
Optimization of the GeneBLAzer® PTGER2-NFAT-bla CHO-K1 Cell Line

GeneBLAzer® PTGER2 CHO-K1 DA Assay Kit**GeneBLAzer® PTGER2 CRE-*bla* CHO-K1 Cells**

Catalog Numbers – K1339 and K1723

Cell Line Descriptions

GeneBLAzer® PTGER2 CHO-K1 DA (Division Arrested) cells and GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 cells contain the human prostaglandin E receptor 2 (PTGER2) receptor (Accession #[NM_000956.2](#)) stably integrated into the CellSensor® CRE-*bla* CHO-K1 cell line. CellSensor® CRE-*bla* CHO-K1 cells (Cat. no. K1535) contain a beta-lactamase (*bla*) reporter gene under control of the Cyclic AMP Response Element (CRE). Division Arrested (DA) cells are available as an Assay Kit, which includes cells and sufficient substrate to analyze 1 x 384-well plate.

DA cells are irreversibly division arrested using a low-dose treatment of Mitomycin-C, and have no apparent toxicity or change in cellular signal transduction. Both GeneBLAzer® PTGER2 CHO-K1 DA cells and GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 cells are functionally validated for Z'-factor and EC₅₀ concentrations of Prostaglandin E2 (Figure 1). In addition, GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 cells have been tested for assay performance under variable conditions, including DMSO concentration, cell number, stimulation time, and substrate loading time. Additional testing data using alternate stimuli are also included.

Target Description

Prostanoids are the metabolites of arachidonic acid by cyclooxygenase and include prostaglandins (PGs) and thromboxanes (TXs). There are five primary types of prostanoids including PGE₂, PGD₂, PGF_{2α}, PGI₂, and TxA₂. Following cell stimulation, prostanoids are synthesized, released, and exert their action on cells in the vicinity of their release. The receptors that mediate the actions of the prostanoids belong to the G protein-coupled receptor (GPCR) superfamily and are termed P receptors with a preceding letter to designate the natural prostanoid to which each receptor is most sensitive: EP, FP, IP, TP, and DP (1-2). The EP class has four receptors named EP1, EP2, EP3, and EP4 based upon the sensitivity to various agonist and antagonists. Each of these receptors belongs to the rhodopsin-type sub-family of GPCRs.

The prostaglandin E2 receptor (PTGER2) also known as EP2 is a G_s coupled receptor which activates adenylate cyclase (3-4). PTGER2 receptors are widely distributed throughout smooth muscle including the myometrium where they mediate relaxation (5). In addition, the Prostaglandin E2 receptor may be involved in bronchodilation (6). Prostaglandin E2 receptor deficient mice developed profound hypertension when on a high salt diet and had impaired ovulation and fertility (7-8).

Validation Summary

Performance of this assay was validated under optimized conditions in 384-well format using LiveBLAzer™-FRET B/G Substrate.

1. Prostaglandin E₂ agonist dose response under optimized conditions

	DA cells	Dividing Cells
EC ₅₀	2 nM	2 nM
Z'-factor	0.89	0.75

Optimum cell no.	= 10K cells/well
Optimum [DMSO]	= up to 1%
Stimulation Time	= 5 hours
Max. [Stimulation]	= 5 μM

2. Alternate agonist dose response

Butaprost EC ₅₀	= 100 nM
Prostaglandin D ₂ EC ₅₀	= 148 μM
AH13205 EC ₅₀	= 64 nM

3. Antagonist dose response

See Antagonist Section

4. Agonist 2nd messenger dose response

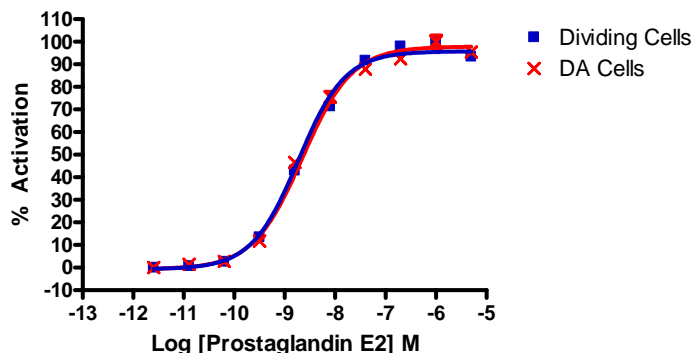
Prostaglandin E ₂ EC ₅₀	= 451 pM
---	----------

