

## Instruction Manual

<b>Catalog Number</b>	PK-RP577-K233
<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-4, 25 units, Orange cap Active Recombinant Caspase-5, 25 units, Clear 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-10, 25 units, Amber cap
<b>Quantity</b>	10 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, WEHD-pNA for caspase-4 and -5, VEID-pNA for caspase-6, DEVD-pNA for caspase-7, IETD-pNA for caspase-8 and -10 and LEHD-pNA for caspase-9) 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 $\mu$ l 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>	Active recombinant caspases are useful in screening caspase inhibitors, studying enzyme kinetics, 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.