
Caspase3 Polyclonal Antibody

(Catalog #A71051)

Background

This gene encodes a protein which is a member of the cysteine-aspartic acid protease (caspase) family. Sequential activation of caspases plays a central role in the execution-phase of cell apoptosis. Caspases exist as inactive proenzymes which undergo proteolytic processing at conserved aspartic residues to produce two subunits, large and small, that dimerize to form the active enzyme. This protein cleaves and activates caspases 6, 7 and 9, and the protein itself is processed by caspases 8, 9 and 10. It is the predominant caspase involved in the cleavage of amyloid-beta 4A precursor protein, which is associated with neuronal death in Alzheimer's disease. Alternative splicing of this gene results in two transcript variants that encode the same protein.

Description

Caspase3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.05% proclin300, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P42574

Purification

Affinity Purification

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 6-105 of human Caspase-3 (NP_004337.2).

Storage

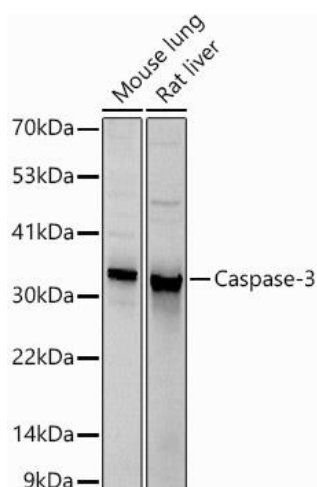
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

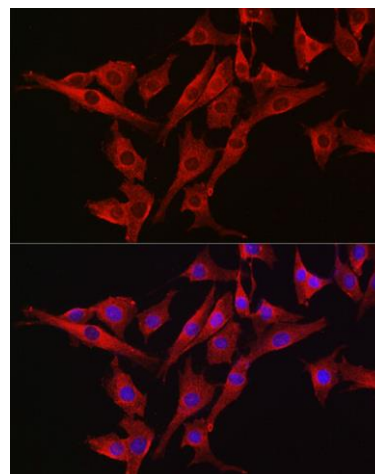
CASP3; CPP32; CPP32B; SCA-1; caspase-3

Application

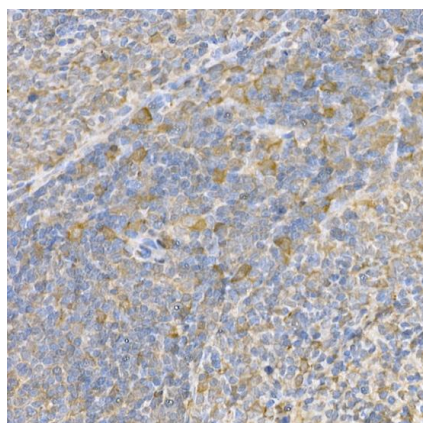
WB, IHC, IF; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:200, IF 1:100 - 1:500



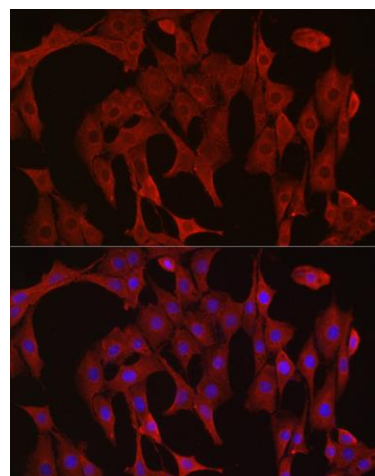
Western blot analysis of extracts of various cell lines, using Caspase3 Polyclonal Antibody at 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25µg per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 5s.



Immunofluorescence analysis of NIH/3T3 cells using Caspase3 Polyclonal Antibody at dilution of 1:300 (40x lens). Blue: DAPI for nuclear staining.



Immunohistochemistry analysis of paraffin-embedded rat spleen using Caspase3 Polyclonal Antibody at dilution of 1:100 (40x lens). Perform high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunofluorescence analysis of PC-12 cells using Caspase3 Polyclonal Antibody at dilution of 1:300 (40x lens). Blue: DAPI for nuclear staining.