

Instruction Manual

Catalog Number	PK-CA577-K131
Description	Contains ready-to-use caspase-1/-2/-3/-5/-6/-8/-9 AFC-labeled substrates. All substrates are provided in liquid ready-to-use form.
Quantity	7 x 25 assays
Sequence / Molecular Weight / Molecular Formula	NA
Purity	see data sheets of individual caspase substrates
Appearance / Formulation / Solubility	Solution in DMSO
Storage & Stability	Store at -20°C. Stable for 1 year 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). 4. Incubate cells on ice for 10 minutes. 5. Add 50 μl of 2X Reaction Buffer (Cat.# PK-CA577-1068-20) containing 10 mM DTT (Cat.# PK-CA577-1201-1) to each sample. 6. Add 5 μl of the 1 mM AFC conjugated substrates (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 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.