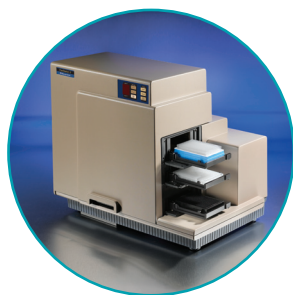


# G<sub>q</sub>-coupled receptor assays using the FlexStation System: M1 muscarinic receptor

APPLICATION BULLETIN NO. 1



**THE FLEXSTATION SYSTEM IS A  
VERSATILE TOOL FOR ASSAY  
DEVELOPMENT AND BASIC  
RESEARCH**

- SIMPLE, RAPID EXPERIMENTAL  
SET-UP AND DATA ANALYSIS
- DETERMINE EC<sub>50</sub> AND IC<sub>50</sub>  
VALUES
- EXCELLENT FOR ASSAY  
DEVELOPMENT AS  
DEMONSTRATED BY Z'  
FACTOR

## INTRODUCTION

G<sub>q</sub>-coupled receptors are transmembrane proteins that are responsible for transmitting extracellular signals to the inside of the cell, which ultimately leads to changes in gene expression. Some G<sub>q</sub>-coupled receptors can be activated through the specific binding of ligands or drugs to the extracellular domains of the protein. This interaction leads to downstream events such as mobilization of calcium from intracellular storage vesicles into the cytoplasm of the cell. Improper G<sub>q</sub>-coupled receptor function is implicated in cancer, neurodegenerative and cardiovascular diseases. Therefore, it is important to compare receptor function in normal and disease states.

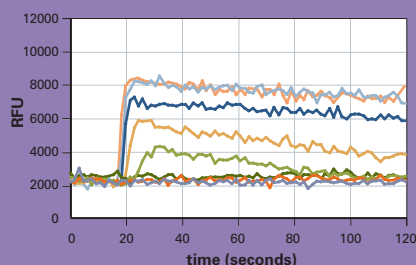
The cell line used in these studies, M1 CHO, is stably transfected with the M1 muscarinic G<sub>q</sub>-coupled receptor<sup>1</sup>. This receptor is stimulated with carbachol, which leads to the release of calcium from the endoplasmic reticulum into the cytosol. Activation of the M1 G<sub>q</sub>-coupled receptor can be monitored using the

FlexStation® Microplate Reader in conjunction with either the calcium-sensitive fluorescent dye, Fluo-3, or the FlexStation Calcium Assay Kit from Molecular Devices. The results demonstrate that a similar range of EC<sub>50</sub> values is obtained for carbachol using both detection methods.

## METHODS

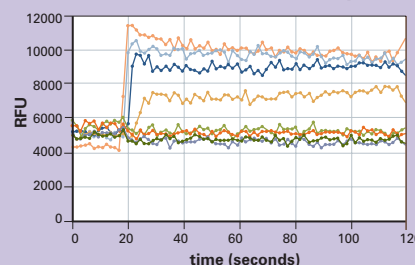
Cells were seeded the night before the experiment at a concentration of 30,000 cells/well in a volume of 100 µL per well of black walled, clear bottomed, 96-well microplates (E&K distributors, cat# 655090). Cells were incubated either for one hour with a buffer containing 4.2 µM Fluo-3 or for one-half hour with 1X Loading Buffer from the FlexStation Calcium Assay Kit (Molecular Devices, cat#R8041) before using FlexStation to dispense carbachol and monitor the mobilization of calcium into the cell cytosol. All experiments were performed at a final concentration of 2.5 mM probenecid in order to inhibit endogenous efflux pumps present in M1 CHO cells.

**carbachol-stimulated M1 CHO  
cells: FLUO-3 (figure 1)**



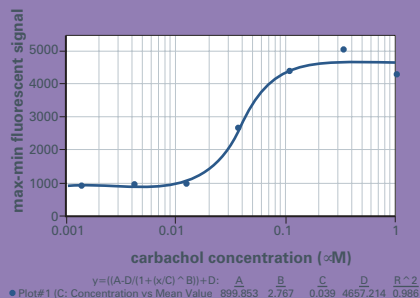
The response of M1 CHO cells to carbachol —  
representative raw data using Fluo-3.

**carbachol-stimulated M1 CHO cells:  
FLEXstation Calcium Assay Kit (figure 2)**



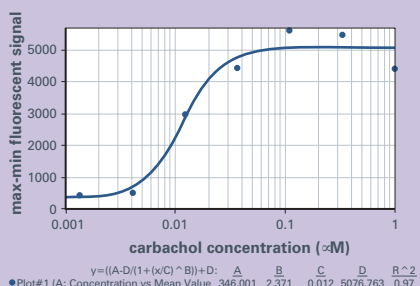
The response of M1 CHO cells to carbachol —  
representative raw data using the FlexStation  
Calcium Assay Kit.

