

PRODUCT INFORMATION



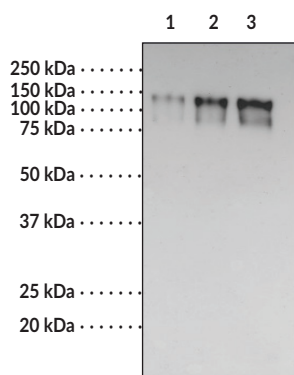
LDL Receptor Polyclonal Antibody - Biotinylated

Item No. 22727

Overview and Properties

Contents:	This vial contains 50 µg of biotinylated IgG.
Synonyms:	LDLR, Low Density Lipoprotein Receptor
Immunogen:	Synthetic peptide from the C-terminal region of mouse LDLR
Species Reactivity:	(+) Human, mouse, and rat
Cross Reactivity:	(+) LDLR
Uniprot No.:	P35951 (murine)
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, 50% glycerol, 0.1% BSA, and 0.02% sodium azide
Host:	Rabbit
MW:	~100-160 kDa
Applications:	Immunocytochemistry (ICC) and Western blot (WB); the recommended starting dilution is 1:250-500. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: rLDLR (50 ng)
Lane 2: rLDLR (100 ng)
Lane 3: rLDLR (150 ng)

WB detection of recombinant LDLR by
LDL Receptor Polyclonal Antibody - Biotinylated (0.5 µg/ml)

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, 07/06/2022

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.¹ 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.² It mediates the endocytosis of LDL by binding to apolipoprotein E (ApoE) or ApoB on the LDL surface, thereby supplying cholesterol to cells.¹ Protein levels of LDLR are decreased in HepG2 cells expressing proprotein convertase subtilisin kexin 9 (PCSK9).³ Knockout of *Ldlr* increases plasma levels of cholesterol and triglycerides and induces the formation of atherosclerotic lesions in mice.⁵ Mutations in *LDLR* are associated with familial hypercholesterolemia.⁴ Cayman's LDL Receptor Polyclonal Antibody - Biotinylated is composed of an LDLR polyclonal antibody conjugated to biotin and can be used for immunocytochemistry (ICC) and Western blot (WB) applications. This antibody recognizes LDLR at approximately 100 to 160 kDa from human, mouse, and rat samples.

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**, 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. 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).
5. 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).

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