

Cell Based Assays on the FLIPR[®]TETRA[®] System: Comparison of the FLIPR[®] Calcium 5 Assay Kit to Other Fluorescence Based Calcium Flux Assays

By Carole Crittenden¹, Jan Dolzer, Ph.D., BingFang Huan¹ Ph.D., Ming Yu¹ Ph.D., Simon Lydford², and Cerys Huggins² Ph.D.
MDS Analytical Technologies, ¹Sunnyvale, California, USA and ²Wokingham, Berkshire, UK

Abstract

Cell-based calcium flux assays on FLIPR[®] fluorometric imaging plate readers are widely used in high throughput screening for identification of GPCR agonists and antagonists within the pharmaceutical industry. Masking technology enables no-wash fluorescence-based detection of changes in intracellular calcium concentration, and remains one of the most popular HTS methods for identifying potential drug candidates. Masking technology significantly lowers background fluorescence and increases the signal-to-noise ratio without the need to wash cells. In this study, we used the FLIPR[®]TETRA[®] instrument to compare the performance of FLIPR[®] Calcium 4 and Calcium 5 Assay Kits to a competitor kit. Several endogenous GPCRs as well as transfected Muscarinic M1 receptor were studied in four cell lines. This evaluation demonstrates that the superior performing FLIPR[®] Calcium 5 Assay Kit provides a larger signal window, increased signal to noise ratio, consistent pharmacology, and equivalent or improved Z factors.

Introduction

The FLIPR[®] Calcium 5 Assay Kit contains a new, superior performance calcium sensitive indicator that does not require media removal and utilizes the same proven quench technology as the FLIPR[®] Calcium 4 Assay Kit. This masking dye technology is licensed to Molecular Devices, Inc. from Bayer AG (patent nos. US 6,420,183, EP 0906572). The masking technology combined with the novel indicator significantly lowers background fluorescence and improves signal-to-noise without washing. Throughput is increased by eliminating labor intensive wash steps, reducing preparation time for wash buffers and wash calibrations. After incubation, the assay functions at room temperature making it applicable to automation using stackers and robots.

FLIPR[®] Calcium 5 Assay Kit Response to Intracellular Calcium Mobilization

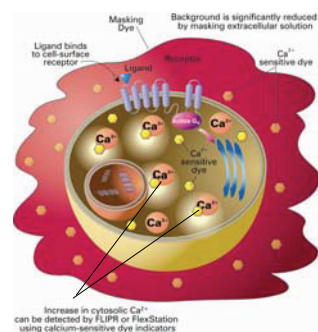


Figure 1. During cellular response to GPCR signaling, fluorescent signal increases as released intracellular calcium molecules bind to the Calcium 5 Assay Kit indicator. Background signal is decreased by extracellular masking technology.

Materials and Methods

Cell lines, receptors, and ligands:

M1WT3 CHO cells (CHO M1) (ATCC Cat# CRL 1985) were plated at 10,000 cells per well in culture media in black-wall clear-bottom 384-well TC coated plates (Corning Cat# 3712). Carbachol (Sigma-Aldrich Cat# C4382) was used to stimulate the Muscarinic M1 receptor and Atropine (Sigma-Aldrich Cat# A0132) was tested as antagonist.

HeLa cells (ATCC Cat# CCL2) were plated at 10,000 cells per well as listed above. Histamine (Sigma-Aldrich Cat# H7125) was used to stimulate the endogenous Histamine H1 agonist and Pyrilamine maleate (Sigma-Aldrich Cat# P5514) was tested as antagonist. Dye for CHO M1 and HeLa cell lines was formulated in each case with 2.5 mM final concentration of water soluble probenecid (Invitrogen Cat# P36400). All cell plates were incubated overnight at 37°C in 5% CO₂.

HEK 293 cells (ATCC Cat# CRL 1573) cells were plated at 15,000 cells per well in culture media in black-wall clear-bottom 384-well poly-D-lysine coated plates (Cat# 356663, BD Life Sciences). Carbachol was used to stimulate the endogenous Muscarinic M3 receptor. Atropine was tested as the antagonist.

Jurkat immortalized T-cells (ATCC Cat# TIB-151) were plated at 75,000/well in 50 µL 1X dye loading media containing 0.1% BSA and centrifuged to insure that they were on the bottom of the well. Cells were incubated at 37 °C in 5% CO₂ for one hour and allowed to come to room temperature prior to the assay on the FLIPR[®]TETRA instrument.

Adenosine 5' triphosphate (Sigma-Aldrich Cat# A2383) was used to stimulate the endogenous P2Y receptor in all cell lines except Jurkat.

