Caspase-4 monoclonal antibody (4B9)

The caspases are a family of cysteine proteases that cleave after certain aspartate residues, and are primarily recognized as mediators of apoptosis. caspases are synthesized as inactive zymogens that can be cleaved to form active enzymes following the induction of apoptosis by stress or death receptors. Initiator caspases (e.g. caspase-8 and -10) are activated by dimerization of the zymogen on a dedicated adaptor protein. These activated initiator caspases in-turn cleave downstream effector or executioner caspases (e.g. caspase-3, -6, and -7) in a cascade-like manner, which cleave key cellular proteins that lead to the morphological changes associated with apoptotic cell death.

This antibody is covered by our Worry-Free Guarantee.

Citations: 9

View Online »

Ordering Information

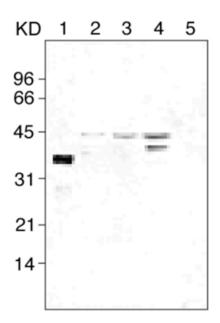
Order Online »

ADI-AAM-114-E

100µg

Manuals, SDS & CofA

View Online »



Western blot analysis of caspase-4 probed with Caspase-4, mAb (4B9). Lane 1: TXp20 (recombinant); Lane 2: U937 cell extract; Lane 3: HL60 cell extract; Lane 4: HUC-Fm cell extract; Lane 5: PC12 cell extract.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application WB

Application NotesDetects a band of ~45kDa by Western blot.

Clone 4B9

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol.

GenBank ID HSU28014

Host Mouse

Immunogen Recombinant human caspase-4 (N-terminal fragment).

lgG1

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000, ECL)Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

Species Reactivity Human

UniProt ID P49662

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

ENZO LIFE SCIENCES,

May 29, 2024

Phone: 800.942.0430

infousa@enzolifesciences.com

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com