Assay Testing Summary

- Assay performance with variable cell number
- Assay performance with variable stimulation time
- Assay performance with variable substrate loading time
- Assay performance with variable DMSO concentration

Primary Agonist Dose Response

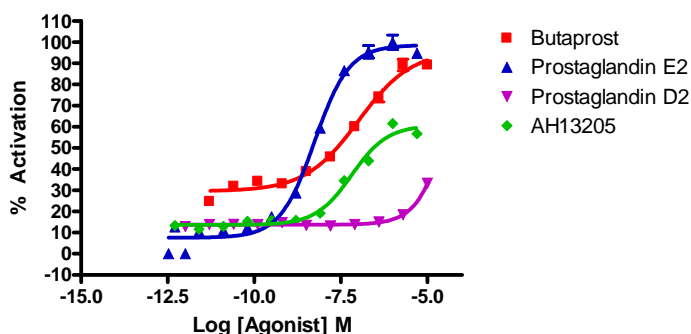
Figure 1 — GeneBLAzer® PTGER2 CHO-K1 DA and GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 dose response to Prostaglandin E₂ under optimized conditions



GeneBLAzer® PTGER2 CHO-K1 DA cells and GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 cells (10,000 cells/well) were plated in a 384-well format and incubated for 16-20 hours. Cells were stimulated with a dilution series of Prostaglandin E₂ in the presence of 0.5% DMSO for 5 hours. Cells were then loaded with LiveBLAzer™-FRET B/G Substrate for 2 hours. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and % Activation plotted against the concentrations of Prostaglandin E₂ (n=6 for each data point).

Alternate Agonist Dose Response

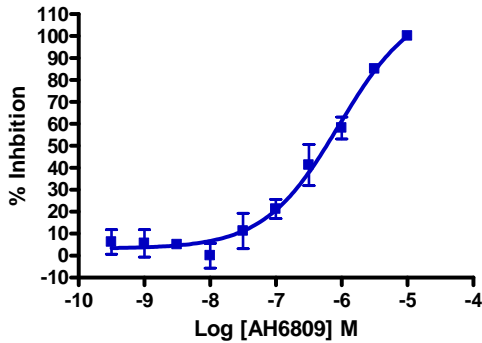
Figure 2 — GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 dose response to Prostaglandin D₂, BW 245C, and Prostaglandin E₂



GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 cells (10,000 cells/well) were plated the day before the assay in a 384-well format. Cells were stimulated with either Prostaglandin D₂ (Sigma #P5172), Butaprost (Sigma #B6309), AH13205 (Sigma #A9102), or Prostaglandin E₂ (Sigma #P5640) over the indicated concentration range in the presence of 0.5% DMSO for 5 hours. Cells were then loaded with LiveBLAzer™-FRET B/G Substrate for 2 hours. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the % Activation plotted against the indicated concentrations of the agonists (n= 8 for each data point). The data shows the correct rank order potency for these agonists.

Antagonist Dose Response

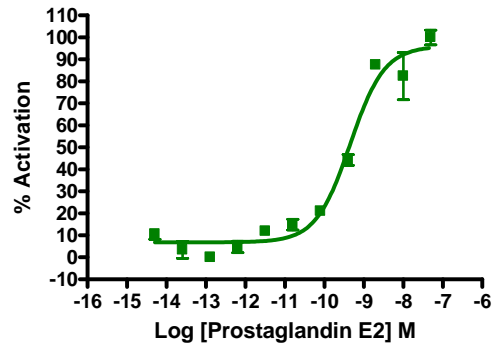
Figure 3 — GeneBLazer® PTGER2-CRE-*bla* CHO-K1 dose response to AH6809



GeneBLazer® PTGER2-CRE-*bla* CHO-K1 cells (10,000 cells/well) were plated the day before the assay in a 384-well black-walled tissue culture assay plate. Cells were treated with AH6809 (Sigma #A1221) and incubated at 37 degrees C for 30 min., followed by 20 nM Prostaglandin E₂ agonist stimulation for 5 hours in 0.5% DMSO. Cells were then loaded for 2 hours with LiveBLazer™-FRET B/G Substrate. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the % Inhibition is shown plotted against the indicated concentrations of the antagonist. AH6809 is known to be most active at DP receptors and weakly active at EP2 receptors.

