
DLK1 Polyclonal Antibody

(Catalog # A71423)

Background

This gene encodes a transmembrane protein that contains multiple epidermal growth factor repeats that functions as a regulator of cell growth. The encoded protein is involved in the differentiation of several cell types including adipocytes. This gene is located in a region of chromosome 14 frequently showing uniparental disomy, and is imprinted and expressed from the paternal allele. A single nucleotide variant in this gene is associated with child and adolescent obesity and shows polar overdominance, where heterozygotes carrying an active paternal allele express the phenotype, while mutant homozygotes are normal.

Description

DLK1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P80370

Purification

Affinity Purification

Immunogen

Recombinant protein of human DLK1

Storage

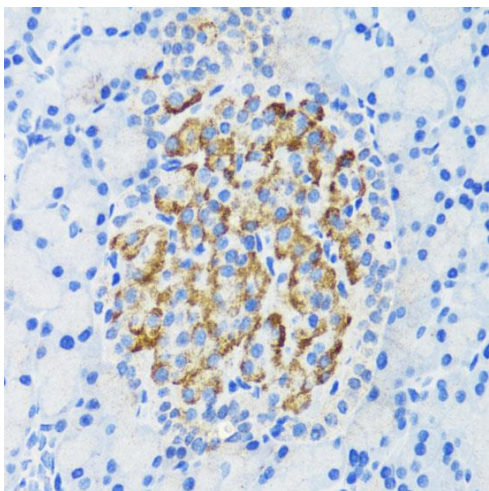
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

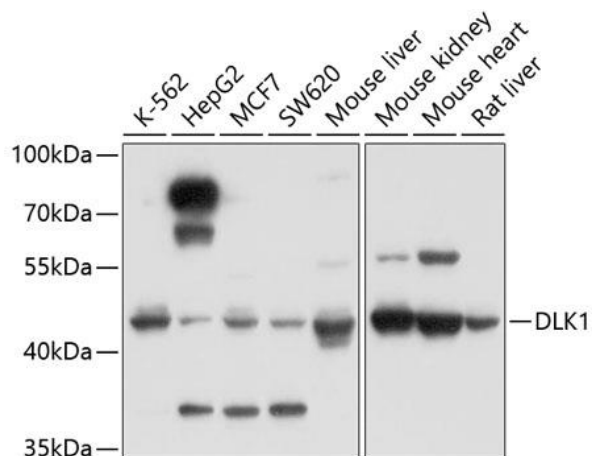
DLK1; DLK; DLK-1; Delta1; FA1; PREF1; Pref-1; ZOG; pG2; protein delta homolog 1

Application

WB, IHC, IF; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:100, IF 1:50 - 1:100



Immunohistochemistry of paraffin-embedded rat pancreatic islet using DLK1 Polyclonal Antibody at dilution of 1:200 (40x lens).



Western blot analysis of extracts of various cell lines, using DLK1 Polyclonal Antibody at 1:1000 dilution.
 Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.
 Exposure time: 30s.