

Instruction Manual

Catalog Number	PK-CA577-1125-20C
Description	
Quantity	20 µl (10 mM)
Sequence / Molecular Weight / Molecular Formula	Z-Leu-Leu-Tyr-FMK (FMK, fluoromethyl ketone); 557 Da
Purity	>95% by HPLC analysis
Appearance / Formulation / Solubility	10 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 1 year under proper storage conditions.
Applications	The effective final concentrations of the Calpain inhibitor is estimated to be 20-50 µM.
References	NA
Caution	NA

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