

Screen Quest™ Calbryte-590 Probenecid-Free and Wash-Free Calcium Assay Kit

 Catalog number: 36200, 36201, 36202
 Unit size: 1 Plate, 10 Plates, 100 Plates

Component	Storage	Amount (Cat No. 36200)	Amount (Cat No. 36201)	Amount (Cat No. 36202)
Component A: Calbryte™ 590 AM	Freeze (< -15 °C), Minimize light exposure	1 vial	1 vial	10 vials
Component B: 10X Pluronic® F127 Plus	Freeze (< -15 °C), Minimize light exposure	1 bottle (1 mL)		
Component B: 10X Pluronic® F127 Plus	Freeze (< -15 °C), Minimize light exposure		1 bottle (10 mL)	10 bottles (10 mL/bottle)
Component C: HHBS (Hanks' buffer with 20 mM Hepes)	Freeze (< -15 °C), Minimize light exposure	1 bottle (9 mL)	1 bottle (100 mL)	Not provided

OVERVIEW

Screen Quest™ Calbryte™ 590 Probenecid-Free and Wash-Free Calcium Assay Kit is designed for red fluorescence-based detection of intracellular calcium mobilization in live cells. Calcium flux assays are widely used in drug discovery for screening G protein-coupled receptors (GPCRs) and calcium channels. The kit uses Calbryte™ 590 AM, a cell-permeant calcium indicator that enters cells and is converted by intracellular esterases into a calcium-responsive fluorescent dye. Calbryte™ 590 AM is designed for improved cellular retention, enabling calcium flux detection without probenecid and without wash steps. Upon cell stimulation, receptor-mediated calcium release or calcium influx produces an increase in red fluorescence that can be measured in a microplate format. Compared with green fluorescence-based calcium assays, the red fluorescence readout can help reduce interference from colored compounds in screening libraries and is compatible with GFP-expressing cell lines for high-content analysis. Screen Quest™ Calbryte™ 590 Probenecid-Free and Wash-Free Calcium Assay Kit is suitable for GPCR calcium signaling studies, calcium channel assays, GFP-compatible calcium imaging, high-throughput screening, and calcium flux analysis in fragile or difficult cell lines.

AT A GLANCE
Protocol Summary

1. Prepare cells in growth medium
2. Add Calbryte™ 590 AM dye-loading solution (100 µL/well for 96-well plate or 25 µL/well for 384-well plate)
3. Incubate at room temperature or 37°C for 45-60 minutes
4. Monitor fluorescence at Ex/Em = 540/590 nm

Important Note

Thaw all the kit components at room temperature before use.

KEY PARAMETERS
Fluorescence microplate reader

Excitation	540 nm
Emission	590 nm
Cutoff	570 nm
Recommended plate	Black wall/clear bottom
Instrument specification(s)	Bottom read mode/programmable liquid handling

Other Instruments: FDSS, NOVOSTar, FlexStation, ViewLux, IN Cell Analyzer, ArrayScan

CELL PREPARATION

For guidelines on cell sample preparation, please visit <https://www.aatbio.com/resources/guides/cell-sample-preparation.html>

PREPARATION OF STOCK SOLUTIONS

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles

Calbryte™ 590 AM stock solution

Add 20 µL (**Cat. # 36200**) or 200 µL (**Cat. # 36201 and # 36202**) of DMSO into the vial of Calbryte™ 590 AM (Component A) and mix them well.

Note: 20 µL of Calbryte™ 590 AM stock solution is enough for one plate. Unused Calbryte™ 590 AM stock solution can be aliquoted and stored at ≤ -20 °C for more than one month if the tubes are sealed tightly.

Note: Protect from light and avoid repeated freeze-thaw cycles.

Assay buffer (1X)

Mix 9 mL of HHBS (Component C, not included in the kit **Cat. # 36202**) with 1 mL of 10X Pluronic® F127 Plus (10X) (Component B) and mix them well.

PREPARATION OF WORKING SOLUTION
Calbryte™ 590 AM dye-loading solution

Add 20 µL of Calbryte™ 590 AM stock solution into 10 mL of Assay Buffer (1X) and mix them well.

Note: This working solution is stable for at least 2 hours at room temperature.

Note: 10 mL dye-loading solution is enough for one 96-wells plate.

SAMPLE EXPERIMENTAL PROTOCOL

1. Add 100 µL/well (96-well plate) or 25 µL/well (384-well plate) of Calbryte™ 590 AM dye-loading solution into the cell plate.
2. Incubate the dye-loading plate in a cell incubator for 60 minutes, and then incubate the plate at room temperature for another 15 - 30 minutes.

Note: If the assay requires 37°C, perform the experiment

immediately without further room temperature incubation. If the cells can function well at room temperature for longer time, incubate the cell plate at room temperature for 1 hour (It is recommended that the incubation time be no longer than 2 hours.)

3. Prepare the compound plate with HHBS or your desired buffer.
4. Run the calcium flux assay by monitoring the fluorescence intensity at Ex/Em = 540/590 nm.

EXAMPLE DATA ANALYSIS AND FIGURES

The reading (SNR x 100%) obtained from the blank standard well is used as a negative control. Subtract this value from the other standards' readings to obtain the base-line corrected values. Then, plot the standards' readings to obtain a standard curve and equation. This equation can be used to calculate ATP samples. We recommend using the Online Four Parameter Logistics Calculator which can be found at:

<https://www.aatbio.com/tools/four-parameter-logistic-4pl-curve-regression-online-calculator>

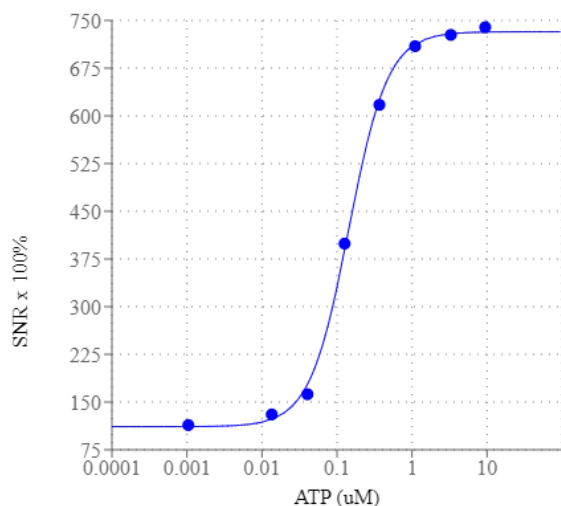


Figure 1. Graph illustrates signal-to-noise ratio (SNR) x 100%. ATP dose response was measured in CHO-K1 cells with Screen Quest™ Calbryte 590 Probenecid-Free and Wash-Free Calcium Assay Kit. CHO-K1 cells were seeded overnight at 50,000 cells/100 μ L/well in a 96-well black wall/clear bottom costar plate. 100 μ L dye loading solution was added and incubated for 45 min at 37°C and 15 min at RT. ATP (50 μ L/well) was added by FlexStation 3 to achieve the final indicated concentrations.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.