

PRODUCT INFORMATION



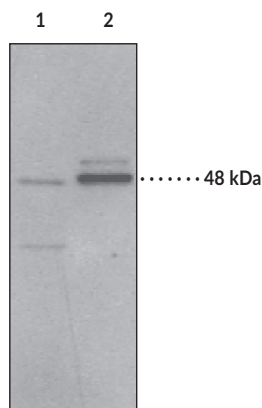
PAF Receptor (human) Polyclonal Antibody

Item No. 160602

Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
Synonym:	Platelet-activating Factor Receptor
Immunogen:	Synthetic peptide from the N-terminal region of human protein PAF receptor
Species Reactivity:	(+) Human and green monkey PAF receptor; other species not tested
Uniprot No.:	P25105
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2 with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Flow cytometry (FC) and Western blot (WB); the recommended starting dilution for FC and WB is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: U937 cell lysate
Lane 2: Raji cell lysate

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

PAF is a potent phospholipid mediator which exerts diverse biological actions by interaction with a G protein-coupled PAF receptor. The PAF receptor has been cloned from a number of species including human, rat, and guinea pig and is characterized as a 7-transmembrane receptor which induces phosphoinositol turnover through G-protein coupling.¹⁻⁵ Northern blot analysis reveals that the receptor is expressed in leukocytes, placenta, lung, spleen, small intestine, kidney, liver, and brain.^{3,4} In leukocyte cell populations the receptor is found on platelets, monocytes, neutrophils, and B-cells, whereas resting T-cells and natural killer cell lines do not express the PAF receptor.⁶ Human monocytes treated with INF- γ have a 2-6 fold increase in PAF receptor expression compared to untreated cells.⁷ PAF receptor is detected on immunoblot at 48 kDa.⁸

References

1. Nakamura, M., Honda, Z., Izumi, T., *et al.* Molecular cloning and expression of platelet-activating factor receptor from human leukocytes. *J. Biol. Chem.* **266**, 20400-20405 (1991).
2. Kunz, D., Gerard, N.P., and Gerard, C. The human leukocyte platelet-activating factor receptor. cDNA cloning, cell surface expression, and construction of a novel epitope-bearing analog. *J. Biol. Chem.* **267**, 9101-9106 (1992).
3. Ye, R.D., Prossnitz, E.R., Zou, A., *et al.* Characterization of a human cDNA that encodes a functional receptor for platelet activating factor. *Biochem. Biophys. Res. Commun.* **180**, 105-111 (1991).
4. Bito, H., Honda, Z., Nakamura, M., *et al.* Cloning, expression and tissue distribution of rat platelet-activating-factor-receptor cDNA. *Eur. J. Biochem.* **227**, 211-218 (1994).
5. Honda, Z., Nakamura, M., Miki, I., *et al.* Cloning by functional expression of platelet-activating factor receptor from guinea-pig lung. *Nature* **349**, 342-346 (1991).
6. Müller, E., Dagenais, P., Alami, N., *et al.* Identification and functional characterization of platelet-activating factor receptors in human leukocyte populations using polyclonal anti-peptide antibody. *Proc. Natl. Acad. Sci. USA* **90**, 5818-5822 (1993).
7. Quellet, S., Müller, E., and Rola-Pleszczynski, M. IFN- γ up-regulates platelet-activating factor receptor gene expression in human monocytes. *J. Immunol.* **152**, 5092-5099 (1994).
8. Marrache, A.M., Gobeil, F.Jr., Bernier, S.G., *et al.* Proinflammatory gene induction by platelet-activating factor mediated via its cognate nuclear receptor. *J. Immunol.* **169(11)**, 6474-6481 (2002).

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