

## Instruction Manual

Catalog Number	PK-CA577-1010-20C
Description	A synthetic peptide that irreversibly inhibits activity of caspase family proteases and blocks apoptosis. The inhibitor is designed as a methyl ester to facilitate cell permeability. The product can be used in 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	20 µl (10 mM)
Sequence / Molecular Weight / Molecular Formula	Z-Val-Ala-Asp-FMK; 468 Da
Purity	One spot by TLC analysis
Appearance / Formulation / Solubility	10 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 1 year under proper storage conditions.
Applications	We recommend using 1000-5000X dilutions for inhibiting caspase activity in Jurkat cell culture (e.g., add 1 µl to 1-5 ml of culture medium). 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.

---