FLIPR[®] Calcium 4 and Calcium 5 Assay Kit Protocol:

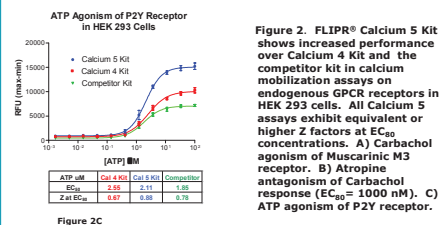
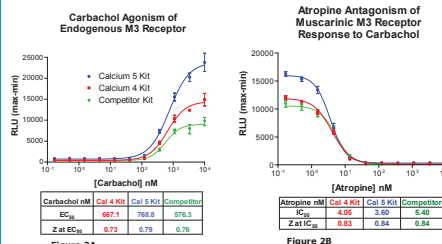
FLIPR[®] Calcium 5 Explorer Kit (Molecular Devices, Inc.) includes 10 vials of dye (Component A) and 1 bottle of Hanks Balanced Salt Solution (HBSS) and 20 mM HEPES adjusted to pH 7.4 (Component B) sufficient for 1 plate each. Bulk kits contain 10 vials of dye sufficient for 10 plates each. Competitor kits contain the same type of components

Dye loading buffer for 1 plate was prepared by dissolving contents of one vial of Component A completely with a final volume of 10 mL Component B loading buffer. Cell plates were removed from the incubator and 25 µL Calcium 4 Kit, Calcium 5 Kit, or competitor kit dye loading buffer was added to each well. Plates were not washed after dye addition. Dye loaded plates were incubated 45 minutes at 37 °C, 5% CO₂ and allowed to come to room temperature 15 minutes prior to reading on the FLIPR[®]TETRA instrument.

Calcium Mobilization Assay on FLIPR[®]TETRA[®]:

A 5X volume of CRC ligand was prepared in HBSS buffer + 20 mM HEPES in 384-well polypropylene plates. Agonist was added during detection on the FLIPR[®]TETRA[®] instrument at optimized parameters. Antagonist was prepared at 5X concentration and added 15 minutes prior to addition of a 6X volume of EC₅₀ concentration of challenge agonist. Relative Fluorescence Units (RFU) were measured for each response for signal maximum minus minimum during approximately 90 seconds after addition. Graphs and EC₅₀/IC₅₀ concentrations were calculated using GraphPad Prism. Z-factor calculations were performed using the method described by Zhang, et al.

Calcium Mobilization in HEK-293 Cells



Calcium Mobilization in CHO M1 Cells

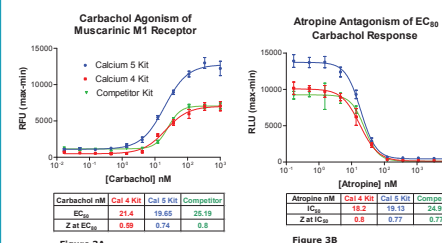
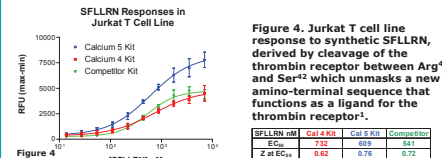
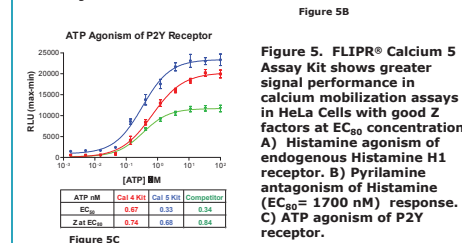
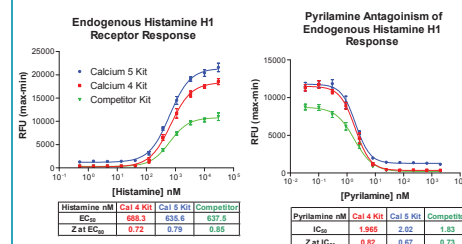


Figure 3. FLIPR[®] Calcium 5 Kit shows superior performance in calcium mobilization assays in CHO M1 Cells with good Z factors at EC₅₀ concentrations. Calcium 4 Kit and the competitor kit exhibit equivalent performance. A) Carbachol agonism of Muscarinic M1 receptor. B) Atropine antagonism of Carbachol (EC₅₀=70 nM) response.

Calcium Mobilization in Jurkat Suspension Cells



Calcium Mobilization in HeLa Cells



ATP Agonism of P2Y Receptor

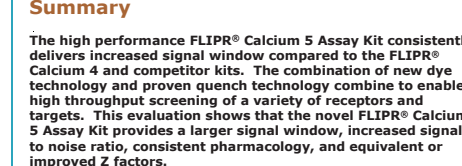


Figure 5. FLIPR[®] Calcium 5 Assay Kit shows greater signal performance in calcium mobilization assays in HeLa Cells with good Z factors at EC₅₀ concentrations. A) Histamine agonism of endogenous Histamine H1 receptor. B) Pyrilamine antagonism of Histamine (EC₅₀=1700 nM) response. C) ATP agonism of P2Y receptor.

Summary

The high performance FLIPR[®] Calcium 5 Assay Kit consistently delivers increased signal window compared to the FLIPR[®] Calcium 4 and competitor kits. The combination of new dye technology and proven quench technology combine to enable high throughput screening of a variety of receptors and targets. This evaluation shows that the novel FLIPR[®] Calcium 5 Assay Kit provides a larger signal window, increased signal to noise ratio, consistent pharmacology, and equivalent or improved Z factors.

Reference

Mari, et al., Thrombin and trypsin-induced Ca²⁺ mobilization in human T cell lines through interaction with different protease-activated receptors, FASEB Journal, February 1996, (10) 309-316.