

# Calcium Detection Upon GPCR Stimulation Using the NOVOstar and Screen Quest™ Fluo-8 NW Calcium Kit



Jinfang Liao<sup>1</sup> and E.J. Dell<sup>2</sup>

<sup>1</sup>ABD Bioquest, Sunnyvale, CA, USA; <sup>2</sup>BMG LABTECH Inc, Durham, NC, USA

Application Note 157

Rev. 06/2007



- Fluo-8 NW is the brightest calcium indicator currently available; >2 times brighter than Fluo-4 NW and >4 times brighter than Fluo-3 AM
- NOVOstar's unique pipetting system allows for instantaneous detection after addition of agonist
- Fluo-8 and Fluo-4 have similar EC<sub>50</sub> values

## Introduction

Analyzing second messengers in cells upon agonist stimulation is an ideal way to measure the activation of G protein coupled receptors (GPCRs)<sup>1</sup>. A calcium flux assay, which measures intracellular calcium, is such a method that is commonly used in drug discovery for screening GPCRs.

The Screen Quest™ Fluo-8 NW Calcium Assay kit from ABD Bioquest provides a homogeneous fluorescence-based assay for detecting intracellular calcium mobilization<sup>2</sup>. Cells expressing a GPCR of interest that signals through calcium are pre-loaded with Fluo-8 NW, which crosses the cell membrane. Once inside the cell, the lipophilic blocking groups of Fluo-8 NW are cleaved by esterases resulting in a negatively charged fluorescent dye that stays inside the cell. Subsequent binding to calcium causes the fluorescence of Fluo-8 NW to be greatly enhanced. Its characteristics of a long wavelength, high sensitivity, and >100 times fluorescence enhancement (when it forms a complex with calcium) make Fluo-8 NW an ideal indicator for measurement of intracellular calcium. The Screen Quest™ Fluo-8 NW Calcium Assay Kits provide an optimized assay method for monitoring G-protein-coupled receptors (GPCRs) and calcium channels. The assay can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation.

The NOVOstar from BMG LABTECH is a multifunctional microplate reader with four measurement modes: fluorescence intensity, luminescence, absorbance, and fluorescence polarization. It is an ideal platform to perform cell-based, fast kinetic assays such as intracellular calcium measurement, because it has a unique pipetting system that can add small amounts of agonist or antagonist (<1.0 µL) to cells followed by instantaneous reading. The NOVOstar also has two injectors that can be used for reagent and/or substrate addition.

## Materials and Methods

- Screen Quest™ Fluo-8 NW Calcium Assay Kit \*1% FBS Medium\* from ABD Bioquest (#36315 (10 plates), #36316 (100 plates))
- Screen Quest™ Fluo-8 NW Calcium Assay Kit \*Medium Removed\* from ABD Bioquest (#36308 (10 plates), #36309 (100 plates))
- NOVOstar Microplate Reader from BMG LABTECH, Inc.
- Black, clear bottom 96-well Costar® microplate
- Other reagents were acquired through normal channels

Detailed protocols for the Screen Quest™ Calcium No Wash Assay, in the presence or absence of medium, are provided with the kits.

HEK293 cells were loaded with either Fluo-8 NW, Fluo-4 NW or Fluo-3 AM dyes. The cells were then stimulated with carbachol using the NOVOstar pipettor, thereby activating endogenous muscarinic receptors and subsequently causing the release of intracellular calcium.

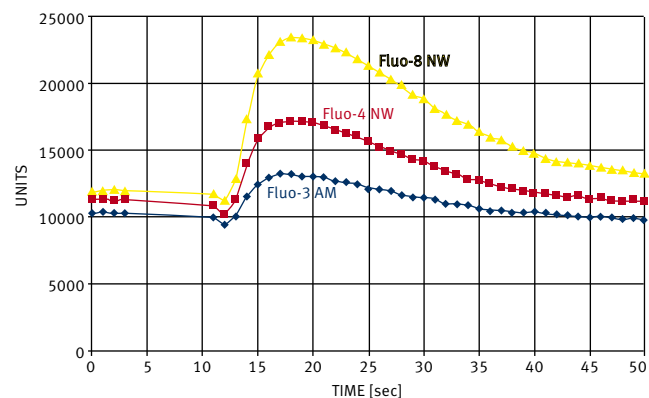
For all assays, HEK293 cells were seeded overnight at 40,000 cells per 100 µl per well in a 96-well black wall/clear bottom Costar® plate. Growth medium was removed in figures 1 and 2, but for figure 3 FBS (fetal bovine serum 1%) was kept in the plate.

Cells were incubated with 100 µl of 4µM Fluo-8 NW, Fluo-4 NW, or Fluo-3 AM for 30 minutes at 37°C, and then another 30 minutes at room temperature. Using the NOVOstar pipettor, either 30 µM of carbachol (25 µl/well) was added (figure 1) or different concentrations of carbachol (25 µl/well for figure 2, and 50 µl/well for figure 3) were added to achieve the final indicated concentrations. Two microplates are used with the NOVOstar, a reagent plate with set dilutions of carbachol and an experimental plate with cells.

Carbachol was added at 12 secs and readings were taken up to 50 secs at 1 sec intervals. The fluorescence intensity mode was used along with a 485/10 excitation filter and a 520/10 emission filter.

