

**ab171868**

**Cal-520, AM**

## Instructions for Use

For detecting calcium in cells using a proprietary fluorescence probe

View kit datasheet: [www.abcam.com/ab171868](http://www.abcam.com/ab171868)

(use [www.abcam.cn/ab171868](http://www.abcam.cn/ab171868) for China, or  
[www.abcam.co.jp/ab171868](http://www.abcam.co.jp/ab171868) for Japan)

This product is for research use only and is not intended for diagnostic use.

# Table of Contents

---

1.	Introduction	2
2.	Kit Contents	3
3.	Storage and Handling	3
5.	Assay Protocol	4
6.	Data Analysis	6
7.	Troubleshooting	8

# 1. Introduction

---

Cal-520, AM (ab171868) provides a robust homogeneous fluorescence-based assay tool for detecting intracellular calcium mobilization. Cal-520 AM is a new fluorogenic calcium-sensitive dye with a significantly improved signal to noise ratio and intracellular retention compared to the existing green calcium indicators (such as Fluo-3 AM and Fluo-4 AM). Cells expressing a GPCR or calcium channel of interest that signals through calcium can be preloaded with Cal-520, AM which can cross cell membrane. Once inside the cell, the lipophilic blocking groups of Cal 520, AM are cleaved by esterases, resulting in a negatively charged fluorescent dye that stays inside cells. Its fluorescence is greatly enhanced upon binding to calcium. When cells stimulated with agonists, the receptor signals the release of intracellular calcium, which significantly increase the fluorescence of Cal-520, AM. The characteristics of its long wavelength, high sensitivity, and >100 times fluorescence enhancement, make Cal-520, AM an ideal indicator for the measurement of cellular calcium. The high S/N ratio and better intracellular retention make the Cal-520, AM calcium assay a robust tool for evaluating GPCR and calcium channel targets as well as for screening their agonists and antagonists.

## **Features**

**Organic anion transporter resistance:** Enable Ca<sup>2+</sup> assays with probenecid-interference GPCRs and Ca<sup>2+</sup> channels.

**Robust:** Significantly higher S/N ratio than those of Fluo-4 AM and any other commercially available fluorescent Ca<sup>2+</sup> assays.

**Convenient:** maximum excitation 492 nm; maximum emission 514 nm.

## **2. Kit Contents**

---

---

<b>Components</b>	<b>Amount</b>
Cal-520, AM	1 mg

---

## **3. Storage and Handling**

---

Keep at -20°C. Protect from light.

## 4. Additional Materials Required

---

- A 96 or 384-well microplate: Tissue culture microplate with black wall and clear bottom.
- A fluorescence microplate reader with a filter set of Ex/Em = 490/525 nm.
- HHBS (1X Hank's with 20 mM Hepes Buffer, pH 7.0).
- 100% DMSO anhydrous.
- Pluronic F127
- Anion transporter inhibitor (optional)

## 5. Assay Protocol

---

This protocol for loading Cal-520, AM esters into live cells. This protocol only provides a guideline, and should be modified according to your specific needs. please read the whole protocol before starting

1. Prepare a 2 to 5 mM stock solution of Cal-520, AM esters in high-quality, anhydrous DMSO.
2. On the day of the experiment, either dissolve Cal-520 in DMSO or thaw an aliquot of the indicator stock solution to room temperature. Prepare a Cal-520 dye working solution of 5 to 10  $\mu$ M in Hanks and Hepes buffer

(HHBS) or the buffer of your choice with 0.02% Pluronic® F-127. The exact concentration of the indicator required for cell loading must be determined empirically.

*Note: The nonionic detergent Pluronic® F-127 is sometimes used to increase the aqueous solubility of Cal-520, AM esters.*

3. If your cells (such as CHO cells) contain the organic anion-transporters, probenecid (0.5–1 mM) may be added to the cell medium to reduce leakage of the de-esterified indicators
4. Incubate the dye-loading plate in a cell incubator for 60 to 90 minutes, and then incubate the plate at room temperature for a further 30 minutes.  
*Note: Incubation with the dye for longer than 2 hours gives a better signal intensity for some cell lines.*
5. Replace the Cal-520, AM dye working solution with HHBS or buffer of your choice (containing an anion transporter inhibitor, such as 1 mM probenecid, if applicable) to remove excess probes.
6. Run the experiments at  $\text{Ex/Em} = 490/525 \text{ nm}$  at room temperature.

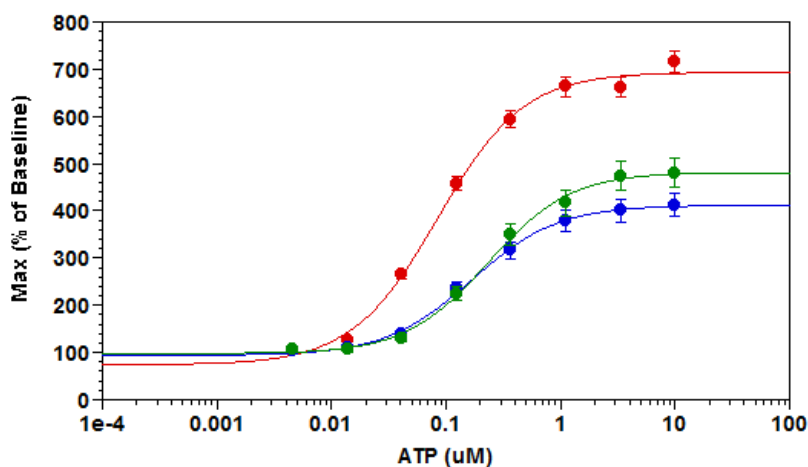
## 6. Data Analysis

---

To determine either the free calcium concentration of a solution or the  $K_d$  of a single-wavelength calcium indicator, the following equation is used:

$$[Ca]_{\text{free}} = K_d[F - F_{\text{min}}]/F_{\text{max}} - F]$$

Where  $F$  is the fluorescence of the indicator at experimental calcium levels,  $F_{\text{min}}$  is the fluorescence in the absence of calcium and  $F_{\text{max}}$  is the fluorescence of the calcium-saturated probe.



**Figure 1. ATP dose response comparison in CHO-K1 cells measured with Cal-520, AM No Wash, FLIPR Calcium 4 or Fluo-4 Direct Calcium Assay Kit.** CHO-K1 cells were seeded overnight in 60,000 cells per 100  $\mu$ L per well in a 96-well black wall/clear bottom costar plate. The cells were incubated with 100  $\mu$ L of the Cal-520, AM No Wash Calcium Assay Kit(Red), FLIPR Calcium 4 Kit(Green) or Fluo-4 Direct Calcium Assay Kit(Blue) for 1.5 hours at 37°C. ATP (50 $\mu$ L/well) was added to achieve the final indicated concentrations.

## 7. 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 samples 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 <b>Deproteinizing sample preparation kit (ab93299)</b>
	Cell/ tissue samples not sufficiently homogenized	Increase sonication time/ number of strokes with the Dounce homogenizer
	Too many freeze-thaw cycles	Aliquot samples to reduce the number of freeze-thaw cycles
	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
Lower/ Higher readings in samples and standards	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
	Incorrect amounts used	Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume)

<b>Problem</b>	<b>Reason</b>	<b>Solution</b>
Standard curve is not linear	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)