
POLG2 Polyclonal Antibody

(Catalog # A70661)

Background

This gene encodes the processivity subunit of the mitochondrial DNA polymerase gamma. The encoded protein forms a heterotrimer containing one catalytic subunit and two processivity subunits. This protein enhances DNA binding and promotes processive DNA synthesis. Mutations in this gene result in autosomal dominant progressive external ophthalmoplegia with mitochondrial DNA deletions.

Description

POLG2 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

Q9UHN1

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 216-485 of human POLG2 (NP_009146.2).

Storage

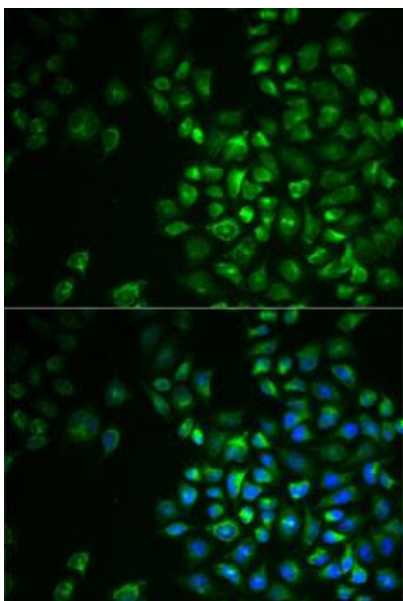
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

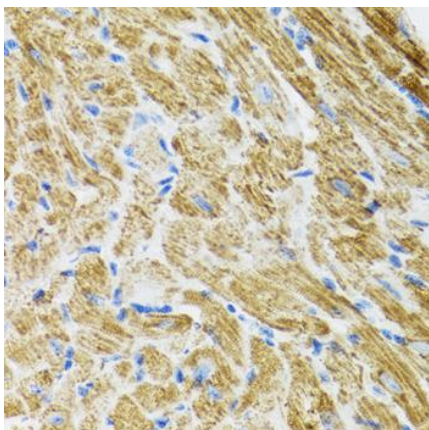
POLG2; HP55; MTPOLB; PEOA4; POLB; POLG-BETA; POLGB; DNA polymerase gamma 2, accessory subunit

Application

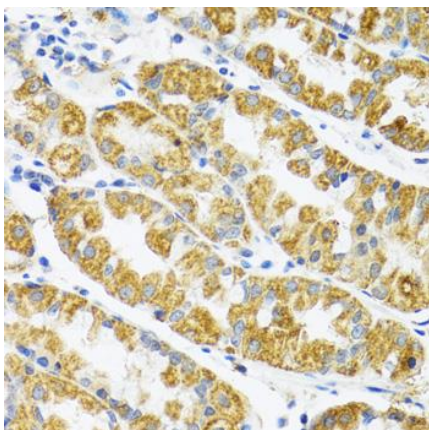
WB, IHC, IF; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:10 - 1:100



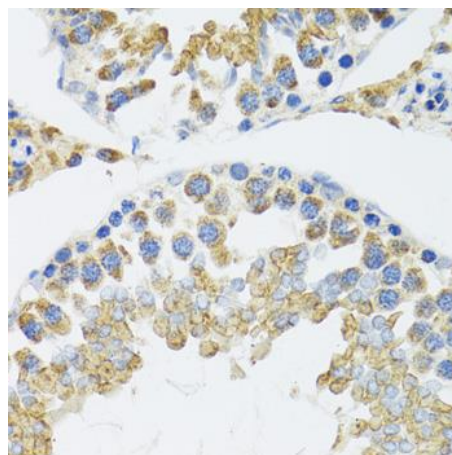
Immunofluorescence analysis of MCF7 cells using POLG2 Polyclonal Antibody. Blue: DAPI for nuclear staining.



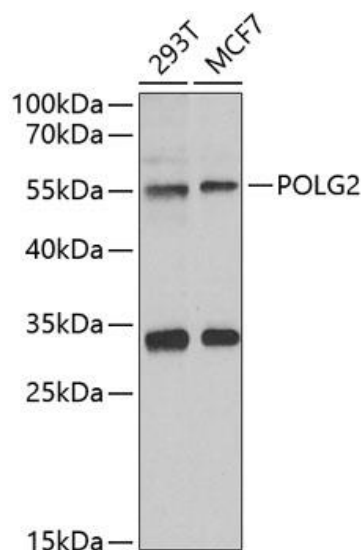
Immunohistochemistry of paraffin-embedded rat heart using POLG2 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human stomach using POLG2 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded mouse testis using POLG2 Polyclonal Antibody at dilution of 1:100 (40x lens).



Western blot analysis of extracts of various cell lines, using POLG2 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.