

## Screen Quest™ Fluo-4 No Wash Calcium Assay Kit

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 36325 (10 plates), 36326 (100 plates)	Keep in freezer and avoid light	FLIPR, FDSS, NOVOSTar, FlexStation, ViewLux, IN Cell Analyzer, ArrayScan

### Introduction

Screen Quest™ Fluo-4 No Wash Calcium Assay Kits provide homogeneous fluorescence-based assays for detecting intracellular calcium mobilization. Cells expressing a GPCR of interest that signals through calcium are pre-loaded with Fluo-4 AM which can cross cell membrane. Once inside the cell, the lipophilic blocking groups of Fluo-4 AM are cleaved by esterases, resulting in a negatively charged fluorescent dye that stays inside cells. Its fluorescence is greatly enhanced upon binding to calcium. When cells stimulated with agonists, the receptor signals the release of intracellular calcium, which significantly increase the fluorescence of Fluo-4. The Screen Quest™ Fluo-4 No Wash Calcium Assay Kits provide an optimized assay method for monitoring the G-protein-coupled receptors and calcium channels. The assay can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation.

### Kit Components

Components	Cat. # 36325 (10 plates)	Cat. # 36326 (100 plates)
Component A: Fluo-4 AM	1 vial, lyophilized	10 vials, lyophilized
Component B: 10X Pluronic® F127 Plus	1 bottle (10 mL/bottle)	10 bottles (10 mL/bottle)
Component C: HHBS (Hanks' Buffer with 20 mM Hepes)	1 bottle (100 mL)	Not included

### Assay Protocol for One Plate

#### Brief Summary

**Prepare cells in growth medium → Add Fluo-4 AM loading solution (100 µL/well/96-well plate or 25 µL/well/384-well plate) → Incubate at 37°C for 1 hour → Monitor fluorescence intensity at Ex/Em = 490/525 nm**

*Warning: Do not add additional probenecid.*

#### 1. Prepare cells:

- 1.1 For adherent cells: Plate cells overnight in growth medium at 40,000 to 80,000 cells/well/100 µL for a 96-well plate or 10,000 to 20,000 cells/well/25 µL for a 384-well plate.
- 1.2 For non-adherent cells: Centrifuge the cells from the culture medium and then suspend the cell pellet in cell growth medium or HHBS at 125,000 to 250,000 cells/well/100 µL for a 96-well poly-D lysine plate or 30,000 to 60,000 cells/well/25 µL for a 384-well poly-D lysine plate. Centrifuge the plate at 800 rpm for 2 minutes with brake off prior to the experiments.

*Note: Each cell line should be evaluated on the individual basis to determine the optimal cell density for the intracellular calcium mobilization.*

#### 2. Prepare Fluo-4 AM dye-loading solution:

- 2.1 Thaw all the kit components at room temperature before use.
- 2.2 Make Fluo-4 AM stock solution: Add 200 µL of DMSO into the vial of Fluo-4 AM (Component A), and mix them well.  
*Note: 20 µL of Fluo-4 NW stock solution is enough for one plate. Unused Fluo-4 AM stock solution can be aliquoted and stored at ≤ -20 °C for more than one month if the tubes are sealed tightly. Protect from light and avoid repeated freeze-thaw cycles.*
- 2.3 Make 1X assay buffer:

Make 1X assay buffer by adding 1 mL of 10 X Pluronic<sup>®</sup> F127 Plus, (10 mL, Component B) into **9 mL** of HHBS buffer (Component C), and mix them well.

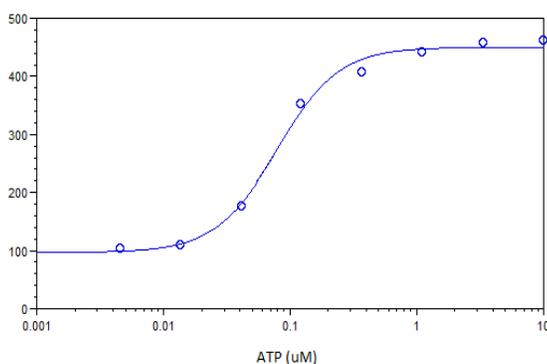
*Note: 10 mL of 1X assay buffer is enough for one plate. Aliquot and store un-used 10X Pluronic<sup>®</sup> F127 Plus (Component B) at  $\leq -20^{\circ}\text{C}$ . Protect from light and avoid repeated freeze-thaw cycles.*

- 2.4 **Make Fluo-4 AM dye-loading solution for one cell plate:** Add 20  $\mu\text{L}$  of Fluo-4 NW stock solution (from Step 2.2) into 10 mL of 1X assay buffer (from Step 2.3), and mix them well. This working solution is stable for at least 2 hours at room temperature.

### 3. Run calcium assay:

- 3.1 Add 100  $\mu\text{L}$ /well (96-well plate) or 25  $\mu\text{L}$ /well (384-well plate) of Fluo-4 AM dye-loading solution (from Step 2.4) into the cell plate.
- 3.2 Incubate the dye-loading plate in a cell incubator for 1 hour, and then incubate the plate at room temperature for another 15 to 30 minutes.
- Note 1: If the assay requires 37°C, perform the experiment immediately without further room temperature incubation.*
- 3.3 Prepare the compound plate with HHBS or your desired buffer.
- 3.4 Run the calcium flux assay by monitoring the fluorescence intensity at Ex/Em = 490/ 525 nm.

## Data Analysis



**Figure 1.** ATP Dose Response was measured in CHO-K1 cells with Screen Quest™ Fluo-4 No Wash Calcium Assay Kit. CHO-K1 cells were seeded overnight at 60,000 cells/100  $\mu\text{L}$ /well in a Costar black wall/clear bottom 96-well plate. The cells were incubated with 100  $\mu\text{L}$  of dye-loading solution using the Screen Quest™ Fluo-4 No Wash calcium assay kit for 1 hour at 37°C, and then at room temperature for another 15 minutes. ATP (50 $\mu\text{L}$ /well) was added by Flexstation 3 to achieve the final indicated concentrations.

## References

1. Falk, S and Rekling, J. C (2009) Neurons in the preBötzing complex and VRG are located in proximity to arterioles in newborn mice. *Neuroscience Letters*. Volume 450, Issue 3, 229-234.
2. Ghoneum, M. Matsuura, M. and Gollapudi, S. (2009) An iron-based beverage, HydroFerrate fluid (MRN-100), alleviates oxidative stress in murine lymphocytes *in vitro*. *Nutrition Journal* 8:18-23.
3. Satoru, T. Kobayashi, K. Takahashi, M. Katahira, K. Goryo, K. Matsushita, N. Yasumoto, K. Fujii-Kuriyama, Y. and Sogawa, K. (2009) Magnesium Deficiency Causes Loss of Response to Intermittent Hypoxia in Paraganglion Cells. *J. Biol. Chem.*, Vol. 284, Issue 28, 19077-19089.

**Warning: This kit is only sold to our authorized distributors and end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of the kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.**