

Instruction Manual

Catalog Number	PK-CA577-1064-100
Description	A synthetic peptide that irreversibly inhibits FLICE and related protease/caspase activity and blocks apoptosis. IETD is the specific recognition sequence found in CPP32 proenzyme. 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-Ile-Glu-Thr-Asp-FMK (FMK, fluoromethyl ketone); 654
Purity	>99% 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-8 activity in Jurkat 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. 2.Nicholson, D.W., et al. (1995) Nature 376:37-40.
Caution	NA

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