

## Instruction Manual

Catalog Number	PK-CA577-1074-100
Description	A synthetic peptide that irreversibly inhibits caspase-9 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-Leu-Glu-His-Asp-FMK (FMK, fluoromethyl ketone); 690.7 Da
Purity	>98% by HPLC analysis
Appearance / Formulation / Solubility	2 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 6 months under proper storage conditions.
Applications	We recommend using 1000X dilutions for inhibiting Caspase-9 activity in cell culture (e.g., add 1 µl to 1 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.
Caution	NA

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