

## Instruction Manual

Catalog Number	PK-CA577-1100-20C
Description	A synthetic peptide that irreversibly inhibits caspase-5 and related caspase activity. 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	20 µl (10 mM)
Sequence / Molecular Weight / Molecular Formula	Z-Trp-Glu-His-Asp-FMK (FMK, fluoromethyl ketone); 763.78 Da
Purity	>99% by HPLC 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-5 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	1.Thornberry, N.A., et al. (1997) J. Biol. Chem. 272:17907-17911. 2.Garcia-Calvo, M., et al. (1998) J. Biol. Chem. 273:32608-32613.
Caution	NA

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