

Active Recombinant Caspase Set IV

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

PACKAGE CONTENT:

Description	Amount	Cap Color	Part Number
Recombinant Caspase-1	25 units	Yellow	1081-25
Recombinant Caspase-2	25 units	Orange	1082-25
Recombinant Caspase-3	25 units	Red	1083-25
Recombinant Caspase-4	25 units	Orange	1084-25
Recombinant Caspase-5	25 units	Clear	1085-25
Recombinant Caspase-6	25 units	Violet	1086-25
Recombinant Caspase-7	25 units	Red	1087-25
Recombinant Caspase-8	25 units	Green	1088-25
Recombinant Caspase-9	25 units	Blue	1089-25
Recombinant Caspase-10	25 units	Amber	1090A-25

UNIT DEFINITION:

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 humans.

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 Minute DNA Ligation Kit
 - 20 Minute Gel Staining/Destaining Kit
- Growth Factors and Cytokines
- Antibodies to Apoptosis and Cell Signaling Molecules