

PRODUCT INFORMATION



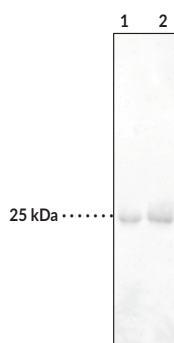
ApoA1 Polyclonal Antibody

Item No. 10008463

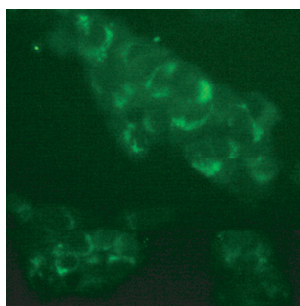
Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
Synonym:	Apolipoprotein A1
Immunogen:	Synthetic peptide from an internal region of human ApoA1
Species Reactivity:	(+) Human, mouse, and rat; other species not tested
Uniprot No.:	P02647
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Immunofluorescence (IF), immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution is 1:75, 1:120, and 1:200, respectively. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

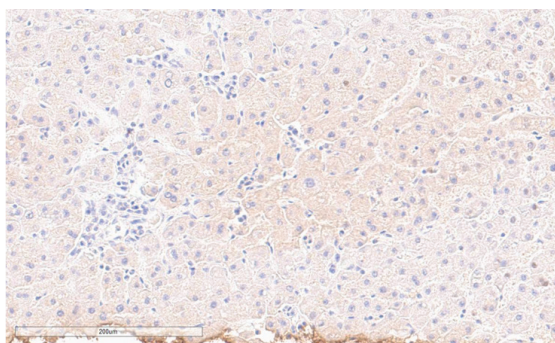
Images



Lane 1: ApoA1 western ready control (5 µl)
Lane 2: ApoA1 western ready control (10 µl)



Immunofluorescent staining of HepG2 cells with Cayman's ApoA1 Polyclonal Antibody at 4 mg/ml. The positive cytoplasm staining was visualized in green with Cayman's Goat Anti-Rabbit IgG FITC (Item No. 10006588).



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human liver tissue after heat-induced antigen retrieval in pH 6.0 citrate buffer. After incubation with ApoA1 Polyclonal Antibody (Item No. 10008463) at a 1:120 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, 10/17/2023

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

ApoA1 is a major protein component in high-density lipoproteins (HDLs). It acts as an acceptor for sequential transfers of phospholipids and free cholesterol from peripheral tissues as well as transporting cholesterol to the liver and other tissues for excretion and steroidogenesis.¹ Serum ApoA1 levels are inversely related to the risk of developing atherosclerosis.² Loss-of-function mutations in ApoA1 are causes of diseases such as HDL deficiency type 1 (or Tangier disease) and type 2 (familial hypoalphalipoproteinemia) and systemic non-neuropathic amyloidosis.^{3,4} Liver and small intestine are two main sources of the protein. Cayman's ApoA1 Polyclonal Antibody detects the protein by WB analysis at 25 kDa in tissue and cell samples such as liver, intestine, and HepG2 cells.

References

1. Ajees, A.A., Anantharamaiah, G.M., Mishra, V.K., *et al.* Crystal structure of human apolipoprotein A-1: Insights into its protective effect against cardiovascular diseases. *Proc. Natl. Acad. Sci. USA* **103(7)**, 2126-2131 (2006).
2. Haas, M.J., Reinacher, D., Li, J.P., *et al.* Regulation of ApoA1 gene expression with acidosis: Requirement for a transcriptional repressor. *J. Mol. Endocrinol.* **27(1)**, 43-57 (2001).
3. Tall, A.R. and Wang, N. Tangier disease as a test of the reverse cholesterol transport hypothesis. *J. Clin. Invest.* **106(10)**, 1205-1207 (2000).
4. Cheung, M.C., Mendez, A.J., Wolf, A.C., *et al.* Characterization of apolipoprotein A-I- and A-II-containing lipoproteins in a new case of high density lipoprotein deficiency resembling tangier disease and their effects on intracellular cholesterol efflux. *J. Clin. Invest.* **91(2)**, 522-529 (1993).

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