

# Active Recombinant Caspase Set IV

**CATALOG #:** K233-10-25  
**LOT #:** \_\_\_\_\_  
**SOURCE:** *E. coli*  
**FORM:** Lyophilized powder  
**STORAGE:** Store at -70°C. Stable for 6-12 months.

**PACKAGE CONTENT:**

Description	Amount	Cap color
Active Recombinant Caspase-1	25 units	Yellow
Active Recombinant Caspase-2	25 units	Orange
Active Recombinant Caspase-3	25 units	Red
Active Recombinant Caspase-4	25 units	Orange
Active Recombinant Caspase-5	25 units	Clear
Active Recombinant Caspase-6	25 units	Violet
Active Recombinant Caspase-7	25 units	Red
Active Recombinant Caspase-8	25 units	Green
Active Recombinant Caspase-9	25 units	Blue
Active Recombinant Caspase-10	25 units	Amber

**UNIT DEFINATION:**

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, VDAD-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.

**RECONSTITUTION:**

The active recombinant caspases can be reconstituted to 0.1-1 unit per µl in PBS or the Reaction Buffer described above (available from BioVision, Cat.# 1068-20) for longer stability. We recommend using 1 unit per assay for analyzing caspase activity.

**APPLICATIONS AND USAGE:**

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 BioVision's Fluorometric or Colorimetric Caspase Activity Assay Kits.

**FOR RESEARCH USE ONLY! Not to be used in human.**

**RELATED PRODUCTS:**

- Apoptosis Detection Kits & Reagents
  - Annexin V Kits & Bulk Reagents
  - Caspase Assay Kits & Reagents
  - Mitochondrial Apoptosis Kits & Reagents
  - Nuclear Apoptosis Kits & Reagents
  - Apoptosis Inducers and Set
  - Apoptotic Cell Isolation Kit
- Cell Fractionation System
  - Mitochondria/Cytosol Fractionation Kit
  - Nuclear/Cytosol Fractionation Kit
  - Membrane Protein Extraction Kit
  - Cytosol/Particulate Rapid Separation Kit
  - Mammalian Cell Extraction Kit
  - FractionPREP Fractionation System
- Cell Proliferation & Senescence
  - Quick Cell Proliferation Assay Kit
  - Senescence Detection Kit
  - High Throughput Apoptosis/Cell Viability Assay Kits
  - LDH-Cytotoxicity Assay Kit
  - Bioluminescence Cytotoxicity Assay Kit
  - Live/Dead Cell Staining Kit
- Cell Damage & Repair
  - HDAC Fluorometric & Colorimetric Assays & Drug Discovery Kits
  - HAT Colorimetric Assay Kit & Reagents
  - DNA Damage Quantification Kit
  - Glutathione Fluorometric & Colorimetric Assay Kits
  - Nitric Oxide Fluorometric & Colorimetric Assay Kits
- Signal Transduction
  - cAMP & cGMP Assay Kits
  - Akt & JNK Activity Assay Kits
  - Beta-Secretase Activity Assay Kit
- Adipocyte & Lipid Transfer
  - Recombinant Adiponectin, Survivin, & Leptin
  - CETP Activity Assay & Drug Discovery Kits
  - PLTP Activity Assay & Drug Discovery Kits
  - Total Cholesterol Quantification Kit
- Molecular Biology & Reporter Assays
  - siRNA Apoptosis and Cloning Vectors
  - Cloning Insert Quick Screening Kit
  - Mitochondrial & Genomic DNA Isolation Kits
  - 5 Minutes DNA Ligation Kit
  - 20 Minutes Gel Staining/Destaining Kit
- Growth Factors and Cytokines
- Antibodies to Apoptosis and Cell Signaling Molecules