

Anti-ALK7

Rabbit Polyclonal Antibody

Catalog # A08-63R

Lot # O2121-15

Cited Applications

WB, ELISA

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

Specificity

Recognizes the ALK7 protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

ALK7 antibody was raised against a 15 amino acid synthetic peptide near the amino terminus of the human ALK7

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

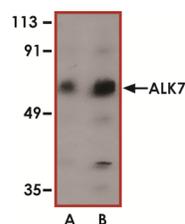
Scientific Background

ALK7 is a type I receptor for the TGFβ family of signaling molecules in which type I receptors phosphorylate cytoplasmic SMAD transcription factors, which then translocate to the nucleus and interact directly with DNA or in complex with other transcription factors (1). Alk7 was expressed in all insulin, glucagon, and somatostatin -positive cells of the pancreas (2). GDF3 regulates adipose tissue homeostasis and energy balance under nutrient overload, in part, by signaling through ALK7 and Alk7 showed reduced fat accumulation and partial resistance to diet-induced obesity, similar to Gdf3.

References

1. Bondestam, J. et.al: cDNA cloning, expression studies and chromosome mapping of human type I serine/threonine kinase receptor ALK7 (ACVR1C). *Cytogenet. Cell Genet.* 95: 157-162, 2001.
2. Bertolino, P. et.al: Activin B receptor ALK7 is a negative regulator of pancreatic beta-cell function. *Proc. Nat. Acad. Sci.* 105: 7246-7251, 2008.

Sample Data



Western blot analysis of ALK7 in human placenta tissue lysate with ALK7 antibody at (A) 1 and (B) 2 µg/ml.

Anti-ALK7

Rabbit Polyclonal Antibody

Catalog Number

A08-63R

Specific Lot Number

O2121-15

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.