

# PRODUCT INFORMATION



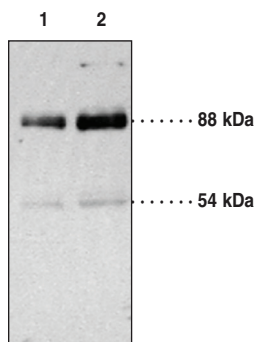
## CD36 Polyclonal Antibody

Item No. 100011

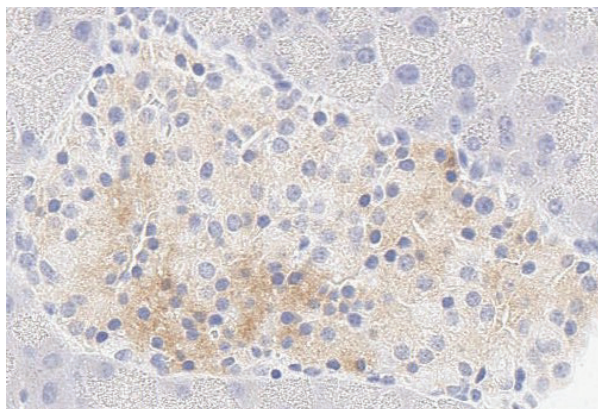
### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl of peptide affinity-purified antibody.
<b>Synonyms:</b>	GPIIb, GPIV, Hexarelin Receptor <sup>1-3</sup> , oxLDL Receptor, Thrombospondin Receptor
<b>Immunogen:</b>	Synthetic peptide from an internal region of human CD36
<b>Species Reactivity:</b>	(+) Human, mouse, and rat CD36
<b>Uniprot No.:</b>	P16671
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	As supplied, 1 year from the QC date provided on the Certificate of Analysis, when stored properly
<b>Storage Buffer:</b>	TBS, pH 7.4, containing 50% glycerol, 0.1% BSA and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Application:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Lane 1: Human platelet lysate (15 µg)  
Lane 2: Human platelet lysate (30 µg)



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) mouse small intestine tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with CD36 Polyclonal Antibody (Item No. 100011) at a 1:200 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/29/2019

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

CD36 is a type-B scavenger receptor that is necessary for the formation of foam cells in atherosclerotic lesions.<sup>4-7</sup> This membrane glycoprotein can internalize fatty acids which activate PPAR $\gamma$  and stimulate further expression of CD36.<sup>8,9</sup> This positive feedback loop combined with the mouse CD36 knock-out studies reinforces the importance of CD36 in lipid metabolism. Additionally CD36 is needed for the phagocytosis of *P. falciparum* infected erythrocytes, retinal pigment epithelial cell photoreceptor fragments, and post-apoptotic monocytes and neutrophils.<sup>1,10-12</sup> CD36 expression has been monitored during hematopoietic cell differentiation and may be an indicator of tumor spreading in lymphocytic leukemia.<sup>13</sup> This transmembrane glycoprotein is detected by immunoblotting at 88 or 54 kDa depending on the degree of glycosylation of the receptor in your sample. Positive controls include adipose and heart tissue, platelets, and macrophages.

## References

---

1. Greenwalt, D.E., Lipsky, R.H., Ockenhouse, C.F., et al. *Blood* **80**(5), 1105-1115 (1992).
2. Bodart, V., Febbraio, M., Demers, A., et al. *Circ. Res.* **90**, 844-849 (2002).
3. Pearce, S.F.A., Roy, P., Nicholson, A.C., et al. *J. Biol. Chem.* **273**(52), 34875-34881 (1998).
4. Febbraio, M., Abumrad, N.A., Hajjar, D.P., et al. *J. Biol. Chem.* **274**, 19055-19062 (1999).
5. Podrez, E.A., Febbraio, M., Sheibani, N., et al. *J. Clin. Invest.* **105**(8), 1095-1108 (2000).
6. Febbraio, M., Podrez, E.A., Smith, J.D., et al. *J. Clin. Invest.* **105**(8), 1049-1056 (2000).
7. Febbraio, M., Hajjar, D.P., and Silverstein, R.L. *J. Clin. Invest.* **108**(6), 785-791 (2001).
8. Han, J., Hajjar, D.P., Febbraio, M., et al. *J. Biol. Chem.* **272**(34), 21654-21659 (1997).
9. Tontonoz, P., Nagy, L., Alvarez, J.G.A., et al. *Cell* **93**, 241-252 (1998).
10. Finnemann, S.C. and Silverstein, R.L. *J. Exp. Med.* **194**(9), 1289-1298 (2001).
11. Navazo, M.D.P., Daviet, L., Savill, J., et al. *J. Biol. Chem.* **271**(26), 15381-15385 (1996).
12. Jiménez, B., Volpert, O.V., Crawford, S.E., et al. *Nature Med.* **6**(1), 41-48 (2000).
13. Rutella, S., Rumi, C., Puggioni, P., et al. *Haematologica* **84**, 419-424 (1999).

## CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM