

## Instruction Manual

Catalog Number	PK-CA577-1103-200
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	200 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"> <li>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).</li> <li>2. Pellet 2-5 x 10<sup>8</sup> cells or use 100-300 µg cell lysates if protein concentration has been measured.</li> <li>3. Resuspend cells in 50 µl of chilled Cell Lysis Buffer (Cat.# PK-CA577-1067-100, PK-CA577-1067-400).</li> <li>4. Incubate cells on ice for 10 minutes.</li> <li>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.</li> <li>6. Add 5 µl of the 1 mM YVAD-AFC substrate (50 µM final concentration) and incubate at 37°C for 1-2 hour.</li> <li>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.</li> </ol>
References	NA
Caution	NA

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