

**GeneBLAzer® M4-Gqo5 CHO-K1 DA Assay Kit****GeneBLAzer® M4-Gqo5 NFAT-*bla* CHO-K1 Cells**

Catalog Numbers – K1377 and K1741

**Cell Line Descriptions**

GeneBLAzer® M4-Gqo5 CHO-K1 DA (Division Arrested) cells and GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 cells contain the human Acetylcholine (muscarinic) subtype 4 receptor (M4), (Accession #[NM\\_000741](#)) stably integrated into the Gqo5-NFAT-*bla* CHO-K1 cell line. Gqo5-NFAT-*bla* CHO-K1 cells (Cat. No. K1536) contain a beta-lactamase (*bla*) reporter gene under control of the Nuclear Factor of Activated T Cells (NFAT) response element and the chimeric G protein, Gqo5. 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® M4-Gqo5 CHO-K1 DA cells and GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 cells are functionally validated for Z'-factor and EC<sub>50</sub> concentrations of carbachol, (Figure 1). In addition, GeneBLAzer® M4-Gqo5-NFAT-*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**

Muscarinic acetylcholine receptors are members of the G protein-coupled receptor (GPCR) superfamily. Muscarinic receptors are widely distributed and mediate the actions of acetylcholine in both the CNS and peripheral tissues. Five muscarinic receptor subtypes have been identified and are referred to as M<sub>1</sub>-M<sub>5</sub> (1-5). The five genes that encode the muscarinic receptors all belong to the rhodopsin-like family (Family A) and share strong sequence homology but have unique regions located at the amino terminus (extracellular) and in the third intracellular loop.

The M<sub>1</sub>, M<sub>3</sub>, and M<sub>5</sub> receptor subtypes couple through the G<sub>q/11</sub> class of G-proteins and activate the phospholipase C pathway. One downstream target activated by the phospholipase C pathway is the transcription factor, NFAT. The M<sub>2</sub> and M<sub>4</sub> receptor subtypes couple through the G<sub>i/o</sub> class of G proteins and inhibit adenylyl cyclase activity.

The clinical implications of the M<sub>4</sub> receptor are unknown; however, studies with knockout mice suggest that M<sub>4</sub> may have implications in dopaminergic function, keratinocyte migration and wound healing, anxiolysis, and analgesia (6-11). In addition, M<sub>4</sub> may be important in memory circuits and have implications in Alzheimer's Disease (12). Additional information on the muscarinic receptors can be found in reviews (13-16).

## Validation Summary

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

### 1. Carbachol agonist dose response under optimized conditions

	DA cells	Dividing Cells
EC <sub>50</sub>	139 nM	176 nM
Z'-factor	0.64	0.60

Recommended cell no.	= 10K cells/well
Recommended [DMSO]	= up to 1%
Recommended Stim. Time	= 5 hours
Max. [Stimulation]	= 3 μM

### 2. Alternate agonist dose response

Oxotremorine M EC <sub>50</sub>	= 31 nM
MCN-A-343 EC <sub>50</sub>	= 844 nM
Bethanecol EC <sub>50</sub>	= 25 μM
Pilocarpine EC <sub>50</sub>	= 66 μM

### 3. Antagonist dose response

Scopolamine IC <sub>50</sub>	= 374 pM
Telenzipine IC <sub>50</sub>	= 3.3 nM
DAMP IC <sub>50</sub>	= 5.1 nM
Methoctramine IC <sub>50</sub>	= 769 nM

### 4. Agonist Dose Response Using Fluo-4NW

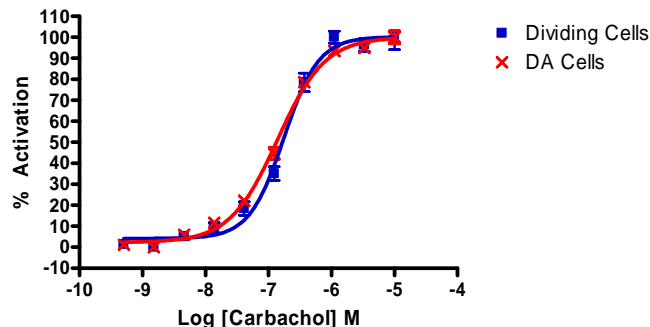
Carbachol EC <sub>50</sub>	= 406 nM
----------------------------	----------

## 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

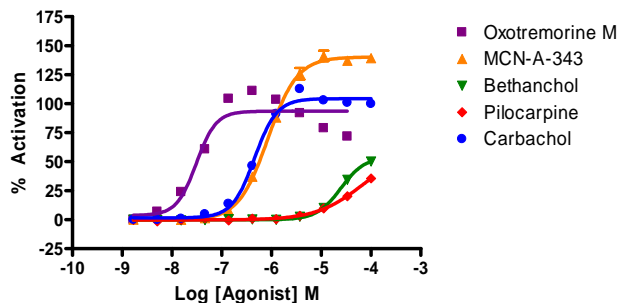
**Figure 1 — GeneBLAzer® M4-Gqo5 CHO-K1 DA and M4-Gqo5-NFAT-*bla* CHO-K1 dose response to carbachol under optimized conditions**



GeneBLAzer® M4-Gqo5 CHO-K1 DA cells and GeneBLAzer® M4-Gqo5-NFAT-*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 carbachol in the presence of 0.5% DMSO for 5 hours. Cells were then loaded with LiveBLAzer™-FRET B/G Substrate (2μM final concentration of CCF4-AM + 1mM Solution D) 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 carbachol (n=6 for each data point).

## Alternate Agonist Dose Response

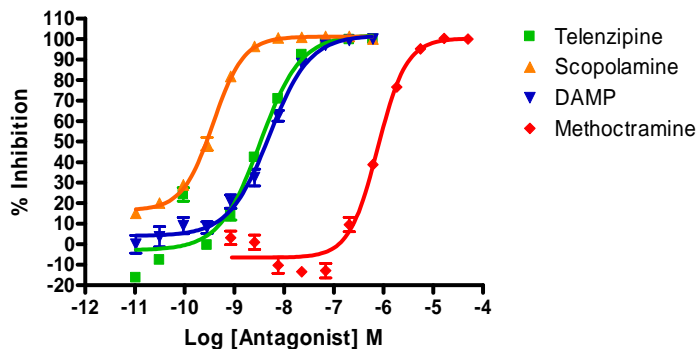
**Figure 2 — GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 dose response to Bethanecol, Oxotremorine M, MCN-A-343, Pilocarpine and Carbachol**



GeneBLAzer® M4-Gqo5-NFAT-*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 Carbachol (Sigma #21760), Bethanecol (Sigma #C5259), Oxotremorine (Sigma #O-100), MCN-A-343 (Sigma #C7041), or Pilocarpine (Sigma #P6503) 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 is shown plotted against the indicated concentrations of the agonists.

### Antagonist Dose Response

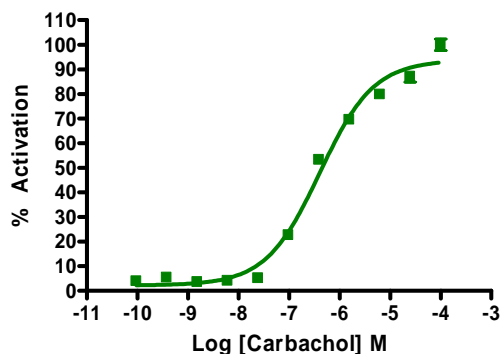
Figure 3 — GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 dose response to Telenzipine, Scopolamine, DAMP and Methoctramine



GeneBLAzer® M5-NFAT-*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 the antagonists Telenzipine (T-122), Scopolamine (S1875), DAMP (D104) and Methoctramine (M-105) and incubated at 37 degrees C for 20 min., followed by 1.5  $\mu$ M Carbachol 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 was plotted against the indicated concentrations of the antagonist. The data shows the correct rank order potency.

