

Product Information

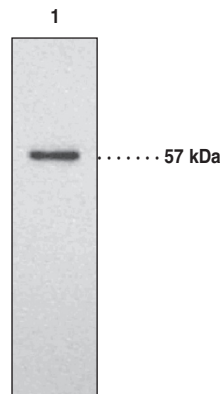


Endothelial Lipase Polyclonal Antibody

Item No. 100030 • Lot. No. XXXXX

Synonyms:	EDL; EL
Contents:	This vial contains (100-500 µg of peptide affinity-purified IgG, <i>lot specific</i>) in 500 µl TBS, pH 7.4, containing 50% glycerol, <i>lot specific</i> mg/ml BSA, and 0.02% sodium azide.
Host:	Rabbit
Antigen:	Human endothelial lipase amino acids 19-32. The antigen alignment with other known species sequences: Human AGSPVPFGPEGRLE Mouse AGS i t t l r P q G s L r
Cross-reactivity:	(+) Human, rat, mouse, porcine, and ovine endothelial lipase
Storage:	≥1 year at -20°C
Applications:	Western blotting and immunohistochemistry (IHC) recommended starting dilutions: western blotting: <i>lot specific</i> µg/ml; IHC (formalin-fixed paraffin-embedded sections): 2 µg/ml. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.
Concentration:	Varies by lot, from 0.2-1.0 mg/ml (100-500 µg/vial). Always 100 µl final working volume for western blotting.

Endothelial lipase (EL) is a member of the triglyceride lipase gene family. It functions primarily as a phospholipase and has low triglyceride lipase activity. It has been shown to be a major genetic determinant for the concentration, structure, and metabolism of high-density lipoprotein, which protects against atherosclerosis.^{1,2} It was originally cloned from endothelial cells and was found to be expressed in a distinct and complementary tissue-restricted fashion, with high-level expression in the liver, placenta, lung, ovary, and macrophage.³ The wide spread distribution of this protein suggests that it plays a general role in lipid metabolism. Immunohistochemical studies demonstrate that EL is expressed in infiltrating cells such as macrophages within atheromatous plaques, in addition to endothelial and smooth muscle cells in non-atherosclerotic coronary arteries. Furthermore, EL expression is detected in the neovasculature within atheromatous plaques in atherosclerotic coronary arteries, indicating that EL may have unique functional roles in atherosclerosis.⁴ Human endothelial lipase has an estimated molecular weight of 57 kDa.



Lane 1: HepG2 cell lysate (~30 µg)

References

1. Ishida, T., Choi, S., Kundu, R.K., *et al.* Endothelial lipase is a major determinant of HDL level. *J. Clin. Invest.* **111**(3), 347-355 (2003).
2. Ma, K., Cilingiroglu, M., Otvos, J.D., *et al.* Endothelial lipase is a major genetic determinant for high-density lipoprotein concentration, structure, and metabolism. *Proc. Natl. Acad. Sci. USA* **100**(5), 2748-2753 (2003).
3. Hirata, K-i., Dichek, H.L., Cioffi, J.A., *et al.* Cloning of a unique lipase from endothelial cells extends the lipase gene family. *J. Biol. Chem.* **274**(20), 14170-14175 (2003).
4. Azumi, H., Hirata, K-i., Ishida, T., *et al.* Immunohistochemical localization of endothelial cell-derived lipase in atherosclerotic human coronary arteries. *Cardiovascular Res.* **58**, 647-654 (2003).

Related Products

Oxidized Lipid HPLC Mixture - Item No. 34004 • 9(R)-HODE cholesteryl ester - Item No. 38406 • 9(S)-HODE cholesteryl ester - Item No. 38411 • (±)13-HODE cholesteryl ester - Item No. 38601 • 13(R)-HODE cholesteryl ester - Item No. 38606 • 13(S)-HODE cholesteryl ester - Item No. 38611 • Cholesteryl Linoleate Hydroperoxides - Item No. 48001 • Azelaoyl PAF - Item No. 60924 • CD36 Polyclonal Antibody - Item No. 100011 • Monoglyceride Lipase Polyclonal Antibody - Item No. 100035 • CD36 Monoclonal Antibody - Item No. 188150 • CD36 Blocking Peptide - Item No. 300011 • Endothelial Lipase (human) Blocking Peptide - Item No. 10004111

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.

Copyright Cayman Chemical Company, 10/05/2011

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