

Instruction Manual

Catalog Number	PK-CA577-1103-1000
Description	Ready-to-use fluorometric substrate for caspase-1/ICE and related caspases that recognize the amino acid sequence YVAD. Caspase-1 and related caspase activity can be quantified by fluorescent detection of free AFC after cleaved from the peptide substrate YVAD-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. The ready-to-use caspase substrate provides an economic alternative for researchers who perform large volume caspase assays. Cell Lysis Buffer (Cat. # PK-CA577-1067-100, PK-CA577-1067-400) and 2X Reaction Buffer (Cat. # PK-CA577-1068-20, PK-CA577-1068-80), and DTT (Cat. # PK-CA577-1201-1) used for caspase assays are also available separately.
Quantity	1000 assays
Sequence / Molecular Weight / Molecular Formula	Ac-Tyr-Val-Ala-Asp-AFC (AFC, 7-amino-4-trifluoromethyl coumarin); 719,7 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 or treat cells by desired method. Concurrently incubate a control culture without treatment. Note: Active recombinant human Caspase-1 is available to use as a positive control (Cat.# PK-RP577-1081-25, PK-RP577-1081-100). 2. Pellet 2-5 x 10⁸ cells or use 100-300 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.# 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 YVAD-AFC substrate (50 µM final concentration) and incubate at 37°C for 1-2 hour. 7. Read samples in a fluorometer equipped with a 400-nm excitation and 505-nm emission filters. For a plate-reading set-up, transfer the samples to a 96-well plate. You may also perform the entire assay directly in a 96-well plate. Fold-increase in YVAD-dependent caspase activity can be determined by comparing these results with the level of the untreated control.
References	NA
Caution	NA

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