

Ion Channel Assay Development using Invitrogen's FRET-based Voltage Sensor Probes

Randall L. Hoffman, Steve Duff and Naveeda Qadir
Invitrogen Corporation, Madison, WI 53719, USA

Application Note 123

Rev. 10/2005



- POLARstar OPTIMA and NOVOstar used for cost effective ion channel assay development
- Plate-to-plate transfer for higher throughput with the NOVOstar
- Injection and FRET measurement at the same time

Introduction

Ion channels are important drug targets because of their critical role in nerve, cardiac, endocrine and skeletal muscle tissues. The lack of sufficiently sensitive screening systems has hampered research in this area. This application note focuses on Voltage Sensor Probes (VSP), a Fluorescence Resonance Energy Transfer (FRET)-based voltage-sensing assay technology from Invitrogen for measuring changes in cellular membrane electrical potential.

This technology enables detection and measurement of rapid changes in membrane voltage and quickly reports them as fluorescence signals from living cells. This technology can be used with any target that changes membrane potential, and is therefore well suited for sodium, potassium, calcium, chloride and ligand-gated ion channel research. The ratiometric method used to detect and quantify changes in cellular membrane potential significantly reduces errors arising from well-to-well variations in cell number, dye loading and signal intensities, plate inconsistencies and temperature fluctuations. These combined features make VSP technology highly amenable for high-throughput screening (HTS) applications.

In order to develop and validate an ion channel assay prior to HTS, it would be helpful for assay development and therapeutic groups to have access to less expensive equipment amenable to VSP technology. This application note demonstrates the use of the POLARstar OPTIMA and NOVOstar plate readers from BMG LABTECH as suitable platforms for development of VSP ion channel assays. These plate readers are much less expensive than HTS ion channel readers and can be used for multiple detection technologies.

Assay Principle

Voltage Sensor Probes (VSP) is a Fluorescence Resonance Energy Transfer (FRET)-based assay technology used for high-throughput ion channel drug discovery. The FRET donor is a membrane-bound, coumarin-phospholipid (CC2-DMPE), which binds only to the exterior of the cell membrane. The FRET acceptor is a mobile, negatively charged, hydrophobic oxonol [either DiSBAC₂(3) or DiSBAC₄(3)], which will bind to either side of the plasma membrane in response to changes in membrane potential. Resting cells have a relatively negative potential, so the two probes associate with the exterior of the cell membrane, resulting in efficient FRET (Figure 1). Exciting the CC2-DMPE donor probe (filter: 400-10 nm) generates a strong red fluorescence signal (filter: 570-20 nm) from the oxonol acceptor probe. When the membrane potential becomes more positive, as occurs with cell depolarization, the oxonol probe rapidly translocates (on a subsecond time scale) to the other face of the membrane. Thus, each oxonol probe “senses” and responds to voltage changes in the cell. This translocation separates the FRET pair, so exciting the CC2-DMPE donor probe now generates a strong blue fluorescence signal (filter: 460-10 nm) from the CC2-DMPE probe.

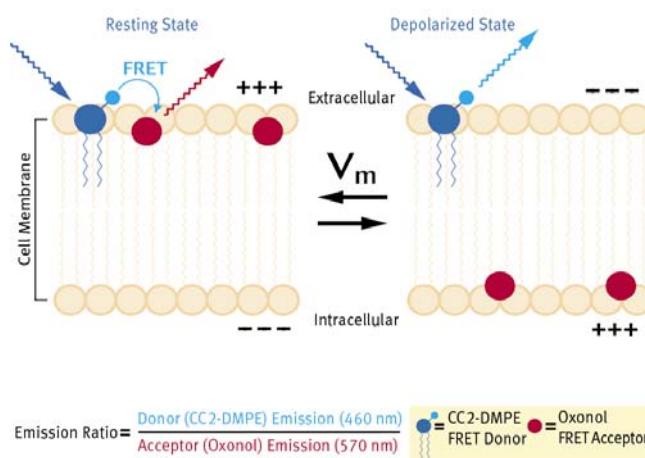


Fig. 1: Schematic illustration of the mechanism of voltage-dependent, FRET-based VSPs

During high-throughput ion channel assay development, a simple model is often first tested for proof of principle for both the assay technology and the instrumentation. In this application note, we demonstrate the compatibility of two instruments from BMG LABTECH for developing ion channel assays using VSP technology from Invitrogen. The ion channel model we have used is endogenously expressed inward rectifying potassium (Kir) channels present in rat basophilic leukemia cells. These channels are opened when high potassium is added to the cell's medium. This causes a change in the cell's membrane potential which can be modulated by barium chloride. In the assay, the cells are manually loaded with the dyes as described below. Barium chloride is then immediately added at various concentrations. After incubation, the plated cells are placed in the reader and in the case of the NOVOstar, the barium chloride is added by the pipettor inside the reader. The reader begins to take baseline readings at the two wavelengths, injects the high potassium buffer (VSP-2 buffer, below), and continues to take readings for a specified time.

