FLIPR® Potassium Assay Kit

The FLIPR[®] Potassium Assay Kit from Molecular Devices[®] provides a fast, simple, and robust fluorescence-based, high-throughput assay for measuring potassium channel activity. This kit, combining a highly sensitive thallium indicator dye with a Molecular Devices proprietary masking dye technology, provides an homogeneous assay with an expanded signal window and an improved signal-to-noise ratio.

Table 1-1: Available Kits

Assay Kit	Evaluation Kit	Explorer Kit	Bulk Kit
FLIPR [®] Potassium Assay Kit	R8330	R8222	R8223
Chapter 1: About the FLIPR Potass Assay Principle	ium Assay Kit		
Advantages			4 4
Chapter 2: Materials and Equipme Kit Components Materials Required But Not R Storage and Handling Compatible Molecular Device	Provided		5 5 5 6 6
Chapter 3: Experimental Protocol Quick Start Protocol Cell Handling Preparing the Loading Buffer Loading Cells Using Loading Preparing the Thallium Sulfa Running the FLIPR Potassium Running the FLIPR Potassium	Buffer te Source Plate Assay Assay on a FlexSt	ation 3 Instrument	7 7 7 8 9 9 9 10
Chapter 4: Data Analysis Example Inhibition of hERG Channel			
Chapter 5: Optimization Guideline Data Analysis Determination of Potassium	and Thallium Conc	entrations	14 14
Obtaining Support			16

This document is provided to customers who have purchased Molecular Devices equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Software that may be described in this document is furnished under a non-transferrable license. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Each product is shipped with documentation stating specifications and other technical information. Molecular Devices products are warranted to meet the stated specifications. Molecular Devices makes no other warranties or representations express or implied, including but not limited to, the fitness of this product for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom. The sole obligation of Molecular Devices and the customer's sole remedy are limited to repair or replacement of the product in the event that the product fails to do as warranted.

For research use only. Not for use in diagnostic procedures.

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: http://www.moleculardevices.com/patents

Product manufactured by Molecular Devices, LLC. 3860 N. First Street, San Jose, California, 95134, United States of America. Molecular Devices, LLC is ISO 9001 registered. ©2019 Molecular Devices, LLC. All rights reserved. ISO 9001

Chapter 1: About the FLIPR Potassium Assay Kit

Assay Principle

The assay exploits the permeation of thallium ions (TI^+) through both voltage-gated and ligand-gated potassium (K^+) channels. In this assay, a novel, highly-sensitive TI^+ indicator dye is used. This dye produces a bright fluorescent signal upon the binding to TI^+ conducted through potassium channels. The intensity of the TI^+ signal is proportional to the number of potassium channels in the open state. Therefore, it provides a functional indication of the potassium channel activities. In addition, one of the Molecular Devices proprietary masking dyes is employed to further reduce background fluorescence for improved signal-to-noise ratio.

During the initial dye-loading step, the acetoxymethyl (AM) ester tagged Tl⁺ indicator dye enters the cells through passive diffusion. Cytoplasm esterases cleave the AM ester and relieve its active thallium-sensitive form. To activate the potassium channels, the cells are then stimulated with either a mixture of K⁺ and Tl⁺, or a ligand in the presence of Tl⁺. The increase of fluorescence in the assay represents the influx of Tl⁺ into the cell specifically through the potassium channel, providing a functional measurement of potassium channel activity.



Increase in cytosolic Tl[⁺] is detected by FLIPR or FlexStation systems using thallium-sensitive dye indicators

Figure 1-1: Assay principle of the FLIPR Potassium Assay Kit

Advantages

Currently, thallium-based potassium assay methodologies are slow and laborious, and can require multiple replacements of assay buffers or medium. In addition, the assay quality is often challenged by well-to-well variations and a small signal window. Therefore, a fast, easy-to-use, and sensitive cell-based assay specific to potassium channels is highly desirable for high-throughput screening. The FLIPR Potassium Assay Kit provides a true homogeneous assay using a no-wash format, with a number of key features including:

- Direct, functional measurement of potassium channel activity
- Homogeneous, no-wash protocol
- Enhanced signal window and higher signal-to-noise ratio
- Rapid procedure with less hands-on time
- Reduced well-to-well variation
- Ease of use with both adherent and non-adherent cells
- Adaptable for use in 96-well, 384-well, or 1536-well formats

Applications

The FLIPR Potassium Assay Kit provides a homogeneous assay for measuring potassium channel activities in a no-wash format. It is designed to work for the majority of ligand-gated or voltage-gated potassium channels.

