

**MULTISCREEN™ MEMBRANE PREPARATION
HUMAN RECOMBINANT IP1 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: MC1206-1

Lot Number: MC1206-1-031314

Quantity: 11.7 mg/ml, 0.5 mg

Packaging Buffer: 20 mM Gly-Gly, 1 mM MgCl₂, 25 mM Sucrose (pH 7.2)

Host cell: CHO-K1

Transfection: Expression vector containing full-length human PTGIR cDNA (GenBank Accession Number: NM_000960) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Background: The human prostaglandin I₂ receptor IP₁ (or PTGIR) mediates the actions of prostaglandin I₂ (PGI₂ or prostacyclin), which is a labile metabolite of arachidonic acid produced in concert with the bis-enoic prostaglandins via the cyclooxygenase pathway. PGI₂ plays a major physiologic role as a strong mediator of vasodilation and inhibitor of platelet activation, and is an antithrombotic agent *in vivo* and mediates inflammation and pain. PGI₂ derived from COX2 plays a critical role in regulating the release of rennin and consequently in renovascular hypertension.

Figure 1

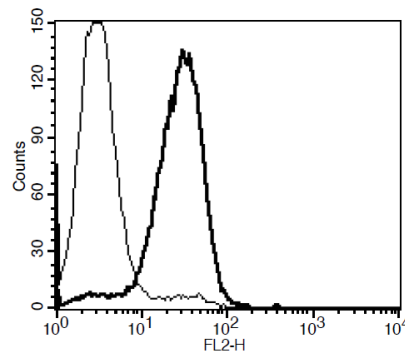


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Boie *et al.* (1994) Cloning and expression of a cDNA for the human prostanoid IP receptor. *J Biol Chem* 269:12173-12178.

Coleman *et al.* (1994) VIII International union of pharmacology classification of prostanoid receptors: properties, distribution, and structure of the receptors and their subtypes. *Pharm Rev* 46:205-229.

FOR RESEARCH USE ONLY.

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com
sales@multispaninc.com
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817
Fax: +1 (510) 887-0863
26219 Eden Landing Road
Hayward, CA 94545-3718
U.S.A.