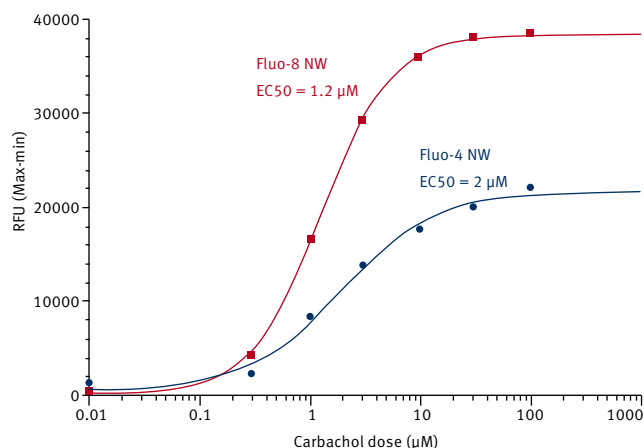
## Results and Discussion

Using the NOVOstar microplate reader with its unique pipetting system, calcium was measured in HEK293 cells after stimulation with carbachol (30µM). Figure 1 shows that Fluo-8 NW has a 2-fold brighter relative fluorescence than Fluo-4 NW and a 4-fold brighter relative fluorescence than Fluo-3 AM.

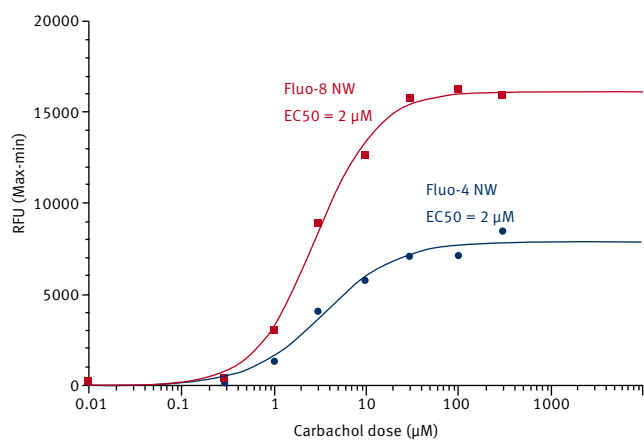


**Fig. 1:** Comparisons of Screen Quest™ Fluo-8 NW Assay Kit, Fluo-4 NW, and Fluo-3 AM on HEK-293 cells after carbachol stimulation (30 µM) with the growth media removed.

This shows that Fluo-8 NW is the brightest fluorescent calcium dye currently available. Further experimentation shows that Fluo-8 NW has a similar binding affinity for calcium as Fluo-4 NW, as determined by their EC<sub>50</sub> values (Figures 2 and 3). In figure 2, the medium was removed from the cells; whereas in figure 3, the medium (1% FBS) was not removed. Note that the EC<sub>50</sub> values derived from the two figures are similar, but that the relative fluorescence units (RFU) are more than twice as high when there is no FBS.



**Fig. 2:** Carbachol Dose Response in HEK-293 cells measured with Screen Quest™ Fluo-8 NW Assay kit and Fluo-4 NW Assay Kit with the growth medium removed. Carbachol (25µl/well) was added by NOVOstar pipette to achieve the final indicated concentrations.



**Fig. 3:** Carbachol Dose Response in HEK-293 cells measured with Screen Quest™ Fluo-8 NW Assay kit and Fluo-4 NW Assay Kit with 1% FBS. Carbachol (50µl/well) was added by NOVOstar pipette to achieve the final indicated concentrations.

## Conclusion

Intracellular calcium measurements can be used to detect the activation of a receptor (i.e. GPCR), which in turn can be used to screen specific agonists or antagonists.

There are two keys aspects to a good intracellular calcium assay:

1. The ability to deliver an agonist/antagonist to cells followed by instantaneous detection of calcium, and
2. A membrane permeable dye that greatly increases in intensity upon binding to calcium.

The NOVOstar (figure 4) from BMG LABTECH fulfils criteria 1 by providing the ideal system to perform all of your fast kinetic, cell-based assays. With its unique pipetting system, the NOVOstar can deliver minute amounts (<1 µL) of agonist or antagonist from the reagent plate to the experimental plate. Detection, using one of its four measurement modes (fluorescence intensity, luminescence, absorbance, or fluorescence polarization), can be initiated prior to, during, and after injection.

The Screen Quest™ Fluo-8 NW calcium kit (in the presence or absence of FBS) from ABD Bioquest fulfils criteria 2. Fluo-8 NW is a membrane permeable dye that increases in fluorescence by 100 fold upon binding to calcium. It is more than twice as bright as Fluo-4 NW and it has a similar affinity for calcium as Fluo-4 NW.

Thus the NOVOstar (figure 4) and Fluo-8 NW, when combined, create the perfect platform for which to measure intercellular calcium.



**Fig. 4:** BMG LABTECH's NOVOstar multifunctional microplate reader with unique pipetting system allows for instantaneous reading upon agonist stimulation.

## References

1. Kiselyov K, Shin DM, Muallem S. (2003) Signalling specificity in GPCR-dependent Ca<sup>2+</sup> signalling. *Cell Signal.* **15**(3), 243-253.
2. For more information on Screen Quest™, Fluo-8 NW and other ABD Bioquest products go to [www.abdbioquest.com](http://www.abdbioquest.com).

<b>Germany:</b>	<b>BMG LABTECH GmbH</b>	<b>Tel: +49 781 96968-0</b>
<b>Australia:</b>	BMG LABTECH Pty. Ltd.	Tel: +61 3 59734744
<b>France:</b>	BMG LABTECH SARL	Tel: +33 1 48 86 20 20
<b>Japan:</b>	BMG LABTECH JAPAN Ltd.	Tel: +81 48 647 7217
<b>UK:</b>	BMG LABTECH Ltd.	Tel: +44 1296 336650
<b>USA:</b>	BMG LABTECH Inc.	Tel: +1 919 806 1735
<b>Internet:</b>	<a href="http://www.bmglabtech.com">www.bmglabtech.com</a>	<a href="mailto:info@bmglabtech.com">info@bmglabtech.com</a>