

Instruction Manual

Catalog Number	PK-CA577-K108
Description	Synthetic peptide inhibitors that irreversibly inhibit the activity of each of the caspase-family proteases. YVAD is the recognition sequence for caspase-1/ICE, VDVAD is the recognition sequence for caspase-2, DEVD is the recognition sequence for caspase-3/CPP32, VEID is the recognition sequence for caspase-6/Mch2, IETD is the recognition sequence for caspase-8/FLICE, LEHD is the recognition sequence for caspase-9, and VAD is the recognition sequence for a caspase(s) that functions upstream of most other caspases.
Quantity	7 x 25 µl
Sequence / Molecular Weight / Molecular Formula	NA
Purity	see data sheets of individual caspase inhibitors
Appearance / Formulation / Solubility	2 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 6-12 months under proper storage conditions.
Applications	We recommend using 1000X dilutions for inhibiting caspase 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	NA
Caution	NA

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