

**Calcein, AM \*UltraPure grade\* \*CAS 890090-35-4\***

 Catalog number: 22003, 22004  
 Unit size: 1 mg, 20x50 ug

Component	Storage	Amount (Cat No. 22003)	Amount (Cat No. 22004)
Calcein, AM *UltraPure grade* *CAS 890090-35-4*	Freeze (< -15 °C), Minimize light exposure	1 vial (1 mg)	20x50 ug

**OVERVIEW**

Calcein-AM is a green fluorogenic dye that stains live cells for cell viability in flow cytometry and fluorescence microscopy. It is the acetoxymethyl (AM) ester form of the fluorescent probe, calcein.

**KEY PARAMETERS**
**Flow cytometer**

Emission	530/30 nm filter
Excitation	488 nm laser
Instrument specification(s)	FITC channel

**Fluorescence microscope**

Emission	FITC filter set
Excitation	FITC filter set
Recommended plate	Black wall/clear bottom

**Fluorescence microplate reader**

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

**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

**Calcein, AM \*UltraPure grade\* stock solution**

1. Prepare a 2 to 5 mM stock solution of Calcein AM in high-quality, anhydrous DMSO.

**Note:** When reconstituted in DMSO, Calcein AM is a clear, colorless solution.

**PREPARATION OF WORKING SOLUTION**
**Calcein, AM \*UltraPure grade\* working solution**

1. Prepare a Calcein AM working solution of 1 to 10 µM in the buffer of your choice (e.g., Hanks and Hepes buffer). For most cell lines, Calcein AM at the final concentration of 4 to 5 µM is recommended. The exact concentration of indicators required for cell loading must be determined empirically.

**Note:** The nonionic detergent Pluronic® F-127 can be used to increase the aqueous solubility of AM esters. In the staining buffer, the final Pluronic® F-127 concentration should be approximately 0.02%. A variety of [Pluronic® F-127 products](#) can be purchased from AAT Bioquest. Avoid long-term storage of AM esters in the

presence of Pluronic® F-127.

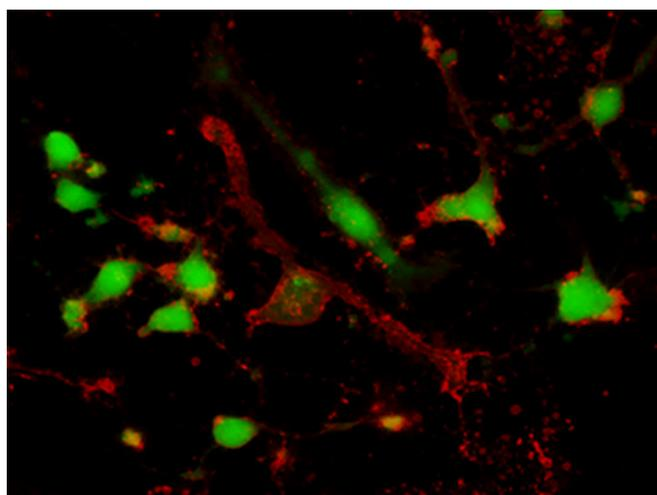
**Note:** If your cells contain organic anion-transporters, probenecid (1–2.5 mM) or sulfapyrazone (0.1–0.25 mM) may be added to the working solution to reduce leakage of the de-esterified indicators. A variety of [ReadiUse™ Probenecid products](#), including water-soluble, sodium salt, and stabilized solutions, can be purchased from AAT Bioquest.

**SAMPLE EXPERIMENTAL PROTOCOL**

1. Prepare cells for imaging.
2. Remove the cell culture medium and wash cells once with serum-free buffer to remove any remaining media.

**Note:** Serum in cell culture media may contain esterase activity, which can increase background interference.

3. Add Calcein AM working solution to the culture.
4. Incubate cells at 37 °C for 30 to 60 minutes.
5. 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.
6. Measure the fluorescence intensity using either a fluorescence microscope equipped with a FITC filter set, a flow cytometer equipped with a blue laser and a 530/30 nm filter (FITC channel), or a fluorescence plate reader at Ex/Em = 490/525 nm cutoff 515 nm.

**EXAMPLE DATA ANALYSIS AND FIGURES**


**Figure 1.** Simultaneous imaging of live and apoptotic HeLa cells labeled using calcein AM (Cat No. 22003) and Annexin V-iFluor® 647 conjugate (Cat No. 20074)

**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](mailto:info@aatbio.com) if you have any questions.