

PRODUCT INFORMATION



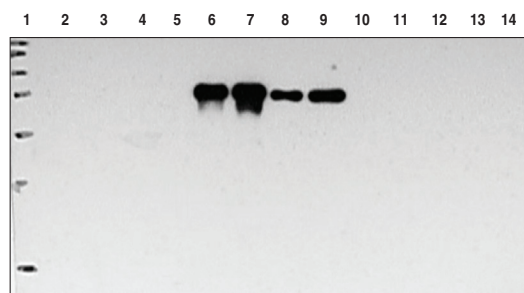
15-Lipoxygenase-2 Polyclonal Antibody

Item No. 10004454

Overview and Properties

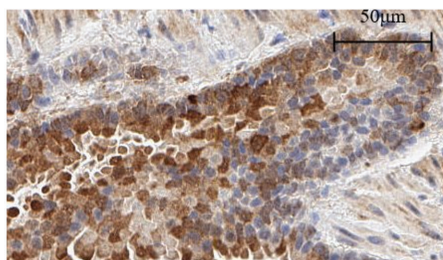
Contents:	This vial contains 50-500 µg of peptide affinity-purified antibody.
Synonyms:	ALOX15B, Arachidonate 15-lipoxygenase B, Arachidonate 15-Lipoxygenase type II, 15-LO-2, 15-LOX-2, Linoleate 13-lipoxygenase 15-Lob
Immunogen:	Synthetic peptide from the internal region of human 15-lipoxygenase-2
Cross Reactivity:	(-) Rabbit reticulocyte 15-lipoxygenase-1, porcine leukocytes 12-lipoxygenase-1, and 5-lipoxygenase (human recombinant)
Species Reactivity:	(+) Human; other species not tested
Uniprot No.:	O15296
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	TBS, pH 7.4, when reconstituted with 500 µl double distilled water
Host:	Rabbit
Applications:	Immunocytochemistry (ICC), immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution for IHC and WB is 1:40 and 1:200, respectively. ICC and other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: Precision Plus Protein Standard
Lane 2: 15-LO1 Western Ready Control (5 µl)
Lane 3: 15-LO1 Western Ready Control (10 µl)
Lane 4: 15-LO1 Recombinant (pure) (0.01 µg)
Lane 5: 15-LO1 Recombinant (pure) (0.02 µg)
Lane 6: 15-LO2 Western Ready Control (5 µl)
Lane 7: 15-LO2 Western Ready Control (10 µl)
Lane 8: 15-LO2 Recombinant (pure) (0.01 µg)
Lane 9: 15-LO2 Recombinant (pure) (0.02 µg)
Lane 10: 5-LO Western Ready Control (1 µl)
Lane 11: 5-LO Western Ready Control (5 µl)
Lane 12: 5-LO Western Ready Control (10 µl)
Lane 13: Porcine Leukocyte lysate (10 µg)
Lane 14: Porcine Leukocyte lysate (20 µg)

Figure 1: 15-Lipoxygenase 2 Polyclonal Antibody used at 1:200 dilution detected 15-LO2 but not 15-LO1, 5-LO or any lipoxygenases present in porcine leukocytes (i.e. 12-LO).



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human adrenal tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with 15-Lipoxygenase-2 Polyclonal Antibody (Item No. 10004454) 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, 06/10/2019

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

Two types of 15-lipoxygenase (15-LO) have been discovered and characterized, both of which metabolize arachidonic acid (AA) to produce 15(S)-hydroxyeicosatetraenoic acid (15(S)-HETE). 15-LO-1 oxygenates AA at both C15 and C12,¹ whereas 15-LO-2 exclusively oxygenates C15 of AA.² Human 15-LO-2 has a molecular mass of approximately 76 kDa and exhibits approximately 40% identity to the reticulocyte 15-LO-1.^{2,3} Expression of 15-LO-2 appears to be restricted to prostate, lung, skin, and cornea and may play a role in the normal development of these tissues.⁴ The protein levels and enzymatic activity of 15-LO-2 are both down-regulated in prostate cancer compared with normal and benign prostate tissues, implicating a possible protective role for 15-LO-2 against tumor formation.⁴⁻⁶

References

1. Kuhn, H., Barnett, J., Grunberger, D., *et al.* Overexpression, purification and characterization of human recombinant 15-lipoxygenase. *Biochem. Biophys. Acta* **1169**(1), 80-89 (1993).
2. Brash, A.R., Chang, M.S., and Boeglin, W.E. Discovery of a second 15S-lipoxygenase in humans. *Proc. Natl. Acad. Sci. USA* **94**, 6148-6152 (1997).
3. Kilty, I., Logan, A., and Vickers, P.J. Differential characteristics of human 15-lipoxygenase isozymes and novel splice variant of 15S-lipoxygenase. *Eur. J. Biochem.* **266**, 83-93 (1999).
4. Tang, S., Bhatia, B., Maldonado, C.J., *et al.* Evidence that arachidonate 15-lipoxygenase 2 is a negative cell cycle regulator in normal prostate epithelial cells. *J. Biol. Chem.* **277**(18), 16189-16201 (2002).
5. Shappell, S.B., Boeglin, W.E., Olson, S.J., *et al.* 15-lipoxygenase-2 (15-LOX-2) is expressed in benign prostatic epithelium and reduced in prostate adenocarcinoma. *Am. J. Pathol.* **155**, 235-245 (1999).
6. Jack, G.S., Brash, A.R., Olson, S.J., *et al.* Reduced 15-lipoxygenase-2 immunostaining in prostate adenocarcinoma: Correlation with grade and expression in high-grade prostatic intraepithelial neoplasia. *Human Pathology* **31**(9), 1146-1154 (2000).

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