

# PRODUCT INFORMATION



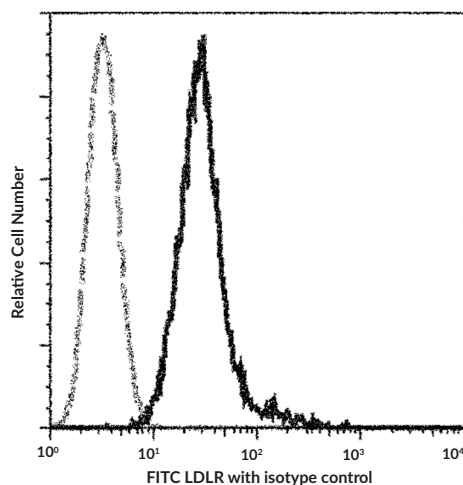
## LDL Receptor Rabbit Monoclonal Antibody (FITC) (Clone 032)

Item No. 37077

### Overview and Properties

<b>Contents:</b>	This vial contains protein A-affinity purified monoclonal antibody.
<b>Synonyms:</b>	LDLR, Low Density Lipoprotein Receptor
<b>Immunogen:</b>	Recombinant mouse LDLR
<b>Cross Reactivity:</b>	(+) LDLR
<b>Species Reactivity:</b>	(+) Mouse
<b>Uniprot No.:</b>	P35951
<b>Form:</b>	Liquid
<b>Storage:</b>	2-8°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS with 0.5% BSA and 0.09% sodium azide
<b>Concentration:</b>	0.1 mg/ml
<b>Clone:</b>	032
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Application:</b>	Flow cytometry (FC); the optimal working concentration/dilution should be determined empirically.

### Image



Profile of anti-LDLR reactivity on Raw264.7 cells analyzed by flow cytometry. Cells should be Fc-blocked by treatment with Mouse BD Fc Block™ purified LDL Receptor Rabbit Monoclonal Antibody (FITC) (Clone 032) prior to staining, washed, then stained with FITC Rabbit anti-LDLR.

**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, 02/07/2024

**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

---

LDL receptor (LDLR) is a cell surface glycoprotein that scavenges LDL from the blood and regulates plasma LDL levels.<sup>1</sup> It is composed of an N-terminal signal sequence, a ligand-binding domain, an EGF precursor homology domain, an O-linked glycosylation domain, a transmembrane region, and a C-terminal cytoplasmic tail. LDLR is primarily expressed in the liver but is also found in the adrenal cortex.<sup>2</sup> It mediates the endocytosis of LDL by binding to apolipoprotein E (ApoE) or ApoB on the LDL surface, thereby supplying cholesterol to cells.<sup>1</sup> Protein levels of LDLR are decreased in HepG2 cells expressing proprotein convertase subtilisin kexin 9 (PCSK9).<sup>3</sup> Knockout of *Ldlr* increases plasma levels of cholesterol and triglycerides and induces the formation of atherosclerotic lesions in mice.<sup>4</sup> Mutations in *LDLR* are associated with familial hypercholesterolemia.<sup>5</sup> Cayman's LDL Receptor Rabbit Monoclonal Antibody (FITC) is composed of an LDLR monoclonal antibody conjugated to fluorescein isothiocyanate (FITC) (Clone 032) and can be used for flow cytometry (FC).

## References

---

1. Goldstein, J.L., Brown, M.S., Anderson, R.G.W., *et al.* Receptor-mediated endocytosis: Concepts emerging from the LDL receptor system. *Annu. Rev. Cell Biol.* **1**, 1-39 (1985).
2. Rudling, M.J., Reihner, E., Einarsson, K., *et al.* Low density lipoprotein receptor-binding activity in human tissues: Quantitative importance of hepatic receptors and evidence for regulation of their expression *in vivo*. *Proc. Natl. Acad. Sci. USA* **87**(9), 3469-3473 (1990).
3. Benjannet, S., Rhainds, D., Essalmani, R., *et al.* NARC-1/PCSK9 and its natural mutants. Zymogen cleavage and effects on the low density lipoprotein (LDL) receptor and LDL cholesterol. *J. Biol. Chem.* **279**(47), 48865-48875 (2004).
4. Praticò, D., Tillmann, C., Zhang, Z.B., *et al.* Acceleration of atherogenesis by COX-1-dependent prostanoid formation in low density lipoprotein receptor knockout mice. *Proc. Natl. Acad. Sci. USA* **98**(6), 3358-3363 (2001).
5. Austin, M.A., Hutter, C.M., Zimmern, R.L., *et al.* Genetic causes of monogenic heterozygous familial hypercholesterolemia: A HuGE prevalence review. *Am. J. Epidemiol.* **160**(5), 407-420 (2004).

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