### Agonist Dose Response Using Fluo-4NW

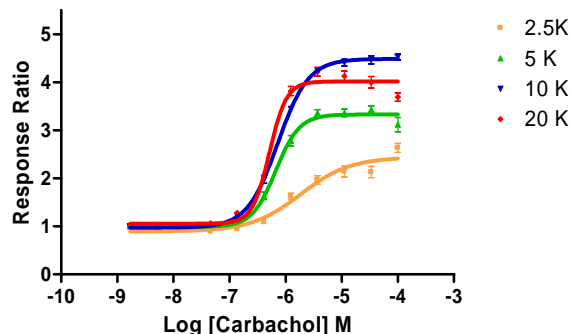
Figure 4 — GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 dose response to carbachol using Fluo-4NW



GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 cells (10,000 cells/well) were plated the day before the assay in a 384-well format. Cells were incubated with Fluo-4NW for 30 min. at 37°C., followed by 30 min. at room temperature. Cells were then stimulated with a dilution series of Carbachol (Sigma #21760) in the presence of 0.5% DMSO. Fluorescence emission values at 516 nm were obtained and the % Activation plotted against the indicated concentrations of carbachol (n=16 for each data point).

### Assay Performance with Variable Cell Number

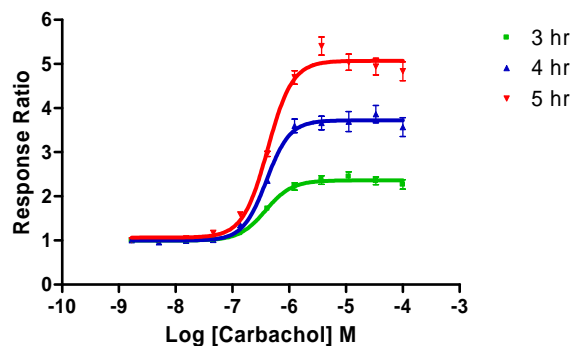
Figure 5— GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 dose response to Carbachol with 2.5, 5, 10, and 20K cells/well



GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 cells were plated one day before the assay at 2,500, 5000, 10,000 or 20,000 cells/well in a 384-well format. On the day of the assay, cells were stimulated with Carbachol (Sigma #21760) in the presence of 0.5% DMSO for 5 hours. Cells were then loaded with LiveBLAzer™-FRET B/G Substrate (2µM final concentration of CCF4-AM + 1mM Solution D) 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

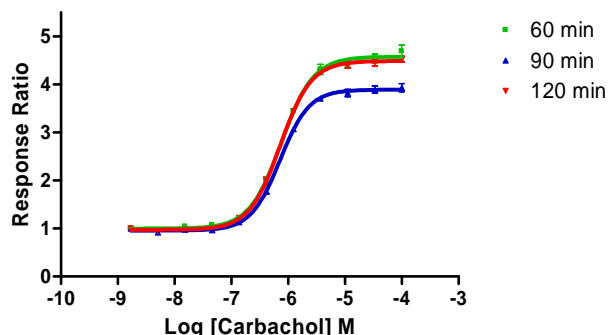
Figure 6 – GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 dose response to Carbachol with 3, 4, and 5 hour stimulation times



GeneBLAzer® M4-Gqo5-NFAT-*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. Carbachol (Sigma #21760) 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 (2µM final concentration of CCF4-AM + 1mM Solution D). Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the Response Ratios for each substrate loading time were plotted against the indicated concentrations of carbachol (n=8 for each data point)

### Assay performance with Variable Substrate Loading Time

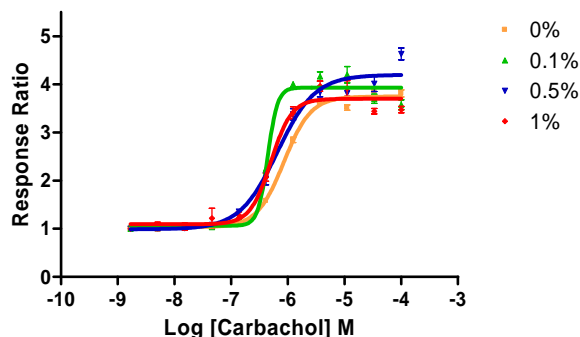
Figure 7 – GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 dose response to Carbachol with 1, 1.5, and 2 hr substrate loading times



GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 cells were plated at 10,000 cells/well in a 384-well format one day before the assay. On the day of the assay, cells were stimulated with Carbachol (Sigma #21760) in the presence of 0.5% DMSO for 5 hours. Cells were then loaded with LiveBLAzer™-FRET B/G Substrate (2µM final concentration of CCF4-AM + 1mM Solution D) for either 1, 1.5, or 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 DMSO concentration

Figure 8 – GeneBLAzer® M4-Gqo5-NFAT-*bla* CHO-K1 dose response to Carbachol with 0, 0.1, 0.5 and 1% DMSO.



