

Instruction Manual

Catalog Number	PK-CA577-K132
Description	Contains ready-to-use caspase-1/-2/-3/-5/-6/-8/-9 pNA-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 $1-5 \times 10^8$ cells. 3. Resuspend cells in 50 μl of chilled Cell Lysis Buffer (Cat.# PK-CA577-1067-100) and incubate cells on ice for 10 minutes. 4. Centrifuge for 1 minute in a microcentrifuge (10,000 x g). 5. Transfer supernatant to a fresh tube and assay protein concentration. 6. Dilute 50-200 μg protein to 50 μl Cell Lysis Buffer for each assay. 7. Add 50 μl of 2X Reaction Buffer (Cat.# PK-CA577-1068-20) containing 10 mM DTT (Cat.# PK-CA577-1201-1) to each sample. 8. Add 5 μl of the 4 mM pNA conjugated substrates (200 μM final conc.) into each tube individually and incubate at 37°C for 1-2 hour. 9. Read samples at 400- or 405-nm in a microtiter plate reader, or spectrophotometer using a 100-μl micro quartz cuvet (Sigma), or dilute sample to 1 ml with Dilution Buffer (Cat.# PK-CA577-1066-100, PK-CA577-1066-500) and using regular cuvet (note: Dilution of the samples proportionally decreases the reading). <p>Fold-increase in caspase activity can be determined by comparing these results with the level of the uninduced control.</p> <p>Note: Background reading from cell lysates and buffers should be subtracted from the readings of both induced and the uninduced samples before calculating fold increase in caspase activity.</p>
References	NA
Caution	NA

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