

---

**Optimization of the GeneBLAzer® M1-NFAT-*bla* CHO-K1 Cell Line**

---

**GeneBLAzer® M1 CHO-K1 DA Cells****GeneBLAzer® M1 NFAT-*bla* CHO-K1 Cells**

Catalog Numbers – K1365 and K1735

**Cell Line Descriptions**

GeneBLAzer® M1 CHO-K1 DA (Division Arrested) cells and GeneBLAzer® M1-NFAT-*bla* CHO-K1 cells contain the human Acetylcholine (muscarinic) subtype 1 receptor (M1), (Accession #NM\_000738) stably integrated into the CellSensor® NFAT-*bla* CHO-K1 cell line. CellSensor® NFAT-*bla* CHO-K1 cells (Cat. no. K1534) contain a beta-lactamase (*bla*) reporter gene under control of the Nuclear Factor of Activated T-cells (NFAT) response element. Division Arrested (DA) cells are available in 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® M1 CHO-K1 DA cells and GeneBLAzer® M1-NFAT-*bla* CHO-K1 cells are functionally validated for Z'-factor and EC<sub>50</sub> concentrations of carbachol (Figure 1). In addition, M1-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-line 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. Activation of this pathway in turn leads to increases in free intracellular calcium levels as inositol triphosphate mediates release of calcium from the endoplasmic reticulum. In addition, protein kinase C is activated via diacylglycerol. 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.

In the brain, M<sub>1</sub> activation mediates "slow" neuronal excitability. Cortical and hippocampal muscarinic receptors are thought to be important in the attentional aspects of cognition. The predominant receptor subtypes in these brain areas are M<sub>1</sub>, M<sub>3</sub>, and M<sub>4</sub>. Therefore, M<sub>1</sub> is a potential target for cognition, Alzheimer's, dementia, and schizophrenia (6). Studies on knock-out mouse models of M<sub>1</sub> are also beginning to reveal potential functions of the receptor (7-9). Additional information on the muscarinic receptors can be found in reviews (10-13).

## 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>	165 nM	105 nM
Z'-factor	0.92	0.87

Optimum cell no.	= 20K cells/well
Optimum [DMSO]	= up to 1%
Optimum Stim. Time	= 5 hours
Max. [Stimulation]	= 10 μM

### 2. Alternate agonist dose response

Bethanecol EC <sub>50</sub>	= 3 μM
Oxotrmorine M EC <sub>50</sub>	= 3.8 nM
MCN -A-343 EC <sub>50</sub>	= 162 nM
Pilocarpine EC <sub>50</sub>	= 366 nM

### 3. Antagonist dose response

Telenzipine IC <sub>50</sub>	= 10 nM
Scopolamine IC <sub>50</sub>	= 2.2 nM
DAMP IC <sub>50</sub>	= 22 nM
Methoctramine IC <sub>50</sub>	= 3.1 μM

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

Carbachol EC <sub>50</sub>	= 55 nM
----------------------------	---------

### 5. Agonist Dose Response with ERK2 Phosphorylation

Methacholine EC <sub>50</sub>	= 74.3 nM
Oxotrmorine M EC <sub>50</sub>	= 9.8 nM
MCN -A-343 EC <sub>50</sub>	= 208 nM
Carbachol EC <sub>50</sub>	= 20.9 nM

### 6. [<sup>3</sup>H] NMS saturation binding analysis

K <sub>D</sub> [ <sup>3</sup> H] NMS	= 0.1 nM
B <sub>max</sub> (pmol/mg)	= 2.0

### 7. Allosteric Modulation

## Assay Testing Summary

### 8. Assay performance with variable cell number

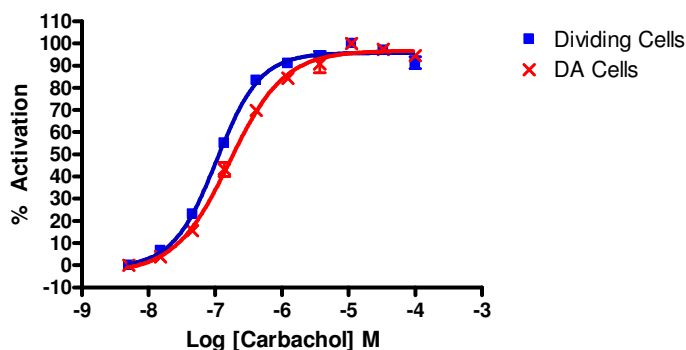
### 9. Assay performance with variable stimulation time

