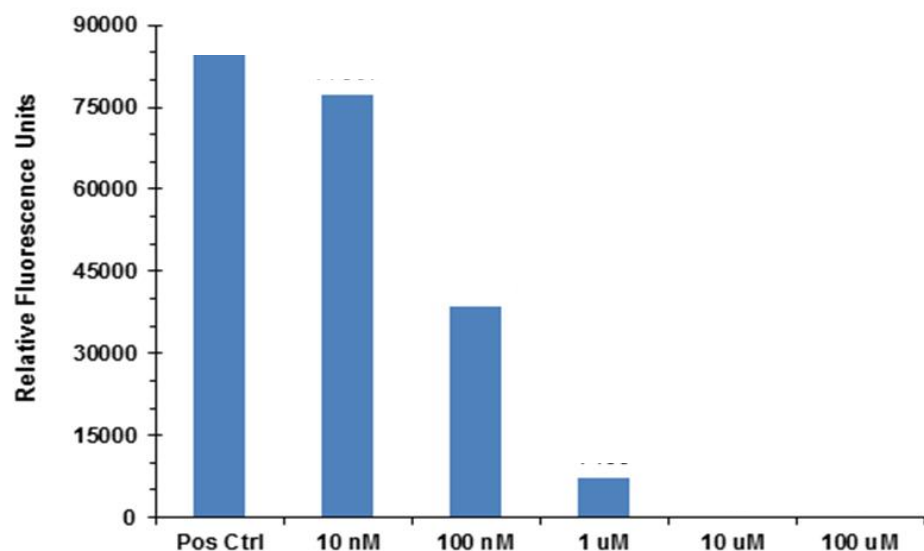


Instruction Manual

Catalog Number	PK-CA577-1176-1
Description	A second-generation caspase-8 inhibitor that is potent, cell permeable, irreversible and nontoxic. (Caution: If the intended use is on purified or recombinant enzymes, esterase should be added to generate free carboxyl groups.)
Quantity	1 mg
Sequence / Molecular Weight / Molecular Formula	785.78 Da; C ₃₈ H ₄₅ F ₂ N ₅ O ₁₁
Chemical Structure	
Purity	NA
Appearance / Formulation / Solubility	Off-white solid; Soluble in DMSO.
Storage & Stability	Store at -20°C. Protect from moisture.
Applications	Inhibition of caspases in <i>in vivo</i> assays.

Q-IETD-OPh Inhibition of Caspase-8 Activity



References	NA
Caution	NA

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