

FMK Negative Control (for caspase inhibitors)

Instruction Manual

Catalog Number	PK-CA577-1122-100
Description	The reagent can be used as a negative control for FMK-labeled caspase inhibitors. (Caution: If the intended use is on purified or recombinant enzymes, esterase should be added to generate free carboxyl groups.)
Quantity	100 µl (2 mM)
Sequence / Molecular Weight / Molecular Formula	Z-Phe-Ala-FMK (FMK, fluoromethyl ketone); 386 Da
Purity	>95% by HPLC analysis
Appearance / Formulation / Solubility	2 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 1 year under proper storage conditions.
Applications	The reagent can be used as a negative control for FMK-labeled caspase inhibitors.
References	NA
Caution	NA

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