

---

## CD59 Polyclonal Antibody

(Catalog # A71164)

---

### Background

This gene encodes a cell surface glycoprotein that regulates complement-mediated cell lysis, and it is involved in lymphocyte signal transduction. This protein is a potent inhibitor of the complement membrane attack complex, whereby it binds complement C8 and/or C9 during the assembly of this complex, thereby inhibiting the incorporation of multiple copies of C9 into the complex, which is necessary for osmolytic pore formation. This protein also plays a role in signal transduction pathways in the activation of T cells. Mutations in this gene cause CD59 deficiency, a disease resulting in hemolytic anemia and thrombosis, and which causes cerebral infarction. Multiple alternatively spliced transcript variants, which encode the same protein, have been identified for this gene.

### Description

CD59 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

P13987

### Purification

Affinity Purification

### Immunogen

Recombinant protein of human CD59

### Storage

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

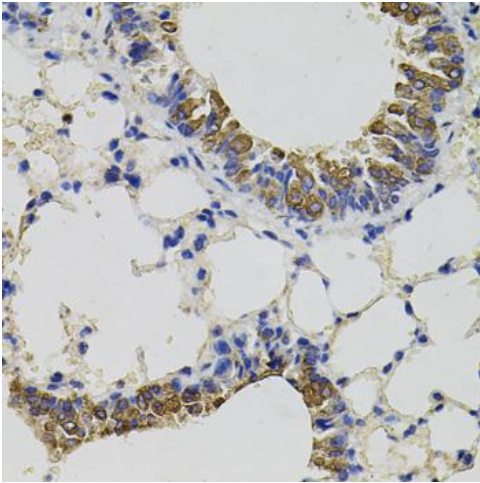
### Alternative Names

CD59; 16.3A5; 1F5; EJ16; EJ30; EL32; G344; HRF-20; HRF20; MAC-IP; MACIF; MEM43; MIC11; MIN1; MIN2; MIN3; MIRL; MSK21; p18-20; CD59 glycoprotein

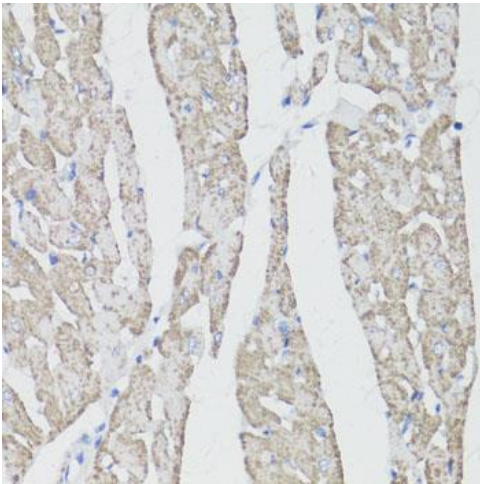
### Application

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

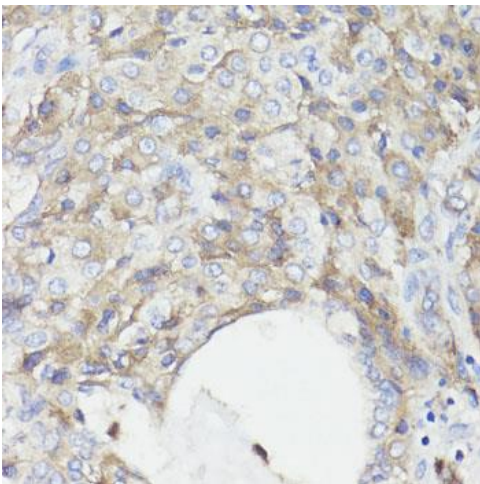
---



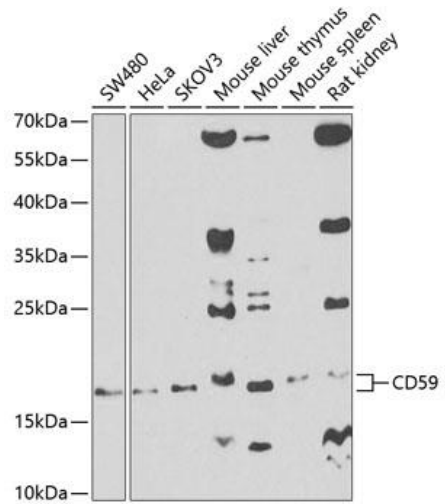
Immunohistochemistry of paraffin-embedded mouse lung using CD59 Polyclonal Antibody (40x lens).



Immunohistochemistry of paraffin-embedded rat heart using CD59 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human lung cancer using CD59 Polyclonal Antibody at dilution of 1:100 (40x lens).



Western blot analysis of extracts of various cell lines, using CD59 Polyclonal Antibody at 1:500 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: 60s.