

---

## SELENBP1 Polyclonal Antibody

(Catalog # A73262)

---

### Background

This gene encodes a member of the selenium-binding protein family. Selenium is an essential nutrient that exhibits potent anticarcinogenic properties, and deficiency of selenium may cause certain neurologic diseases. The effects of selenium in preventing cancer and neurologic diseases may be mediated by selenium-binding proteins, and decreased expression of this gene may be associated with several types of cancer. The encoded protein may play a selenium-dependent role in ubiquitination/de-ubiquitination-mediated protein degradation. Alternatively spliced transcript variants encoding multiple isoforms have been observed for this gene.

### Description

SELENBP1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

Q13228

### Purification

Affinity Purified

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 243-472 of human SELENBP1 (NP\_003935.2).

### Storage

Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

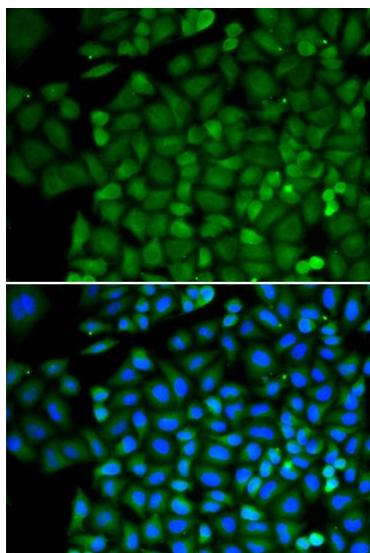
### Alternative Names

SELENBP1; HEL-S-134P; LPSB; SBP56; SP56; hSBP

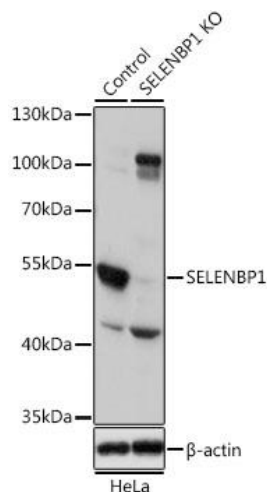
### Application

WB, IF; Recommended dilution: WB 1:500 - 1:2000, IF/ICC 1:50 - 1:200

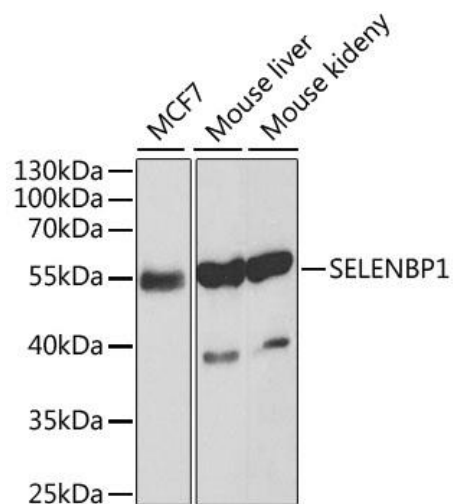
---



Immunofluorescence analysis of HeLa cells using SELENBP1 Polyclonal Antibody. Blue: DAPI for nuclear staining.



Western blot analysis of extracts from normal (control) and SELENBP1 knockout (KO) HeLa cells, using SELENBP1 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 of various cell lines, using SELENBP1 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.