

Anti-DAPK2

Rabbit Polyclonal Antibody

Catalog # D02-63R

Lot # O2109-15

Cited Applications

WB, ELISA, IHC

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

Specificity

Recognizes the DAPK2 protein and has no cross response with DAPK1

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

DAPK2 antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human DAPK2. This antigenic peptide is identical to the corresponding amino acids of mouse origin.

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

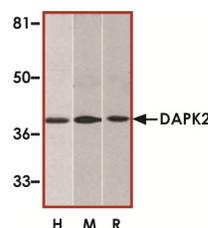
Scientific Background

DAPK2 or death-associated protein kinase 2 belongs to a family of proapoptotic Ca(2+)/calmodulin-regulated serine/threonine kinases. Overexpression of DAPK2 induces cell apoptosis. DAPK2 has been shown to be a novel Sp1-dependent target gene for E2F1 and Krüppel-like factor 6 (KLF6) in cell death response (1). Both E2F1 and KLF6 strongly activate the DAPK2 promoter. DAPK2 plays a role in granulopoiesis where it is highly expressed. β-catenin can block anoikis of malignant kidney and intestinal epithelial cells and promote their anchorage-independent growth by down-regulating DAPK2 (2). β-catenin-induced down-regulation of DAPK2 requires the presence of the transcription factor TCF-4.

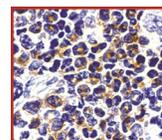
References

1. Britschgi A. et al: DAPK2 is a novel E2F1/KLF6 target gene involved in their proapoptotic function. *Oncogene*. 2008 Sep 25;27(43):5706-16.
2. Li, H. et al: Down-regulation of death-associated protein kinase-2 is required for beta-catenin-induced anoikis resistance of malignant epithelial cells. *J Biol Chem*. 2009 Jan 23;284(4):2012-22.

Sample Data



Western blot analysis of DAPK2 in A431 (H), mouse spleen (M), and rat kidney (R) lysates with DAPK2 antibody at 1 µg/ml.



Immunohistochemistry of DAPK2 in mouse spleen cells with DAPK2 antibody at 2 µg/ml.

Anti-DAPK2

Rabbit Polyclonal Antibody

Catalog Number

D02-63R

Specific Lot Number

O2109-15

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.