

Cal Green[™] 1, AM [Equivalent to Calcium Green-1, AM]

Catalog number: 20501, 20502 Unit size: 10x50 ug, 1 mg

Component	Storage	Amount (Cat No. 20501)	Amount (Cat No. 20502)
Cal Green™ 1, AM [Equivalent to Calcium Green	Freeze (< -15 °C), Minimize light exposure	10x50 ug	1 vial (1 mg)
-1, AM]			

OVERVIEW

Cal Green[™] 1 (AAT Bioquest) is the same molecule to Calcium Green-1 (Invitrogen). It exhibits an increase in fluorescence intensity upon binding calcium ion. The cell-permeant dye, Cal Green 1 AM (Calcium Green-1 AM) is a 488 nm-excitable calcium indicator. Compared to Fluo-3 AM, Cal Green-1 AM is more fluorescent at low calcium concentrations in cells, facilitating the determination of baseline calcium levels and increasing the visibility of resting cells. It has been used in many calcium signaling investigations, including measuring intracellular calcium, following calcium influx and release, and multiphoton excitation imaging of calcium in living tissues. Cells can be loaded with Cal Green-1 AM by adding the dissolved indicator directly to cultured cells in medium. The fluorescence signal from these cells is generally measured using fluorescence microscopy, fluorescence microplate assays, or flow cytometry.

KEY PARAMETERS

Fluorescence microscope

Excitation	FITC
Emission	FITC
Recommended plate	Black wall/clear bottom

Fluorescence microplate reader

Excitation	490
Emission	525
Cutoff	515
Recommended plate	Black wall/clear bottom
Instrument specification(s)	Bottom read mode/Programmable liquid
	handling

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.

Cal Green[™] 1 AM Stock Solution

Prepare a 2 to 5 mM stock solution of Cal Green[™] 1 AM in high-quality, anhydrous DMSO.

PREPARATION OF WORKING SOLUTION

Cal Green[™] 1 AM Working Solution

On the day of the experiment, either dissolve Cal GreenTM 1 AM in DMSO or thaw an aliquot of the indicator stock solution to room temperature. Prepare a dye working solution of 2 to 20 μ M in a buffer of your choice (e.g., Hanks and Hepes buffer) with 0.04% Pluronic® F-127. For most cell lines, Cal GreenTM 1 AM at a final concentration of 4-5 μ M is recommended. The exact concentration of indicators 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 Green[™] 1 AM. A variety of Pluronic® F-127 solutions can be purchased from AAT Bioquest.

Note If your cells contain organic anion-transporters, probenecid (1-2 mM) may be added to the dye working solution (final in well concentration will be 0.5-1 mM) to reduce leakage of the de-esterified indicators. A variety of ReadiUse™

probenecid products, including water-soluble, sodium salt, and stabilized solution, can be purchased from AAT Bioquest.

SAMPLE EXPERIMENTAL PROTOCOL

Following is our recommended protocol for loading AM esters into live cells. This protocol only provides a guideline and should be modified according to your specific needs.

- 1. Prepare cells in growth medium overnight.
- On the next day, add 1X Cal Green[™] 1 AM working solution into your cell plate.

Note If your compound(s) interfere with the serum, replace the growth medium with fresh HHBS buffer before dye-loading.

 Incubate the dye-loaded plate in a cell incubator at 37 °C for 30 to 60 minutes.

Note Incubating the dye for longer than 1 hour can improve signal intensities in certain cell lines.

- Replace the dye working solution with HHBS or buffer of your choice (containing an anion transporter inhibitor, such as 1 mM probenecid, if applicable) to remove any excess probes.
- Add the stimulant as desired and simultaneously measure fluorescence using either a fluorescence microscope equipped with a FITC filter set or a fluorescence plate reader containing a programmable liquid handling system such as an FDSS, FLIPR, or FlexStation, at Ex/Em = 490/525 nm cutoff 515 nm.

EXAMPLE DATA ANALYSIS AND FIGURES



Figure 1. Chemical structure for Cal Green[™] 1, AM [Equivalent to Calcium Green-1, AM]

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.