
CD14 Polyclonal Antibody

(Catalog #A72619)

Background

The protein encoded by this gene is a surface antigen that is preferentially expressed on monocytes/macrophages. It cooperates with other proteins to mediate the innate immune response to bacterial lipopolysaccharide. Alternative splicing results in multiple transcript variants encoding the same protein.

Description

CD14 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.05% proclin300, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P08571

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 20-330 of human CD14 (NP_000582.1).

Storage

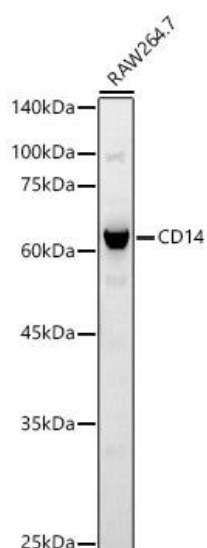
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

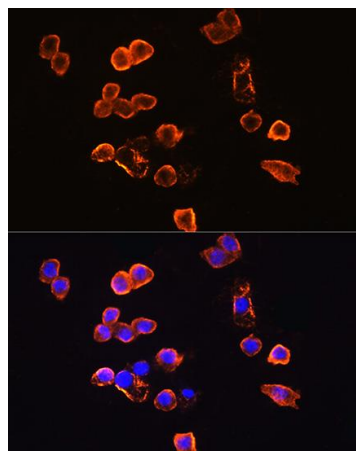
CD14; CD14 molecule

Application

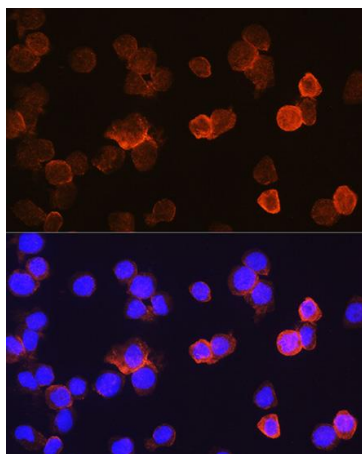
ELISA, WB, IF; Recommended dilution: WB: 1:500-1:1000, IF:1:50-1:200



Western blot analysis of extracts of RAW264.7 cells, using CD14 Polyclonal Antibody at 1:600 dilution. Secondary antibody: HRP 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: 30s.



Immunofluorescence analysis of RAW264.7 cells using CD14 Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.



Immunofluorescence analysis of THP-1 cells using CD14 Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.