

# Product Information



## IP Receptor (mouse) Blocking Peptide

Item No. 360070

Prostaglandin I<sub>2</sub> (PGI<sub>2</sub>) is a potent vasorelaxant and inhibitor of human platelet aggregation acting *via* the G-protein coupled PGI<sub>2</sub> receptor (IP receptor).<sup>1</sup> The IP receptor also participates in the signal transduction of the pain response, cardioprotection, and inflammation.<sup>2-6</sup> Cloning of the mouse and human IP receptors revealed a significant amino-terminal truncation of the human IP receptor compared to the mouse receptor.<sup>7,8</sup> Accordingly, the Cayman IP receptor (mouse) polyclonal antibody is useful only for immunochemical analysis of mouse and rat samples.

### Laboratory Procedures

This vial contains 200 µg of lyophilized peptide. This blocking peptide can be used in conjunction with Cayman's IP Receptor (mouse) Polyclonal Antibody (Item No. 160070) to block protein-antibody complex formation during immunochemical detection of the IP receptor.

Reconstitute the lyophilized peptide with 200 µl of PBS or distilled water. Store this peptide solution at -20°C. It should be stable for at least three years. To block antibody/protein complex formation, the following procedure is recommended:

1. Mix the IP Receptor (mouse) Polyclonal Antibody (Item No. 160070) and blocking peptide together in a 1:4 (v/v) ratio in a microfuge tube. For example, mix 20 µl of antibody and 80 µl of peptide.\*
2. Incubate for one hour at room temperature with occasional mixing prior to further dilution and application of the mixture to the immunoblot.
3. Dilute the mixture to the final working antibody concentration and apply to the slide or membrane as usual.

\*This is a recommended mixture. The minimum amount of peptide needed for complete blocking has not been precisely determined and may vary depending on the sample being analyzed. The amount of peptide required may need to be increased if sufficient blocking does not occur.

### 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, EP<sub>3</sub>C and EP<sub>4</sub>, mediate the prostaglandin E<sub>2</sub>-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 A<sub>2</sub>. *Science* **296**, 539-541 (2002).
5. 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**, 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).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/360070](http://www.caymanchem.com/catalog/360070)

### 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](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)

**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 or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 09/06/2012