

Product Information



IP Receptor (human) Polyclonal Antibody

Item No. 10005518 • Lot No. XXXXXX

Synonyms:	PGI ₂ Receptor; Prostacyclin Receptor; Prostaglandin I ₂ Receptor
Contents:	This vial contains XXX µg of affinity-purified IgG in 500 µl TBS, pH 7.4, containing 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide
Host:	Rabbit
Antigen:	Human IP receptor N-terminal amino acids 1-16 (MADSCRNLTYVRGSVG). Homology of peptide antigen with the IP receptor from other species is as follows: Human MADSCRNLTYVRGSVG Bovine MADSCRNLTYVR d SVG Mouse and rat MA g SCwN iTYV q d SVG
Cross-reactivity:	(+) Human, mouse, and rat IP receptor; other species not tested but expected to detect bovine IP receptor
Stability:	≥1 year at -20°C
Application:	Western blotting at X µg/ml (X:XXX) is the recommended starting dilution. Other applications have not been tested, therefore optimal working dilutions should be empirically determined

Prostaglandin I₂ (PGI₂) is a potent vasorelaxant and inhibitor of human platelet aggregation.¹ The PGI₂ receptor (IP receptor) also participates in signal transduction of the pain response, cardioprotection, and inflammation.²⁻⁶ Cloning of the murine and human IP receptors revealed a significant amino-terminal truncation of the human IP receptor compared to the murine receptor.^{7,8} Cayman Chemical's IP receptor antibody detects a 67 kDa band on immunoblot. The predicted migration of the deglycosylated protein is 40 kDa and bands at either 67 or 40 kDa may be detected depending on the degree of post-translational modification of the sample.⁹

References

1. Aristoff, P.A., Johnson, P.D., and Harrison, A.W. Synthesis of 9-substituted carbacyclin analogues. *J. Org. Chem.* **48**, 5341-5348 (1983).
2. Southall, M.D. and Vasko, M.R. Prostaglandin receptor subtypes, EP₃C and EP₄, mediate the prostaglandin E₂-induced cAMP production and sensitization of sensory neurons. *J. Biol. Chem.* **276(19)**, 16083-16091 (2001).
3. Murata, T., Ushikubi, F., Matsuoka, T., *et al.* Altered pain perception and inflammatory response in mice lacking prostacyclin receptor. *Nature* **388**, 678-682 (1997).
4. Cheng, Y., Austin, S.C., Rocca, B., *et al.* Role of prostacyclin in the cardiovascular response to thromboxane A₂. *Science* **296**, 539-541 (2002).
5. 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**, 301-304 (1999).
6. McLaughlin, V.V., Genthner, D.E., Panella, M.M., *et al.* Reduction in pulmonary vascular resistance with long-term epoprostenol (prostacyclin) therapy in primary pulmonary hypertension. *N. Engl. J. Med.* **338**, 273-277 (1998).
7. 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).
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).
9. Smyth, E.M., Nestor, P.V., and Fitzgerald, G.A. Agonist-dependent phosphorylation of an epitope-tagged human prostacyclin receptor. *J. Biol. Chem.* **271(52)**, 33698-33704 (1996).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10005518

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 of any kind, expressed or implied, including, but not limited to, the warranties of fitness for a particular purpose and merchantability, which extends beyond the description of the chemicals on the face hereof, except that the material will meet our specifications at the time of delivery. Buyer's exclusive remedy and Cayman Chemical Company's sole liability hereunder shall be limited to refund of the purchase price of, or at Cayman Chemical Company's option the replacement of, all material that does not meet our specifications. Cayman Chemical Company shall not be liable otherwise or for incidental or consequential damages, including, but not limited to, the costs of handling. Said refund or replacement is conditioned on Buyer giving written notice to Cayman Chemical Company within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within said thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

Copyright Cayman Chemical Company, 12/12/2011

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