



# Examining the FP receptor-triggered CREB activation in cell-based assays

P1554

C.S. Ho<sup>†</sup>, C. Patel<sup>‡</sup>, B. Saepoo<sup>†</sup>, and I. Morano<sup>‡</sup>

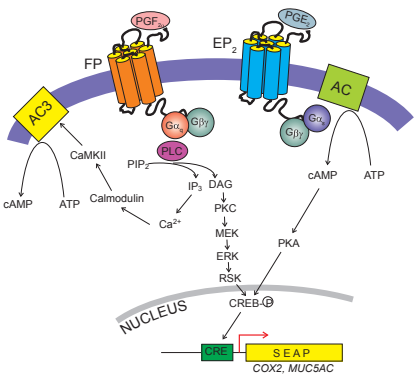
<sup>†</sup>Assay Services Division and <sup>‡</sup>Drug Discovery Division, Cayman Chemical, Ann Arbor, Michigan

## Abstract

Prostaglandins (PGs) play a crucial role in mediating a wide range of cellular processes and progression in inflammation, cancer, cardiovascular, and other diseases. Therefore, enzymes involved in the production of prostanoids, such as cyclooxygenases and prostaglandin synthases, as well as the receptors that mediate their biological responses, are important therapeutic targets. Several families of G protein-coupled receptors have been identified for prostanoids with different agonist selectivity and subsequent signaling pathways. The PGF<sub>2α</sub> receptor (FP) was initially characterized as a G<sub>q</sub>-coupled receptor, which leads to an increase in inositol triphosphate/diacylglycerol, intracellular calcium mobilization, and protein kinase C (PKC) activation. Interestingly, PGF<sub>2α</sub> has also been suggested to induce CREB-regulated genes in human amniotic fibroblasts and tracheobronchial epithelial cells through PKC activity. The activation of CREB by the FP receptor has been reproduced in this study by co-transfecting the FP receptor with a CRE-SEAP reporter into HEK293T cells using reverse transfection. Replicate transfection plates were used to examine receptor activations by agonists in a calcium mobilization assay and an ELISA for cAMP. In the calcium assay, cloprostenol, as well as the free acids of bimatoprost and latanoprost, exhibited similar potencies and efficacies of PGF<sub>2α</sub> with EC<sub>50</sub> values around 50 nM. In contrast, agonists induced only subtle increases of cAMP unless a G<sub>i</sub>/G<sub>o</sub> chimeric G protein was co-expressed to redirect the receptor activation signal to the adenylate cyclase pathway. This verified that CREB activation by the FP receptor is not mediated by the adenylate cyclase pathway. In the presence of the G<sub>i</sub>/G<sub>o</sub> chimeric G protein, all four agonists that were tested led to the same maximal increase in cAMP levels after a 45-minute incubation. Latanoprost free acid and PGF<sub>2α</sub> exhibited comparable EC<sub>50</sub> values in the cAMP assay performed in parallel with the calcium assay on replicate plates, whereas bimatoprost and cloprostenol were one order more potent in the G<sub>i</sub>/G<sub>o</sub>-facilitated cAMP assay than in the calcium mobilization assay. This suggests the FP receptor may preferentially interact with the chimeric G protein upon certain agonist binding. On the other hand, the expression of G<sub>i</sub>/G<sub>o</sub> is not required in the CRE-SEAP reporter assay for FP receptor activity, although the addition of G<sub>i</sub>/G<sub>o</sub> produced a left shift in the dose-response curves. This direct stimulation of CRE-SEAP reporter using the same assay platform was not observed with other G<sub>q</sub>-coupled receptors, such as orexin 1 and 2 receptors, that can also activate the PKC pathway. In other words, the activation of PKC alone may not be sufficient to mediate the effect of the FP receptor on CREB.

