

## Anti-ATG13

Rabbit Polyclonal Antibody

**Catalog # A113-30G**

Lot # O2361-12

### Cited Applications

E, WB, IHC, IF

*Ideal working dilutions for each application should be empirically determined by the investigator.*

### Specificity

Recognizes the ATG13 protein

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

ATG13 antibody was raised against a 15 amino acid synthetic peptide near the carboxy terminus of human ATG13

### Formulation

PBS + 0.02% sodium azide.

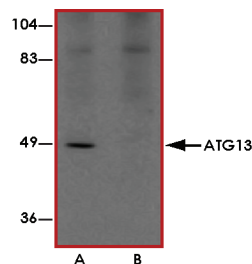
### Stability

1yr at -20°C from date of shipment.

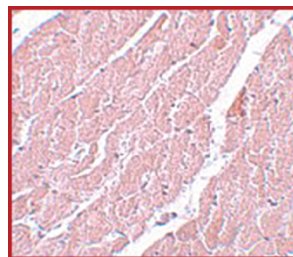
### References

1. Gozuacik D et al. Autophagy as a cell death and tumor suppressor mechanism. *Oncogene*2004; 23:2891-906.
2. Kisen GO, et al: Reduced autophagic activity in primary rat hepatocellular carcinoma and ascites hepatoma cells. *Carcinogenesis*1993; 14:2501-5.
3. Kamada Y et al: Tor-mediated induction of autophagy via Apg1 protein kinase complex. *J. Cell. Biol.*2000; 150:1507-13.
4. Jung CH et al: ULK-Atg13-FIP200 complexes mediate mTOR signaling to the autophagy machinery. *Mol. Biol. Cell* 2009; 20:1992-2003.

### Sample Data



Western blot analysis of ATG13 in rat heart tissue lysate with ATG13 antibody at 1 µg/mL in (A) the absence and (B) the presence of blocking peptide.



Immunohistochemistry of ATG13 in mouse heart with ATG13 antibody at 5 µg/mL.

## Anti-ATG13

Rabbit Polyclonal Antibody

Catalog Number

A113-30G

Specific Lot Number

O2361-12

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**