

# PRODUCT INFORMATION



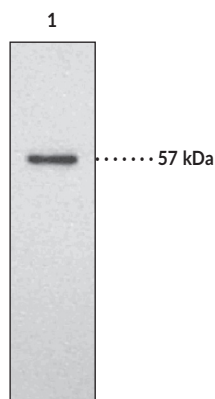
## Endothelial Lipase Polyclonal Antibody

Item No. 100030

### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
<b>Synonyms:</b>	EDL, EL
<b>Immunogen:</b>	Synthetic peptide from the N-terminal region of human endothelial lipase
<b>Species Reactivity:</b>	(+) Human, mouse, ovine, porcine, and rat
<b>Uniprot No.:</b>	Q9Y5X9
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



Lane 1: HepG2 cell lysate (~30 µg)

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, 10/10/2023

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

---

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.<sup>1,2</sup> 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.<sup>3</sup> 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.<sup>4</sup> Human endothelial lipase has an estimated molecular weight of 57 kDa.

## 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).

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