

Instruction Manual

Catalog Number	PK-CA577-1009-100
Description	A synthetic peptide inhibitor that irreversibly inhibits caspase-3 and related protease/caspase activity and blocks apoptosis. DEVD is the specific recognition sequence found in poly (ADP-ribose) polymerase (PARP). 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-Asp-Glu-Val-Asp-FMK (FMK, fluoromethyl ketone); 668 Da
Purity	>98% by HPLC analysis
Appearance / Formulation / Solubility	2 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 1 year under proper storage conditions.
Applications	We recommend using 1000X dilutions for inhibiting Caspase-3 activity in Jurkat cell culture (e.g., Add 1 µl to 1 ml of culture medium). However, the optimal doses may vary for different cells and culture conditions.
References	Xiang, J., et al. (1996) Proc. Natl. Acad. Sci. USA 93:14559. Nicholson, D. W., et al. (1995) Nature 376:37. Lazebnic, Y. A., et al. (1994) Nature 371:346.
Caution	NA

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