

Optimization of the GeneBLAzer® ADRB3 CRE-*bla* CHO-K1 Cell Line**GeneBLAzer® ADRB3 CHO-K1 DA Cells****GeneBLAzer® ADRB3-CRE-*bla* CHO-K1 Cells**

Catalog Numbers – K1473 and K1745

Cell Line Descriptions

GeneBLAzer® ADRB3 CHO-K1 DA (Division Arrested) cells and GeneBLAzer® ADRB3-CRE-*bla* CHO-K1 cells contain the human Adrenrgic Beta 3 Receptor (ADRB3) (Accession # [NM_000025.1](#)) 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).

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® ADRB3 CHO-K1 DA cells and GeneBLAzer® ADRB3-CRE-*bla* CHO-K1 cells are functionally validated for Z'-factor and EC₅₀ concentrations of Isoproterenol, (Figure 1). In addition, GeneBLAzer® ADRB3-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

Adrenoceptors are a class of GPCRs that mediate the actions of the neurotransmitter noradrenalin and the hormone/neurotransmitter adrenaline. The adrenoceptors were separated into two major types, based on the rank order potencies of their agonist's, into α and β subtypes (1). The α subtype has been further broken down into 1α , 2α , and 3α subtypes based on molecular and pharmacological evidence (2). Initially the β -adrenoceptors were divided into two subtypes. The β_1 receptor is dominant in the heart and adipose tissue and is equally activated by noradrenaline and adrenaline. The β_2 receptor is responsible for relaxation of uterine, vascular and airway smooth muscle and is less sensitive to noradrenaline than adrenaline (3). The β_3 receptor is atypical as it is insensitive to common β -antagonists (4, 5).

Adrenoceptor B₃ (ADRB3) is a major factor in lipolysis and thermogenesis (6). These receptors are also involved in relaxation of the colon (7, 8). ADRB3 targeted drugs may have a negative inotropic effect (9), meaning that these drugs affect the force in which the heart contracts.

ADRB3 receptors distributed widely throughout the body, however they are found most frequently in adipose tissue and the gall bladder. Other locations where ADRB3 receptors are found is the small intestine, stomach, prostate, left atrium and bladder (10, 11).

Validation Summary

Testing and validation of this assay was evaluated in a 384-well format using LiveBLAzer™-FRET B/G Substrate.

1. Isoproterenol dose response under optimized conditions

	<u>DA cells</u>	<u>Dividing Cells</u>
EC ₅₀	5.5 nM	4 nM
Z'-factor	0.75	0.66

Recommended cell no.	= 10K cells/well
Recommended [DMSO]	= up to 1%
Recommended Stim. Time	= 5 hours
Max. [Stimulation]	= 600 nM

2. Alternate agonist dose response

(-)-Epinephrine EC ₅₀	= 0.84 nM
BRL 37344 EC ₅₀	= 0.74 nM

3. Antagonist dose response

ICI118551 IC ₅₀	= 19 μM
----------------------------	---------

4. Agonist 2nd Messenger Dose Response

Isoproterenol EC ₅₀	= 1.1 nM
--------------------------------	----------

Assay Testing Summary

5. Assay performance with variable cell number

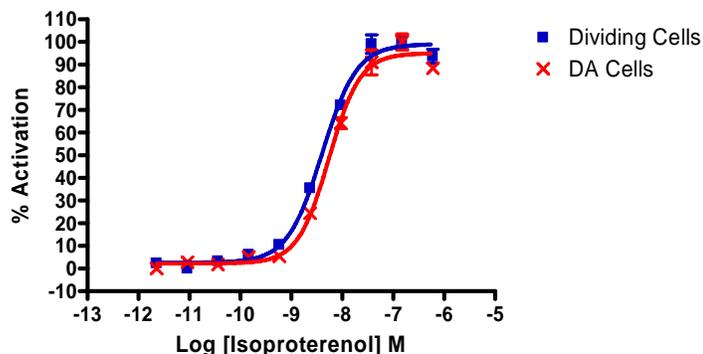
6. Assay performance with variable stimulation time

7. Assay performance with variable substrate loading time

8. Assay performance with variable DMSO concentration

Primary Agonist Dose Response

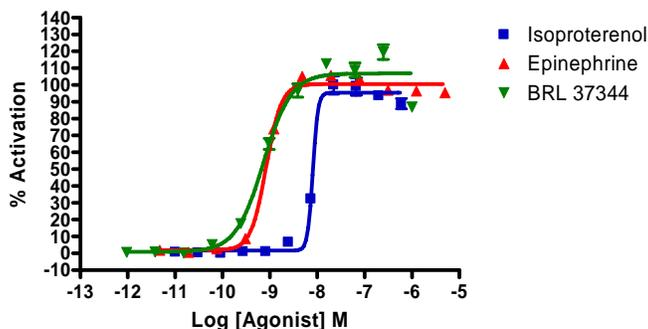
Figure 1 — GeneBLAzer® ADRB3-CRE-*bla* CHO-K1 and GeneBLAzer® ADRB3 CHO-K1 DA Isoproterenol dose response under optimized conditions



GeneBLAzer® ADRB3 CHO-K1 DA and GeneBLAzer® ADRB3-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 Isoproterenol (Sigma #I5627) 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 for each replicate against the concentrations of Isoproterenol (n=16 for each data point).

