

Instruction Manual

Catalog Number	PK-CA577-1079-100
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	100 µl (2 mM)
Sequence / Molecular Weight / Molecular Formula	Z-Ala-Thr-Ala-Asp(OMe)-FMK (FMK, fluoromethyl ketone); 540 Da
Purity	>95% by HPLC analysis
Appearance / Formulation / Solubility	2 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 6-12 months under proper storage conditions.
Applications	We recommend using 1000X dilutions for inhibiting Caspase-12 activity in tissue and cell lysate (add 1 µl to 1 ml of tissue and cell lysate). The optimal doses may vary for different tissue and cell preparations.
References	NA
Caution	NA

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