

# PRODUCT INFORMATION



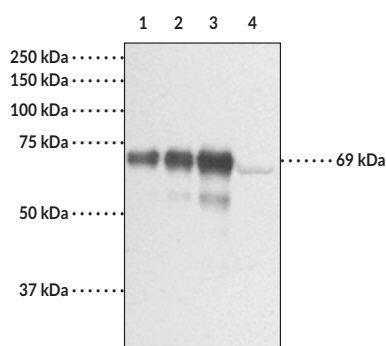
## Leukotriene A<sub>4</sub> Hydrolase Polyclonal Antibody

Item No. 160250

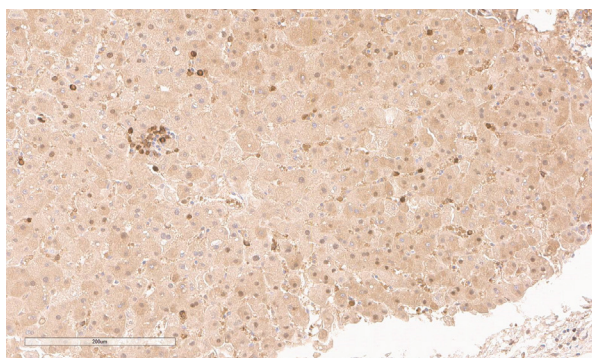
### Overview and Properties

<b>Contents:</b>	This vial contains 100 µg of protein A-purified polyclonal antibody.
<b>Synonym:</b>	LTA <sub>4</sub> Hydrolase
<b>Immunogen:</b>	Human recombinant leukotriene A <sub>4</sub> hydrolase (LTA <sub>4</sub> hydrolase) <sup>1,2</sup>
<b>Species Reactivity:</b>	(+) Human; other species not tested
<b>Uniprot No.:</b>	P09960
<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 1:40 and 1:200, respectively. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

### Images



Lane 1: LTA<sub>4</sub> Hydrolase Standard (0.005 µg)  
Lane 2: LTA<sub>4</sub> Hydrolase Standard (0.01 µg)  
Lane 3: LTA<sub>4</sub> Hydrolase Standard (0.02 µg)  
Lane 4: Raji cell lysate (50 µg)



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 Leukotriene A<sub>4</sub> hydrolase Polyclonal Antibody (Item No. 160250) at a 1:40 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, 11/16/2023

**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

---

Leukotriene A<sub>4</sub> (LTA<sub>4</sub>) hydrolase has been cloned from a variety of species including human, rat, and guinea pig.<sup>1,3-5</sup> These species exhibit ~90% homology at the amino acid level. Human LTA<sub>4</sub> hydrolase has a calculated molecular mass of 69 kDa based on the deduced amino acid sequence.<sup>1,3</sup>

## References

---

1. Funk, C.D., Rådmark, O., Fu, J.Y., *et al.* Molecular cloning and amino acid sequence of leukotriene A<sub>4</sub> hydrolase. *Proc. Natl. Acad. Sci. USA* **84**, 6677-6681 (1987).
2. Mancini, J.A. and Evans, J.F. Cloning and characterization of the human leukotriene A<sub>4</sub> hydrolase gene. *Eur. J. Biochem.* **231**, 65-71 (1995).
3. Minami, M., Ohno, S., Kawasaki, H., *et al.* Molecular cloning of a cDNA coding for human leukotriene A<sub>4</sub> hydrolase. Complete primary structure of an enzyme involved in eicosanoid synthesis. *J. Biol. Chem.* **262**, 13873-13876 (1987).
4. Makita, N., Funk, C.D., Imai, E., *et al.* Molecular cloning and functional expression of rat leukotriene A<sub>4</sub> hydrolase using the polymerase chain reaction. *FEBS Lett.* **299**, 273-277 (1992).
5. Minami, M., Mutoh, H., Ohishi, N., *et al.* Amino-acid sequence and tissue distribution of guinea-pig leukotriene A<sub>4</sub> hydrolase. *Gene* **161**, 249-251 (1995).

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