

## Calcein Deep Red™ acetate

Catalog number: 22010  
Unit size: 1 mg

Component	Storage	Amount
Calcein Deep Red™ acetate	Freeze (< -15 °C), Minimize light exposure	1 vial (1 mg)

### OVERVIEW

This product is now replaced by #22011. Calcein AM is one of the most popular fluorescent probes used for labeling and monitoring cellular functions of live cells. However, the single color of Calcein AM makes it impossible to use this valuable reagent in the multicolor applications. For example, it is impossible to use Calcein AM in combination of GFP-transfected cells due to the same color to GFP. To address this color limitation of Calcein AM, we have developed Calcein Orange™, Calcein Red™ and Calcein Deep Red™. These new Calcein AM analogs enable the multicolor labeling and functional analysis of live cells in combination with Calcein AM. Non-fluorescent Calcein Deep Red™ acetate can easily get into live cells and hydrolyzes to generate strongly fluorescent Calcein Deep Red™ (Cat#: 21902) dye. Calcein Deep Red™ dye can be monitored with the common Cy5 filter set.

### KEY PARAMETERS

#### Flow cytometer

Excitation	640 nm laser
Emission	660/20 nm filter
Instrument specification(s)	APC channel

#### Fluorescence microscope

Excitation	Cy5 filter set
Emission	Cy5 filter set
Recommended plate	Black wall/clear bottom

#### Fluorescence microplate reader

Excitation	620
Emission	660
Cutoff	630
Recommended plate	Solid black

### 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 Deep Red™ Acetate Stock Solution

Prepare a 2 to 5 mM stock solution of Calcein Deep Red™ acetate in high-quality, anhydrous DMSO.

**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.

### PREPARATION OF WORKING SOLUTION

#### Calcein Deep Red™ Acetate Working Solution

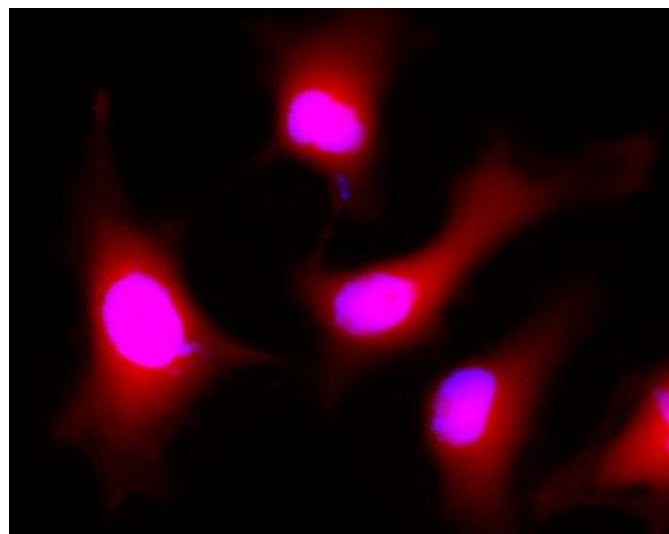
Prepare a Calcein Deep Red™ acetate working solution of 1 to 10 µM in the buffer of your choice (e.g., Hanks and Hepes buffer). For most cell lines, Calcein Deep Red™ acetate at the final concentration of 4 to 5 µM is recommended. The exact concentration of indicators required for cell loading must be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

**Note** 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.

### 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 Deep Red™ acetate 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 Cy5 filter set, a flow cytometer equipped with a 660/20 nm filter (APC channel), or a fluorescence plate reader at Ex/Em = 620/660 nm cutoff 630 nm.

### EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.** Images of Live HeLa cells stained with Calcein Deep Red™ acetate (Cat.22010). Cell nuclei were stained with Hoechst 33342 (Blue, Cat#17535).

### 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.