

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 thrombospondins, 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.

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com