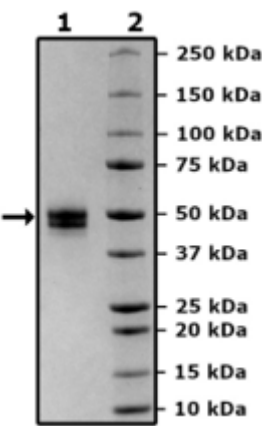


Product Information

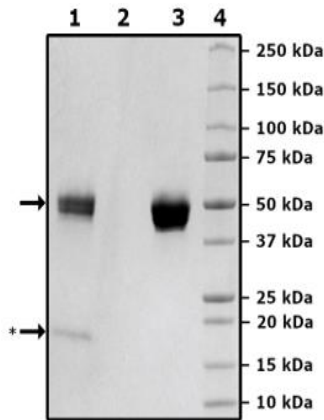
Description:	Recombinant human CD14 molecule, encompassing amino acids 20-336. This construct contains a C-terminal Avi-Tag™ followed by a His-tag (6xHis). The protein was affinity purified.
Background:	CD14 (Cluster of Differentiation-14) is a surface antigen expressed predominantly by monocytes and macrophages. It is considered a patterned recognition receptor (PRR), which recognizes pathogen markers. Specifically, CD14 acts as a co-receptor for lipopolysaccharides (LPS) found in the outer membrane of gram-negative bacteria. Through this interaction, CD14 supports initiation of the innate immune response during bacterial infection. CD14 also contributes toward immune responses to viral pathogens such as human respiratory syncytial virus (RSV) and may amplify the inflammatory response observed in severe cases of SARS-CoV-2 infection (COVID-19). Monoclonal antibodies against CD14 are currently undergoing clinical trial as a candidate treatment to limit severe inflammatory responses common in patients hospitalized with COVID-19.
Species:	Human
Construct:	CD14 (20-336-Avi-His)-(Biotin)
Concentration:	2.67 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	37 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_000591
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown



- 1. Beads
- 2. Flow thru
- 3. Control
- 4. Standards

* Avidin from beads.