

# PRODUCT INFORMATION

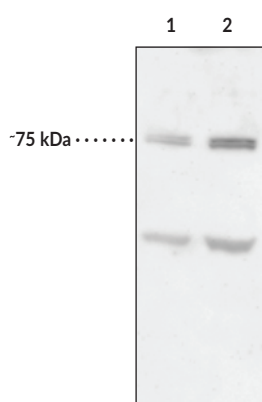


## 12-Lipoxygenase (leukocyte-type; mouse) Polyclonal Antibody Item No. 32822

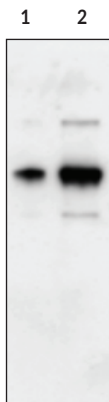
### Overview and Properties

<b>Contents:</b>	This vial contains 500 µg of protein A-purified polyclonal antibody.
<b>Synonyms:</b>	Alox12, Arachidonate 12-lipoxygenase, 12-LO, 12-LOX
<b>Immunogen:</b>	Full-length recombinant mouse 12-lipoxygenase (leukocyte-type)
<b>Species Reactivity:</b>	(+) Mouse, porcine
<b>Uniprot No.:</b>	P39654
<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 and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	ELISA and Western blot (WB); the recommended starting dilution is 1:200 for ELISA and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Lane 1: Porcine leukocyte lysate (10 µg)  
Lane 2: Porcine leukocyte lysate (25 µg)



Lane 1: Recombinant 12-Lipoxygenase (leukocyte-type; mouse) protein (1 ng)  
Lane 2: Recombinant 12-Lipoxygenase (leukocyte-type; mouse) protein (5 ng)

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, 12/13/2023

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

12-Lipoxygenase (12-LO) is one of several LOs that catalyzes the peroxidation of polyunsaturated fatty acids in a position-specific manner.<sup>1</sup> There are four 12-LO subtypes, encoded by distinct genes, that can be categorized according to their tissue expression in mice.<sup>2</sup> Leukocyte-type 12-LO is encoded by *Alox15* in mice and shares 74% sequence identity with human *ALOX15* but, in contrast to its human ortholog, encodes a 12-lipoxygenating *ALOX15* isoform.<sup>3-5</sup> Leukocyte-type 12-LO exists as a single polypeptide chain composed of an N-terminal regulatory domain and a C-terminal catalytic domain.<sup>1</sup> It catalyzes the oxidation of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) to 12(S)-HpETE (Item No. 44570) and, to a lesser extent, 15(S)-HpETE (Item No. 44720), which are further reduced to 12(S)-HETE (Item No. 34570) and 15(S)-HETE (Item No. 34720), respectively.<sup>1,6,7</sup> Other substrates of leukocyte-type 12-LO include eicosapentaenoic acid, docosahexaenoic acid, linoleic acid, and  $\gamma$ -linolenic acid, which are oxygenated at various positions in a substrate-dependent manner.<sup>8</sup> Leukocyte-type 12-LO is constitutively expressed at high levels in resident peritoneal macrophages, and its activity is increased by the Th2 cytokines IL-4 and IL-13 *in vitro*.<sup>5</sup> *Alox15*<sup>-/-</sup> mice are protected from increases in systolic blood pressure in a model of hypertension induced by L-NAME (Item No. 80210).<sup>9</sup> Cayman's 12-Lipoxygenase (leukocyte-type; mouse) Polyclonal Antibody can be used for ELISA and Western blot (WB) applications. The antibody recognizes 12-LO (leukocyte-type) at ~73 kDa from porcine and mouse samples.

## References

1. Kühn, H., Banthiya, S., and van Leyen, K. Mammalian lipoxygenases and their biological relevance. *Biochim. Biophys. Acta* **1851**(4), 308-330 (2015).
2. Kühn, H. and O'Donnell, V.B. Inflammation and immune regulation by 12/15-lipoxygenases. *Prog. Lipid Res.* **45**(4), 334-356 (2006).
3. Ivanov, I., Kühn, H., and Heydeck, D. Structural and functional biology of arachidonic acid 15-lipoxygenase-1 (*ALOX15*). *Gene* **573**(1), 1-32 (2015).
4. Adel, S., Karst, F., González-Lafont, À., *et al.* Evolutionary alteration of *ALOX15* specificity optimizes the biosynthesis of antiinflammatory and proresolving lipoxins. *Proc. Nat. Acad. Sci. USA* **113**(30), E4266-E4275 (2016).
5. Kühn, H., Gehring, T., Schröter, A., *et al.* Cytokine-dependent expression regulation of *ALOX15*. *J. Cytokine Biol.* **1**(2), 1000106 (2016).
6. Mashima, R. and Okuyama, T. The role of lipoxygenases in pathophysiology; new insights and future perspectives. *Redox Biol.* **6**, 297-310 (2015).
7. Chen, X.-S., Kurre, U., Jenkins, N.A., *et al.* cDNA cloning, expression, mutagenesis of C-terminal isoleucine, genomic structure, and chromosomal localizations of murine 12-lipoxygenases. *J. Biol. Chem.* **269**(19), 13979-13987 (1994).
8. Kutzner, L., Goloshchapova, K., Heydeck, D., *et al.* Mammalian *ALOX15* orthologs exhibit pronounced dual positional specificity with docosahexaenoic acid. *Biochim. Biophys. Acta Mol. Cell Biol. Lipids* **1862**(7), 666-675 (2017).
9. Kriska, T., Cepura, C., Magier, D., *et al.* Mice lacking macrophage 12/15-lipoxygenase are resistant to experimental hypertension. *Am. J. Physiol. Heart Circ. Physiol.* **302**(11), H2428-H2438 (2012).

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