

## Active Recombinant Mouse Caspase-3

**CATALOG #:** 1183-25 25 units  
1183-100 100 units

**LOT #:** \_\_\_\_\_

### UNIT DEFINITION:

One unit of the recombinant caspase-3 is the enzyme activity that cleaves 1 nmol of the caspase substrate DEVD-pNA (pNA: *p*-nitroaniline) 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.

**FORM:** Lyophilized powder

**RECONSTITUTION:** Reconstitute to 1 unit per  $\mu$ l in PBS containing 15% glycerol.

**SPECIFIC ACTIVITY:** 300,000 units/mg

### STORAGE CONDITIONS:

The lyophilized caspase-3 is stable for 1 year at  $-70^{\circ}\text{C}$ . Following reconstitution in PBS, the enzyme should be aliquoted and immediately stored at  $-70^{\circ}\text{C}$ . Avoid multiple freeze/thaw cycles as activity might decrease.

### DESCRIPTION:

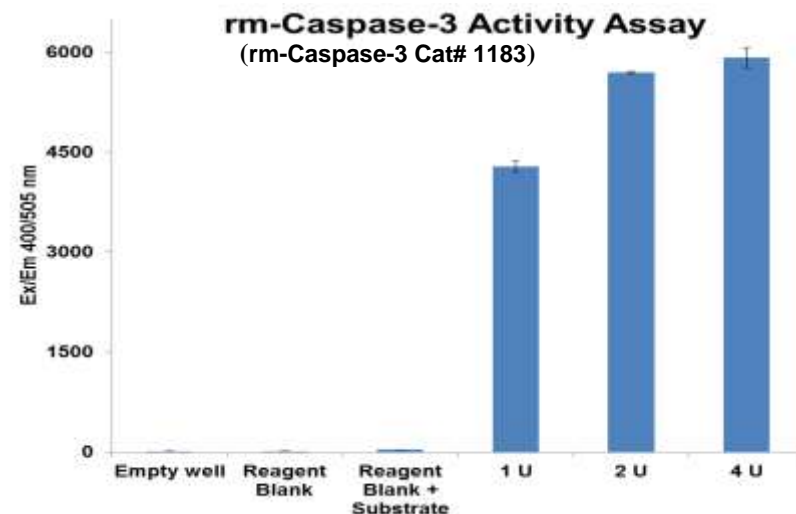
Caspase-3 (also known as CPP32, Yama and apopain) is a major member of the caspase-family of cysteine proteases. Caspase-3 exists in cells as an inactive 32 kDa proenzyme. During apoptosis procaspase-3 is processed at aspartate residues by self-proteolysis and/or cleavage by upstream caspases, such as caspase-6 (Mch2), caspase-8 (Flice) and granzyme B. The processed form of caspase-3 consists of large (17 kD) and small (11 kD) subunits which associate to form the active enzyme. The active caspase-3 has been shown involving in the proteolysis of several important molecules, such as poly (ADP-ribose) polymerase (PARP), the sterol regulatory element binding proteins (SREBPs), focal adhesion kinase (FAK), and others. The recombinant active human caspase-3 expressed in *E. coli* spontaneously undergoes autoprocessing to yield subunits characteristic of the native enzyme. The active caspase-3 preferentially cleaves caspase-3 substrates (e.g., DEVD-AFC or DEVD-pNA) and is routinely tested at BioVision for its ability to enzymatically cleave these two substrates Ac-DEVD-pNA (Cat. #1008-200) or Ac-DEVD-AFC (1007-200).

### APPLICATIONS AND USAGE:

Active caspase-3 is useful in studying enzyme regulation, determining target substrates, screening caspase inhibitors, or as a positive control in caspase activity assays. We recommend using 1 unit/assay for analyzing caspase activity.

For a complete caspase-3 assay protocol, please refer to BioVision's Caspase-3/ CPP32 Fluorometric or Colorimetric Assay Kits (Cat.#: K105 and K106).

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



**Recombinant Mouse Caspase-3 Activity Profile Using the Caspase-3 Fluorometric Assay Kit (K105-25)**

### 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
- Apoptosis siRNA Vectors

#### 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