

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT IP1 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1206-1

Lot Number: 2/26/13

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

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

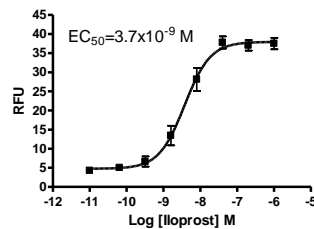
Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS

Stability: Stable for 1-2 days after thawing

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.

Application: Functional assays



Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

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.

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