## Introduction

The prostaglandin F<sub>2α</sub> (PGF<sub>2α</sub>) receptor, PTGFR or FP, is a G protein-coupled receptor that has been shown to activate the G<sub>q</sub> signaling pathway leading to elevation of intracellular calcium levels.<sup>1</sup> HEK293 cells stably transfected with FP have been tested with various agonists in calcium mobilization assays and exhibited comparable potencies in these cell lines compared to endogenous expression of the FP receptor.<sup>2</sup> Nevertheless, the stimulation of FP receptors has also been linked to the activation of CREB and upregulation of COX2 and MUC5AC, suggesting a cross-talk between signaling pathways through PKC.<sup>3-5</sup> On the other hand, a different model has been proposed that the FP receptor-triggered calcium signal may potentiate the EP<sub>2</sub>/G<sub>q</sub> pathway-mediated cAMP increase through CaMKII and isoform 3 of adenylate cyclase.<sup>6</sup> In contrary, a recent study reported the opposite effects of EP<sub>2</sub>/cAMP/PKA and FP/calcium/PKC pathways in the expression of OVGP1 in bovine oviductal epithelial cells.<sup>7</sup> In this study, a transfection complex with an optimized ratio of FP expression plasmid and CRE-SEAP reporter construct was coated on 96-well microplates for reverse transfection of HEK293T cells. The FP receptor was stimulated with several known agonists the day after seeding cells, and the receptor activation and downstream gene regulation were monitored by calcium mobilization, ELISA for cAMP, and reporter gene assays. By comparing the relative potencies of these agonists in different assays, with and without the supplement of chimeric G protein, and the use of kinase inhibitors, the signaling mechanism models of the FP receptor were examined on this reverse transfection platform.



## Materials and Methods

### DNA constructs

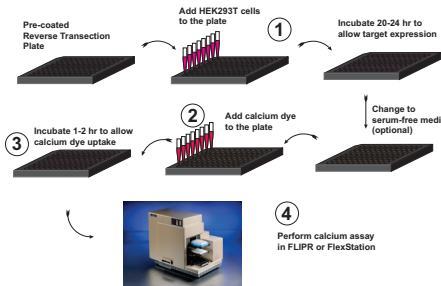
An FP2A cDNA fragment with optimized codon usage and Kozak sequence was designed and assembled into pCMV6-XL5 vector. A G<sub>i</sub>/G<sub>o</sub> chimera expression construct was produced by replacing the coding sequence for the C-terminal 5 amino acid of G<sub>i</sub> cDNA with corresponding sequence from G<sub>o</sub> as described by Komatsuzaki, et al.<sup>8</sup> CRE-SEAP reporter plasmid contains the cDNA of a secreted form of the human placental alkaline phosphatase (SEAP) under the regulation of five cyclic AMP response elements as previously described.<sup>3,10</sup>

### Cell culture and transfection

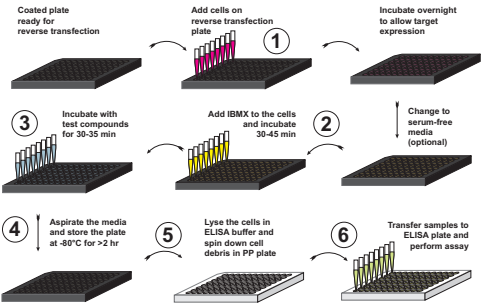
Transfection complex containing lipids for transfection and optimized amount of FP2A expression construct plasmid and CRE-SEAP reporter DNA was immobilized on tissue culture treated 96-well microplates using a proprietary protocol. Expression construct for G<sub>i</sub>/G<sub>o</sub> was added to the DNA mixture in the experiment that required the chimeric G protein. HEK293T/17 cells were grown in complete medium (Dulbecco's modified essential medium supplemented with 10% fetal calf serum). Actively dividing HEK293T cells were plated on the microplates with transfection complex at 50,000 cells per well 18-24 hours before assay in complete medium, Opti-MEM with 0.5% FBS, or Opti-MEM alone depending on the assay.

### Cell-based assays

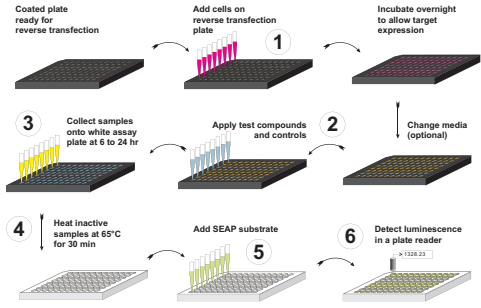
For the calcium mobilization assay, cells were plated in complete medium, which was changed to 100 µl/well serum-free DMEM at 20-24 hours after plating cells. Calcium 6 dye (Molecular Devices) was dissolved in assay buffer with probenecid according to manufacturer's instruction. The dye was dispensed at 100 µl/well to the plate and incubated at 37°C for 90 minutes. Test compounds were dissolved in DMSO at 10 mM stock concentration and diluted in assay buffer to 5x of final concentration on a polypropylene deep well plate. Assay was performed in FlexStation 3 after equilibrating the cell plate to ambient temperature in the machine for 30 minutes. Fluorescence at ex/em 485/525 nm was monitored every 3 seconds. Dilutions of testing compounds were delivered to the well in 50 µl volume at 20 seconds.