Agonist 2nd Messenger Response

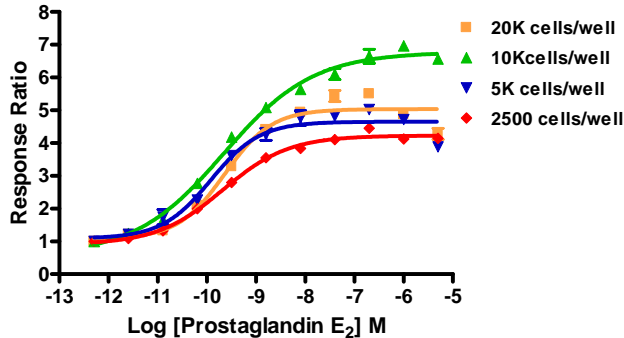
Figure 4— GeneBLazer® PTGER2-CRE-*bla* CHO-k1 2nd messenger dose response to Prostaglandin E2 under optimized conditions



GeneBLazer® PTGER2-CRE-*bla* CHO-K1 cells were tested for a response to Prostaglandin E2 with a TR-FRET cAMP assay.

Assay Performance with Variable Cell Number

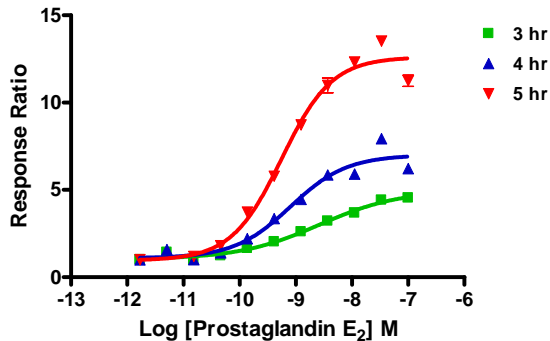
Figure 5— GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 dose response to Prostaglandin E₂ with 2.5, 5, 10 and 20K cells/well



GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 cells were plated in assay medium the day before the assay at 2,500, 5,000, 10,000 or 20,000 cells/well in a 384-well format. On the day of the assay, cells were stimulated with Prostaglandin E₂ (Sigma #P5640) in the presence of 0.5% DMSO for 5 hours. Cells were then loaded with LiveBLAzer™-FRET B/G Substrate for 2 hours. Fluorescence emission values at 460 nm and 530 nm for the various cell numbers were obtained using a standard fluorescence plate reader and the Response Ratios plotted against the indicated concentrations of Carbachol (n=8 for each data point).

Assay performance with Variable Stimulation Time

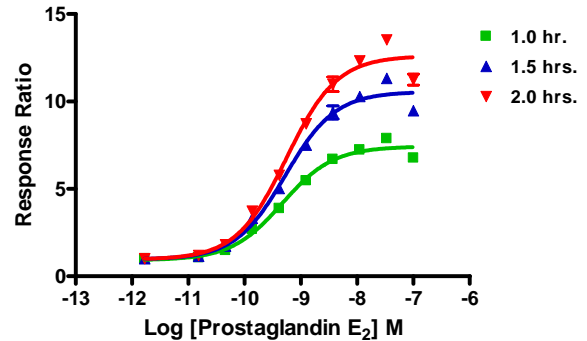
Figure 6 – GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 dose response to Prostaglandin E₂ with 3, 4 and 5 hour stimulation times



GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 cells (10,000 cells/well) were plated the day before the assay in a 384-well black-walled tissue culture assay plate. Prostaglandin E₂ (Sigma #P5640) was then added to the plate over the indicated concentration range for 3, 4, or 5 hrs in 0.5% DMSO and then loaded for 2 hours with LiveBLAzer™-FRET B/G Substrate. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the Response Ratios plotted against the indicated concentrations of carbachol (n=16 for each data point).