### 10. Assay performance with variable substrate loading time

### 11. Assay performance with variable DMSO concentration

## Primary Agonist Dose Response

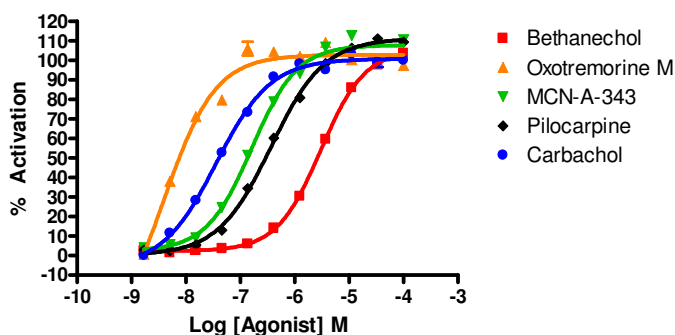
**Figure 1 — GeneBLazer® M1-NFAT-*bla* CHO-K1 DA and GeneBLazer® M1-NFAT-*bla* CHO-K1 dose response to carbachol under optimized conditions**



GeneBLazer® M1-NFAT-*bla* CHO-K1 DA cells and GeneBLazer® M1-NFAT-*bla* CHO-K1 cells (20,000 cells/well) were plated the day of the assay in a 384-well format and stimulated with Carbachol (Sigma #21760) 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 carbachol (n= 4 for each data point).

## Alternate Agonist Dose Response

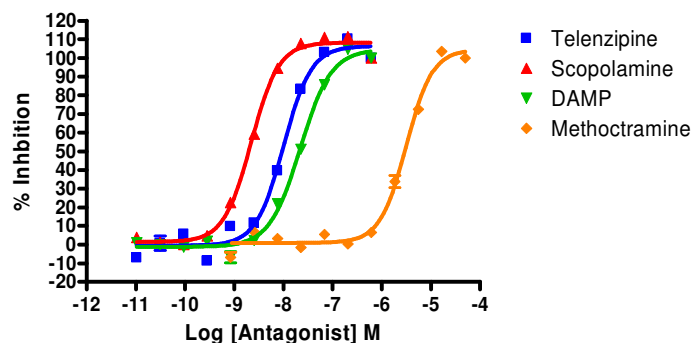
**Figure 2 — GeneBLazer® M1-NFAT-*bla* CHO-K1 dose response to Bethanechol, Oxotremorine M, MCN-A-343, Pilocarpine and Carbachol**



GeneBLazer® M1-NFAT-*bla* CHO-K1 cells (20,000 cells/well) were plated the day of the assay in a 384-well format. Cells were stimulated with either Carbachol (Sigma #21760), Bethanechol (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 plotted against the indicated concentrations of the agonists (n= 8 for each data point).

## Antagonist Dose Response

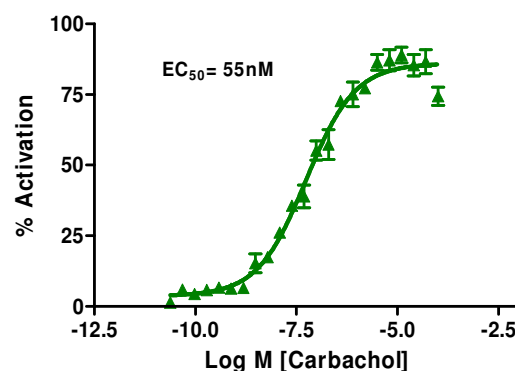
**Figure 3 — GeneBLazer® M1-NFAT-*bla* CHO-K1 dose response to Telenzipine, Scopolamine, DAMP and Methoctramine**



GeneBLazer® M1-NFAT-*bla* CHO-K1 cells (20,000 cells/well) were plated the day of the assay in a 384-well black-walled tissue culture assay plate. Cells were treated with telenzipine (Sigma #T-122), Scopolamine (Sigma #S1875), DAMP (Sigma #D104), or Methoctramine (Sigma #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 is shown plotted against the indicated concentrations of the antagonists. The data shows the correct rank order potency.

