

Active Recombinant Caspase Set III

CATALOG #: ALX-850-243

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	ALX-850-243-1
Recombinant Caspase-2	25 units	Orange	ALX-850-243-2
Recombinant Caspase-3	25 units	Red	ALX-850-243-3
Recombinant Caspase-6	25 units	Violet	ALX-850-243-6
Recombinant Caspase-7	25 units	Red	ALX-850-243-7
Recombinant Caspase-8	25 units	Green	ALX-850-243-8
Recombinant Caspase-9	25 units	Blue	ALX-850-243-9
Recombinant Caspase-10	25 units	Amber	ALX-850-243-10

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, 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.

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:

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 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 Minutes DNA Ligation Kit
- 20 Minute Gel Staining/Destaining Kit

Growth Factors and Cytokines

Antibodies to Apoptosis and Cell Signaling Molecules