

# PRODUCT INFORMATION



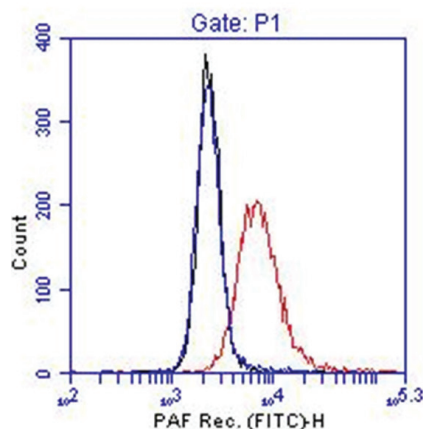
## PAF Receptor (human) Monoclonal Antibody (11A4, Clone 21)

Item No. 160600

### Overview and Properties

|                            |                                                                                                                                                                                                                                                                                        |
|----------------------------|----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Contents:</b>           | This vial contains 100 µg of protein A-purified monoclonal antibody.                                                                                                                                                                                                                   |
| <b>Synonym:</b>            | Platelet-activating Factor Receptor                                                                                                                                                                                                                                                    |
| <b>Immunogen:</b>          | Synthetic peptide from the internal region of human PAF receptor                                                                                                                                                                                                                       |
| <b>Species Reactivity:</b> | (+) Human, bovine, and porcine; other species not tested                                                                                                                                                                                                                               |
| <b>Uniprot No.:</b>        | P25105                                                                                                                                                                                                                                                                                 |
| <b>Form:</b>               | Lyophilized                                                                                                                                                                                                                                                                            |
| <b>Storage:</b>            | -20°C (as supplied)                                                                                                                                                                                                                                                                    |
| <b>Stability:</b>          | ≥3 years                                                                                                                                                                                                                                                                               |
| <b>Storage Buffer:</b>     | TBS, pH 7.4, when reconstituted in 500 µl of deionized water                                                                                                                                                                                                                           |
| <b>Clone:</b>              | 11A4, Clone 21                                                                                                                                                                                                                                                                         |
| <b>Host:</b>               | Mouse                                                                                                                                                                                                                                                                                  |
| <b>Isotype:</b>            | IgG2a                                                                                                                                                                                                                                                                                  |
| <b>Applications:</b>       | ELISA, Flow Cytometry (FC), and Immunocytochemistry (ICC); the recommended starting dilution for all applications is 1:1,000. Not recommended for Western blot. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

### Image



5000 THP-1 cells per sample

**Black:** IgG-FITC control

**Blue:** Secondary Antibody Alone (1:2000)

**Red:** 160600 (2µg), then second Ab (1:2000)

**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.<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- $\gamma$  have a 2-6 fold increase in PAF receptor expression compared to untreated cells.<sup>7</sup>

## References

---

1. Nakamura, M., Honda, Z., Izumi, T., *et al.* *J. Biol. Chem.* **266**, 20400-20405 (1991).
2. Kunz, D., Gerard, N.P., and Gerard, C. *J. Biol. Chem.* **267**, 9101-9106 (1992).
3. Ye, R.D., Prossnitz, E.R., Zou, A., *et al.* *Biochem. Biophys. Res. Commun.* **180**, 105-111 (1991).
4. Bito, H., Honda, Z., Nakamura, M., *et al.* *Eur. J. Biochem.* **227**, 211-218 (1994).
5. Honda, Z., Nakamura, M., Miki, I., *et al.* *Nature* **349**, 342-346 (1991).
6. Müller, E., Dagenais, P., Alami, N., *et al.* *Proc. Natl. Acad. Sci. USA* **90**, 5818-5822 (1993).
7. Quellet, S., Müller, E., and Rola-Pleszczynski, M. *J. Immunol.* **152**, 5092-5099 (1994).

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