GeneBLAzer® M4-Gqo5-NFAT-*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. Carbachol (Sigma #21760) 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 (2µM final concentration of CCF4-AM + 1mM Solution D). Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the Response Ratios are for each DMSO concentration were plotted against the indicated concentrations of carbachol (n=8 for each data point).

## References

1. Bonner, T. I., *et al* (1987) **Identification of a family of muscarinic acetylcholine receptor genes.** *Science*, **237**, 527 - 532.
2. Bonner, T. I., *et al* . (1988) **Cloning and expression of the human and rat m5 muscarinic acetylcholine receptor genes.** *Neuron*, **1**, 403 - 410.
3. Kubo, T. *et al.* (1986) **Cloning, sequencing and expression of complementary DNA encoding the muscarinic acetylcholine receptor.** *Nature*, **323**, 411 - 416.
4. Kubo, T. *et al.* (1986) **Primary structure of porcine cardiac muscarinic acetylcholine receptor deduced from the cDNA sequence.** *FEBS Lett.*, **209**, 367 - 372.
5. Peralta, E.G., *et al.* (1987) **Distinct primary structures, ligand binding properties and tissue-specific expression of four human muscarinic acetylcholine receptors.** *EMBO J.*, **6**, 3923 – 3929.
6. Bymaster, F.P., Carter, P.A., Zhang, L., Falcone, J.F., Stengel, P.W., Cohen, M.L., Shannon, H.E., Gomeza, J., Wess, J. and Felder, C.C. (2001) **Investigations into the physiological role of muscarinic M2 and M4 muscarinic and M4 receptor subtypes using receptor knockout mice.** *Life Sci.*, **68**, 2473-2479.
7. Bymaster, F.P., McKinzie, D.L., Felder, C.C. and Wess, J. (2003) **Use of M1-M5 muscarinic receptor knockout mice as novel tools to delineate the physiological roles of the muscarinic cholinergic system.** *Neurochem Res.*, **28**, 437-442.
8. Tzavara, E.T., Bymaster, F.P., Davis, R.J., Wade, M.R., Perry, K.W., Wess, J., McKinzie, D.L., Felder, C., and Nomikos, G.G. (2004) **M4 muscarinic receptors regulate the dynamics of cholinergic and dopaminergic neurotransmission: relevance to the pathophysiology and treatment of related CNS Pathologies.** *FASEB J.*, **18**, 1410-1412.
9. Chernvavsky, A.I., Arredondo, J., Wess, J., Karlsson, E. and Grando, S.A. (2004) **Novel signaling pathways mediating reciprocal control of keratinocyte migration and wound epithelialization through M3 and M4 muscarinic receptors.** *J. Cell Biol.*, **19**, 261-272.
10. Degroot, A. and Nomikos, G.G. (2006) **Genetic deletion of muscarinic M4 receptors is anxiolytic in the shock-probe burying model.** *Eur. J. Pharmacol.*, **531**, 183-186.
11. Chen, S.R., Wess, J., and Pan, H.L. (2005) **Functional activity of the M2 and M4 receptor subtypes in the spinal cord studied with muscarinic acetylcholine receptor knockout mice.** *J. Pharmacol. Exp. Ther.*, **313**, 765-770.
12. Levey, A. (1996) **Muscarinic acetylcholine receptor expression in memory circuits: Implications for treatment of Alzheimer disease.** *Proc. Natl. Acad. Sci. USA*, **93**, 13541-13546.
13. Caulfield, M. P. and Birdsall, N. J. M. (1998) **International Union of Pharmacology. XVII. Classification of muscarinic acetylcholine receptors.** *Pharmacol. Rev.*, **50**, 279 - 290.
14. Caulfield, M.P. (1993) **Muscarinic receptors characterization, coupling and function.** *Pharmacol. Ther.*, **58**, 319 - 379.
15. Eglén, R. M., Hegde, S. S. and Watson, N. (1996) **Muscarinic receptor subtypes and smooth muscle function.** *Pharmacol. Rev.*, **48**, 531 - 565.
16. Hulme, E. C., Birdsall, N. J. M. and Buckley, N. J. (1990) **Muscarinic receptor subtypes.** *Annu. Rev. Pharmacol. Toxicol.*, **30**, 633 - 673.