

Instruction Manual

Catalog Number	PK-CA577-1128-20C
Description	A synthetic peptide that irreversibly inhibits granzyme B activity. The inhibitor is designed as a methyl ester to facilitate cell permeability.
Quantity	20 µl (10 mM)
Sequence / Molecular Weight / Molecular Formula	Z-Ala-Ala-Asp(Ome)-CMK; 455.89 Da
Purity	Single spot by TLC
Appearance / Formulation / Solubility	10 mM in DMSO
Storage & Stability	Store at -20°C. Protect from light and moisture.
Applications	see Description
References	Gong, B., et al. (1999) Cell Growth Diff. 10:491-496. Caputo, et al. (1993) J. Biol. Chem. 268:17672-17675. Shi, L., et al. (1992) J. Exp. Med. 176:1521-1528. Otake, S., et al. (1991) Biochemistry 30:2217-2220.
Caution	NA

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