

## Active Caspase-8 Polyclonal Antibody

<b>CATALOG #:</b>	3259-100BFGFAF
<b>AMOUNT:</b>	100 µg
<b>LOT #:</b>	_____
<b>HOST:</b>	Rabbit
<b>IMMUNOGEN:</b>	Synthetic peptide corresponding to residues surrounding amino acid 392 of mouse caspase-8 (ID# BV-9).
<b>SPECIES REACTIVITY:</b>	Mouse, rat
<b>ANTIBODY FORMULATION:</b>	100 µg (0.5 mg/ml) affinity purified rabbit polyclonal antibody in PBS (No azide, No glycerol, no BSA).

### STORAGE CONDITIONS:

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

### DESCRIPTION:

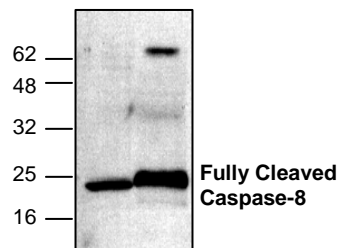
Caspase-8, a member of the caspase-family of proteases, plays a key role in mediating Fas (CD95) and TNF induced apoptosis. Caspase-8 is synthesized as inactive pro-enzyme and activation of the enzyme involves proteolytic cleavage that leads to the release of the active p18 and p10 subunits. Activated caspase-8 is able to cleave and activate downstream caspases, such as caspase-3, -6, -7 and a death agonist member of the Bcl-2/Bcl-xL family, Bid, leading to apoptosis.

### SPECIFICITY:

The antibody detects the 25 kDa large subunit of caspase-8. It does not detect the full-length of caspase-8.

### APPLICATIONS AND USAGE:

The antibody can be used for Western blotting (0.5-4 µg/ml) in samples from mouse and rat origins. However, the optimal conditions should be determined individually.



### Cat.# 3259-100

Western blot analysis of caspase-8 expression in mouse intestine and rat kidney tissue lysate.

### 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