

---

## C5 Polyclonal Antibody

(Catalog #A71922)

---

### Background

This gene encodes a component of the complement system, a part of the innate immune system that plays an important role in inflammation, host homeostasis, and host defense against pathogens. The encoded preproprotein is proteolytically processed to generate multiple protein products, including the C5 alpha chain, C5 beta chain, C5a anaphylatoxin and C5b. The C5 protein is comprised of the C5 alpha and beta chains, which are linked by a disulfide bridge. Cleavage of the alpha chain by a convertase enzyme results in the formation of the C5a anaphylatoxin, which possesses potent spasmogenic and chemotactic activity, and the C5b macromolecular cleavage product, a subunit of the membrane attack complex (MAC). Mutations in this gene cause complement component 5 deficiency, a disease characterized by recurrent bacterial infections. Alternative splicing results in multiple transcript variants.

### Description

C5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: PBS with 0.09% sodium azide, 50% glycerol, pH7.3.

### Specificity

Human, Mouse, Rat

### Isotype

IgG

### Uniprot ID

P01031

### Purification

Affinity Purified

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1567-1676 of human C5 (NP\_001726.2)

### Storage

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

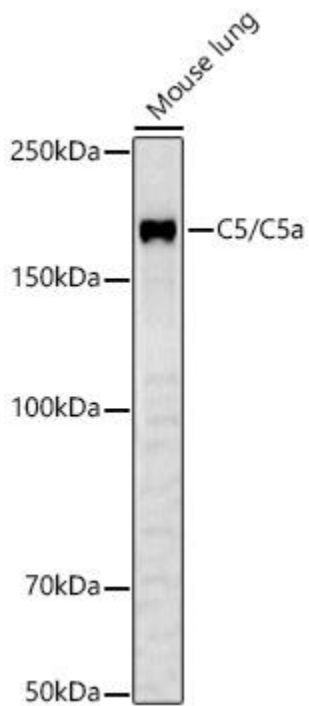
### Alternative Names

C5; C5D; C5a; C5b; CPAMD4; ECLZB; complement C5

### Application

WB, IF, ELISA; Recommended dilution: WB:1:500-1:1000, ELISA - recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.

---



Western blot analysis of lysates from Mouse lung, using C5 Polyclonal Antibody at 1:600 dilution.  
Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution.  
Lysates/proteins: 25µg per lane.  
Blocking buffer: 3% nonfat dry milk in TBST.  
Exposure time: 90s.