

---

## CD14 Polyclonal Antibody

(Catalog #A64104)

---

### Background

Coreceptor for bacterial lipopolysaccharide (PubMed:1698311, PubMed:23264655). In concert with LBP, binds to monomeric lipopolysaccharide and delivers it to the LY96/TLR4 complex, thereby mediating the innate immune response to bacterial lipopolysaccharide (LPS) (PubMed:20133493, PubMed:23264655). Acts via MyD88, TIRAP and TRAF6, leading to NF-kappa-B activation, cytokine secretion and the inflammatory response. (PubMed:8612135).

### Description

CD14 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

P08571

### Purification

Protein G purified

### Immunogen

Recombinant Human Monocyte differentiation antigen CD14 protein (20-345AA)

### Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

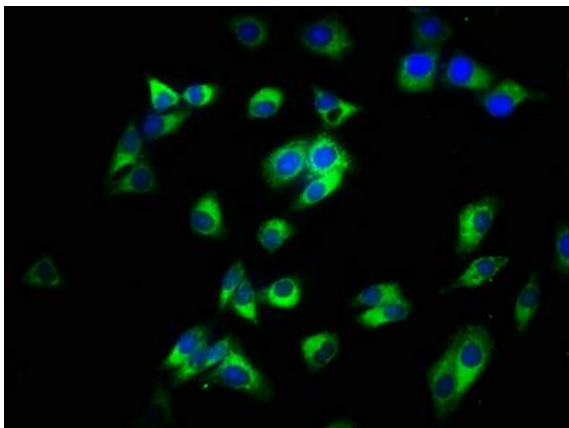
### Alternative Names

Monocyte differentiation antigen CD141, Myeloid cell-specific leucine-rich glycoprotein, CD14, CD14

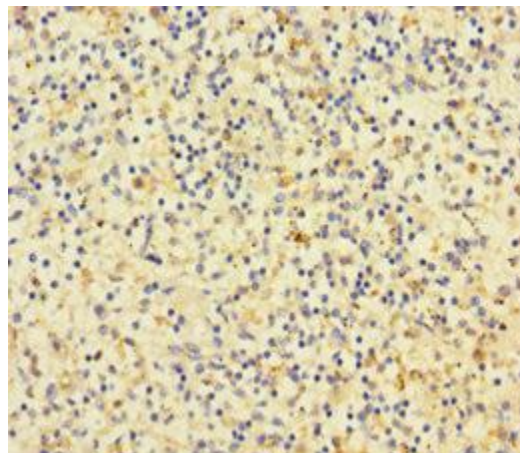
### Application

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

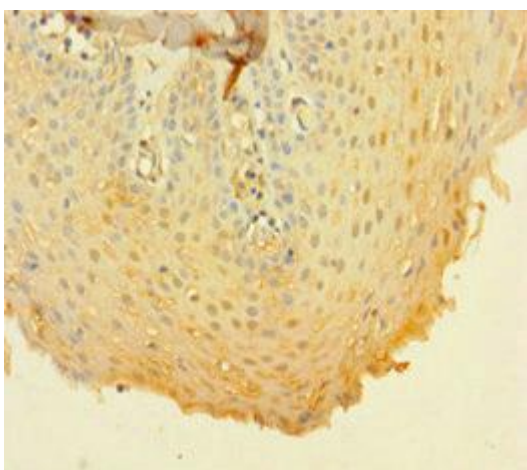
---



Immunofluorescence staining of HepG2 cells with CD14 Antibody at 1:200, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).



Immunohistochemistry of paraffin-embedded human spleen tissue using CD14 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human tonsil tissue using CD14 Antibody at dilution of 1:100