

Instruction Manual

Catalog Number	PK-CA577-1115-20C
Description	A synthetic peptide inhibitor that inhibits caspase-13 and related caspase activity. FMK inhibitors are irreversible and non-toxic, and can be used in both in vivo and in vitro studies. 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-Leu-Glu-Glu-Asp-FMK (FMK, fluoromethyl ketone); 696 Da
Purity	>95% 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-13 and related caspase activity (e.g., add 1 µl to 1-5 ml of culture medium). The optimal doses may vary among different cells and culture conditions. The inhibitor is designed as a methyl ester to facilitate cell permeability.
References	Humke, E.W., et al. (1998) J. Biol. Chem. 273:15702-15707.
Caution	NA

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