

# PRODUCT INFORMATION



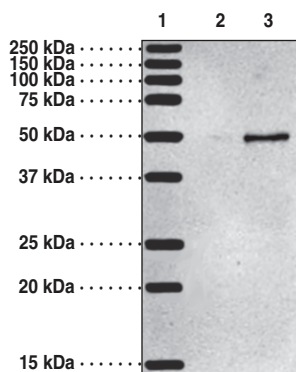
## 5-OxoETE Receptor Polyclonal Antibody

Item No. 100025

### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
<b>Synonyms:</b>	GPR170, OXER1, OXE Receptor, Oxoeicosanoid Receptor 1, R527, TG1019
<b>Immunogen:</b>	Synthetic peptide from the C-terminal region of human 5-OxoETE receptor
<b>Species Reactivity:</b>	(+) Human, Cos-7 (African green monkey), mouse, porcine, and rat; other species not tested
<b>Uniprot No.:</b>	Q8TDS5
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol, and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	Immunocytochemistry and Western blot; the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



Lane 1: Precision Plus Protein Standard  
Lane 2: Cos1 cell lysate (30 µg)  
Lane 3: Cos1 cell lysate (60 µg)

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, 10/10/2023

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

---

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

## References

---

1. 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).
2. 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).

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