

# Product Information

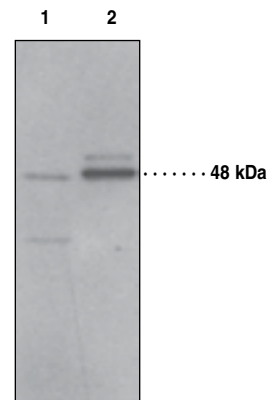


## PAF Receptor (human) Polyclonal Antibody

Item No. 160602 • Lot. No. XXXXXX

- Contents:** This vial contains (100-500 µg, *lot specific*) of peptide affinity-purified IgG in 500 µl of TBS, pH 7.4, containing 0.5 mg/ml BSA, 50% glycerol, and 0.02% sodium azide.
- Host:** Rabbit
- Antigen:** Synthetic peptide from the human PAF receptor (amino acids 1-17; MEPHDSHMDSEFRYTL) conjugated to KLH<sup>1-3</sup>
- Cross-reactivity:** (+) Human, mouse, rat, and porcine PAF receptor  
Homology of human PAF receptor peptide antigen with other species:  
Human            M E P H D S S H M D S E F R Y T L  
Rat                M E q n g S f r v D S E F R Y T L  
Guinea Pig      M E l n s S S r v D S E F R Y T L
- Stability:** ≥1 year at -20°C
- Applications:** Flow cytometry, immunocytochemistry, and western blot. For western blot use at approximately 1:200.
- Concentration:** Varies by lot; from 0.2-1.0 mg/ml (100-500 µg/vial). Always 100 µl final working volume for western blotting.

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.<sup>1-5</sup> Northern blot analysis reveals that the receptor is expressed in leukocytes, placenta, lung, spleen, small intestine, kidney, liver, and brain.<sup>3,4</sup> 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.<sup>6</sup> Human monocytes treated with INF-γ have a 2-6 fold increase in PAF receptor expression compared to untreated cells.<sup>7</sup> PAF receptor is detected on immunoblot at 48 kDa.<sup>8</sup>



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

### 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. Quillet, S., Müller, E., and Rola-Pleszczynski, M. INF-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).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/160602](http://www.caymanchem.com/catalog/160602)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 02/03/2012

## Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
[www.caymanchem.com](http://www.caymanchem.com)