For the cAMP ELISA, cells were plated in Opti-MEM with 0.5% FBS, which was replaced with 160 µl/well Opti-MEM containing 0.5 mM IBMX at 20-24 hours after plating cells on the transfection plate. After 45 minutes of incubation, 40 µl of serial dilutions of test compounds in Opti-MEM+IBMX were added to the well. After 30-45 minutes of incubation at 37°C, media were aspirated, and the plates were put on dry ice before transferring to -80°C. After storing the plates at -80°C for at least two hours, the plates were removed from the freezer and thawed on bench. Cells were lysed in ELISA buffer containing 0.1% Tween-20, and the cell lysates were centrifuged to pellet the cell debris. The supernatants were examined with a cAMP ELISA kit (Cayman Item No. 501040).



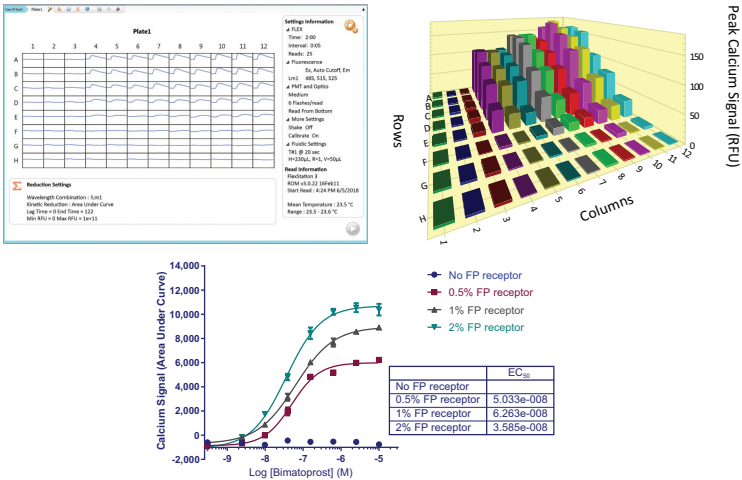
For the reporter assay, cells were plated in serum-free Opti-MEM and serial dilutions of test compounds were applied to the cells after they were seeded on the transfection plate for 20-24 hours before stimulation. Media samples were collected at 6-24 hours after stimulation and assayed for SEAP reporter activity using a luminescence substrate.



## Results

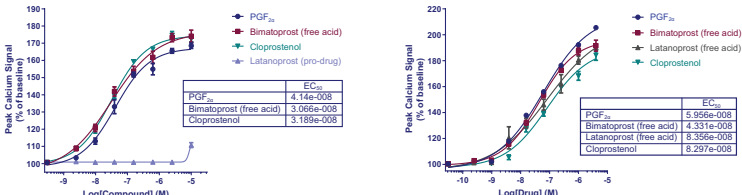
### Titration of FP cDNA in a Reverse Transfection Complex by Calcium Assay

The expression level of FP receptors was titrated by diluting the expression plasmid in empty vector DNA for transfection. Transfection complexes with a fixed amount of total DNA, but different percentages of FP2A, were coated on TC-treated 96-well plates in triplicated columns using a proprietary protocol. Cells were added to the plates and stimulated with serial dilutions of a control agonist, bimatoprost (free acid), according to the calcium assay protocol. An optimal transfection complex was based on the robustness of assay and EC<sub>50</sub> values.



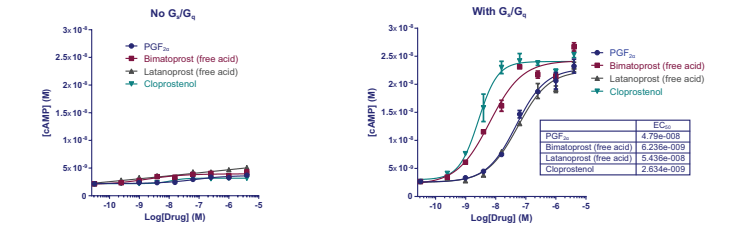
### Validation of the Assay with Selected Agonists

Using the same transfection complex recipe, two batches of plates were produced separately. One plate from each batch was plated with different passages of HEK293T/17 cells on different days and was tested in the calcium mobilization assay with several known agonists serially diluted in slightly different schemes. The EC<sub>50</sub> values obtained for PGF<sub>2α</sub>, bimatoprost (free acid), and cloprostenol are comparable between the two experiments.



