
TNRC6A Polyclonal Antibody

(Catalog # A72897)

Background

This gene encodes a member of the trinucleotide repeat containing 6 protein family. The protein functions in post-transcriptional gene silencing through the RNA interference (RNAi) and microRNA pathways. The protein associates with messenger RNAs and Argonaute proteins in cytoplasmic bodies known as GW-bodies or P-bodies. Inhibiting expression of this gene delocalizes other GW-body proteins and impairs RNAi and microRNA-induced gene silencing.

Description

TNRC6A 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

Q8NDV7

Purification

Affinity Purification

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 950-1050 of human TNRC6A (NP_055309.2).

Storage

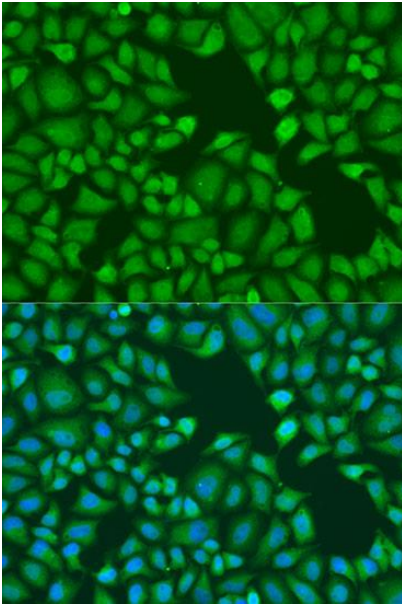
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

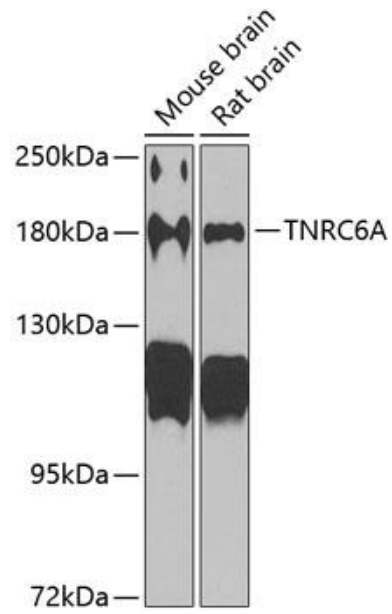
TNRC6A; CAGH26; GW1; GW182; TNRC6; trinucleotide repeat containing 6A

Application

WB, IF; Recommended dilution: WB 1:500 - 1:1000, IF 1:50 - 1:100



Immunofluorescence analysis of U2OS cells using TNRC6A Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.



Western blot analysis of extracts of various cell lines, using TNRC6A 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.