

## Active Caspase-3 Polyclonal Antibody

<b>CATALOG #:</b>	3015-100
<b>AMOUNT:</b>	100 µg
<b>LOT #:</b>	
<b>HOST:</b>	Rabbit
<b>IMMUNOGEN:</b>	Synthetic peptide mapping to the N-terminus adjacent to the cleavage site of human caspase-3 (Internal ID# BV-2306)
<b>SPECIES REACTIVITY:</b>	Human, Mouse, Rat

**FORMULATION:**

500 µl (0.2 mg/ml) antigen affinity purified rabbit anti-Active Caspase-3 antibody in phosphate-buffered saline (PBS) containing 30% glycerol, 0.5% BSA, and 0.01% thimerosal.

**STORAGE CONDITIONS:**

Store at -20°C. For long term storage, aliquot and freeze at -70°C. Avoid repeated freeze/thaw cycles.

**DESCRIPTION:**

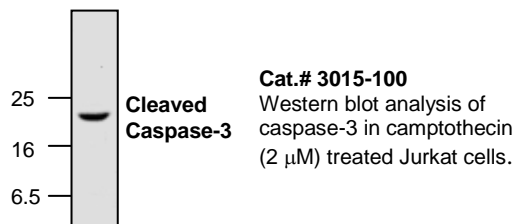
Caspase-3 has been extensively studied and implicated to play an important role in apoptosis. Active caspase-3 proteolytically cleaves and activates other caspases, as well as relevant targets in the cells (e.g., PARP). The antigen affinity purified antibody recognizing the active form of caspase-3 provides a new tool for identifying apoptotic cell populations in both tissue sections and cultured cells.

**SPECIFICITY:**

The antibody preferentially recognizes the p17 fragment of the active caspase-3 in samples from human, mouse, and rat origins.

**APPLICATION AND USAGE:**

The antibody can be used for Western blotting (1:50 -1:500 dilutions) and immunohistochemistry (1:10-1:20 dilutions). However, the optimal concentrations should be determined individually.



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

**RELATED PRODUCTS:**

## Apoptosis Detection Kits &amp; 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 &amp; 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 &amp; Repair

- HDAC Fluorometric & Colorimetric Assays & Drug Discovery Kits
- HAT Colorimetric Assay Kit & Reagents
- DNA Damage Quantification Kit
- Glutathione & Nitric Oxide Fluorometric & Colorimetric Assay Kits

## Signal Transduction

- cAMP & cGMP Assay Kits
- Akt & JNK Activity Assay Kits
- Beta-Secretase Activity Assay Kit

## Adipocyte &amp; Lipid Transfer

- Recombinant Adiponectin, Survivin, & Leptin
- CETP Activity Assay & Drug Discovery Kits
- PLTP Activity Assay & Drug Discovery Kits
- Total Cholesterol Quantification Kit

## Molecular Biology &amp; Reporter Assays

- siRNA Vectors
- Cloning Insert Quick Screening Kit
- Mitochondrial & Genomic DNA Isolation Kits
- 5 Minutes DNA Ligation Kit
- 20 Minutes Gel Staining/Destaining Kit
- β-Galactosidase Staining Kit & Luciferase Reporter Assay Kit

## Growth Factors and Cytokines

## Monoclonal and Polyclonal Antibodies