

Anti-DAPK3

Rabbit Polyclonal Antibody

Catalog # D03-63R

Lot # O2109-7

Cited Applications

WB, ELISA, ICC, IF

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the DAPK3 protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

DAPK3 antibody was raised against a peptide corresponding to amino acids near the center of human DAPK3.

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

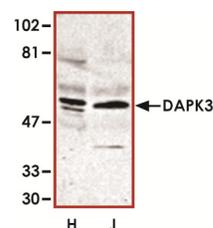
Scientific Background

DAPK3 or Death-associated protein kinase 3 (also known as ZIP) plays a role in apoptosis (1). DAPK3 is a nuclear serine/threonine-specific kinase that phosphorylates core histones H3 and H4, and myosin light chain in vitro. DAPK3 interacts with transcription and splicing factors as well as with pro-apoptotic protein Par-4 suggesting that it participates in multiple cellular processes. DAPK3 contains a leucine zipper structure at its C terminus and this region is responsible for binding to ATF4. The leucine zipper domain is necessary for the homodimerization of DAPK3 as well as for the activation of the kinase (2).

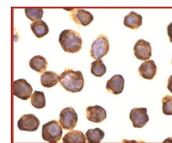
References

1. Kawai, T. et al: ZIP kinase, a novel serine/threonine kinase which mediates apoptosis. *Mol. Cell Biol.* 1998;18(3):1642-51.
2. Preuss, U. et al. Novel mitosis-specific phosphorylation of histone H3 at Thr11 mediated by Dlk/ZIP kinase. *Nucleic Acids Res.* 2003; 31(3):878-85.

Sample Data



Western blot analysis of DAPK3 in HeLa (H) and Jurkat (J) whole cell lysates with DAPK antibody at 1 µg/ml.



Immunocytochemistry of DAPK3 in Jurkat cells with DAPK3 antibody at 10 µg/ml.

Anti-DAPK3

Rabbit Polyclonal Antibody

Catalog Number

D03-63R

Specific Lot Number

O2109-7

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.