

Instruction Manual

Catalog Number	PK-CA577-1007-1000
Description	Ready-to-use fluorometric substrate for caspase-3/CPP32 ($K_m = 9.7 \mu\text{M}$) and related caspases that recognize the amino acid sequence DEVD. The sequence DEVD is based on caspase-3 cleavage site in poly (ADP-ribose) polymerase (PARP). CPP32 and related caspase activity can be quantified by fluorescent detection of free AFC after cleaved from the peptide substrate DEVD-AFC at Ex. = 400 nm and Em. = 505 nm, using a fluorometer or a multi-well fluorescence plate reader. Alternatively, a shift in fluorescence from blue to green upon cleavage can be visualized using a hand-held long-UV lamp.
Quantity	1000 assays
Sequence / Molecular Weight / Molecular Formula	Ac-Asp-Glu-Val-Asp-AFC (AFC, 7-amino-4-trifluoromethyl coumarin); 729 Da
Purity	>98% by HPLC analysis.
Appearance / Formulation / Solubility	1 mM in DMSO
Storage & Stability	Store at -20°C , protected from light. Stable for 6 months under proper storage conditions.
Applications	<ol style="list-style-type: none"> 1. Induce apoptosis in cells by desired method. Concurrently incubate a control culture without induction. 2. Count cells and pellet $1-5 \times 10^6$ cells or use 50-200 μg cell lysates if protein concentration has been measured. 3. Resuspend cells in 50 μl of chilled Cell Lysis Buffer (PK-CA577-1067-100, PK-CA577-1067-400) 4. Incubate cells on ice for 10 minutes. 5. Add 50 μl of 2X Reaction Buffer (Cat. # PK-CA577-1068-20, PK-CA577-1068-80) containing 10 μM DTT (Cat. # PK-CA577-1201-1) to each sample. 6. Add 5 μl of the 1 mM DEVD-AFC (50 μM final conc.) into each tube individually and incubate at 37°C for 1-2 hour. 7. Read samples in a fluorometer equipped with a 400-nm excitation filter and 505-nm emission filter. For a plate-reading set-up, transfer the samples to a 96-well plate. You may perform the entire assay directly in a 96-well plate. 8. Fold-increase in Caspase-3 activity can be determined by comparing these results with the level of the uninduced control.
References	NA
Caution	NA

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