

## ab112129 Calcium Flux Assay Kit (Fluo-8, No Wash)

For detecting calcium in cells by using our proprietary fluorescence probe  
This product is for research use only and is not intended for diagnostic use.

### Storage and Stability

On receipt entire assay kit should be stored at -20°C, protected from light.

### Materials Supplied

Item	Amount
Component A: Fluo-8	1 vial
Component B: 10X Pluronic® F127 Plus	1 mL
Component C: HHBS	9 mL

### Materials Required, Not Supplied

- 96 or 384-well microplate: Tissue culture microplate with black wall and clear bottom.
- HTS fluorescence microplate reader with a filter set of Ex/Em = 490/525 nm. We recommend the following instruments: FLIPR™, FDSS, BMG NOVOstar™, FlexStation, ViewLux, IN Cell Analyzer or Arrayscan.
- HHBS (1X Hank's with 20 mM Hepes Buffer, pH 7.0).
- 100% DMSO
- (Optional) Carbachol (ab141354) or other GPCR agonists to induce agonist-mediated calcium release in cells.

### Assay Procedure

- *Warning: Do not add additional probenecid.*

### Prepare Cells

1. For adherent cells: Plate cells overnight in growth medium with 5-10% FBS at 40,000 to 80,000 cells/well/100 µL for a 96-well plate or 10,000 to 20,000 cells/well/25 µL for a 384-well plate.
2. For non-adherent cells: Centrifuge the cells from the culture medium and then suspend the cell pellet in HHBS at 125,000 to 250,000 cells/well/100 µL for a 96-well poly-D-lysine plate or 30,000 to 60,000 cells/well/25 µL for a 384-well poly-D-lysine plate. Centrifuge the plate at 800 rpm for 2 minutes with brake off prior to the experiments.

Δ **Note:** Each cell line should be evaluated on an individual basis to determine the optimal cell density for the intracellular calcium mobilization.

### Preparation of Fluo-8 dye-loading solution

1. Thaw all the kit components at room temperature before use.
2. Make Fluo-8 stock solution: Add 20 µL of DMSO into Fluo-8 (Component A) and mix them well.

Δ **Note:** 20 µL of Fluo-8 stock solution is enough for one plate. Unused Fluo-8 stock solution can be aliquoted and stored at -20°C for more than one month if the tubes are sealed tightly. Keep protected from light and avoid repeated freeze-thaw cycles.

3. Make 1X assay buffer: Add 9 mL of HHBS (Component C) into 1 mL of 10X Pluronic® F127 Plus (Component B) and mix them well.

Δ **Note:** 10 mL of 1X assay buffer is suitable for one plate. Aliquot and store unused 1X assay buffer at -20°C. Keep from light and avoid repeated freeze-thaw cycles.

4. Make Fluo-8 dye-loading solution for one cell plate: Add 20 µL of Fluo-8 stock solution (from Step B.2) into 10 mL of 1X assay buffer (from Step B.3) and mix them well. This working solution is stable for at least 2 hours at room temperature.

### Run Calcium Assay

1. Add 100 µL/well (96-well plate) or 25 µL/well (384-well plate) of Fluo-8 dye-loading solution (from Step B.4) into the cell plate.

Δ **Note:** Alternatively, grow the cells in growth medium with 5 - 10% FBS to improve cell growth. In this case, it is important to replace the growth medium with HHBS buffer in order to minimize background fluorescence, and compound interference with serum.

2. Incubate the dye-loading plate in a cell incubator for 30 minutes, and then incubate the plate at room temperature for another 30 minutes.

Δ **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 1-2 hours.

3. Prepare the compound plates with HHBS or the desired buffer.
4. Using your instrument that contains a pipettor (such as Flexstation from Molecular Devices), dispense compounds directly onto the cell plate while collecting the data simultaneously (see next step for Ex/Em instructions).

