

Anti-ALK4

Rabbit Polyclonal Antibody

Catalog # A07-63R

Lot # O2121-14

Cited Applications

WB, ELISA

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

Specificity

Recognizes the ALK4 protein

Cross Reactivity

Human and Mouse

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

ALK4 antibody was raised against a 14 amino acid synthetic peptide near the carboxy terminus of the human ALK4

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

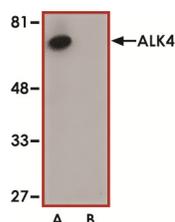
Scientific Background

ALK4 is a member of the subfamily of receptor ser/thr kinases that mediates signaling by the Activins. ALK4 is expressed in many human tissues, including kidney, pancreas, brain, lung, and liver. Truncated ALK4, predominantly expressed in human pituitary adenomas, function as dominant negative receptors to interfere with wild-type receptor function and blocks the antiproliferative effect of activin possibly contributing to development of human pituitary tumors (1). ALK4 is able to mediate Nodal signaling in the presence of Cripto during vertebrate development (2).

References

1. Zhou, Y. et al: Truncated activin type I receptor Alk4 isoforms are dominant negative receptors inhibiting activin signaling. Mol Endocrinol. 2000 Dec;14(12):2066-75.
2. Reissmann, E. et al: The orphan receptor ALK7 and the Activin receptor ALK4 mediate signaling by Nodal proteins during vertebrate development. Genes Dev. 2001 Aug 1; 15(15):2010-22.

Sample Data



Western blot analysis of ALK4 in human kidney tissue lysate with ALK4 antibody at 1 µg/ml in (A) the absence and (B) the presence of blocking peptide.

Anti-ALK4

Rabbit Polyclonal Antibody

Catalog Number

A07-63R

Specific Lot Number

O2121-14

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.