

PRODUCT INFORMATION



CD36 Monoclonal Antibody (Clone JC63.1)

Item No. 188150

Overview and Properties

Contents:	This vial contains 100 µg of protein L purified IgA.
Synonyms:	GPIIb, GPIV, Hexarelin Receptor, oxLDL Receptor, Thrombospondin Receptor ¹⁻³
Immunogen:	Adenovirus expressing full-length recombinant mouse CD36
Species Reactivity:	(+) Human, mouse, and rat CD36
Uniprot No.:	Q08857
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, with 50% glycerol, and 0.02% sodium azide
Clone:	JC63.1
Host:	CD36 null mouse
Isotype:	IgA
Application:	Flow cytometry; the recommended starting concentration is 0.5 µg/ml. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image

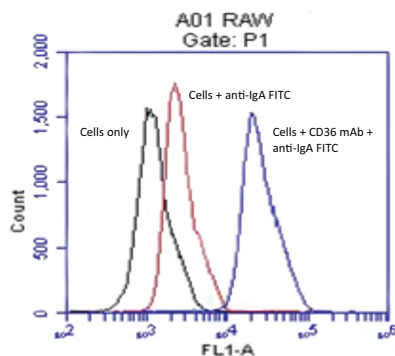


Figure 1: Flow cytometry with RAW 264.7 cells

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, 10/23/2019

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 and thereby atherosclerotic lesions.¹⁻⁴ This membrane glycoprotein can internalize fatty acids which activate PPAR γ and stimulate further expression of CD36.^{5,6} This positive feedback loop combined with the murine 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.⁷⁻¹⁰ CD36 expression has been monitored during hematopoietic cell differentiation and may be an indicator of tumor spreading in lymphocytic leukemia.¹¹ Positive controls include cells from adipose and heart tissue, platelets, and macrophages. Cayman Chemical's CD36 Monoclonal Antibody (Clone JC63.1) is recommended for flow cytometry on CD36-containing cells of mouse, rat, and human origin. The antibody can also be used to antagonize CD36-mediated cellular activity. Functional blocking inhibits CD36 binding of oxLDL (macrophages) and prevents foam cell formation of macrophages.

References

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