

Caspase-9 Polyclonal Antibody

CATALOG #:	3409-100
AMOUNT:	100 µg
LOT #:	_____
HOST:	Rabbit
IMMUNOGEN:	Synthetic peptide surrounding amino acid 345 of mouse caspase-9 (ID# BV-13-14)
SPECIES REACTIVITY:	mouse, rat

FORMULATION:

100 µg (0.2 mg/ml) protein A 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:

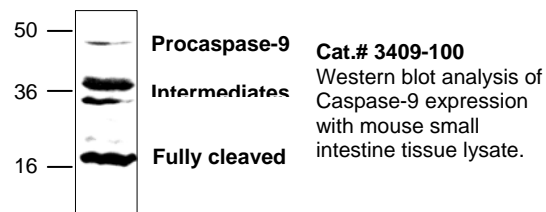
Caspase-9 is one of the most important caspases among the caspase family members. Upon induction of apoptosis, cytochrome c released from mitochondria associates with procaspase-9/Apaf-1. The complex processes procaspase-9 into a large subunit (35 kDa or 17 kDa) and small (10 kDa) by self-cleavage at D315. Activated caspase-9 further cleaves other caspase members including caspase-3, one of the proteases responsible for the proteolytic cleavage of many key proteins in apoptosis. In addition to self-cleavage, procaspase-9 can also be cleaved in vivo by caspase-3 at D330. The process served as a positive feedback to amplify the apoptotic signal in caspase activation pathway.

APPLICATION AND USAGE:

The antibody can be used for Western blotting (0.5-4 µg/ml). However, the optimal conditions should be determined individually. Mouse small intestine and rat kidney tissue lysates can be used as positive controls. Blocking peptide (Cat.# 3409BP-50) is available separately.

SPECIFICITY:

The antibody detects 45 kDa procaspase-9, 35/37 kDa cleaved large fragment and the 25 kDa fully cleaved 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