

---

## GAPDH Polyclonal Antibody

(Catalog # A70921)

---

### Background

This gene encodes a member of the glyceraldehyde-3-phosphate dehydrogenase protein family. The encoded protein has been identified as a moonlighting protein based on its ability to perform mechanistically distinct functions. The product of this gene catalyzes an important energy-yielding step in carbohydrate metabolism, the reversible oxidative phosphorylation of glyceraldehyde-3-phosphate in the presence of inorganic phosphate and nicotinamide adenine dinucleotide (NAD). The encoded protein has additionally been identified to have uracil DNA glycosylase activity in the nucleus. Also, this protein contains a peptide that has antimicrobial activity against *E. coli*, *P. aeruginosa*, and *C. albicans*. Studies of a similar protein in mouse have assigned a variety of additional functions including nitrosylation of nuclear proteins, the regulation of mRNA stability, and acting as a transferrin receptor on the cell surface of macrophage. Many pseudogenes similar to this locus are present in the human genome. Alternative splicing results in multiple transcript variants.

### Description

GAPDH Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: PBS with 0.01% thiomersal, 50% glycerol, pH7.3

### Specificity

Human, Mouse, Rat

### Isotype

IgG

### Uniprot ID

P04406

### Purification

Affinity Purification

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-335 of human GAPDH (NP\_002037.2).

### Storage

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

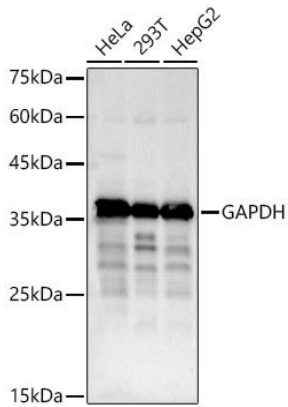
### Alternative Names

GAPDH; G3PD; GAPD; HEL-S-162eP; glyceraldehyde-3-phosphate dehydrogenase

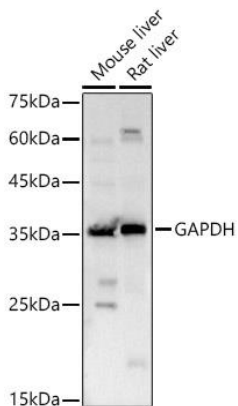
### Application

WB, IHC, IF, IP; Recommended dilution: WB 1:10000 - 1:30000, IHC-P 1:50 - 1:200, IF/ICC 1:50 - 1:200, IP 1:2000 - 1:20000

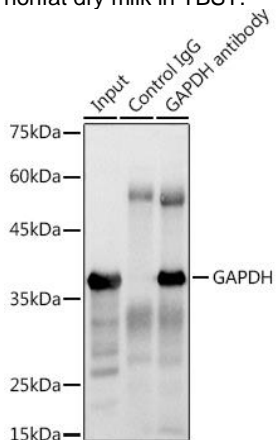
---



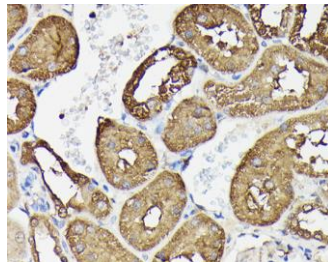
WB analysis of extracts of various cell lines, using GAPDH pAb at 1:30000 dilution. Secondary Ab: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug/lane. Blocking buffer: 3% nonfat dry milk in TBST.



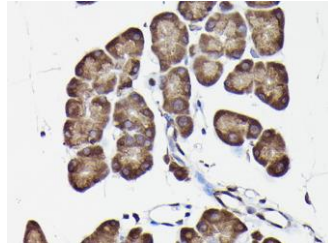
WB of extracts of various cell lines, using GAPDH pAb at 1:30000 dilution. Secondary Ab: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug/lane. Blocking buffer: 3% nonfat dry milk in TBST.



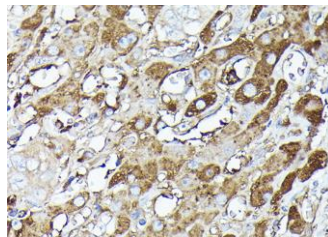
IP analysis of 300ug extracts of HeLa cells using 3ug pAb. WB dilution of 1:20000.



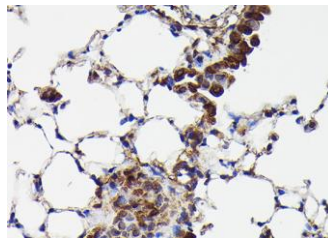
IHC of paraffin-embedded rat kidney using GAPDH pAb at dilution of 1:100 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



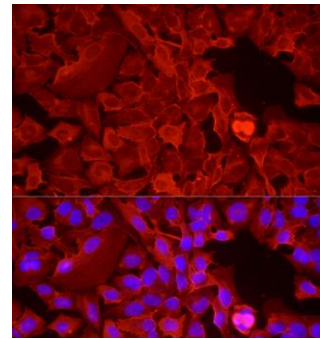
IHC of paraffin-embedded rat pancreas using GAPDH pAb at dilution of 1:100 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



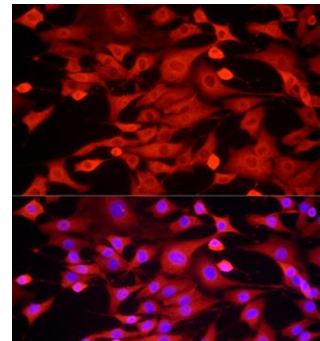
IHC of paraffin-embedded human liver cancer using GAPDH PAb at dilution of 1:100 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



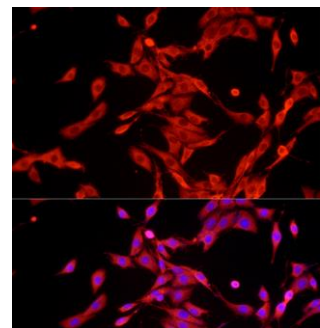
IHC of paraffin-embedded mouse lung using GAPDH pAb at dilution of 1:100 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



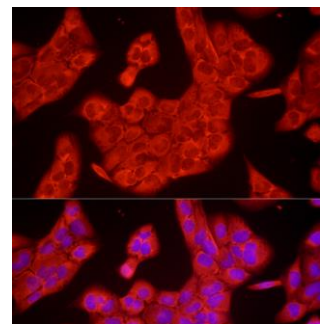
IF analysis of HeLa cells using GAPDH pAb at dilution of 1:20 (40x lens). Blue: DAPI for nuclear staining.



IF analysis of NIH/3T3 cells using GAPDH pAb at dilution of 1:20 (40x lens). Blue: DAPI for nuclear staining.



IF analysis of PC-12 cells using GAPDH pAb at dilution of 1:20 (40x lens). Blue: DAPI for nuclear staining.



IF analysis of U2OS cells using GAPDH pAb at dilution of 1:20 (40x lens). Blue: DAPI for nuclear staining.