

## Instruction Manual

Catalog Number	PK-CA577-1012-100
Description	A synthetic peptide that irreversibly inhibits ICE (interleukin-1 $\beta$ converting enzyme, $K_i = 0.76$ nM) activity. YVAD is the specific recognition sequence for ICE/caspase-1. The inhibitor is designed as a methyl ester to facilitate cell permeability. (Caution: If the intended use is on purified or recombinant enzymes, esterase should be added to generate free carboxyl groups.)
Quantity	100 $\mu$ l (2 mM)
Sequence / Molecular Weight / Molecular Formula	Z-Tyr-Val-Ala-Asp-FMK (FMK, fluoromethyl ketone); 630 Da
Purity	>99% by HPLC analysis
Appearance / Formulation / Solubility	2 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-1 activity in Jurkat cell culture (e.g., add 1 $\mu$ l to 1-5 ml of culture medium). The optimal doses may vary for different cells and culture conditions.
References	Schlegel, J., et al. (1996) J. Biol. Chem. 271:1841-1485. Nicholson, D.W., et al. (1995) Nature 376:37-40. Walker, N.P.C., et al. (1994) Cell 78:343-349. Thornberry, N., et al. (1992) Nature 356:768-771.
Caution	NA

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