

PRODUCT INFORMATION



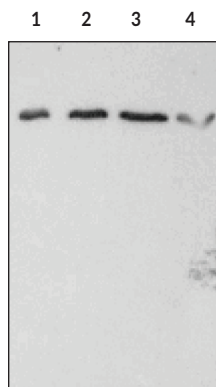
FP Receptor Polyclonal Antibody

Item No. 101802

Overview and Properties

Contents:	This vial contains peptide-affinity purified antibody lyophilized from 500 µl.
Synonyms:	PGF _{2α} Receptor, Prostaglandin F _{2α} Receptor
Immunogen:	Synthetic peptide from the N-terminal cytoplasmic region of mouse protein FP receptor
Species Reactivity:	(+) Human, bovine, mouse, ovine, and rat; other species not tested
Uniprot No.:	P43117
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	TBS, pH 7.4, with 0.5% BSA when reconstituted in 500 µl deionized water
Host:	Rabbit
Applications:	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: Bovine cornea (20 µg)
Lane 2: Bovine cornea (30 µg)
Lane 3: Bovine cornea (40 µg)
Lane 4: Ram seminal vesicle microsomes (30 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, 11/02/2023

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

The prostaglandin F_{2α} receptor (FP receptor) is a G protein-coupled receptor known to elicit smooth muscle contraction in a variety of tissues.⁶ FP receptor mRNA is found in reproductive, gastric, neural, and ocular tissues as well as specialized cells of the kidney.^{2,3,7,8} Two FP receptor isoforms, FP_A and FP_B have been identified in sheep.⁵ FP_B receptor is truncated at the carboxy-terminus thereby preventing desensitization (which normally occurs by protein kinase C phosphorylation of FP_A) and produces aberrant signalling in FP_B transfectants.^{9,10} The existence of a truncated human FP isoform responsible for signal-amplification contributing to luteolysis, cell growth, or malignant transformation is currently unknown.^{5,9-11} Cayman's FP Receptor Polyclonal Antibody can be used for Western blot applications. The calculated size of FP Receptor is 41 kDa and this antibody recognizes multiple bands from 50-67 kDa from human, bovine, mouse, ovine, and rat samples.

References

1. Sugimoto, Y., Hasumoto, K., Namba, T., *et al.* *J. Biol. Chem.* **269**, 1356-1360 (1994).
2. Abramovitz, M., Boie, Y., Nguyen, T., *et al.* *J. Biol. Chem.* **269**, 2632-2636 (1994).
3. Lake, S., Gullberg, H., Wahlqvist, J., *et al.* *FEBS Lett.* **355**, 317-325 (1994).
4. Sakamoto, K., Ezashi, T., Miwa, K., *et al.* *J. Biol. Chem.* **269**, 3881-3886 (1994).
5. Pierce, K.L., Bailey, T.J., Hoyer, P.B., *et al.* *J. Biol. Chem.* **272**, 883-887 (1997).
6. Narumiya, S., Sugimoto, Y., and Ushikubi, F. *Physiol. Rev.* **79**, 1193-1226 (1999).
7. Kitanaka, J., Hasimoto, H., Sugimoto, Y., *et al.* *Prostaglandins* **48**, 31-41 (1994).
8. Mukhopadhyay, P., Bian, L., Yin, H., *et al.* *Invest. Ophthalmol. Vis. Sci.* **42**(2), 424-428 (2001).
9. Fujino, H., Srinivasan, D., Pierce, K.L., *et al.* *Mol. Pharmacol.* **57**, 353-358 (2000).
10. Fujino, H. and Regan, J.W. *J. Biol. Chem.* **276**(16), 12489-12492 (2001).
11. Sugimoto, Y., Yamasaki, A., Segi, E., *et al.* *Science* **277**, 681-683 (1997).

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