Alternate Agonist Dose Response

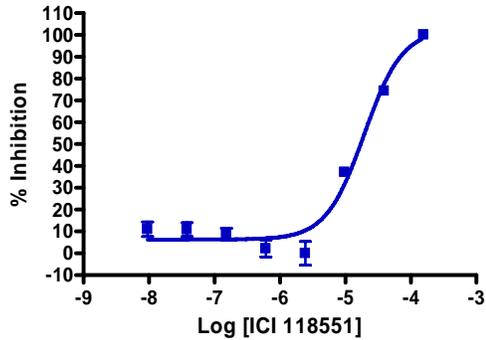
Figure 2 — GeneBLAzer® Isoproterenol, (-)-Epinephrine and BRL 37344 Dose Response



GeneBLAzer® ADRB3-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 dilution series of Isoproterenol (Sigma #I5627), BRL 37344 (Sigma #B169) and (-)-Epinephrine (Sigma #E4250) 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 shown plotted against the concentrations of Isoproterenol, BRL 37344, and (-)-Epinephrine (n=8 for each data point). The cell lines show the correct rank order potency for these compounds.

Antagonist Dose Response

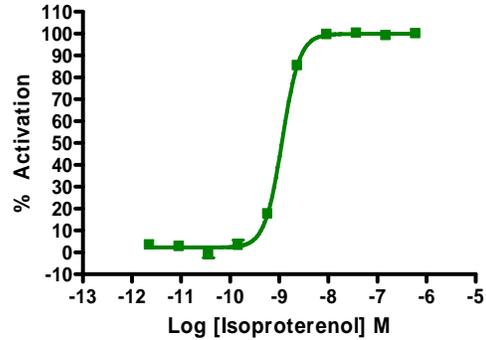
Figure 3 — GeneBLAzer® ADRB3-CRE-*bla* CHO-K1 ICI 118,551 dose response



GeneBLAzer® ADRB3-CRE-*bla* CHO-K1 cells were plated 16-20 hours prior to assay at 10,000 cells per well in a 384-well format. Cells were treated with a dilution series of ICI 118,551 (Sigma #I1127) in the presence of 0.25% DMSO. Cells were then incubated at 37°C & 5% CO₂ for 30 min. Isoproterenol (Sigma #I5627) was then added to the plate at the EC₈₀ concentration of 26 nM along with 0.25% DMSO (0.5% final concentration). Cells were incubated for 5 hours 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 % Inhibition shown plotted against the concentrations of the antagonists. This data shows the proper functioning of the assay in antagonist mode. (n=16 for each data point).

Agonist 2nd Messenger Dose Response

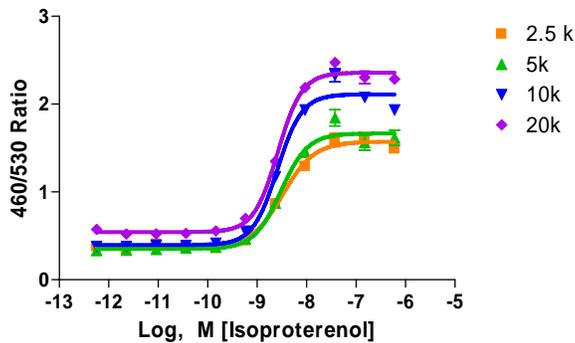
Figure 4— GeneBLAzer® ADRB3-CRE-*bla* CHO-k1 2nd messenger dose response to Isoproterenol under optimized conditions.



GeneBLAzer® ADRB3-CRE-*bla* CHO-K1 cells were tested for a response to isoproterenol with a TR-FRET cAMP assay

Assay Performance with Variable Cell Number

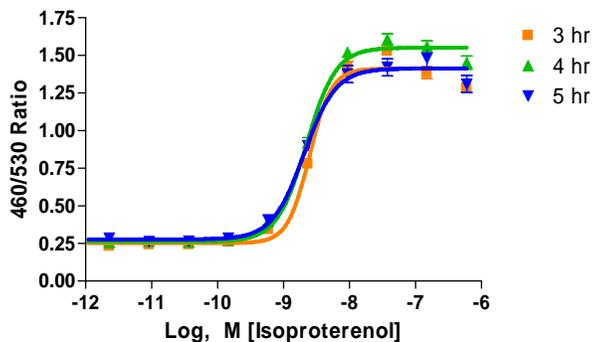
Figure 4 – Isoproterenol dose response with 2.5, 5, 10, and 20K cells/well



GeneBLAzer[®] ADRB3-CRE-*bla* CHO-K1 cells were plated at 2,500 5,000 10,000 or 20,000 cells/well in a 384-well format and incubated for 16-20 hours. Cells were stimulated with a dilution series of Isoproterenol (Sigma #I5627) 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 and plotted for each cell number against the concentrations of Isoproterenol (n=8 for each data point).