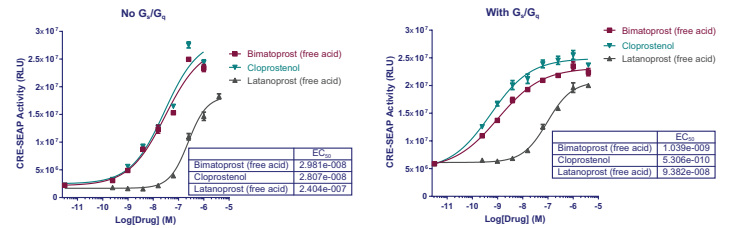
### Examining the Role of cAMP in FP Receptor Signaling Pathway

The same serial dilutions of FP receptor agonists were tested on an identical reverse transfection plate from the calcium mobilization assay above in a cAMP assay. With the endogenous G<sub>q</sub>/G<sub>o</sub> signaling pathway, there are only marginal increases in cAMP levels upon activation of the FP receptor by these agonists. This indicates the activation of a CREB-regulated gene by the FP receptor cannot be mediated by the cAMP pathway directly. With the supplement of an expression plasmid for the G<sub>i</sub>/G<sub>o</sub> chimeric G protein α-subunit in the transfection complex, the originally G<sub>q</sub>-coupled receptor activation was redirected to a G<sub>q</sub> signaling pathway, leading to drastic increases in cAMP. PGF<sub>2α</sub> and latanoprost (free acid) exhibited EC<sub>50</sub> values comparable to the calcium mobilization assays above, whereas cloprostenol and bimatoprost (free acid) shifted the dose-response curves to the left by around 1 to 1.5 orders of magnitude. This suggests the chimeric G protein may mimic the effect of EP<sub>2</sub>/G<sub>q</sub> pathway coactivation when the FP receptor is activated by certain agonists. In other words, the activation of the FP receptor may trigger both the G<sub>q</sub> and G<sub>i</sub> (through G<sub>i</sub>/G<sub>o</sub>) pathways and exhibits a synergistic effect in cAMP increase depending on the agonist.



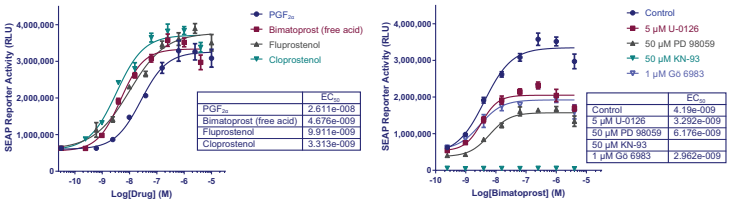
### Downstream Signaling Activity Revealed by Reporter Assay

The same batch of plates examined in the calcium mobilization and cAMP assays above was used to show the downstream effect of FP receptor activity by SEAP reporter assay after overnight incubation with test compounds. In the reverse transfection plate without G<sub>i</sub>/G<sub>o</sub> chimeric G protein, both cloprostenol and bimatoprost (free acid) exhibited comparable EC<sub>50</sub> values as in the calcium mobilization assay while latanoprost (free acid) showed a few-fold higher EC<sub>50</sub>. This confirmed the activation of CREB even without a significant increase in cAMP level by FP receptor activation. With the supplement of G<sub>i</sub>/G<sub>o</sub> chimeric G protein, the FP receptor activation can be coupled to the cAMP/PKA/CREB pathway directly. Under this condition, the basal reporter activity was elevated presumably due to a constitutive activity by the expression of G<sub>i</sub>/G<sub>o</sub> chimeric G protein. In addition, the dose-response curves of bimatoprost (free acid) and cloprostenol shifted to the left by 1.5 orders of magnitude, while the latanoprost (free acid) exhibited a much smaller shift. These findings are consistent with the kinetic changes by G<sub>i</sub>/G<sub>o</sub> observed in the assays for secondary messengers in the upstream signaling pathways upon FP receptor activation above.



