

PRODUCT INFORMATION



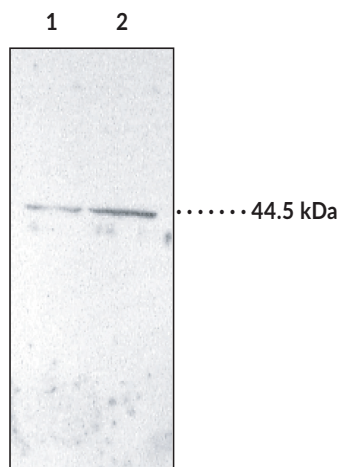
IP Receptor (mouse) Polyclonal Antibody

Item No. 160070

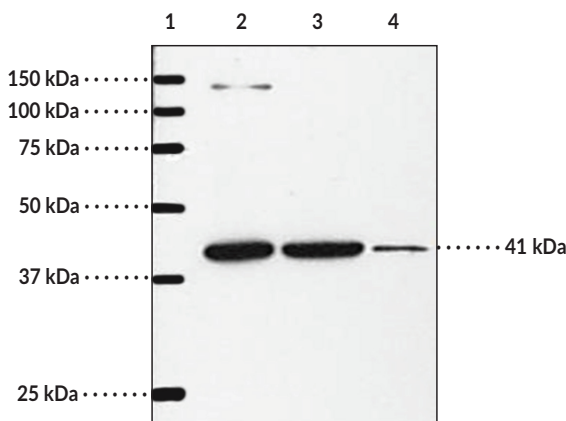
Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
Synonyms:	PGI ₂ Receptor, Prostacyclin Receptor, Prostaglandin I ₂ Receptor
Immunogen:	Peptide from the N-terminal region of mouse IP receptor
Cross Reactivity:	(+) IP Receptor
Species Reactivity:	(+) Human, mouse; other species not tested
Uniprot No.:	P43252
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Application:	Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: Mouse brain homogenate (40 µg)
Lane 2: Mouse brain homogenate (80 µg)



Lane 1: Precision Plus Protein Standard
Lane 2: Hek293T cell lysate (50 µg)
Lane 3: DLD1 cell lysate (35 µg)
Lane 4: A549 cell lysate (50 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, 11/13/2023

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 IP Receptor is a class A rhodopsin-like G protein-coupled receptor that mediates the actions of prostaglandin I₂ (PGI₂).¹ The C-terminal intracellular tail of the IP receptor undergoes isoprenylation and palmitoylation that results in anchoring of the tail to the plasma membrane. The IP receptor is expressed in platelets and vascular smooth muscle cells and in the aorta, lungs, heart, and kidneys.^{1,2} It signals through G proteins in a cell type- and expression-dependent manner and is involved in cardiovascular, inflammatory, and immune functions, as well as the pain response.^{1,3-6} An arginine-to-cysteine mutation at position 212 in the IP receptor inhibits its ability to activate adenylyl cyclase, which leads to increased platelet aggregation *ex vivo* and increases disease severity and the incidence of cardiovascular events in patients with a high risk of cardiovascular disease.⁷ The mouse IP receptor contains 28 additional N-terminal amino acids compared with the human receptor.^{2,8} Cayman's IP Receptor (mouse) Polyclonal Antibody can be used for Western blot. The antibody recognizes the IP receptor at approximately 45 kDa from human and mouse samples.

References

1. Stitham, J., Arehart, E.J., Gleim, S.R., *et al.* Human prostacyclin receptor structure and function from naturally-occurring and synthetic mutations. *Prostaglandins Other Lipid Mediat.* **82(1-4)**, 95-108 (2007).
2. Nakagawa, O., Tanaka, I., Usui, T., *et al.* Molecular cloning of human prostacyclin receptor cDNA and its gene expression in the cardiovascular system. *Circulation* **90**, 1643-1647 (1994).
3. Cheng, Y., Austin, S.C., Rocca, B., *et al.* Role of prostacyclin in the cardiovascular response to thromboxane A₂. *Science* **296(5567)**, 539-541 (2002).
4. Cui, Y., Kataoka, Y., Satoh, T., *et al.* Protective effect of prostaglandin I₂ analogs on ischemic delayed neuronal damage in gerbils. *Biochem. Biophys. Res. Commun.* **265(2)**, 301-304 (1999).
5. Norlander, A.E. and Peebles, R.S.J. Prostaglandin I₂ and T regulatory cell function: Broader impacts. *DNA Cell Biol.* (2021).
6. Murata, T., Ushikubi, F., Matsuoka, T., *et al.* Altered pain perception and inflammatory response in mice lacking prostacyclin receptor. *Nature* **388(6643)**, 678-682 (1997).
7. Arehart, E.J., Stitham, J., Asselbergs, F.W., *et al.* Acceleration of cardiovascular disease by a dysfunctional prostacyclin receptor mutation: Potential implications for cyclooxygenase-2 inhibition. *Circ. Res.* **102(8)**, 986-993 (2008).
8. Namba, T., Oida, H., Sugimoto, Y., *et al.* cDNA cloning of a mouse prostacyclin receptor. Multiple signaling pathways and expression in thymic medulla. *J. Biol. Chem.* **269**, 9986-9992 (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