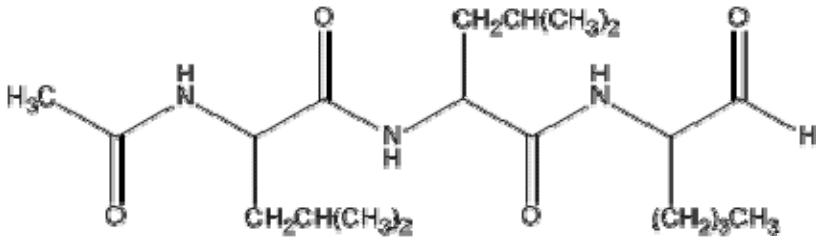


Calpain Inhibitor I

A cell-permeable inhibitor of calpain I, calpain II, cathepsin B, and cathepsin L

Instruction Manual

Catalog Number	PK-CA577-1834-25
Description	A cell-permeable inhibitor of calpain I (K _i = 190 nM), calpain II (K _i = 220 nM), cathepsin B (K _i = 150 nM), and cathepsin L (K _i = 0.5 nM). Inhibits neutral cysteine proteases and the proteasome (K _i = 6 μM). Can initiate apoptosis in HL-60 cells but blocks dexamethasone-induced apoptosis in thymocytes and cycloheximide-induced apoptosis in metamyelocytes. Inhibits cyclin B degradation and arrests the cell cycle at G1/S and at meta-/anaphase.
Quantity	25 mg
Sequence / Molecular Weight / Molecular Formula	383.54 Da; C ₂₀ H ₃₇ N ₃ O ₄
Chemical Structure	 <p>The chemical structure shows a cyclic peptide backbone with four amide bonds. The side chains are: a methyl group (H₃C), an isobutyl group (CH₂CH(CH₃)₂), another isobutyl group (CH₂CH(CH₃)₂), and a tert-butyl group ((CH₃)₃CH₂).</p>
Purity	≥ 95% as determined by HPLC.
Appearance / Formulation / Solubility	White powder; Soluble in DMSO (~25 mg/ml) or EtOH.
Storage & Stability	Store at -20°C.
Applications	see Description
References	NA
Caution	Do not take internally. Wear gloves and mask when handling the product! Avoid contact by all modes of exposure.

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