Chapter 2: Materials and Equipment

Kit Components

Table 2-1: Components of the FLIPR Potassium Assay Kits

Item	Evaluation Kit (R8330)	Explorer Kit (R8222)	Bulk Kit (R8223)
Component A	2 vials	10 vials	10 vials
Component C	2 vials	10 vials	10 vials
Chloride-free Assay Buffer (5X)	1 bottle	1 bottle	1 bottle
Potassium Sulfate solution 200 mM of K ₂ SO4 in water	1 bottle	1 bottle	1 bottle
Thallium Sulfate solution 50 mM of Tl ₂ SO4 in water	1 bottle	1 bottle	1 bottle
Assay Buffer (Component B) 20 mM HEPES buffer + 1X Hank's Balanced Salt solution (HBSS), pH 7.4	1 bottle	1 bottle	None

- The entire Evaluation Kit is sufficient for two (2) 96-well or 384-well microplates. The entire Explorer Kit is sufficient for ten (10) 96-well or 384-well microplates. Each vial is sufficient for one (1) microplate.
- The entire Bulk Kit is sufficient for one-hundred (100) 96-well or 384-well microplates. Each vial is sufficient for ten (10) microplates.

Materials Required But Not Provided

Table 2-2: Reagents and Supplies

Item	Suggested Vendor
Dimethyl Sulfoxide (DMSO) is used for dissolving Component A. It is important that the DMSO used is of high quality and stored properly.	Sigma #D8418, or equivalent
Component B*: 20 mM HEPES buffer + 1X Hank's Balanced Salt solution (HBSS), pH 7.4 *Component B is provided for Evaluation kits and Explorer kits only.	10X Hank's Balanced Salt Solution (Gibco #14065-056, or equivalent) 1M HEPES buffer solution (Irvine Scientific #9319, or equivalent) Water for cell culture (Irvine Scientific #9312, or equivalent)
FLIPR® validated pipette tips, 96, 384 or 1536 gasket	Molecular Devices
Probenecid, an inhibitor for the anion-exchange protein, might be required for some cell lines to ensure that the dye stays inside the cell and is not pumped back out. Prepare a stock solution of 500 mM in 1N NaOH, and then dilute to 250 mM in HH buffer. Prepare Loading Buffer such that the final in-well concentration of probenecid is 2.5 mM when added to cells.	Sigma #P8761, or other chemical suppliers Tip: Use of water-soluble probenecid is also possible following individual manufacturer instructions.

Storage and Handling

On receipt of the FLIPR Potassium Assay Kit, store Component A and Component C at -20° C (-4°F), and store the other components at 4°C. Under these conditions, the reagents are stable for six (6) months in the original packaging.

After reconstitution, the Loading Buffer is stable for up to eight (8) hours at room temperature. Aliquots can be frozen and stored (without probenecid) for up to one (1) week without loss of activity.



WARNING! Reagents can contain chemicals that are harmful. Exercise care when handling reagents as described in the related safety data sheet (SDS or MSDS). The safety data sheet is available in the Knowledge Base on the Molecular Devices support web site: www.moleculardevices.com/service-support



DANGER! Thallium sulfate is harmful if swallowed or touched. Disposal of this product, solutions, and any by-products must comply with the requirements of environmental-protection and waste-disposal legislation, and any regional local authority requirements. Exercise care when handling this product as described in the related safety data sheet (SDS or MSDS). The safety data sheet is available in the Knowledge Base on the Molecular Devices support web site: www.moleculardevices.com/service-support

Compatible Molecular Devices Instruments

The FLIPR Potassium Assay Kit is designed to be used with the following Molecular Devices instruments and their corresponding tips:

- FLIPR Penta High-Throughput Cellular Screening System
- FLIPR[®] Tetra High-Throughput Cellular Screening System
- FlexStation 3 Multi-Mode Microplate Reader

Chapter 3: Experimental Protocol

Quick Start Protocol

To run a potassium assay protocol:

- 1. Plate the cells in microplates and incubate overnight at 37°C, 5% CO₂.
- 2. Prepare the Loading Buffer the following day.
- 3. Remove the cell plates from the incubator and add an equal volume of Loading Buffer to each well. For example, 25 μL of Loading Buffer to 25 μL of cells and media for a 384-well microplate.
- 4. Incubate the prepared plates for one hour at room temperature in the dark.
- 5. Prepare the compound plates.
- 6. Run the experiment on a FLIPR Penta instrument.

Cell Handling

The Molecular Devices FLIPR Potassium Assay Kit is designed to work with many cell types, both adherent and non-adherent. Standard procedures vary across laboratories, and Molecular Devices recognizes that a variety of cell handling conditions can be used at the discretion of the user. This section provides general guidelines for preparing cells for use with the FLIPR Potassium Assay Kit.

