

Anti-VLK

Rabbit Polyclonal Antibody

Catalog # V05-163R

Lot # O2121-26

Cited Applications

WB, ELISA, IF, IHC

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

Specificity

Recognizes the VLK protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

VLK antibody was raised against a 15 amino acid synthetic peptide from near the center of human VLK

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

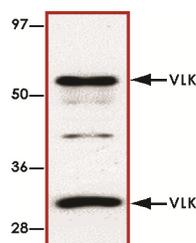
Scientific Background

VLK or vertebrate lonesome kinase is a protein kinase which is first expressed in E-cadherin-positive anterior visceral endoderm and mesendoderm, and later its expression is confined to E-cadherin-negative mesenchyme. VLK regulates the rate of protein export from the Golgi and thereby plays an important role in the formation of functional stroma by mesenchymal cells (1). Targeted disruption of VLK leads to a defect in lung development and delayed ossification of endochondral bone. VLK deficient mice display neonatal lethality due to respiratory failure, with a suckling defect arising from a cleft palate. VLK is required for the appropriate timing of flat proliferative chondrocyte differentiation (2).

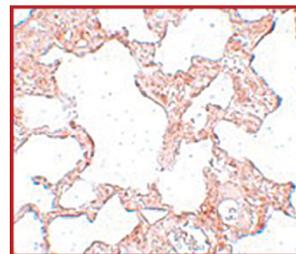
References

1. Kinoshita, et.al: The novel protein kinase Vlk is essential for stromal function of mesenchymal cells. *Development* 136: 2069-2079, 2009.
2. Imuta, Y. et al: Short limbs, cleft palate, and delayed formation of flat proliferative chondrocytes in mice with targeted disruption of a putative protein kinase gene, Pkdc (AW548124). *Dev Dyn.* 2009 Jan; 238(1):210-22.

Sample Data



Western blot analysis of VLK in human lung tissue lysate with VLK antibody at 1 ug/ml.



Immunohistochemistry of VLK in rat lung tissue with VLK antibody at 2.5 ug/ml.

Anti-VLK

Rabbit Polyclonal Antibody

Catalog Number

V05-163R

Specific Lot Number

O2121-26

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.