

APPLICATION NOTE

Development of a cell-based potassium-chloride transporter assay using the FLIPR Potassium Assay Kit

Introduction

Potassium-chloride transporter member 5 (SLC12A5, or KCC2) is one of nine cationchloride cotransporters (CCCs) encoded by the SLC12 family of genes, and is the only CCC transporter preferentially expressed in neurons. KCC2 plays a critical role in the correct functioning of the central nervous system where it is pivotal in maintaining neuronal intracellular CI⁻ concentration ([Cl-]). KCC2 is involved in the control of numerous neuronal processes and its impaired activity has been identified in epilepsy, neuropathic pain and spasticity following spinal cord injury¹. This suggests that positive modulators of the activity or expression of KCC2 may provide effective therapy for neurological conditions arising from defects in KCC2 functionality.

The FLIPR® Potassium Assay Kit contains a novel, highly-sensitive thallium (TI⁺) indicator dye that produces bright fluorescence upon binding to TI+ introduced through potassium channels. The intensity of the signal is proportional to the number of open potassium channels on the cells; therefore, it acts as a surrogate indicator of potassium ion channel activity. The kit also employs Molecular Devices proprietary masking dye to reduce background fluorescence resulting in an improved signal/background ratio. As TI⁺ is also transported by cation-chloride cotransporters, the FLIPR Potassium Assay Kit can be used to easily assess the effect of compounds on KCC2 activity by measuring the initial rate of KCC2-mediated TI⁺ transport/influx.

Assay principle

The assay kit's TI⁺ indicator dye has an expanded signal window compared to non-homogenous assays. BTC-AM is a Ca2+ indicator that has also been used for monitoring potassium channel since TI⁺ ions also enhance the fluorescence of BTC. During the dye-loading step, the TI⁺ indicator dyes enter the cells as acetoxymethyl (AM) esters by passive diffusion across the cell membrane. Cytoplasmic esterases cleave the AM esters and release the active fluorogenic forms. A proprietary extracellular masking dye is included with the assay kit to reduce background fluorescence (Figure 1). To activate the potassium-chloride transporter, the cells are stimulated with a mixture of K⁺ and TI⁺ in the presence or absence of test compounds (e.g. KCC2 inhibitors). Under the assay conditions used, the increase in fluorescent signal represents

Benefits

- Kinetic measurement of K⁺/Cl⁻ cotransporter activity in a cellbased assay provides information rich data
- Homogenous no-wash protocol reduces well-to-well variation, improving assay robustness
- Non-radioactive reagent (unlike traditional Rb⁺ efflux assays)
- Amenable to
 high-throughput screening



Figure 1. The FLIPR Potassium Assay Kit principle. Increase in cytosolic TI⁺ is detected by FLIPR or FlexStation systems using thallium-sensitive dye indicators.

the influx of TI⁺ into the cell specifically through the cotransporter, and therefore represents a functional measurement of the cotransporter activity. Modulation of the cotransporter activity is achieved by the inclusion of KCC2 inhibitors.

The FLIPR Potassium Assay Explorer Kit (Molecular Devices Cat. # R8222) contains a TI⁺-sensitive dye, masking dye for homogenous operation, 200 mM K₂SO₄, 50 mM TI₂SO₄, 5X chloride-free buffer, and HBSS with 20 mM HEPES buffer. The kit is sufficient for ten 96-, 384-, or 1536-well plates. The assay workflow is shown in Figure 2.

Experimental procedures Cell culture and transfection

HEK293T (ECACC Cat. # 12022001) cells were grown to 80–90% confluence in T_{7r} flasks containing growth medium (DMEM supplemented with 10% FBS and 2 mM L-Glutamine). Cultures were maintained at 37°C in the presence of 5% CO₂. Cells were passaged every 3-4 days at a ratio of 1:6. Two days before the assay, cells were transiently transfected with human KCC2 (hKCC2) using Lipofectamine 2000 (Thermo Fisher Cat. # 11668027) according to the manufacturer's instructions. The transfection mixture was mixed with growth medium, plated directly into a 96-well, black-walled, clear bottom plate at 10,000 cells/well and incubated at 37°C and 5% CO₂ for 48 hours.

KCC2 inhibition assay

Growth medium was removed from the plates which were then incubated with either FLIPR Potassium Assay Kit or BTC-AM (2 μ M) for one hour at 37°C in the dark. The loading buffer contained 10 μ M bumetanide to reduce the endogenous NKCC TI⁺ influx/transport signal, and 0.1 mM ouabain, a Na⁺/K⁺-ATPase blocker, to inhibit ATP-dependent sodium-potassium exchange across the cell membrane. Following the dye loading step, cells incubated with BTC-AM were washed to remove excess dye. The KCC2 inhibitors were added and incubated for the last 45 minutes of the dye loading time (FLIPR Potassium Assay Kit only). Optimized stimulus buffer (9 mM K⁺ / 0.9 mM TI⁺) was added to the well during







Figure 3. SoftMax Pro 6.4.2 import workflow.

