

## Instruction Manual

Catalog Number	PK-CA577-1171-1
Description	A synthetic peptide that is cell permeable, nontoxic and used as a negative control to the caspase inhibitor Q-VD-OPH, the new generation of caspase inhibitor. It is stable in aqueous environment and exhibits several folds less inhibitory activity than the corresponding caspase Q-VD-OPH. It is the perfect choice for both in vitro and in vivo studies.
Quantity	1 mg
Sequence / Molecular Weight / Molecular Formula	C <sub>27</sub> H <sub>27</sub> F <sub>2</sub> N <sub>3</sub> O <sub>6</sub> , 527.19
Purity	>95% by HPLC analysis
Appearance / Formulation / Solubility	Pure thin resin. Soluble in DMSO.
Storage & Stability	Store at -20°C. Stable for 1 year under proper storage conditions.
Applications	We recommend using 1-10 µM concentration for in vitro enzyme kinetics studies. For in vivo usage, quantity needs to be determined by researcher. 20 mg/Kg of body weight is a proper place to start.
References	NA
Caution	NA

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