

Product Information



CD36 Monoclonal Antibody (Clone JC63.1)

Item No. 188150

Synonyms:	GPIIb, GPIV, Hexarelin Receptor, oxLDL Receptor, Thrombospondin Receptor ¹⁻³
Contents:	Purified mouse anti-CD36 IgA (1 mg/ml) in PBS, pH 7.2, containing 50% glycerol and 0.02% sodium azide.
Host:	CD36 null mouse, clone JC63.1
Antigen:	Adenovirus expressing full-length recombinant mouse CD36 ⁴
Isotype:	IgA
Cross Reactivity:	(+) Mouse, rat, and human CD36
Stability:	≥1 year at -20°C
Application:	Flow cytometry: 0.5 µg/ml (1:2,000); Functional blocking: 2 µg/ml (1:500) inhibits CD36 binding of ox-LDL (macrophages) and prevents foam cell formation of macrophages.

CD36 is a type-B scavenger receptor that is necessary for the formation of foam cells and thereby atherosclerotic lesions.⁵⁻⁸ This membrane glycoprotein can internalize fatty acids which activate PPAR γ and stimulate further expression of CD36.^{9,10} 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.^{1,4,11,12} CD36 expression has been monitored during hematopoietic cell differentiation and may be an indicator of tumor spreading in lymphocytic leukemia.¹³ Positive controls include cells from adipose and heart tissue, platelets, and macrophages.

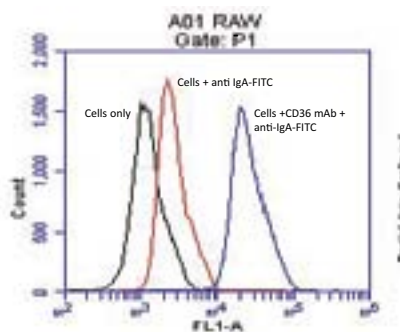


Figure 1: Flow Cytometry with RAW 264.7 cells

References

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

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/188150

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/20/2011

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com