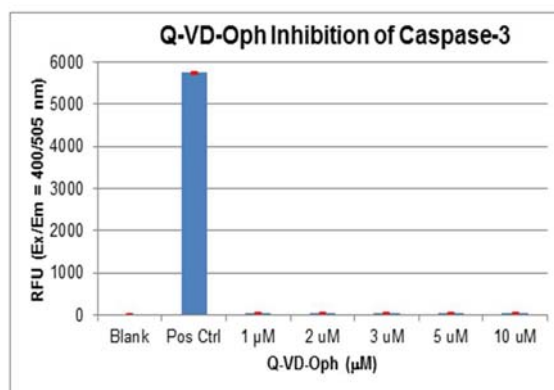
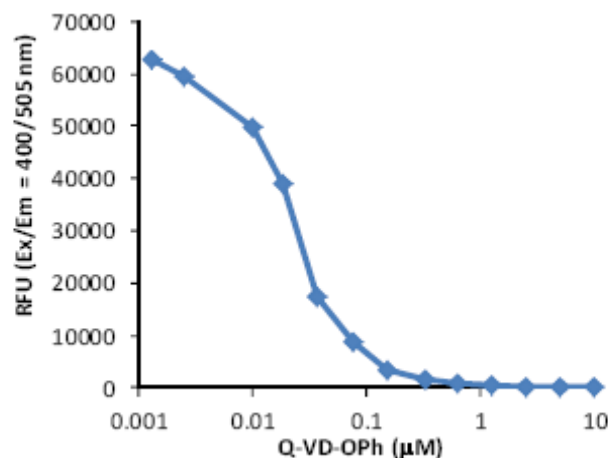


Instruction Manual

Catalog Number	PK-CA577-1170-5
Description	A synthetic peptide that is potent, cell permeable, nontoxic and irreversibly inhibits caspase activity to blocks apoptosis. The new generation of caspase inhibitor is more stable in aqueous environment and exhibits several folds higher activity than the corresponding FMK caspase inhibitors. Q-VD-OPH is the inhibitor of choice for both in vitro and in vivo studies. (Caution: If the intended use is on purified or recombinant enzymes, esterase should be added to generate free carboxyl groups.). IC ₅₀ = 20-40 nM.
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	Quinolyl-Val-Asp-OPH; 514 Da
Purity	>95% by HPLC analysis
Appearance / Formulation / Solubility	White to off-white solid. Soluble in DMSO (200 mg/ml).
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.
References	NA
Caution	NA

Inhibition of Caspase-3 activity by next generation caspase inhibitor, QVD-OPh:
 Different concentrations of inhibitor Q-VD-OPh were tested to check the inhibition of Caspase-3 activity. Active Caspase-3 was incubated with the inhibitor Q-VD-OPh for 7 minutes prior to addition of synthetic peptide substrate DEVD-AFC (AFC, 7-amino-4-trifluoromethyl coumarin). Fluorescence was measured at Ex/Em = 400/505 nm.



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