

Instruction Manual

Catalog Number	PK-CA577-1123-20C
Description	A synthetic peptide that irreversibly inhibits activity of caspase family proteases and blocks apoptosis. The inhibitor is designed as a methyl ester to enhance cell permeability. The conjugation with biotin allows detection of activated caspases in conjunction with avidin-tags. (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	Biotin-Val-Ala-Asp-FMK (FMK, fluoromethyl ketone); 672.81 Da
Purity	>95% by HPLC analysis
Appearance / Formulation / Solubility	10 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 or binding to activated caspases (e.g., add 1 µl to 1 ml of culture medium). However, the optimal doses should be determined individually.
References	Slee, E.A., et al. (1996) <i>Biochem. J.</i> 315:21-24. Armstrong, et al. (1996) <i>J. Biol. Chem.</i> 271:16850-16857. Qi, X.M., et al. (1997) <i>Oncogene</i> 15:1207-1212.
Caution	NA

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