Assay performance with Variable Substrate Loading Time

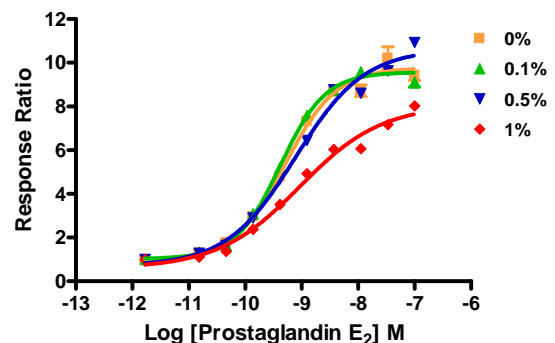
Figure 7 – GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 dose response to Prostaglandin E₂ with 1, 1.5 and 2 hour loading times



GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 cells were plated at 10,000 cells/well in a 384-well format the day before the assay. On the day of the assay, cells were stimulated with Prostaglandin E₂ (Sigma #P5640) in the presence of 0.5% DMSO for 5 hours. Cells were then loaded with LiveBLAzer™-FRET B/G Substrate for either 1, 1.5, and 2 hours. Fluorescence emission values at 460 nm and 530 nm for the various loading times were obtained using a standard fluorescence plate reader and the Response Ratios plotted against the indicated concentrations of Prostaglandin D₂ (n=16 for each data point).

Assay Performance with variable DMSO concentration

Figure 8 – GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 dose response to Prostaglandin D₂ with 0, 0.1, 0.5 and 1% DMSO



GeneBLAzer® PTGER2-CRE-*bla* CHO-K1 cells (10,000 cells/well) were plated the day before the assay in a 384-well black-walled tissue culture assay plate. Prostaglandin D₂ (Sigma #P5640) was then added to the plate over the indicated concentration range. DMSO was added to the assay at concentrations from 0% to 1%. Cells were stimulated for 5 hrs with agonist and loaded for 2 hours with LiveBLAzer™-FRET B/G Substrate. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the Response Ratios are shown plotted for each DMSO concentration against the indicated concentrations of carbachol (n=8 for each data point).

References

1. Coleman, R.A., Smith, W.L. and Narumiya, S. (1994) **VIII. International Union of Pharmacology classification of prostanoid receptors: properties, distribution and structure of the receptors and their subtypes.** *Pharmacol. Rev.*, **46**, 205-229.
2. Narumiya, S., Sugimoto, Y. and Ushikubi, F. (1999) **Prostanoid receptors: structures, properties, and functions.** *Physiol. Rev.*, **79**, 1193-1226.
3. Alvarez, R., Taylor, A., Fazzari, J.J. and Jacobs, J.R. (1981) **Regulation of cyclic AMP metabolism in human platelets. Sequential activation of adenylate cyclase and cyclic AMP phosphodiesterase by prostaglandins.** *Mol. Pharmacol.*, **20**, 302-309.
4. Breyer, R.M., Bagadassarian, C.K., Myers, S.A. and Breyer, M.D. (2001) **Prostanoid receptors: subtypes and signaling.** *Annu. Rev. Pharmacol. Toxicol.*, **41**, 661-690.
5. Senior, J., Marshall, K., Sangha, R., Baxter, G.S. and Clayton, J.K. (1991) **In vitro characterization of prostanoid EP-receptors in the non-pregnant human myometrium.** *Br. J. Pharmacol.* **102**, 747-753.
6. Sheller, J.R., Mitchell, D., Meyrick, B., Oates, J. and Breyer, R. (2000) **EP(2) receptor mediates bronchodilation by PGE(2) in mice.** *J. Appl. Physiol.* **88**, 2214-2218.
7. Kennedy, C.R., Zhang, Y., Brandon, S., Guan, Y., Coffee, K., Funk, C.D., Magnuson, M.A., Oates, J.A., Breyer, M.D. and Breyer, R.M. (1999) **Salt-sensitive hypertension and reduced fertility in mice lacking the prostaglandin EP₂ receptor.** *Nat. Med.* **5**, 217-220.
8. Hizaki, H., Segi, E., Sugimoto, Y., Hirose, M., Saji, T., Ushikubi, F., Matsuoka, T., Noda, Y., Tanaka, T., Yoshida, N., Narumiya, S. and Ichikawa, A. (2000) **Abortive expansion of the cumulus and impaired fertility in mice lacking the prostaglandin E receptor subtype EP₂.** *Proc. Natl. Acad. Sci. U.S.A.* **96**, 10501-10506.