

# PRODUCT INFORMATION



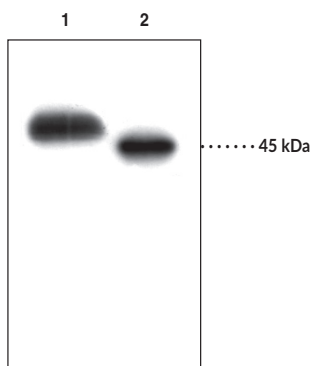
## PAF Acetylhydrolase (human) Polyclonal Antibody

Item No. 160603

### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl of peptide affinity-purified polyclonal antibody
<b>Synonyms:</b>	Lp-PLA <sub>2</sub> , PAF-AH, Platelet-activating Factor Receptor Acetylhydrolase
<b>Immunogen:</b>	Synthetic peptide corresponding to the C-terminal region of human PAF-AH
<b>Cross Reactivity:</b>	(+) PAF-AH
<b>Species Reactivity:</b>	(+) Human; (-) Mouse
<b>Uniprot No.:</b>	Q13093
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	TBS, pH 7.4, with 50% glycerol, 0.1% BSA, and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Application:</b>	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: LDL fraction from human plasma  
Lane 2: Recombinant human PAF-AH (20 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, 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

---

Platelet-activating factor acetylhydrolase (PAF-AH) is a calcium-independent member of the phospholipase A<sub>2</sub> (PLA<sub>2</sub>) family that hydrolyzes the acetyl group at the *sn*-2 position of PAF, a phospholipid mediator involved in a variety of biological activities.<sup>1,2</sup> There are three types of PAF-AH: two intracellular forms, PAF-AH1 and PAF-AH2, which are ubiquitously expressed, and an extracellular form, lipoprotein-associated phospholipase A<sub>2</sub> (Lp-PLA<sub>2</sub>), which is found in the plasma in association with LDL and HDL.<sup>3-5</sup> Unlike PAF-AH1, which is largely restricted to hydrolysis of PAF, PAF-AH2 and Lp-PLA<sub>2</sub> exhibit broader substrate specificity, hydrolyzing phospholipids with oxidized or short *sn*-2 acyl chains in addition to PAF.<sup>2,3</sup> PAF-AH1 has roles in neuronal development, protein trafficking, and spermatogenesis, and PAF-AH2 and Lp-PLA<sub>2</sub> have a protective roles against oxidative stress and atherosclerosis, respectively.<sup>6</sup> Cayman's PAF-acetylhydrolase (human) Polyclonal Antibody can be used for Western blot (WB) applications.

## References

---

1. Kono, N. and Arai, H. Platelet-activating factor acetylhydrolases: An overview and update. *Biochim. Biophys. Acta Mol. Cell Biol. Lipids* **1864**(6), 922-931 (2019).
2. McIntyre, T.M., Prescott, S.M., and Stafforini, D.M. The emerging roles of PAF acetylhydrolase. *J. Lipid Res.* **50**(Suppl), S255-S259 (2009).
3. Kono, N. and Arai, H. Intracellular platelet-activating factor acetylhydrolase, type II: A unique cellular phospholipase A<sub>2</sub> that hydrolyzes oxidatively modified phospholipids. *Enzymes* **38**, 43-54 (2015).
4. Murakami, M. Lipoquality control by phospholipase A2 enzymes. *Proc. Jpn. Acad. Ser. B Phys. Biol. Sci.* **93**(9), 677-702 (2017).
5. Karasawa, K., Harada, A., Satoh, N., et al. Plasma platelet activating factor-acetylhydrolase (PAF-AH). *Prog. Lipid Res.* **42**(2), 93-114 (2003).
6. Ashraf, M.A. and Nookala, V. In: StatPearls [Internet]. Treasure Island (FL): StatPearls Publishing (2021). Available from: <https://www.ncbi.nlm.nih.gov/books/NBK557392/>

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