

Instruction Manual

Catalog Number	PK-CA577-1142-5
Description	A synthetic peptide that irreversibly inhibits activity of caspase-2. The inhibitor is designed as a methyl ester to facilitate cell permeability. The product can be used for both in vitro and in vivo studies. (Caution: If the intended use is on purified or recombinant enzymes, esterase should be added to generate free carboxyl groups.)
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	Z-Val-Asp(OMe)-Val-Ala-Asp(OMe)-FMK; 695 Da
Purity	One spot by TLC
Appearance / Formulation / Solubility	Solid; Reconstitute in DMSO (20 mM).
Storage & Stability	Store at -20°C. Stable for 1 year under proper storage conditions.
Applications	We recommend using 2-10 μ M for inhibiting caspase activity in cell culture. The optimal doses may vary for different cells and culture conditions.
References	NA
Caution	NA

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.