Materials and Methods

Cell Culture

RBL-2H3 (Rat Basophilic Leukemia, ATCC #CRL2256) cells were plated at 50,000 cells/well in Corning® 3603 96-well plates 18-24 hours prior to experimental procedure.

Preparation of VSP Loading Buffers

- 5 μM CC2-DMPE Loading Buffer: Premix 10 μL of 5 mM CC2-DMPE (Invitrogen) and 10 μL of 100 mg/mL Pluronic® F-127 (Sigma). Add 10 mL of VSP Solution 1 (160 mM NaCl, 4.5 mM KCl, 2 mM CaCl₂, 1 mM MgCl₂, 10 mM glucose, 10 mM HEPES, pH 7.4) and vortex vigorously to mix. Protect from light prior to use.
- 10 μM DiSBAC₂(3) Loading Buffer: Premix 8.3 μL of 12 mM DiSBAC₂(3) (Invitrogen) and 12.5 μL of 200 mM VABSC-1 (Invitrogen). Add 10 mL of VSP Solution 1 and vigorously vortex to mix. Protect from light prior to use.

Loading Cells

Media was removed from all wells of the 96-well plates and replaced with 100 μ L VSP Solution 1 (VSP-1). The VSP-1 was immediately removed and replaced with 100 μ L CC2-DMPE Loading Buffer and incubated at room temperature for 30 minutes, covered and protected from light. The CC2-DMPE Loading Buffer was removed after 30 minutes and the plates washed once with 100 μ L VSP-1. The VSP-1 was immediately replaced with 100 μ L DiSBAC₂(3) Loading Buffer and incubated at room temperature for 30 minutes, covered and protected from light.

BaCl₂ (Sigma), when appropriate, was added either immediately following addition of DiSBAC₂(3) Loading Buffer or immediately prior to VSP-2 high K⁺ buffer (164.5 mM KCl, 2 mM CaCl₂, 1 mM MgCl₂, 10 mM glucose, 10 mM HEPES, pH 7.4).

Results and Discussion

Immediately following the DiSBAC₂(3) Loading Buffer 30 minute incubation, the plates were analyzed with either the POLARstar OPTIMA or NOVOstar. Both instruments basically follow the same protocol of initially reading the plate with two “simultaneous” emissions before VSP-2 addition and continuing after VSP-2 addition.

Data is analyzed by comparing the baseline subtracted ratio of donor (at 460 nm) to acceptor (at 570 nm) before (Ro) and after (Rf) addition of VSP-2: the final normalized assay ratio = Rf/Ro.

The POLARstar OPTIMA can inject VSP-2 and collect data at two wavelengths once per second. The software enables real-time visual readout in both plate and well views. Data is automatically exported to the BMG LABTECH evaluation software package for analysis.

A POLARstar OPTIMA plate readout view is shown in Figure 2A) for a “positive control” assay with a 96-well plate with 50,000 RBL cells/well. Column 1 is the no cells control.

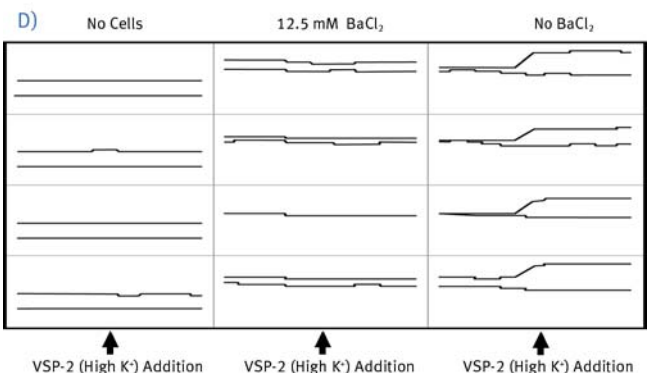
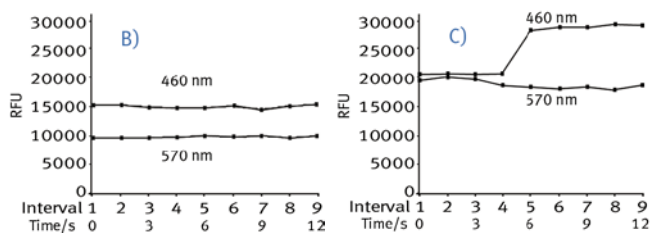
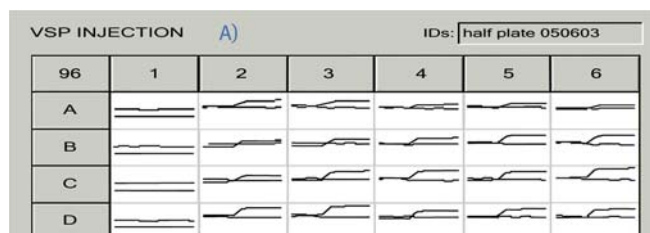


