

Caspase-8 Polyclonal Antibody

CATALOG #:	3158-100
AMOUNT:	100 µg
LOT #:	_____
HOST:	Rabbit
IMMUNOGEN:	Synthetic peptide surrounding the cleavage site of human Caspase-8 (ID# BV-006)
SPECIES REACTIVITY:	human.

FORMULATION:

100 µg (0.5 mg/ml) affinity purified rabbit anti-caspase-8 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 DESCRIPTION:

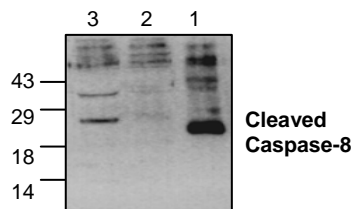
Caspase family of cysteine proteases has been shown to play a key role in apoptosis. Caspase-8 is a 55 kDa cytosolic protein that is synthesized as an inactive pro-enzyme. Activation of caspase-8 involves a two-step proteolysis: the cleavage of caspase-8 to generate a 43 and a 12 kDa fragment which is further processed to 10 kDa. The p43 is then cleaved to yield p26 and the release of the active site containing p18.

SPECIFICITY:

The affinity purified antibody mainly detects the 26 kDa fully cleaved fragment and the 35-43 kDa intermediate of caspase-8 in apoptotic samples.

APPLICATION AND USAGE:

The antibody can be used for Western blotting (0.5-4 µg/ml). Based on researchers feedback, it can also be used in immunohistochemistry (10-20 mg/ml). However, the optimal conditions should be determined individually. Blocking peptide (Cat.# 3158BP-50) is available separately.



Cat.# 3158-100

Western blot analysis of Caspase-8.
Lane 1: Recombinant h-caspase-8;
Lane 2: Untreated Jurkat cells;
Lane 3: Camptothecin induced cells.

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