

Instruction Manual

Catalog Number	PK-CA577-1113-200
Description	Ready-to-use fluorometric substrate for AEVD-dependent caspases. Activity of AEVD-dependent caspase activity can be quantified by fluorescent detection of free AFC after cleavage from the peptide substrate AEVD-AFC at Ex. = 400 nm and Em. = 505 nm, using a fluorometer or 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. Cell Lysis Buffer (Cat. # PK-CA577-1067-100, PK-CA577-1067-400), 2X Reaction Buffer (Cat. # PK-CA577-1068-20, PK-CA577-1068-80), DTT (Cat. # PK-CA577-1201-1) and other reagents used for caspase assays are also available separately.
Quantity	200 assays
Sequence / Molecular Weight / Molecular Formula	Ac-Ala-Glu-Val-Asp-AFC (AFC, 7-amino-4-trifluoromethyl coumarin); 685 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 (Cat.# 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 AEVD-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. <p>Fold-increase in AEVD-dependent caspase activity can be determined by comparing these results with the level of the uninduced control.</p>
References	NA
Caution	NA

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