Adherent cells are the most frequently used cells with the kits. They are generally plated the day before an experiment and then incubated in a 37° C, 5% CO₂ incubator overnight. See Table 3-1 for suggested plating volumes and seeding densities to create an 80% to 90% confluent cell monolayer for the assay in a FLIPR[®] Penta instrument.

Cell Type (cells/well)	96-well microplate (100 μL growth medium)	384-well microplate (25 µL growth medium)
Adherent cells	20000 to 80000	5000 to 20000
Non-adherent cells	40000 to 200000	10000 to 50000

Table	3-1:	Suggested	Plating	Volumes	and	Seeding	Densities
Table	3-T.	Juggesteu	indung	volumes	ana	Jecume	Densities

For non-adherent cells, Molecular Devices recommends centrifuging cells from culture medium and re-suspending the pellet in the culture medium or applicable buffer of choice on the day of the experiment. Cells can be dye-loaded in a tube or while plated. Molecular Devices recommends that after the cells are plated, centrifuge the microplates at 100 x g for up to 4 minutes, with the brake off. As an alternative, non-adherent cells can be treated like adherent cells, plating the day before the assay using the same plating volumes and seeding densities, as long as the cells are seeded onto coated plates (for example, poly-D-lysine or collagen) to make sure that they are well attached to the microplate bottom.

Preparing the Loading Buffer

To prepare the loading buffer:

- 1. Remove one vial each of Component A and Component C from the freezer, and then equilibrate to room temperature.
- Equilibrate Component B to room temperature. For the Bulk Kit, prepare 100 mL of 20 mM HEPES plus 1X HBSS, pH 7.4 as Component B.
- 3. Dissolve the contents of the Component C vial in DMSO, and then mix thoroughly by vortexing.
 - For the Evaluation Kit or Explorer Kit, use 30 μL of DMSO.
 - For the Bulk Kit, use 300 μL of DMSO.
- Combine the vial of Component A with 10 mL of the Component B buffer. For the Bulk Kit, use only 10 mL of the prepared Component B buffer in this step.
- Combine the Component C solution from step 3 to the Component A solution from step 4, and then mix by vortexing for ~1 to 2 minutes until the contents of the vial are dissolved.

It is important that the contents are completely dissolved to ensure reproducibility between experiments.

6. For Bulk kit only, combine the solution from step 5 with the remaining 90 mL of the prepared Component B buffer, and then mix thoroughly.

		1	1		
í		-	-		
I	F			-	
I	H	-	-		
J	-	-	-	-	

Note: If the cells require probenecid (such as CHO or other cells containing an organic anion transporter), then add probenecid in the Loading Buffer to 5 mM so that the final in-well working concentration is 2.5 mM. Adjust Loading Buffer pH to 7.4 after addition of probenecid. See the procedure for making probenecid in Materials Required But Not Provided on page 5.

Ē

Note: Do not store frozen aliquots of Loading Buffer with probenecid, and always prepare new probenecid on the day of the experiment. Use of watersoluble probenecid is also possible following individual manufacturer instructions.

Loading Cells Using Loading Buffer

To load cells using loading buffer:

- 1. Remove the cell plates from the incubator or centrifuge.
- 2. Add an equal volume of Loading Buffer to each well: 100 μL per well for 96-well plates and 25 μL per well for 384-well plates.



Note: Molecular Devices does not recommend washing cells before dye loading. However, growth medium and serum might interfere with certain assays. In this case, the supernatant can be aspirated and replaced with an equal volume of serum-free 20mM HEPES buffer + 1X HBSS, pH 7.4 before adding the Loading Buffer.

3. After adding dye, incubate the cell plates for 1 hour at room temperature. The loading time should be optimized for each cell line and target.



Note: Do NOT wash the cells after dye loading.

Preparing the Thallium Sulfate Source Plate

To introduce TI⁺ into the assay plate, prepare a 5X solution by diluting ligands or K_2SO_4 (for voltage-gated targets only) with TI_2SO_4 in 1X Chloride-free Buffer.

Concentrations of Tl_2SO_4 and K_2SO_4 (for voltage-gated targets only) used for assays should be optimized for each target channel. Molecular Devices suggests a titration of Tl^+ or a combination of Tl^+ and K^+ at 0.5 mM to 3 mM Tl^+ and 5 mM to 30 mM K^+ (final concentration in wells). See Optimization Guidelines on page 14.

Note: Since thallium sulfate and potassium sulfate have two equivalents of cation per mole, consider them as 2X in their respective cation concentrations.



