

## Instruction Manual

Catalog Number	PK-CA577-1009-20C
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	20 µl (10 mM)
Sequence / Molecular Weight / Molecular Formula	Z-Asp-Glu-Val-Asp-FMK (Z-DEVD-FMK)(FMK, fluoromethyl ketone); 668 Da
Purity	>98% 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-3 activity in Jurkat cell culture (e.g., add 1 µl to 1-5 ml of culture medium). The inhibitor can be added to culture medium together with apoptosis inducer or 1 hr before inducer is added. 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.