

## Instruction Manual

<b>Catalog Number</b>	PK-CA577-1117-1000
<b>Description</b>	Ready-to-use fluorometric substrate for caspases that recognize the amino acid sequence ATAD. Caspase activity can be quantified by fluorescent detection of free AFC after cleaved from the peptide substrate ATAD-AFC at Ex/Em = 400/505 nm using a fluorometer or a 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 amount of caspase assays. Cell Lysis Buffer (Cat. 1067-100), 2X Reaction Buffer (PK-CA577-1068-20, -80), DTT (Cat. # PK-CA577-1201-1), and other reagents used for caspase activity assays are also available separately.
<b>Quantity</b>	1000 assays
<b>Sequence / Molecular Weight / Molecular Formula</b>	Ac-Ala-Thr-Ala-Asp-AFC (AFC, 7-amino-4-trifluoromethyl coumarin); 630,56 Da
<b>Purity</b>	>90% by HPLC analysis
<b>Appearance / Formulation / Solubility</b>	4 mM in DMSO
<b>Storage &amp; Stability</b>	Store at -20°C, protected from light. Stable for 6 months
<b>Applications</b>	<ol style="list-style-type: none"> <li>1. Induce apoptosis in cells by desired method. Concurrently incubate a control culture without induction.</li> <li>2. Count cells and pellet <math>1-5 \times 10^6</math> cells or use 50-200 <math>\mu\text{g}</math> cell lysates if protein concentration has been measured.</li> <li>3. Resuspend cells in 50 <math>\mu\text{l}</math> 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 <math>\mu\text{l}</math> of 2X Reaction Buffer (Cat.# PK-CA577-1068-20, PK-CA577-1068-80) containing 10 <math>\mu\text{M}</math> DTT (Cat. # PK-CA577-1201-1) to each sample.</li> <li>6. Add 5 <math>\mu\text{l}</math> of the 1 mM ATAD-AFC (50 <math>\mu\text{M}</math> final conc.) into each tube individually and incubate at 37°C for 1-2 hour.</li> <li>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.</li> </ol> <p>Fold-increase in ATAD-dependent caspase activity can be determined by comparing these results with the level of the uninduced control.</p>
<b>References</b>	NA
<b>Caution</b>	NA

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