
DKK3 Polyclonal Antibody

(Catalog #A71962)

Background

This gene encodes a protein that is a member of the dickkopf family. The secreted protein contains two cysteine rich regions and is involved in embryonic development through its interactions with the Wnt signaling pathway. The expression of this gene is decreased in a variety of cancer cell lines and it may function as a tumor suppressor gene. Alternative splicing results in multiple transcript variants encoding the same protein.

Description

DKK3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.05% proclin300, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q9UBP4

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 22-200 of human DKK3 (NP_056965.3)

Storage

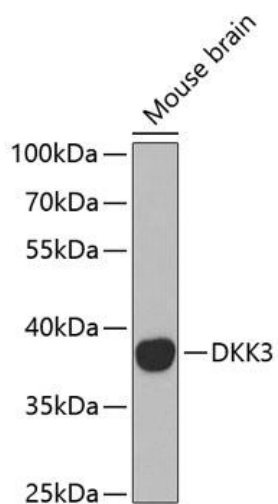
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

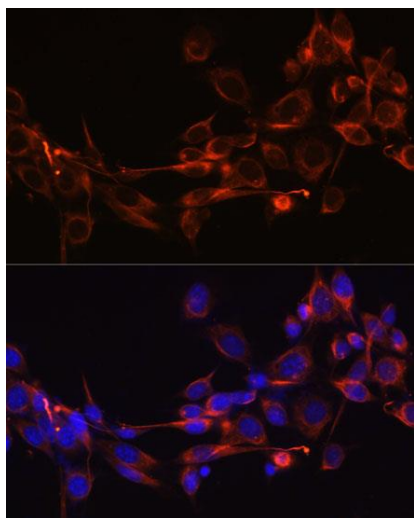
DKK3; REIC; RIG; dickkopf-related protein 3

Application

WB,IF; Recommended dilution: WB 1:500 – 1:1000, IF 1:50 – 1:200



Western blot analysis of extracts of mouse brain, using DKK3 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: 90s.



Immunofluorescence analysis of NIH-3T3 cells using DKK3 Polyclonal Antibody at dilution of 1:100 (40x lens).
Blue: DAPI for nuclear staining.