

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

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1205

Lot Number: 03/06/13

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

Freeze Medium: Sigma Freezing Medium (C-6164)

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

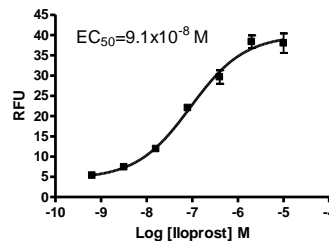
Propagation Medium: DMEM, 10% FBS

Stability: Stable for 1-2 days after thawing

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



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.