

Caspase-10 Polyclonal Antibody

CATALOG #:	3410-200
AMOUNT:	200 µg
LOT #:	_____
HOST:	Rabbit
IMMUNOGEN:	Synthetic peptide surrounding amino acid 404 of human caspase-10 (ID# BV-15)
SPECIES REACTIVITY:	Human, Mouse, Rat

FORMULATION:

200 µg (200 µg/ml) affinity purified rabbit anti-caspase-10 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:

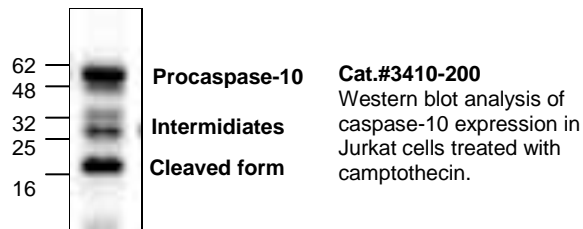
Caspase family of cysteine proteases play a key role in apoptosis. Caspase-10 is expressed in many tissues and cell lines, and has been identified as caspase-10/a (short form) and caspase-10/b (long form), respectively. Similar as other caspases, caspase-10 is also synthesized as an inactive pro-enzyme that is processed in cells undergoing apoptosis. Activated caspase-10 activates caspase-3,-4,-6,-7,-8, and -9, which are then involved in proteolytic cleavage of many cellular proteins. Caspase-10 has been shown involving in the Fas and TNFR1 induced apoptosis.

SPECIFICITY:

The antibody detects the proform (~55 kDa), intermediates (30 kDa and 34 kDa), and cleaved form (~20 kDa) of caspase-10 on SDS-PAGE immunoblots.

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.# 3410BP-50).



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

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