

## ANXA2 Polyclonal Antibody

(Catalog # A71015)

### Background

This gene encodes a member of the annexin family. Members of this calcium-dependent phospholipid-binding protein family play a role in the regulation of cellular growth and in signal transduction pathways. This protein functions as an autocrine factor which heightens osteoclast formation and bone resorption. This gene has three pseudogenes located on chromosomes 4, 9 and 10, respectively. Multiple alternatively spliced transcript variants encoding different isoforms have been found for this gene.

### Description

ANXA2 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

P07355

### Purification

Affinity Purification

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-339 of human ANXA2 (NP\_004030.1).

### Storage

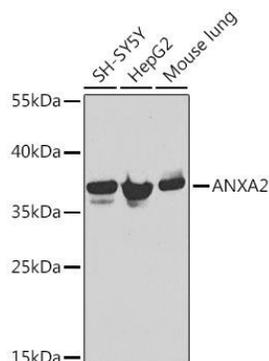
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

### Alternative Names

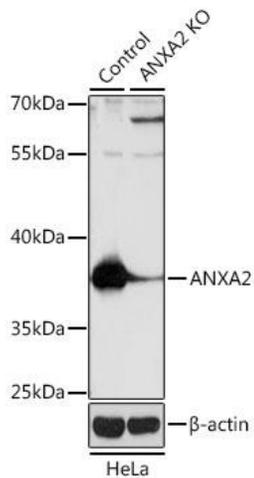
ANXA2; ANX2; ANX2L4; CAL1H; HEL-S-270; LIP2; LPC2; LPC2D; P36; PAP-IV; annexin A2

### Application

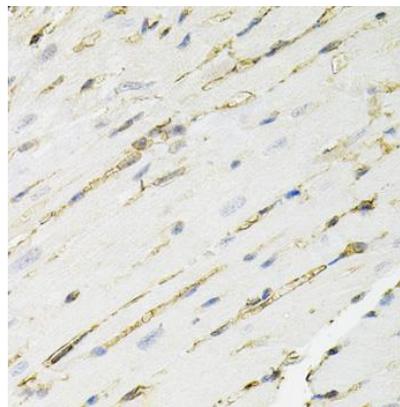
WB, IHC, IF; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:20 - 1:50



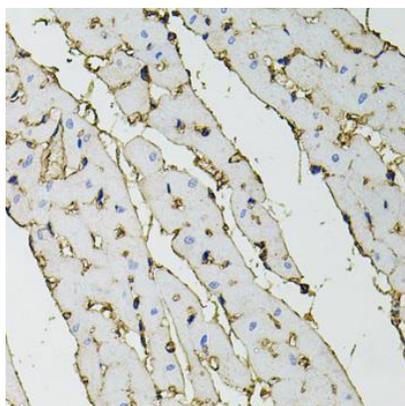
Western blot analysis of extracts of various cell lines, using ANXA2 Polyclonal Antibody at 1:1000 dilution. Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST.



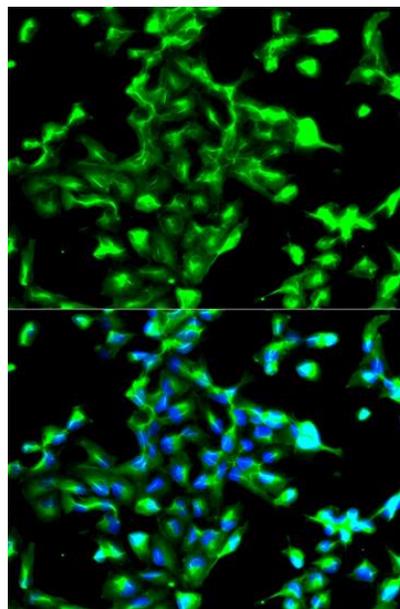
Western blot analysis of extracts from normal (control) and ANXA2 knockout (KO) HeLa cells, using ANXA2 Polyclonal Antibody at 1:1000 dilution.  
 Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.  
 Lysates/proteins: 25ug per lane.  
 Blocking buffer: 3% nonfat dry milk in TBST.  
 Exposure time: 3s.



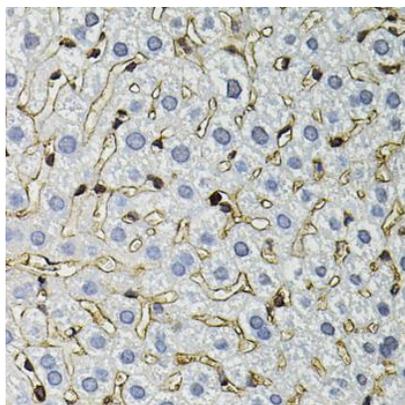
Immunohistochemistry of paraffin-embedded mouse heart using ANXA2 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded rat heart using ANXA2 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunofluorescence analysis of A549 cells using ANXA2 Polyclonal Antibody. Blue: DAPI for nuclear staining.



Immunohistochemistry of paraffin-embedded mouse liver using ANXA2 Polyclonal Antibody at dilution of 1:100 (40x lens).