

Anti-ALK2

Rabbit Polyclonal Antibody

Catalog # A06-63R

Lot # O2121-13

Cited Applications

WB, ELISA

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

Specificity

Recognizes the ALK2 protein

Cross Reactivity

Human and Mouse

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

ALK2 antibody was raised against a 14 amino acid synthetic peptide near the amino terminus of the human ALK2

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

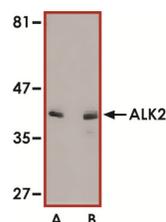
Scientific Background

ALK2 is a receptor serine/threonine kinase that is member of the ALK family and is upstream of signaling pathway involving the SMAD proteins especially SMAD1/5/8. Knockdown of ALK2, but not TGFβR1 (ALK5), abrogates endoglin-mediated decrease in cell motility of prostate cancer cells and constitutively active ALK2 is sufficient to restore a low-motility phenotype in endoglin deficient cells (1). Therefore, endoglin decreases prostate cancer cell motility through activation of the ALK2-Smad1 pathway. ALK2 is the key gene involved in Fibrodysplasia ossificans progressiva (FOP), a rare autosomal dominant congenital disorder characterized by progressive heterotopic bone formation in muscle tissues (2).

References

1. Craft, C.S. et al: Endoglin inhibits prostate cancer motility via activation of the ALK2-Smad1 pathway. *Oncogene*. 2007 Nov 8;26(51):7240-50.
2. Shore, E. M. et al: A recurrent mutation in the BMP type I receptor ACVR1 causes inherited and sporadic fibrodysplasia ossificans progressiva. *Nature Genet.* 38: 525-527, 2006.

Sample Data



Western blot analysis of ALK2 in A549 cell lysate with ALK2 antibody at 1 µg/ml in (A) the absence and (B) the presence of blocking peptide.

Anti-ALK2

Rabbit Polyclonal Antibody

Catalog Number

A06-63R

Specific Lot Number

O2121-13

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.