Fig. 2: VSP data on the BMG LABTECH POLARstar OPTIMA

In Figure 2B) a current state window view of a no cell control shows no change in 460 nm upon VSP-2 addition, whereas in Figure 2C) an increase in the 460 nm signal upon injection of VSP-2 is detected. Figure 2D) shows enlarged views of current state windows from 12 wells of a BaCl₂ dose response. Note the lack of 460 nm emission increase in both the no cell control and 12.5 mM BaCl₂ wells compared to untreated wells.

The NOVOstar was programmed to use its auto pipettor to add BaCl₂ dilutions directly from one 96-well plate via plate-to-plate transfer to the assay plate with RBL cells. It then injected VSP-2 using one of its two additional injectors and collected data at two wavelengths once per second (Figure 3).

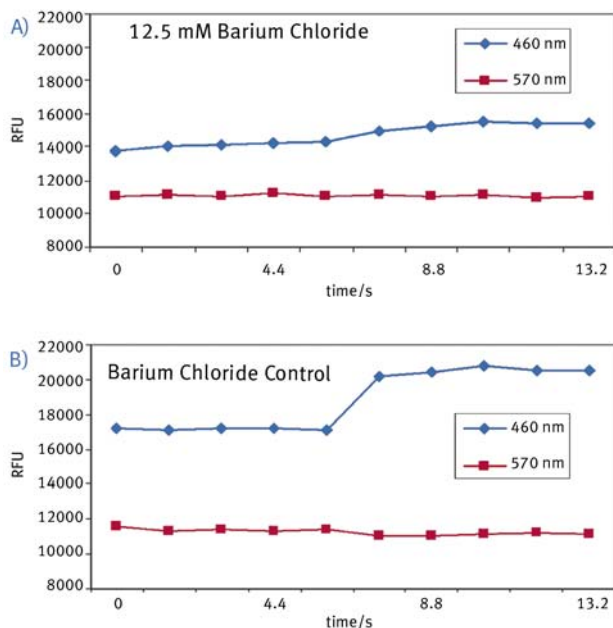


Fig. 3: VSP data on the BMG LABTECH NOVOstar

The raw fluorescent readout was evaluated in Excel and the addition of 12.5 mM BaCl₂ avoids a significant change in 460 nm emission (Figure 3A), whereas an untreated control well shows as expected an increase in 460 nm fluorescence signal upon VSP-2 addition (Figure 3B).

Conclusion

This application note demonstrates the compatibility of Voltage Sensor Probe technology with two plate readers from BMG LABTECH. The key to the technology is having instrumentation that can measure two wavelengths in rapid succession. This takes advantage of the ratiometric FRET readout, which greatly reduces variation common to many cell based assays. Our model for ion channel assay development is endogenously expressed inward rectifying Kir channels present in rat basophilic leukemia cells. With both instruments we showed real-time readouts of the ion channel following K⁺ addition and blockage of the Kir channel with BaCl₂. The NOVOstar injectors were able to inject both the BaCl₂ inhibitor and the K⁺ stimulant. The ability of both the POLARstar OPTIMA and NOVOstar to inject ion channel stimulant, switch filters at approximately 1 Hz, and continue to take readings during the kinetic event make these suitable instruments for therapeutic groups in developing ion channel assays for lower throughput applications and screening.

Germany: **BMG LABTECH GmbH** Tel: +49 781 96968-0
 Australia: BMG LABTECH Pty. Ltd. Tel: +61 3 59734744
 China: BMG LABTECH Co. Ltd. Tel: +86 10 84110632
 France: BMG LABTECH SARL Tel: +33 1 48 86 20 20
 Japan: BMG LABTECH JAPAN Ltd. Tel: +81 48 647 7217
 UK: BMG LABTECH Ltd. Tel: +44 1296 336650
 USA: BMG LABTECH Inc. Tel: +1 919 806 1735

Internet: www.bmglabtech.com info@bmglabtech.com