

PRODUCT INFORMATION



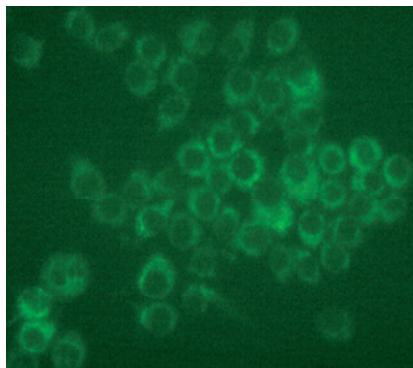
LDL Receptor Polyclonal Antibody

Item No. 10007665

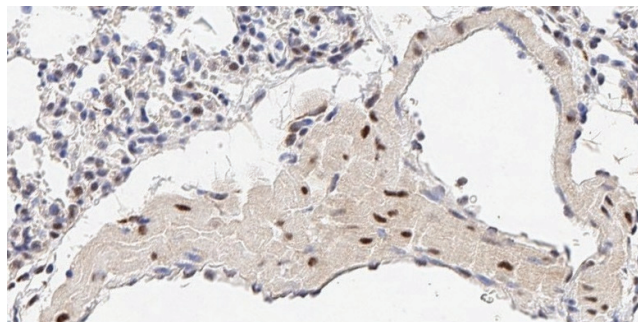
Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified antibody.
Synonyms:	LDLR, Low Density Lipoprotein Receptor
Immunogen:	Synthetic peptide from the C-terminal region of mouse LDL Receptor
Species Reactivity:	(+) Human, mouse, and rat; other species not tested
Uniprot No.:	P35951
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Immunocytochemistry (ICC), Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution for ICC and IHC is 1:100 and 1:200 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Immunofluorescent staining of RAW 264.7 cells with Cayman's LDLR Polyclonal Antibody at 4 µg/ml. The positive cytoplasm staining was visualized in green with Cayman's Goat Anti-Rabbit IgG FITC (Item No. 10006588).



Immunohistochemistry (IHC) analysis of formalin-fixed, paraffin-embedded (FFPE) mouse lung tissue after heat-induced antigen retrieval in pH 6.0 citrate buffer. After incubation with LDL Receptor Polyclonal Antibody (Item No. 10007665), at a 1:100 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, 05/01/2019

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

Cholesterol is an essential element of cell membranes and is carried around the body packaged in lipoproteins, primarily low-density lipoproteins (LDLs). The LDL receptors (LDLRs) are cell surface glycoproteins that scavenge LDL from the blood and regulate plasma LDL cholesterol. LDLRs contain five primary domains: the ligand binding domain, the homology with the EGF precursor domain, the O-linked sugars domain, the membrane-spanning region, and a cytoplasmic tail.¹ In humans, more than 60% of the LDL R3 are found in the liver.² LDLR expression is under hormonal control both *in vivo* and *in vitro*. Mutations in the LDLR gene cause disorders such as familial hypercholesterolemia and atherosclerosis.¹ Murine LDLR is 864 amino acids in length with an estimated molecular weight of 95 kDa. The protein is highly glycosylated through N- and O-linkages and thus migrates at 100 to 160 kDa bands on SDS-PAGE.¹ Cayman's LDL Receptor Polyclonal Antibody detects both glycosylated and unglycosylated proteins in tissue/cell samples such as liver, HepG2, and RAW 264.7 cells.

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).

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