

Product Information

Description

Z-VDVAD-FMK is a synthetic peptide that irreversibly inhibits the activity of caspase-2. It attenuates oxyhemoglobin-induced cleavage of PARP and apoptosis in endothelial cells.

Target:

Caspase-2

Pathway

Apoptosis

Chemical Name:

methyl (3S)-5-fluoro-3-[[[(2S)-2-[[[(2S)-2-[[[(2S)-4-methoxy-2-[[[(2S)-3-methyl-2-(phenylmethoxycarbonylamino)butanoyl]amino]-4-oxobutanoyl]amino]-3-methylbutanoyl]amino]propanoyl]amino]-4-oxopentanoate

Molecular Formula:

C₃₂H₄₆N₅O₁₁F

Supplied As:

Solid

Solubility:

≥34.8 mg/mL in DMSO; insoluble in EtOH; insoluble in H₂O. For obtaining a higher solubility, please warm the tube at 37°C and shake it in the ultrasonic bath for a while. Long term storage in solution is not recommended. Not stable in diluted solutions.

Stability:

At least 6 months at -20°C. Avoid freeze/thaw cycles.

Storage:

-20°C

MW:

695.73 Da

Cas No.

210344-92-6

Purity:

≥98%

Chemical Structure