DANGER! Thallium sulfate is harmful if swallowed or touched. Disposal of this product, solutions, and any by-products must comply with the requirements of environmental-protection and waste-disposal legislation, and any regional local authority requirements. Exercise care when handling this product as described in the related safety data sheet (SDS or MSDS). The safety data sheet is available in the Knowledge Base on the Molecular Devices support web site: www.moleculardevices.com/service-support

Running the FLIPR Potassium Assay

After incubation, transfer the microplates to the FLIPR instrument and start the FLIPR Potassium Assay.

Do the assay signal test before the experiment. The LED intensity, exposure time, and the gate or gain can be adjusted to get the desired RFU range.

- For a FLIPR Tetra System or FLIPR Penta instrument with an EMCCD camera, adjust the typical average baseline counts to a range from 1500 RFU to 2500 RFU.
- For a FLIPR Penta instrument with an HS EMCCD camera, adjust the typical average baseline counts to a range from 2,500 to 5,000 RFU.
- For a FLIPR Tetra System or FLIPR Penta instrument with an ICCD camera, adjust the typical average baseline counts to a range from 8000 RFU to 10000 RFU.

Update the assay signal test to transfer the adjusted settings to the protocol.

Before reading the microplate, set up the FLIPR instrument using the recommended experimental setup parameters listed in Table 3-2. and Table 1.

Faster addition speeds closer to the cell monolayer are recommended to ensure better mixing of compounds and lower signal variance across the plate. Molecular Devices recommends further assay development and adjustment of the volume, height, and speed of dispense to optimize the individual cell response.

Recommended Settings for the FLIPR Instrument

Before reading the microplate, set up your FLIPR system using the following recommended experiment settings. Your settings depend on the camera installed in your instrument.

Parameter	96-well microplate	384-well microplate	
Excitation LED (nm)	470–495	470–495	
Emission Filter (nm)	515–575	515–575	
Camera Mode	EMCCD camera: NA HS EMCCD camera: Normal ICCD camera: NA	EMCCD camera :NA HS EMCCD camera: Normal ICCD camera: NA	
Camera Gain	EMCCD camera: 70 to 150 HS EMCCD camera: 4 ICCD camera: Fixed at 2000	EMCCD camera: 70 to 150 HS EMCCD camera: 4 ICCD camera: Fixed at 2000	
Exposure (sec)	EMCCD camera: 0.1 to 0.5 HS EMCCD camera: 0.2 ICCD camera: 0.4	EMCCD camera: 0.1 to 0.5 HS EMCCD camera: 0.2 ICCD camera: 0.04	
LED Intensity (%)	EMCCD camera: 50 to 100 HS EMCCD camera: 40 ICCD camera: 40	EMCCD camera: 50 to 100 HS EMCCD camera: 40 ICCD camera: 40	
Camera Gate (%) (ICCD camera only)	6	6	
Addition Volume (µL)	50	12.5	

Table 3-2: Experimental setup parameters for the FLIPR Penta instrument

Parameter	96-well microplate	384-well microplate
Compound Concentration (Fold)	бх	6х
Addition Height (µL)	120 to 160	30 to 40
Addition Speed (µL/sec)	Adherent Cells: 60 to 100 Non-Adherent Cells: 20 to 40	Adherent Cells: 30 to 40 Non-Adherent Cells: 10 to 20
Tip-Up Speed (mm/sec)	20	20

Table 3-2: Experimental setup parameters for the FLIPR Penta instrument (continued)

Running the FLIPR Potassium Assay on a FlexStation 3 Instrument

Before reading the microplate, set up the FlexStation 3 instrument using a SoftMax[®] Pro Software protocol. The recommended experimental setup parameters are listed in Table 3-3: Experiment Setup Parameters for the FlexStation 3 Instrument on page 12.

Molecular Devices recommends that parameters be optimized for each cell line and target to deliver the best performance for your assay.

After incubation, transfer the microplates directly to the FlexStation 3 instrument assay plate carriage and run the assay.

In an individual well or column of wells, the peak(s) should be complete within 1 to 3 minutes after addition. For an entire plate however, the protocol will not complete until all chosen columns are finished. The assays are run one column at a time.

Analyze the data using the SoftMax Pro Software.

