
TWNK Polyclonal Antibody

(Catalog #A69095)

Background

Involved in mitochondrial DNA (mtDNA) metabolism. Could function as an adenine nucleotide-dependent DNA helicase. Function inferred to be critical for lifetime maintenance of mtDNA integrity. In vitro, forms in combination with POLG, a processive replication machinery, which can use double-stranded DNA (dsDNA) as template to synthesize single-stranded DNA (ssDNA) molecules. May be a key regulator of mtDNA copy number in mammals.

Description

TWNK Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300. 50% Glycerol, 0.01M PBS, pH 7.4.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q96RR1

Purification

Protein G purified

Immunogen

Recombinant Human Twinkle protein, mitochondrial protein (559-684AA)

Storage

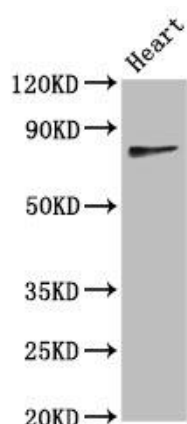
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Twinkle protein, mitochondrial, Progressive external ophthalmoplegia 1 protein, T7 gp4-like protein with intramitochondrial nucleoid localization, T7-like mitochondrial DNA helicase, Twinkle mtDNA helicase, TWNK

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500 - 1:5000, IHC:1:500 - 1:1000



Western Blot

Positive WB detected in: Mouse heart tissue

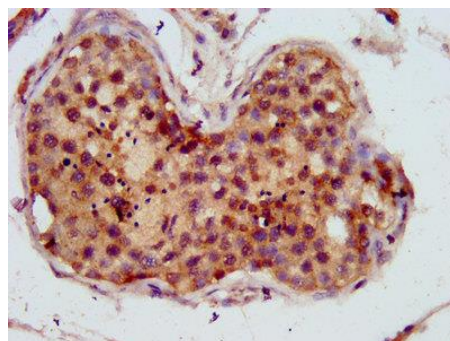
All lanes: TWNK antibody at 4.4ug/ml

Secondary

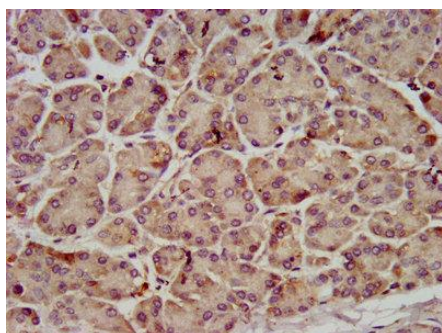
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 78, 67, 61 KDa

Observed band size: 78 KDa



IHC image of TWNK Polyclonal Antibody diluted at 1:800 and staining in paraffin-embedded human testis tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



IHC image of TWNK Polyclonal Antibody diluted at 1:800 and staining in paraffin-embedded human pancreatic tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.