

Instruction Manual

Catalog Number	PK-CA577-1119
Description	A synthetic peptide that inhibits Granzyme B.
Quantity	1 mg
Sequence / Molecular Weight / Molecular Formula	Ac-LEPD-CHO; 498.53 Da; C ₂₂ H ₃₄ N ₄ O ₉
Purity	>95% by HPLC analysis
Appearance / Formulation / Solubility	Lyophilized solid. Soluble in DMSO.
Storage & Stability	Store at -20°C. Protect from light and moisture. Stable for 1 year as supplied under proper storage conditions.
Applications	NA
References	NA
Caution	NA

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