## Agonist Dose Response Using Fluo-4NW

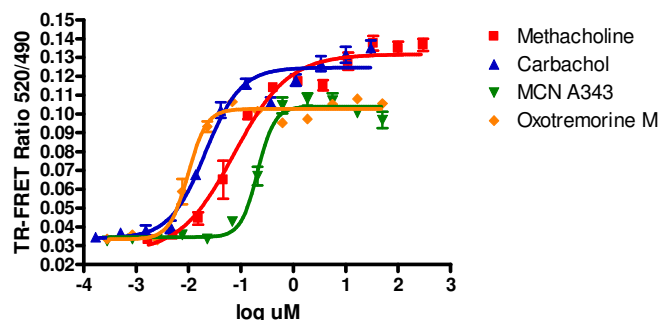
**Figure 4 — GeneBLazer® M1-NFAT-*bla* CHO-K1 dose response to Carbachol using Fluo-4NW**



GeneBLazer® M1-NFAT-*bla* CHO-K1 cells (10,000 cells/well) were plated the day before the assay in a 384-well format. Cells were then 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 plotted as % Activation against the indicated concentrations of carbachol (n=16 for each data point).

## ERK2 Phosphorylation

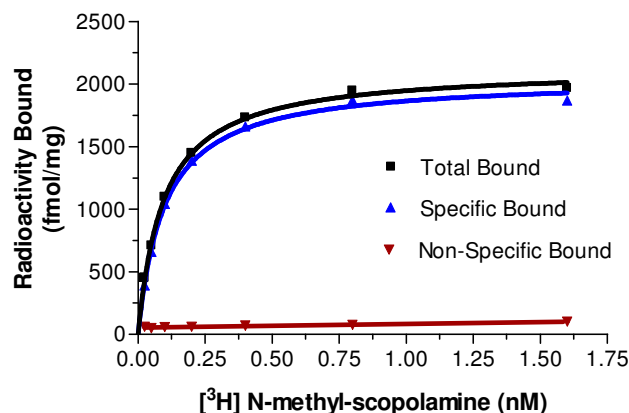
**Figure 5 — GeneBLazer® M1-NFAT-*bla* CHO-K1 agonist dose responses with the BacMam ERK2 [pThr185/pTyr187] Cellular Assay Kit**



GeneBLazer® M1-NFAT-*bla* CHO-K1 cells were assayed with the BacMam ERK2 [pThr185/pTyr187] Cellular Assay Kit. Transduced cells were stimulated for 6 minutes with a dilution series of either Methacholine, Carbachol, MCN-A-343, or Oxitremorine M prior to detection of ERK2 phosphorylation with the TR-FRET kit. Time resolved fluorescence emission values at 520 nm and 490 nm were obtained using a standard fluorescence plate reader and the TR-FRET ratio plotted.

## Radioligand Binding

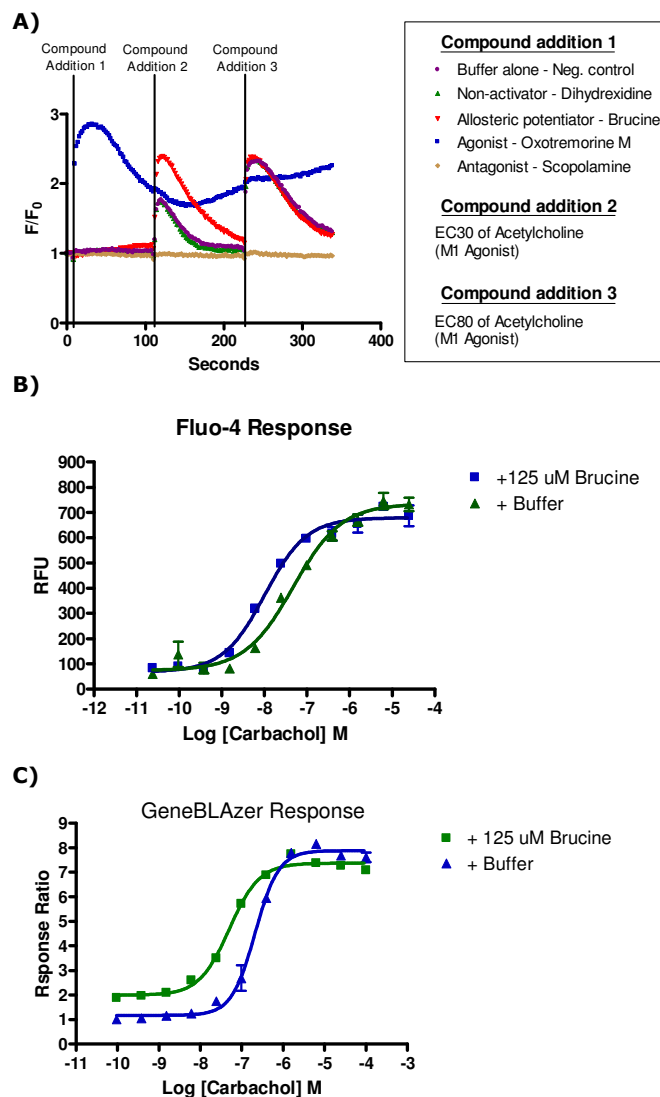
**Figure 6 — [<sup>3</sup>H] NMS saturation binding to GeneBLazer® M1-NFAT-*bla* CHO-K1 Membranes**



