

## Instruction Manual

<b>Catalog Number</b>	PK-RP577-K232
<b>Components</b>	Active Recombinant Caspase-1, 25 units, Yellow cap Active Recombinant Caspase-2, 25 units, Orange cap Active Recombinant Caspase-3, 25 units, Red cap Active Recombinant Caspase-6, 25 units, Violet cap Active Recombinant Caspase-7, 25 units, Red cap Active Recombinant Caspase-8, 25 units, Green cap Active Recombinant Caspase-9, 25 units, Blue cap Active Recombinant Caspase-10a, 25 units, Amber cap
<b>Quantity</b>	8 x 25 units
<b>Unit Definition</b>	One unit of the recombinant human caspases is the enzyme activity that cleaves 1 nmol of the individual caspase substrate (YVAD-pNA for caspase-1, VDVAD-pNA for caspase-2, DEVD-pNA for caspase-3, VEID-pNA for caspase-6, DEVD-pNA for caspase-7, IETD-pNA for caspase-8, LEHD-pNA for caspase-9, and IETD-pNA for caspase-10) per hour at 37°C in a reaction solution containing 50 mM Hepes, pH 7.2, 50 mM NaCl, 0.1% Chaps, 10 mM EDTA, 5% Glycerol, and 10 mM DTT.
<b>Source</b>	E. coli
<b>Formulation</b>	Lyophilized powder
<b>Reconstitution</b>	The active recombinant caspases can be reconstituted to 0.1-1 unit per µl in PBS or - for longer stability - in PBS containing 15% glycerol or the Reaction Buffer described above (also available separately from PromoKine, Cat. No. PK-CA577-1068-20 and PK-CA577-1068-80). We recommend using 1 unit per assay for analyzing caspase activity.
<b>Storage</b>	Store at -70°C. Stable for 6-12 months.
<b>Applications</b>	Highly active recombinant human caspases were expressed in E. coli and purified. Active recombinant caspases are useful in screening caspase inhibitors, studying enzyme kinetics and regulation, determining substrate specificity, and as a positive control in caspase activity assays. For a complete caspase activity assay protocol, please refer to PromoKine's Fluorometric or Colorimetric Caspase Activity Assay Kits.

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