

PAF Acetylhydrolase Assay Kit ab133088

[2 References](#) [1 Image](#)

Overview

Product name PAF Acetylhydrolase Assay Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
PAF-AH	89			3.5%

Inter-assay

Sample	n	Mean	SD	CV%
PAF-AH	89			10%

Sample type Serum, Plasma, Cell Lysate, Tissue Homogenate

Assay type Enzyme activity

Range 0.02 $\mu\text{mol}/\text{min}/\text{ml}$ - 0.2 $\mu\text{mol}/\text{min}/\text{ml}$

Species reactivity **Reacts with:** Mammals, Other species

Product overview Abcam's PAF Acetylhydrolase Assay Kit (ab133088) provides an accurate and convenient method for measurement of PAF acetylhydrolase (PAF-AH) activity (both cytosolic and extracellular). The assay uses 2-thio PAF which serves as a substrate for all PAF-AHs. Upon hydrolysis of the acetyl thioester bond at the *sn*-2 position by PAF-AH, free thiols are detected using 5,5'-dithio-*bis*-(2- nitrobenzoic acid).

Notes Platelet-activating factor (PAF) is a biologically active phospholipid synthesized by a variety of cells upon stimulation. PAF is converted to the biologically inactive lyso-PAF by the enzyme PAF acetylhydrolase (PAF-AH). PAF-AHs are located intra- and extra-cellularly (for instance, in both cytosol and plasma) Plasma PAF-AH is highly selective for phospholipids with very short acyl groups at the *sn*-2 position and is associated with lipoproteins.

Platform Microplate reader

Properties

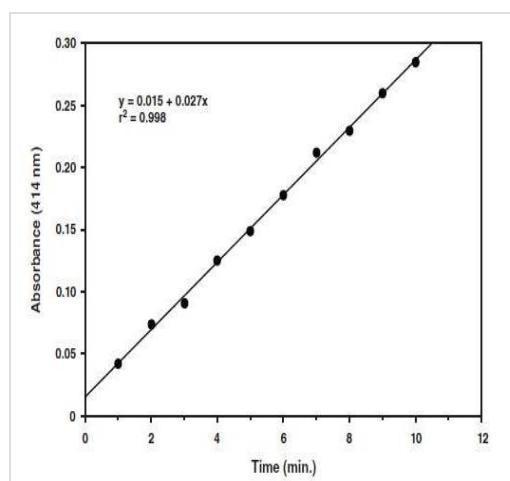
Storage instructions Store at -20°C. Please refer to protocols.

Components	96 tests
2-thio PAF (substrate)	2 vials
96-Well Plate Cover	1 unit
96-Well Solid Plate (Colorimetric Assay)	1 unit
DTNB	4 vials
Human Plasma PAF-AH Standard	1 vial
PAF Acetylhydrolase Assay Buffer 1	1 vial
PAF Acetylhydrolase Assay Buffer 2	1 vial

Function Exhibits lysophospholipase, transacylase, PAF acetylhydrolase and asparaginase activities.

Sequence similarities In the N-terminal section; belongs to the asparaginase 1 family.
Contains 5 ANK repeats.

Images



Representative Standard Curve using ab133088.

Activity of Human PAF-AH

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you

- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors