

Product Information



ApoAI Western Ready Control

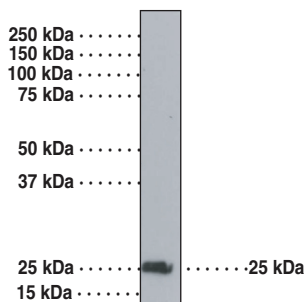
Item No. 10009751

Synonym:	Apolipoprotein AI
Source:	Human recombinant protein
M_r:	~25 kDa
Purity:	Whole cell lysate; a single band is detected on a western blot when probed with Cayman's ApoAI Polyclonal Antibody (Item No. 10008463)
Supplied in:	Laemmli buffer (2% SDS, 10% glycerol, 62.5 mM Tris-HCl pH 6.8, containing 5% 2-mercaptoethanol, bromophenol blue)
Stability:	≥2 years at -20°C
Applications:	Positive control for western blotting

ApoAI is a major protein component in high-density lipoproteins (HDL). 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 ApoAI levels are inversely related to the risk of developing atherosclerosis.² Loss-of-function mutations in ApoAI 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.

Laboratory Procedures

Western ready controls are designed to serve as positive controls on western blots. This ApoAI western ready control is supplied pre-boiled in 55 µl of Laemmli buffer. To use, thaw and load 10 µl of ApoAI western ready control per well and follow a standard western blotting protocol. The amount needed per well will depend on the antibody and detection system used.



Western blot of human recombinant ApoAI probed with Cayman's ApoAI Polyclonal Antibody (Item No. 10008463).

References

1. Ajees, A.A., Anantharamaiah, G.M., Mishra, V.K., *et al.* Crystal structure of human apolipoprotein A-I: 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 ApoAI gene expression with acidosis: Requirement for a transcriptional repressor. *J. Mol. Endocrinol.* **27**, 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**, 522-529 (1993).

Related Products

ACAT-2 Polyclonal Antibody - Item No. 100027 • ACAT-1 Polyclonal Antibody - Item No. 100028 • Hormone Sensitive Lipase Polyclonal Antibody - Item No. 10006371 • SREBP-2 Polyclonal Antibody - Item No. 10007663 • LDL Receptor Polyclonal Antibody - Item No. 10007665

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

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.