### dose-response to carbachol: FLUO-3 (figure 3)



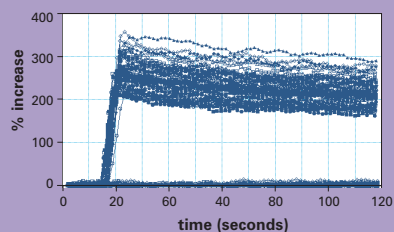
SOFTmax<sup>®</sup> PRO graph window. Representative dose response curve for carbachol is shown using Fluo-3.

### dose-response to carbachol: FLEXstation Calcium Assay Kit (figure 4)



SOFTmax PRO graph window. Representative dose response curve for carbachol is shown using the FlexStation Calcium Assay Kit.

### Z' factor studies using FLEXstation (figure 5)



Z' factor studies using FlexStation. Data is displayed as "Max-Min" and "Percent Increase" by choosing these parameters in SOFTmax PRO software.

### DETERMINATION OF EC<sub>50</sub> VALUES

Cells were exposed to a range of carbachol (0 to 1 μM) using half log dilutions of the agonist. Results are representative of at least four independent experiments performed over a period of at least two different days. The raw data was analyzed as "Max-Min" and "Absolute" RFU values (Figures 1, 2). The average value of replicate samples was plotted on a graph using a curve fit of four parameter, for which the "C" value is the concentration of drug that causes 50% of maximal response (EC<sub>50</sub> value, see Figure 3). The EC<sub>50</sub> values for carbachol using Fluo-3 ranged from 19–39 nM, whereas it ranged from 12–30 nM using the FlexStation Calcium Assay Kit (see Figures 3 and 4).

### DETERMINATION OF Z' FACTOR

Cells in half of the microplate were treated with 0.5 μM carbachol, while cells in the other half of the plate were treated with 1X Reagent Buffer (1X HBSS/20 mM HEPES pH 7.4). The FlexStation Calcium Assay Kit was used to monitor carbachol-mediated calcium mobilization (Figure 5). Z' factor was calculated using the formula<sup>2</sup>:

$$\frac{(3\sigma_{c+} - 3\sigma_{c-})}{|\mu_{c+} - \mu_{c-}|}$$

The terms  $\sigma_{c-}$  and  $\sigma_{c+}$  denote the standard deviation of the negative control and positive control, respectively, whereas  $|\mu_{c+} - \mu_{c-}|$  denotes the absolute value of the difference in the mean of the positive control and the mean of the negative control. The Z' factor obtained for FlexStation using the FlexStation Calcium Assay Kit was 0.65. This indicates a large separation band between the negative and positive controls and an optimized, high quality assay.

### SUMMARY

FlexStation generates highly reproducible results for calcium mobilization assays used to monitor the activity of G<sub>q</sub>-coupled receptors (Fig. 3). First, the EC<sub>50</sub> values obtained for carbachol were consistent between detection methods, Fluo-3 and the FlexStation Calcium Assay Kit. Also, assays analyzed on the same day showed less than two-fold differences, and assays analyzed on different days showed less than four-

fold differences, which further demonstrates the reliable performance of the FlexStation System (data not shown). Lastly, the Z' factor determination indicates that FlexStation provides an excellent tool for assay development for the model system examined. In addition, other benefits include SOFTmax PRO software that makes it simple to set up the instrument and analyze the data. Moreover, the dual monochromators provide increased flexibility for assay development or basic research purposes. In conclusion, FlexStation provides an easy-to-use, versatile tool to help basic research and assay development laboratories reach their goals.

For more detailed information, please refer to the application note "Comparison of FLIPR<sup>®</sup> and FlexStation<sup>®</sup> Systems for Calcium Mobilization Assays", available for download from Molecular Devices' web site at: [www.moleculardevices.com/pages/flex\\_lit1.html](http://www.moleculardevices.com/pages/flex_lit1.html)

### REFERENCES

- 1 Buck, M. A., and C. M. Fraser. 1990. *Biochem. Biophys. Res. Commun.* 173: 666–672.
- 2 Zhang, J.-H., T.D.Y. Thomas, and K. R. Oldenburg. 1999. *J. Biomol. Screening* 4: 67–73.

### SALES OFFICES

- USA & Canada +1-800-635-5577
- Brazil +55-11-3616-6607
- China (Beijing) +86-10-6410-8669
- China (Shanghai) +86-21-6887-8820
- Germany +49-89/96-05-88-0
- Japan (Osaka) +81-6-6399-8211
- Japan (Tokyo) +81-3-5282-5261
- South Korea +82-2-3471-9531
- United Kingdom +44-118-944-8000

Check our web site for a current listing of our worldwide distributors.

[www.moleculardevices.com](http://www.moleculardevices.com)

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

The trademarks used herein are the property of Molecular Devices, Inc. or their respective owners.

Specifications subject to change without notice.