Saturation binding analysis was performed by incubating GeneBLazer® M1 NFAT-*bla* CHO-K1 membranes with increasing concentrations of [<sup>3</sup>H] N-methyl-scopolamine (NMS) in PBS, pH 7.4, with 100 µg/ml BSA. Non-specific binding was determined in the presence of 10 µM atropine. Samples were incubated for 2 hrs prior to filtration through a GF/B 96-well filter plate, which was pre-treated with 0.5% PEI. Filters were washed with cold 50 mM Tris-Cl, pH 7.4.

## Allosteric Modulation

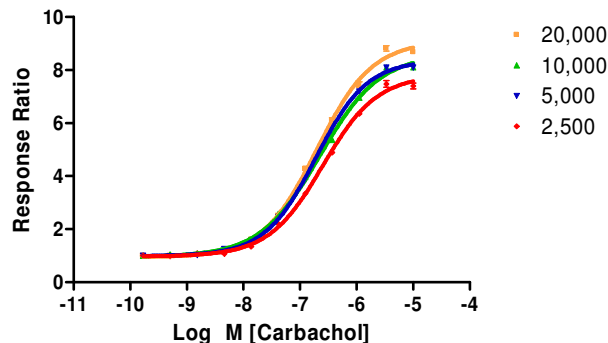
**Figure 7 — Allosteric modulation of GeneBLazer® M1-NFAT-*bla* CHO-K1 detected with Fluo-4 or the GeneBLazer® Assay**



GeneBLazer® M1-NFAT-*bla* CHO-K1 were assayed with either Fluo-4 (Figures A and B) or beta-lactamase (Figure C). Brucine is an allosteric modulator of the M1 muscarinic receptor which when applied allows carbachol to have greater potency at lower doses.

### Assay Performance with Variable Cell Number

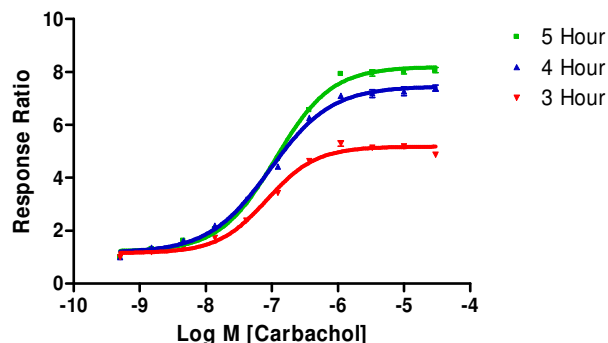
**Figure 8— GeneBLazer® M1-NFAT-*bla* CHO-K1 dose response to carbachol using 2.5, 5, 10, and 20K cells/well**



GeneBLazer® M1-NFAT-*bla* CHO-K1 cells were plated at 2500, 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 for 2.5 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

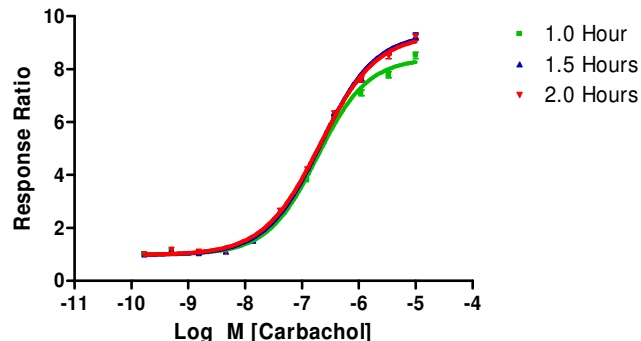
**Figure 9 – GeneBLazer® M1-NFAT-*bla* CHO-K1 dose response to carbachol using 3, 4, and 5 hour stimulation times.**



GeneBLazer® M1-NFAT-*bla* CHO-K1 cells (20,000 cells/well) were plated the day of 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. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the Response Ratios plotted (n=8 for each data point).