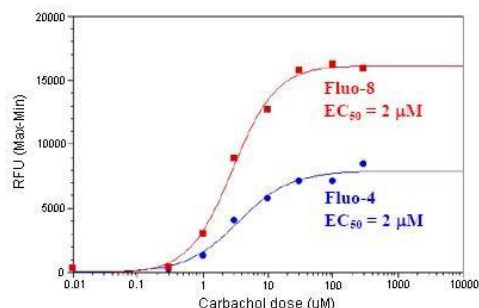
Δ **Note:** Typically the instrument collects the data every second for 100 seconds. Use the max signal to generate the plot. You can use a fluorescence microscope by adding the stimuli while taking the picture simultaneously.

5. Run the calcium flux assay by monitoring the fluorescence intensity at Ex/Em = 490/525 nm.

Δ **Note:** It is important to run the signal test before the experiment. Different instruments have their own range. Adjust the signal test intensity to the level of 10% to 15% of the maximum instrument intensity counts.

### Data Analysis

The reading (RFU (Max-Min)) 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 Carbachol dose samples.



**Figure 1.** Carbachol Dose Response was measured in HEK-293 cells with ab112129. HEK-293 cells were seeded overnight at 40,000 cells/100  $\mu$ L/well in black wall/clear bottom 96-well plate. The cells were incubated with 100  $\mu$ L of dye-loading solution using ab112129 for 1 hour at room temperature. Carbachol (50  $\mu$ L/well) was added to achieve the final indicated concentrations.

### Troubleshooting

Problem	Reason	Solution
Assay not working	Assay buffer at wrong temperature	Assay buffer must not be chilled - needs to be at RT
	Protocol step missed	Re-read and follow the protocol exactly
	Plate read at incorrect wavelength	Ensure you are using appropriate reader and filter settings (refer to datasheet)
	Unsuitable microtiter plate for assay	Fluorescence: Black plates (clear bottoms); Luminescence: White plates; Colorimetry: Clear plates. If critical, datasheet will indicate whether to use flat- or U-shaped wells
Unexpected results	Measured at wrong wavelength	Use appropriate reader and filter settings described in datasheet
	Samples contain impeding substances	Troubleshoot and also consider deproteinizing samples
	Unsuitable sample type	Use recommended sample types as listed on the datasheet
	Sample readings are outside linear range	Concentrate/ dilute samples to be in linear range
Samples with inconsistent readings	Unsuitable sample type	Refer to datasheet for details about incompatible samples
	Samples prepared in the wrong buffer	Use the assay buffer provided (or refer to datasheet for instructions)
	Samples not deproteinized (if indicated on datasheet)	Use the <b>10kDa spin column (ab93349)</b> or appropriate deproteinization protocol
	Too many freeze-thaw cycles	Aliquot samples to reduce the number of freeze-thaw cycles

Lower/ Higher readings in samples and standards	Samples contain impeding substances	Troubleshoot and also consider deproteinizing samples
	Samples are too old or incorrectly stored	Use freshly made samples and store at recommended temperature until use
	Not fully thawed kit components	Wait for components to thaw completely and gently mix prior use
	Out-of-date kit or incorrectly stored reagents	Always check expiry date and store kit components as recommended on the datasheet
	Reagents sitting for extended periods on ice	Try to prepare a fresh reaction mix prior to each use
	Incorrect incubation time/ temperature	Refer to datasheet for recommended incubation time and/ or temperature
Standard curve is not linear	Incorrect amounts used	Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume)
	Not fully thawed kit components	Wait for components to thaw completely and gently mix prior use
	Pipetting errors when setting up the standard curve	Try not to pipette too small volumes
	Incorrect pipetting when preparing the reaction mix	Always prepare a master mix
	Air bubbles in wells	Air bubbles will interfere with readings; try to avoid producing air bubbles and always remove bubbles prior to reading plates
	Concentration of standard stock incorrect	Recheck datasheet for recommended concentrations of standard stocks
	Errors in standard curve calculations	Refer to datasheet and re-check the calculations
	Use of other reagents than those provided with the kit	Use fresh components from the same kit

### Technical Support

Copyright © 2025 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

**For all technical or commercial enquiries please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)