sub> F <sub>2</sub> N <sub>3</sub> O <sub>6</sub>
<b>Supplied As:</b>	Solid
<b>Solubility:</b>	≥26.35 mg/mL in DMSO; insoluble in H <sub>2</sub> O; ≥97.4 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.
<b>Stability:</b>	At least 6 months at -20°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-20°C
<b>MW:</b>	527 Da
<b>Purity:</b>	≥98% by HPLC

## Chemical Structure