### Dissecting the Signaling Pathway Leading to CREB Activation by the FP Receptor

To investigate the signaling pathway linking the FP/G<sub>q</sub> activation to CREB phosphorylation, inhibitors for MEK1/2, PKC, and CaMKII were examined in the reporter assay. To avoid the toxic effect of prolonged incubation with kinase inhibitors, a 6-hour (instead of 16-hour) stimulation protocol was used after longer post-transfection incubation. The new protocol was found to have lower EC<sub>50</sub> values with several control agonists examined, while the relative potencies were not changed. Serial dilutions of bimatoprost (free acid) were added to the cells after pre-incubation with kinase inhibitors for 20 minutes. The CaMKII inhibitor KN-93 at 50 µM revealed severe cytotoxic effects and abolished the activity of the FP receptor. Both MEK1/2 inhibitors U-0126 (5 µM) and PD 98059 (50 µM) showed significant suppression of the FP receptor trigger reporter activity, while the latter also led to a decrease in basal reporter activity. The PKC inhibitor Gö 6983 (1 µM) produced a suppressive effect similar to that of the MEK1/2 inhibitors. With these MEK1/2 and PKC inhibitors, the EC<sub>50</sub> of bimatoprost (free acid) was not affected. These findings confirmed the involvement of PKC and MEK1/2 in the signaling pathway connecting the FP receptor to the phosphorylation of CREB.



## Conclusion and Discussion

1. A robust assay platform has been developed to study FP receptor activity.
2. The same reverse transfection plate can be used in calcium mobilization, cAMP, and reporter assay.
3. G<sub>i</sub>/G<sub>o</sub> chimeric G protein can efficiently couple the FP receptor to the G<sub>q</sub> signaling pathway and may mimic the effect of EP<sub>2</sub> co-stimulation.
4. FP receptor activation leads to phosphorylation of CREB through the PKC/MEK pathway without increasing intracellular cAMP levels.

### References

1. Ito, S., Sakamoto, K., Morizaki, Oda, N., et al. Prostaglandin F<sub>2α</sub> receptor is coupled to G<sub>q</sub> in cDNA-transfected Chinese hamster ovary cells. *Biochem. Biophys. Res. Commun.* **2002**; 275:752-752 (1994).
2. Kelly, C.E., Williams, G.W., and Shatt, N.A. Real-time intracellular Ca<sup>2+</sup> mobilization by bimatoprost, unoprostone, and other analogs via endogenous muscarinic, calc, and cloned human FP prostaglandin receptors. *J. Pharmacol. Exp. Ther.* **304**(1): 238-245 (2003).
3. Chung, W.-C., Ren, S.-H., Sun, H., et al. CREB-mediated prostaglandin F<sub>2α</sub>-induced MUC5AC overexpression. *J. Immunol.* **162**(4): 2349-2354 (2009).
4. Jones, T., and Fagan, K. Novel suppression mechanism operating in early phase of adenosine by positive feedback loop for enhancement of cyclooxygenase-2 expression through prostaglandin F<sub>2α</sub> receptor-mediated activation of MEK/ERK/CREB cascade. *FEBS J.* **278**(14): 2903-2912 (2011).
5. Gao, C.M., Kaurizumi, N., Sun, K., et al. Cross-talk between PKC and CREB in the induction of COX-2 by PGF<sub>2α</sub> in human choroid plexus. *Endocrinology* **153**(5): 4938-4945 (2012).
6. Aben, A.B., Sales, K.J., Catalano, R.J., et al. EP2 receptor-mediated cAMP release is augmented by PGF<sub>2α</sub> activation of the FP receptor via the calcium-calmodulin pathway. *Cell Signal.* **22**(1): 71-79 (2010).
7. Zhang, N., Mao, W., Zhang, Y., et al. The prostaglandin F<sub>2α</sub> receptor (PTGFR) and prostaglandin F<sub>2α</sub> receptor (PTGFR) mediate oviductal glycoprotein 1 expression in bovine oviductal epithelial cells. *J. Reprod. Dev.* **64**(2): 101-108 (2018).
8. Komatsuzaki, K., Matsuyama, Y., Gotoh, M., et al. A novel system that reports the G protein-linked to a given receptor: a study of type 3 cannabinoid receptor. *FEBS Lett.* **406**(1-2): 165-170 (1997).
9. Chen, W., Shultz, T.S., Stok, R.S., et al. A colorimetric assay for measuring activation of G<sub>i</sub> and G<sub>o</sub> coupled signaling pathways. *Anal. Biochem.* **226**(2): 349-354 (1995).
10. Danneberg, Y., Herms, S., Thibaudou, E., et al. A reporter gene assay for high-throughput screening of G-protein-coupled receptors: stably or transiently expressed in HEK293 EBNA cells grown in suspension culture. *Anal. Biochem.* **340**(2): 315-325 (2005).