Recommended Settings for the FlexStation 3 Instrument

Parameter	96-well microplate	384-well microplate
Read Mode	Fluorescence	Fluorescence
Read Position	Bottom Read	Bottom Read
Excitation Wavelength (nm)	485	485
Emission Wavelength (nm)	525	525
Automatic Emission Cut-Off (nm)	515	515
PMT Sensitivity	6	6
Pipette Height (μL)	230	50
Transfer Volume (μL)	50	12.5
Compound Concentration (Fold)	5X	5X
Addition Speed (Rate)	Adherent Cells: 3 Non-Adherent Cells: 1	Adherent Cells: 2 to 3 Non-Adherent Cells: 1

Table 3-3: Experiment Setup Parameters for the FlexStation 3 Instrument

Chapter 4: Data Analysis Example

Given a state in the state of the state of

Inhibition of hERG Channel

Compound	IC ₅₀ (nM) FLIPR Instrument	IC ₅₀ (nM) Electrophysiology
Terfenadine	274	332
Astemizole	76	62
Pimozide	135	147

In this experiment, the concentration-dependent inhibition of hERG channels by three reference compounds was analyzed with the FLIPR Potassium Assay Kit. The hERG channel was stably expressed in the Chinese hamster ovary (CHO) cells, and all compounds were incubated with cells for 30 minutes after 1 hour of dye loading at room temperature. Stimulus buffer containing 1X chloride-free buffer and final concentrations of 10 mM K⁺, and 1 mM TI⁺ was added to the wells while reading on the FLIPR Tetra Instrument. The IC₅₀ values derived from this assay were compared to the electrophysiology results collected with the IonWorks Barracuda Instrument.

Chapter 5: Optimization Guidelines

Note: Performance of the Molecular Devices Reagent Kits on Molecular Devices instruments have been validated for use with Molecular Devices pipette tips.

Data Analysis

Correct data analysis is crucial to correctly assess the functionality of the assay. It is very important to choose the correct time point to calculate final data. In the assay, the assay background trace (optimized Thallium only or maximal inhibition condition) shows a slow upward slope over the course of the assay. Using a late time point results in a reduced assay window impacting the quality of the data and also the IC_{50} determination.

Molecular Devices recommends using a time point that results in the greatest assay window as defined as Max signal vs. Thallium only trace. Identification of the optimal time point can be done by assessing the raw data manually or by using the Negative Control Correction feature in the FLIPR® ScreenWorks® Software. This feature subtracts the assigned background curve, resulting in bell shaped traces. It can be used for identification of the optimal time point or as a general analysis tool. Using Negative Control Correction, the Max-Min analysis automatically uses the data at the optimal time point.

If other analysis methods are preferred, the optimal time point determined can be manually chosen (maximum of bell shaped curve).

Use of the Negative Control Correction feature to determine the optimal time point and signal window for the assay



Figure 5-1: Before and after Negative Control Correction

 $K_\nu 1.3$ channel activity was measured with 2 mM TI⁺ and different K^+ concentrations. The data are displayed with and without Negative Control Correction.

Determination of Potassium and Thallium Concentrations

Potassium channels are grouped into several classes that have very different properties. It is very important to determine the target-specific optimal concentrations of TI⁺, or TI⁺ and K⁺ for each assay. To determine the optimal concentrations to use for the assay, Molecular Devices recommends setting up an initial experiment with a matrix of ligand and TI⁺ or K⁺ and TI⁺ at 5 mM to 30 mM K⁺ and 0.5 mM to 3M mM TI⁺ (final concentration in the wells).

Titration of K⁺ and TI⁺ concentrations to optimize assay performance for different targets



Figure 5-2: Kv1.3 Channel Activity Assay

 $K_\nu 1.3$ channel activity was measured with a titration of different TI^* and K^* concentrations. The data are displayed with Negative Control Correction.



Figure 5-3: hERG Channel Activity Assay

hERG channel activity was measured with a titration of different TI⁺ and K⁺ concentrations. The data are displayed with Negative Control Correction.

Obtaining Support

Molecular Devices is a leading worldwide manufacturer and distributor of analytical instrumentation, software, and reagents. We are committed to the quality of our products and to fully supporting our customers with the highest level of technical service.

Our Support website, www.moleculardevices.com/service-support, has a link to the Knowledge Base, which contains technical notes, software upgrades, safety data sheets, and other resources. If you still need assistance after consulting the Knowledge Base, you can submit a request to Molecular Devices Technical Support.

You can contact your local representative or Molecular Devices Technical Support at 800-635-5577 (North America only) or +1 408-747-1700. In Europe, call +44 (0) 118 944 8000.

Please have the instrument serial number (on the rear of the instrument), or reagent kit lot number, and any related sample data files available when you call.

Contact Us

Phone: +1-800-635-5577 Web: moleculardevices.com info@moldev.com Email:

Visit our website for a current listing of worldwide distributors.

The trademarks used herein are the property of Molecular Devices, LLC or their respective owners Specifications subject to change without notice. Patents: www.moleculardevices.com/productpatents All rights reserved. FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

R3622 D

