

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

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1205

**Lot Number:** DC1205-100623

**Quantity:** 1 vial (4 x 10<sup>6</sup>) frozen cells

**Freeze Medium:** Cell Banker 2

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human PTGFR cDNA (GenBank accession number NM\_000959) with FLAG tag sequence at N-terminus

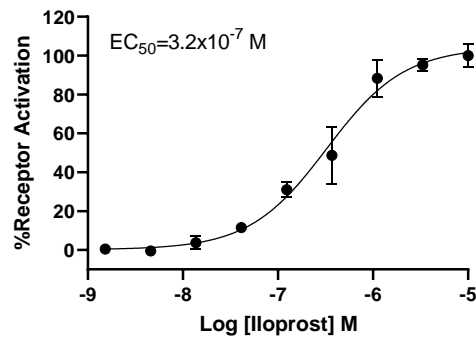
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS

**Background:** Prostaglandin F (2-alpha) is involved in a number of physiologic processes. It serves as a potent luteolytic agent in many species, has been implicated as a modulator of intraocular pressure, smooth muscle contraction in the uterus and elsewhere. Its effects on cells are mediated through specific interaction with the prostaglandin F receptor (FP or PTGFR). Knockout mice lacking the FP receptor are unable to deliver normal fetuses at term due to a lack of response to oxytocin. The mice also failed to show the decline in serum progesterone expected to precede parturition.

**Application:** Functional assays

**Figure 1**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

**References:**

Abramovitz *et al.* (1994) Cloning and expression of a cDNA for the human prostanoid FP receptor. *J Biol Chem* 269(4):2632-6.

Kim *et al.* (2012) Prostaglandin FP receptor inhibitor reduces ischemic brain damage and neurotoxicity. *Neurobiol Dis* 48(1):58-65.

**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.