

Caspase-9 Polyclonal Antibody

CATALOG #:	3016-100
AMOUNT:	100 µg
LOT #:	_____
HOST:	Rabbit
IMMUNOGEN:	Synthetic peptide surrounding amino acid 290 of human caspase-9 (ID# BV-12)
SPECIES REACTIVITY:	Human, Mouse, Rat, Hamster, Rabbit, Pig, Bovine

FORMULATION:

100 µg (0.2 mg/ml) affinity purified rabbit anti-caspase-9 polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 0.01% thimerosal.

STORAGE CONDITIONS:

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

BACKGROUND:

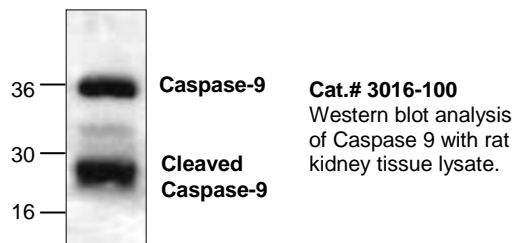
Caspase-9 is one of the most important caspases among the caspase family members. Caspase-9 is synthesized as inactive pro-enzyme that is processed in cells undergoing apoptosis. The processed caspase-9 consists of large (32-35 kD) and small (10 kD) subunits which associate to form an active enzyme. Activated caspase-9 in turn cleaves and activates caspase-3, one of the proteases responsible for the proteolytic cleavage of many key proteins in apoptosis.

APPLICATION AND USAGE:

The antibody can be used for Western blotting (0.5-4 µg/ml). However, the optimal conditions should be determined individually. Blocking peptide is available separately (Cat.# 3016BP-50).

SPECIFICITY:

The antibody detects proform (~46 kDa), intermediate (~33-36 kDa), and the fully cleaved (~20-25 kDa) caspase-9.



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

Cell Damage & 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 & 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 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