

# PRODUCT INFORMATION



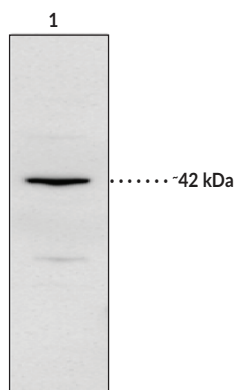
## IP Receptor (human) Polyclonal Antibody

Item No. 10005518

### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
<b>Synonyms:</b>	PGI <sub>2</sub> Receptor; Prostacyclin Receptor; Prostaglandin I <sub>2</sub> Receptor
<b>Immunogen:</b>	Synthetic peptide from the N-terminal region of human IP receptor
<b>Cross Reactivity:</b>	(+) IP Receptor
<b>Species Reactivity:</b>	(+) Human and mouse; other species not tested
<b>Uniprot No.:</b>	P43119
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	TBS, pH 7.4, with 50% glycerol, 0.1% BSA, and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Application:</b>	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: RT4 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, 10/12/2023

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 I<sub>2</sub> (PGI<sub>2</sub>).<sup>1</sup> 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.<sup>1,2</sup> 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.<sup>1,3-6</sup> 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.<sup>7</sup> The human IP receptor is N-terminally truncated by 28 amino acids compared with the murine receptor.<sup>8,9</sup> Cayman's IP Receptor (human) Polyclonal Antibody can be used for Western blot (WB). The antibody recognizes the IP receptor at approximately 42 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(4)**, 1643-1647 (1994).
3. Cheng, Y., Austin, S.C., Rocca, B., *et al.* Role of prostacyclin in the cardiovascular response to thromboxane A<sub>2</sub>. *Science* **296(5567)**, 539-541 (2002).
4. Cui, Y., Kataoka, Y., Satoh, T., *et al.* Protective effect of prostaglandin I<sub>2</sub> 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 I<sub>2</sub> 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. 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(4)**, 1643-1647 (1994).
9. 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(13)**, 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