Assay Performance with Variable Stimulation Time

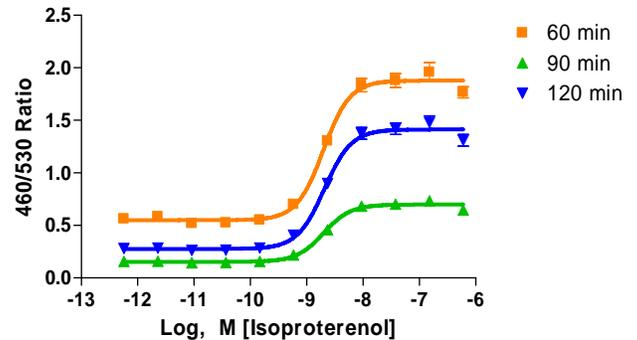
Figure 5 – Isoproterenol dose response with 3, 4 and 5 hr stimulation times



GeneBLAzer[®] ADRB3-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 Isoproterenol (Sigma #I5627) for 3, 4, or 5 hrs in the presence of 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 plotted for each stimulation time against the concentrations of Isoproterenol (n=8 for each data point).

Assay Performance with Variable Substrate Loading Times

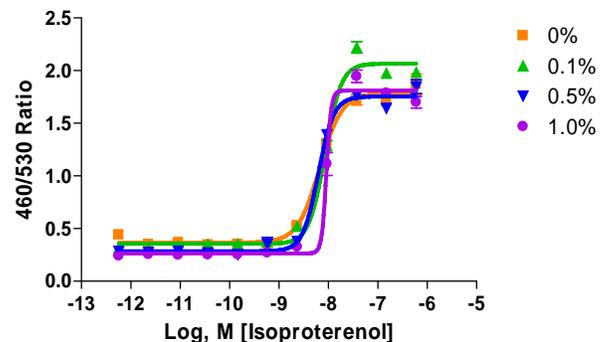
Figure 6 – GeneBLAzer[®] Isoproterenol dose response with 1, 1.5, and 2 hour substrate loading times.



GeneBLAzer[®] ADRB3-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 Isoproterenol (Sigma #I5627) in the presence of 0.5% DMSO for 5 hours. Cells were then loaded for either 1, 1.5 or 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 plotted for each substrate loading time against the concentrations of Isoproterenol (n=8 for each data point).

Assay Performance with Variable DMSO Concentration

Figure 7 – GeneBLAzer[®] Isoproterenol dose response with 0, 0.1, 0.5 and 1% DMSO



GeneBLAzer[®] ADRB3-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 Isoproterenol (Sigma I5627) for 5 hours. DMSO was added to the cells at concentrations from 0% to 1%. 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 plotted for each DMSO concentration against the concentrations of Isoproterenol (n=8 for each data point).

References

1. Ahlquist, R.P. **A study of adrenotropic receptors.** *AM. J Physiol.* **153**, 586-600 (1948).
2. Bylund, D.B. **Subtypes of α_2 - adrenoceptors: pharmacological and molecular biological evidence converge.** *Trends Pharmacol. Sci.* **9**, 356-361 (1988).
3. Lands, A. M., et al. **Differentiation of receptor systems activated by sympathomimetic amines.** *Nature*, **214**, 597 - 598. (1967).
4. Arch, J. R. S. *et al.*, **Atypical β -adrenoceptors on brown adipocytes as a target for anti-obesity drugs.** *Nature*, **309**, 163 - 165. (1984).
5. Bond, R. A. and Clarke, D. E. **Agonist and antagonist characterization of a putative adrenoceptor with distinct pharmacological properties from the alpha- and beta-subtypes.** *Br. J. Pharmacol.*, **95**, 723 - 734. (1988)
6. Jockers, R., et al. **Desensitization of the beta-adrenergic response in human brown adipocytes.** *Endocrinology.*, **139**, 2676 - 2684. (1998).
7. Severi, C., et al. **Beta3-adrenoceptors: relaxant function and mRNA detection in smooth muscle cells isolated from the human colon.** *Can J Physiol Pharmacol.*, **82**, 515 - 522. (2004).
8. Bardou, M., et al. **In vitro inhibition of human colonic motility with SR 59119A and SR 59104A: evidence of a beta3-adrenoceptor-mediated effect.** *Eur J Pharmacol.*, **353**, 281 - 287. (1998)
9. Gauthier, C., et al. **Functional beta3-adrenoceptor in the human heart.** *J Clin Invest.*, **98**, 556 - 562. (1996)
10. Krief, S., et al. **Tissue distribution of beta 3-adrenergic receptor mRNA in man.** *J Clin Invest.*, **91**, 344 - 349. (1993).
11. Berkowitz, D. E., et al. **Distribution of beta 3-adrenoceptor mRNA in human tissues.** *Eur J Pharmacol.*, **289**, 223 - 228. (1995).