

Caspase-6 Polyclonal Antibody

CATALOG #:	3256-100
AMOUNT:	100 µg
LOT #:	_____
HOST:	Rabbit
IMMUNOGEN:	Synthetic peptide mapping to the C-terminus of rat caspase-6 large fragment (ID# BV-24)
SPECIES REACTIVITY:	human, mouse, rat

FORMULATION:

100 µg (0.2 mg/ml) rabbit anti-caspase-6 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:

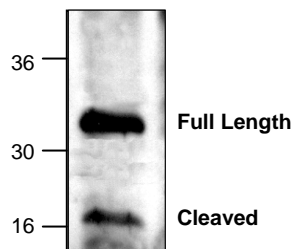
Caspase-6 is one of the major executioner caspases functioning in cellular apoptotic processes. Similar as other caspases, caspase-6 is also synthesized as an inactive proenzyme that is processed in cells undergoing apoptosis. Caspase-6 has also been shown involving in the proteolysis of poly (ADP-ribose) polymerase (PARP) and nuclear lamin A.

SPECIFICITY:

The antibody detects a full length (35 kDa) and the cleaved fragment (18 kDa) of caspase-6. Blocking Peptide is available separately (Cat.# 3256BP-50).

APPLICATION AND USAGE:

The antibody can be used for Western blotting (0.5-4 µg/ml). However, the optimal conditions should be determined individually. Jurkat cell lysate (Cat.# 2401-100), mouse intestine (Cat.# 2601-100), and rat kidney (Cat. 2702-100) tissue lysate can be used as a positive controls.



Cat.# 3256-100
Western blot analysis of Caspase-6 expression rat kidney tissue lysate. Both full length and the cleaved large fragment have been detected.

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