Assay cells with FLIPR Potassium Assay Kit

Acquire data with the

FLIPR Tetra System



Paste raw results into Excel using the SoftMax Pro import template



Import data into SoftMax Pro Software



Figure 4. Representative fluorescence signal traces (\geq 6 wells per treatment) obtained on the FLIPR Tetra System in HEK293 cells loaded with the FLIPR Potassium Assay Kit or BTC-AM. Data show the effect of buffer only addition (–) or TI⁺ addition (–) in hKCC2 transfected cells, and TI⁺ addition (–) in mock transfected cells loaded with the FLIPR Potassium Assay Kit. For comparison, the effect of TI⁺ addition (–) in transfected cells loaded with BTC-AM is also shown.



Figure 5. Representative fluorescence traces obtained on the FLIPR Tetra System in hKCC2-transfected cells loaded with the FLIPR Potassium Assay Kit. Data show the effect of buffer only addition (–) or TI⁺ addition in the absence (–) or presence (–) of 30 μ M R-(+)-DIOA.

Specifications subject to change without notice. Patents: www.moleculardevices.com/productpatents FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

©2015 Molecular Devices, LLC 11/15 1994A Printed in USA

We have shown that the FLIPR Potassium Assay Kit can be used to measure

K⁺/TI⁺-evoked increase in fluorescence

are very consistent from well to well, and

R-(+)-DIOA signals is significant (P < 0.001;

the separation between the control and

In the absence of inhibitor the K^+/TI^+

stimulus produced a Vmax of 53.9 ±

5.1 units per second, whereas in the

presence of 30 µM R-(+)-DIOA this was

reduced to 17.5 \pm 1.9 units per second

(Figure 6) where $n \ge 30$ wells over four

independent experiments.

Student's t-test).

Conclusion

the functional activity of the hKCC2 cation-chloride cotransporter using a homogeneous, no-wash protocol. The assay kit displayed a large assay window and excellent reproducibility, and was able to successfully detect a known modulator of hKCC2 activity. Compared to traditional dyes such as BTC-AM, the simplified protocol and robust assay quality, in combination with the high-throughput capability of the FLIPR Tetra System, provides a powerful solution to screen for modulators of the hKCC2 cotransporter.

Acknowledgements

Andrea Townsend-Nicholson and Stephanie Schorge, University College London, and Simon Lydford, Molecular Devices (UK) Ltd

References

- 1. Kahle KT, Staley KJ, Nahed BV, Gamba G, Hebert SC, Lifton RP, et al. Roles of the cation-chloride cotransporters in neurological disease. Nat Clin Pract Neurol. 2008:4:490-503
- 2. Zhang JH, Chung TD & Oldenburg KR. A Simple Statistical Parameter for Use in Evaluation and Validation of High Throughput Screening Assays. J Biomol Screen. 1999;4:67-73.

detection on the FLIPR® Tetra System using the 470-495 nm excitation LEDs and 515–575 nm emission filter. Data files were imported into SoftMax® Pro Data Acquisition and Analysis Software for subsequent analysis using the Import Feature (Figure 3), available for SoftMax Pro 6.4.1 or higher. The initial rate (Vmax in units per second) was calculated from the first 10 seconds of data post stimulus addition.

Results

After loading hKCC2-transfected or mocktransfected HEK293 cells with the novel FLIPR Potassium Assay Kit or BTC-AM, we measured the increase in fluorescence triggered by the addition of external TI⁺ during detection on the FLIPR Tetra System. The hKCC2-driven FLIPR signal exhibits a rapid increase over the first ten seconds after TI⁺ addition in cells loaded with the FLIPR Potassium Assay Kit, followed by a slower increase and eventually a plateau phase. The increase was significantly higher (P < 0.001; Student's t-test) in hKCC2expressing cells, compared to mocktransfected cells as shown in Figure 4. The average Z factor was calculated as 0.60 \pm 0.05 indicating a robust assay². In contrast, with BTC-AM loaded cells the signal was much slower to develop, showed a large degree of variability and the magnitude of the signal was far lower. For these reasons, no further comparisons between the assay kit and BTC-AM were made.

A series of experiments were then carried out with hKCC2-expressing cells loaded with the FLIPR Potassium Assay Kit to assess whether the TI+ influx assay could be used to detect modulators of hKCC2 activity. In each 96-well plate, cells were treated with either vehicle control or 30 µM R-(+)-DIOA (an alkanoic acid that has been shown to be a potent and selective inhibitor of K⁺/Cl⁻ cotransporters). As seen in Figure 5, the resultant signals from the

Contact Us

Phone: +1-800-635-5577 Web: www.moleculardevices.com Email: info@moldev.com Check our website for a current listing of worldwide distributors.



Figure 6. The effect of different challenges on hKCC2-transfected cells. Data shows the effect of three different treatment groups: 1. buffer only addition (■), and 2. TI⁺/K⁺ addition in the absence (■) or 3. presence (\blacksquare) of 30 μ M R-(+)-DIOA. Bars represent mean V_{max} (units/sec.) ± SEM (n \ge 30).

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



