

Product Information

Description:	Recombinant human CD36 (also known as Platelet glycoprotein 4), encompassing amino acids 30-439. The C-terminus of this construct is fused with the Fc fragment of a human IgG1 followed by an Avi-Tag™. The recombinant protein was affinity purified.
Background:	CD36 (cluster of differentiation 36) is also known as platelet glycoprotein 4, fatty acid translocase (FAT), scavenger receptor class B member 3 (SCARB3), and glycoprotein 88 (GP88). It is expressed in many cell types including erythrocytes, monocytes, differentiated adipocytes, skeletal muscle cells, mammary and intestinal epithelial cells, and endothelial cells. CD36 is one of the major glycoproteins present at the surface of platelets. CD36 has intricate biological functions, depending on the cell type and ligand. CD36 is involved in various diseases that involve the immune and vascular systems and in diseases for which lipid metabolism is important such as diabetes and obesity, inflammation, atherosclerosis, heart disease, and Alzheimer's disease. In addition, since lipids are an important energy source for tumor cells, CD36 is associated with cancer. Upregulated CD36 expression has been observed in multiple types of cancer and is correlated with poor clinical outcomes.
Species:	Human
Construct:	CD36 (30-439-Fc(IgG1)-Avi)
Concentration:	0.45 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:	75 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_000072.3
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

