# CD36 (human), (recombinant)

CD36 binds to collagen, thrombospondin, anionic phospholipids and oxidized low-density lipoprotein (oxLDL). It also may function as a cell adhesion molecule. It directly mediates cytoadherence of Plasmodium falciparum parasitized erythrocytes. It also binds long chain fatty acids and may function in the transport and/or as a regulator of fatty acid transport. It also serves as a receptor for thombospondins, THBS1 AND THBS2, mediating their antiangiogenic effects. As a co-receptor for TLR4-TLR6 heterodimer, it promotes inflammation in monocytes/macrophages. Upon ligand binding, such as oxLDL or amyloid-beta 42, it rapidly induces the formation of a heterodimer of TLR4 and TLR6, which is internalized and triggers inflammatory response, leading to NF-κ-B-dependent production of CXCL1, CXCL2 and CCL9 cytokines, via MyD88 signaling pathway, and CCL5 cytokine, via TICAM1 signaling pathway, as well as IL-1B secretion.

### **Ordering Information**

Order Online »

ENZ-PRT170-0010	10µg
-----------------	------

Manuals, SDS & CofA

**View Online »** 

### **Handling & Storage**

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name SCARB3, GP3B, GP4, Platelet glycoprotein 4

Formulation Lyophilized from 0.22µm filtered solution in PBS, pH 7.4.

Generally 5-8% mannitol or trehalose is added as a

protectant before lyophilization.

**MW** ~47.5 kDa

Purity ≥95% (SDS-PAGE)

**Source** Produced in HEK 293 cells.

UniProt ID P16671

Last modified: May 29, 2024

