

## FITC-Palmitate Conjugate

Catalog number: 22192  
Unit size: 5 mg

Component	Storage	Amount (Cat No. 22192)
FITC-Palmitate Conjugate	Freeze (< -15 °C), Minimize light exposure	5 mg

### OVERVIEW

FITC-Palmitate Conjugate is a highly hydrophobic fatty acid lipid that is tagged with fluorescein. Palmitic acid is a saturated C16 fatty acid that can interact well with the hydrophobic cell membranes or lysosome membranes. Fluorescein-modified palmitic acid has strong green fluorescence that can well be imaged with the standard FITC filter set. AAT Bioquest also offer other dye-labeled lipids, such as phospholipids, sphingolipids, sterols, glycerolipids, and others.

### KEY PARAMETERS

#### Fluorescence microscope

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

### 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.

#### FITC-Palmitate Conjugate stock solution

Prepare a 5 to 10 mM stock solution by dissolving an appropriate amount of FITC-Palmitate Conjugate in DMSO.

**Note:** For example, to prepare a 5 mM DMSO stock solution dissolve 1 mg of the FITC-Palmitate conjugate in 235 µL of DMSO.

### PREPARATION OF WORKING SOLUTION

#### FITC-Palmitate Conjugate working solution

Prepare a 5 to 10 µM FITC-Palmitate conjugate working solution by diluting the stock solution in HHBS.

**Note:** Prepare a fresh working solution before use.

### SAMPLE EXPERIMENTAL PROTOCOL

The following protocol can be used as a guideline. If needed, optimize the protocol to achieve the desired results.

1. Culture cells as desired.
2. Remove cell culture medium.

**Note:** Cell culture medium must be removed since it may interfere with the conjugate staining.

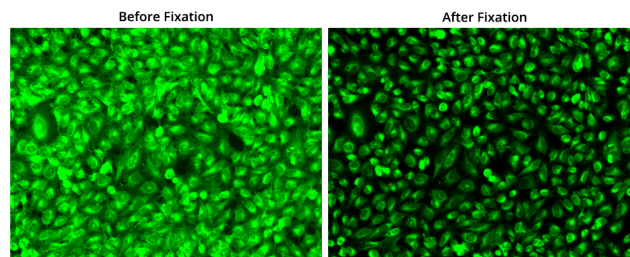
3. Add 100 µL of FITC-Palmitate Conjugate working solution per well.
4. Incubate the cells in the dark for 30 minutes at room temperature.
5. Remove the FITC-Palmitate Conjugate working solution and wash cells twice with HHBS.
6. Add HHBS buffer to the cells and observe the cells using a

fluorescence microscope with a FITC filter set.

7. **Optional:** The cells can be fixed after staining.

### EXAMPLE DATA ANALYSIS AND FIGURES

Placeholder for image details



**Figure 1.** Image of HeLa cells stained with FITC-Palmitate Conjugate and then fixed with 4% formaldehyde. Live HeLa cells were cultured in a 96-well plate overnight, and the medium was removed before staining. 100 µL of the FITC-Palmitate conjugate (10 µM) in HH buffer was added to each well. Cells were stained for 30 minutes at room temperature, washed twice with HH buffer, and then fixed with 4% formaldehyde. The cells were imaged with a FITC filter before and after fixation.

### 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.