

Product Information



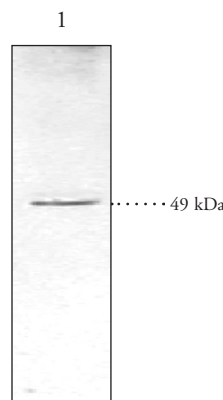
5-OxoETE Receptor Polyclonal Antibody

Item No. 100025 • Lot No. XXXX

- Synonyms:** R527, TG1019
- Contents:** This vial contains 400 µg of peptide affinity-purified IgG in *lot specific* µl TBS, pH 7.4, containing 50% glycerol, *lot specific*% BSA, and 0.02% sodium azide
- Host:** Rabbit
- Antigen:** Human 5-OxoETE receptor C-terminal amino acids 408-423
- Cross-reactivity:** (+) Human, mouse, rat, porcine, and Cos-7 (African green monkey) 5-OxoETE receptor; other species not tested
- Stability:** ≥1 year at -20°C
- Applications:** Recommended starting dilution for western blot - 4 µg/ml and for immunocytochemistry - 4 µg/ml. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

5-OxoETE (5-oxo-6E,8Z,11Z,14Z-eicosatetraenoic acid) is a potent stimulator of chemotaxis for eosinophils and neutrophils that mediates its action *via* a putative G-protein coupled receptor (GPCR).^{1,2} The 5-OxoETE receptor couples to G_{i/o} to inhibit cyclic AMP production and to mobilize intracellular calcium.^{1,2} Northern blot analysis reveals that the receptor is highly expressed on eosinophils and neutrophils, suggesting a key role in the inflammatory response.² In addition, the protein can also be detected in tissues such as kidney and liver.^{1,2}

Cayman Chemical's 5-OxoETE receptor polyclonal antibody can be used in western blot analysis (detection at 48 kDa) and immunocytochemistry. The high expression level of the receptor in leukocytes makes it the best positive control sample. When probing other tissue samples such as liver and kidney, membrane preparations or higher sensitivity detection systems may be required for visualization on immunoblot.



1. Porcine leukocyte lysate (25 µg)

References

- Hosoi, T., Koguchi, Y., Sugikawa, E., *et al.* Identification of a novel human eicosanoid receptor coupled to G_{i/o}. *J. Biol. Chem.* **277**(35), 31459-31465 (2002).
- Jones, C.E., Holden, S., Tenaillon, L., *et al.* Expression and characterization of a 5-oxo-6E,8Z,11Z,14Z-eicosatetraenoic acid receptor highly expressed on human eosinophils and neutrophils. *Mol. Pharmacol.* **63**(3), 471-477 (2003).

Related Products

5-OxoETE - Item No. 34250 • Cyclic AMP EIA Kit - Item No. 581001

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

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, 10/05/2011