### Assay performance with Variable Substrate Loading Time

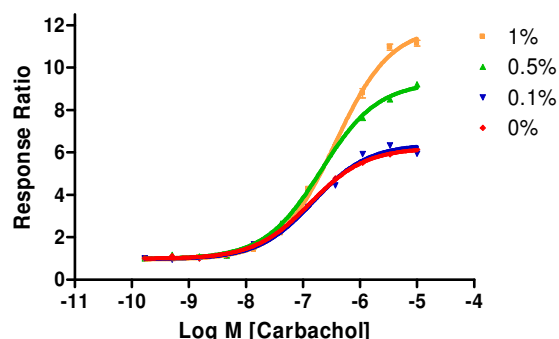
**Figure 10 – GeneBLazer® M1-NFAT-*bla* CHO-K1 dose response to carbachol using 1, 1.5, and 2 hour loading times**



GeneBLazer® M1-NFAT-*bla* CHO-K1 cells were plated at 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 for either 1, 1.5, and 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 11 – GeneBLazer® M1-NFAT-*bla* CHO-K1 dose response to carbachol using 0, 0.1, 0.5 and 1% DMSO.**



GeneBLazer® M1-NFAT-*bla* CHO-K1 cells (10,000 cells/well) were plated the day of 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. 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. Bonner, T. I., Buckley, N. J., Young, A. C. and Brann, M. R. (1987) **Identification of a family of muscarinic acetylcholine receptor genes.** *Science*, **237**, 527 - 532.
2. Bonner, T. I., Young, A. C., Brann, M. R. and Buckley, N. J. (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., Ashkenazi, A., Winslow, J. W., Smith, D. H., Ramachandran, J. and Capon, D. J. (1987) **Distinct primary structures, ligand binding properties and tissue-specific expression of four human muscarinic acetylcholine receptors.** *EMBO J.*, **6**, 3923 - 3929.
6. Levey, A.I. (1996) **Muscarinic acetylcholine receptor expression in memory circuits: Implications for treatment of Alzheimer's disease.** *Proc. Natl. Acad. Sci. U.S.A.*, **93**, 13541-13546.
7. Hamilton, S. E. et al. (1997) **Disruption of the m1 receptor gene ablates muscarinic receptor-dependent M current regulation and seizure activity in mice.** *Proc. Natl. Acad. Sci. U.S.A.*, **94**, 13311 - 13316.
8. Rouse, S. T., Hamilton, S. E., Potter, L. T., Nathanson, N. M. and Conn, P. J. (2000) **Muscarinic-induced modulation of potassium conductances is unchanged in mouse hippocampal pyramidal cells that lack functional M1 receptors.** *Neurosci. Lett.*, **278**, 61 - 64.
9. Shapiro, M. S., Loose, M. D., Hamilton, S. E., Nathanson, N. M., Gomeza, J., Wess, J. and Hille, B. (1999) **Assignment of muscarinic receptor subtypes mediating G-protein modulation of Ca<sup>2+</sup> channels by using knockout mice.** *Proc. Natl. Acad. Sci. U.S.A.*, **96**, 10899 - 10904.
10. Caulfield, M. P. and Birdsall, N. J. M. (1998) International Union of Pharmacology. XVII. **Classification of muscarinic acetylcholine receptors.** *Pharmacol. Rev.*, **50**, 279 - 290.
11. Caulfield, M.P. (1993) **Muscarinic receptors characterization, coupling and function.** *Pharmacol. Ther.*, **58**, 319 - 379.
12. Eglen, R. M., Hegde, S. S. and Watson, N. (1996) **Muscarinic receptor subtypes and smooth muscle function.** *Pharmacol. Rev.*, **48**, 531 - 565.
13. Hulme, E. C., Birdsall, N. J. M. and Buckley, N. J. (1990) **Muscarinic receptor subtypes.** *Annu. Rev. Pharmacol. Toxicol.*, **30**, 633 - 673.