

Instruction Manual

Catalog Number	PK-CA577-1079-20C
Description	A synthetic peptide that irreversibly inhibits caspase-12 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-Ala-Thr-Ala-Asp-FMK I(FMK, fluoromethyl ketone); 540 Da
Purity	>95% by HPLC analysis
Appearance / Formulation / Solubility	10 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 6-12 months under proper storage conditions.
Applications	We recommend using 1000-5000X dilutions for inhibiting Caspase-12 activity in tissue and cell lysate (e.g., add 1 µl to 1 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.