

## Cell Navigator® Fluorimetric Lipid Droplet Assay Kit \*Blue Fluorescence\*

Catalog number: 22731  
Unit size: 200 Tests

| Component                    | Storage                                    | Amount (Cat No. 22731) |
|------------------------------|--|------------------------|
| Component A: Droplite™ Blue  | Freeze (< -15 °C), Minimize light exposure | 1 Vial                 |
| Component B: Staining Buffer | Freeze (< -15 °C), Minimize light exposure | 1 Bottle (20 mL)       |

### OVERVIEW

Droplite™ Blue is a lipophilic stain that can specifically stain lipid droplets/adiposomes present in the cells.

Lipid droplets, also known as lipid bodies or adiposomes, are essential cellular organelles responsible for the storage and hydrolysis of neutral lipids. These lipid-rich structures serve as key reservoirs for fatty acids and cholesterol and play a crucial role in biological processes such as energy production, cellular signaling, and membrane maintenance. Impaired cytoplasmic lipid droplet accumulation is associated with various diseases and pathological conditions including metabolic disorders, liver dysfunction, and cardiovascular diseases.

AAT Bioquest's Cell Navigator® Fluorimetric Lipid Droplet Assay Kit provides a powerful and reliable method for accurately studying lipid droplet accumulation within the cells. This kit features Droplite™ Blue, a lipophilic stain that exhibits bright blue fluorescence in a lipid-rich environment while giving minimal fluorescence in aqueous media. Droplite™ Blue is an ideal vital stain for visualizing intracellular lipid droplets using fluorescence microscopy, flow cytometry, or fluorescence microplate readers. The blue fluorescence can be easily detected using standard DAPI filter sets, offering a convenient and efficient tool for lipid droplet analysis.

### AT A GLANCE

#### Protocol Summary

1. Prepare cells with test compounds.
2. Add Droplite™ Blue working solution.
3. Incubate at room temperature or 37°C for 10 to 30 minutes.
4. Read fluorescence intensity with a fluorescence microscope using a DAPI filter set.

#### Important Note

This protocol is our recommended guideline for live cells, but it can be adjusted to meet your specific requirements. Since Droplite™ Blue exhibits minimal fluorescence in aqueous media, removing the growth medium and staining solution after staining is optional. Stained cells can be fixed with 3–4% formaldehyde. Alternatively, prefixed cells (fixed with 3–4% formaldehyde) can also be stained using the Droplite™ Blue staining solution.

### KEY PARAMETERS

#### Fluorescence microplate reader

|                             |                         |
|-----------------------------|-------------------------|
| Cutoff                      | 420 nm                  |
| Emission                    | 450 nm                  |
| Excitation                  | 350 nm                  |
| Recommended plate           | Black wall/clear bottom |
| Instrument specification(s) | Bottom read mode        |

#### Fluorescence microscope

|                   |                         |
|-------------------|-------------------------|
| Emission          | DAPI filter set         |
| Excitation        | DAPI filter set         |
| Recommended plate | Black wall/clear bottom |

### CELL PREPARATION

For guidelines on cell sample preparation, please visit:

<https://www.aatbio.com/resources/guides/cell-sample-preparation.html>

### PREPARATION OF WORKING SOLUTION

1. To prepare the Droplite™ Blue working solution, dilute 5 µL of the Droplite™ Blue (Component A) in 1 mL of Staining Buffer (Component B).

**Note:** 50 µL of Droplite™ Blue (Component A) is enough for one 96-well plate. Protect the solution from light. The optimal concentration of Droplite™ Blue may vary depending on the application. Adjust staining conditions based on the cell type and the permeability of cells or tissues to the probe.

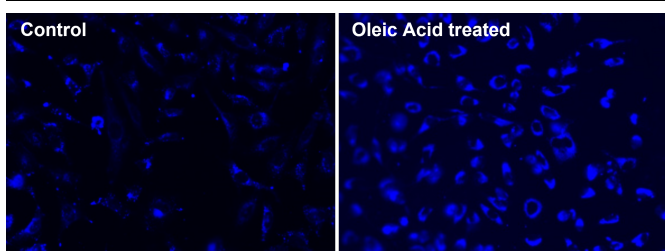
### SAMPLE EXPERIMENTAL PROTOCOL

#### For adherent cells:

1. Grow cells either in a 96-well black wall/clear bottom plate (100 µL/well/96-well) or on cover-slips inside a petri dish filled with the appropriate culture medium.
2. Gently aspirate the culture medium and add equal volume (such as 100 µL/well/96-well plate) of the Droplite™ Blue staining solution.
3. Incubate the cells in a 37°C, 5% CO<sub>2</sub> incubator for 10 - 30 minutes.
4. Remove Droplite™ Blue working solution (Optional).
5. Measure fluorescence at Ex/Em = 350/450 nm with a microplate reader or observe the cells using a fluorescence microscope equipped with a DAPI filter set.

**For suspension cells:**

1. Centrifuge the cells at 1000 rpm for 5 minutes to get  $1 - 5 \times 10^5$  cells per tube.
2. Resuspend cells in 500  $\mu$ L of Droplite™ Blue working solution.
3. Incubate at room temperature or 37°C for 10 to 30 min, protected from light.
4. Centrifuge to remove the Droplite™ Blue working solution, and resuspend cells in 500  $\mu$ L of pre-warmed medium or buffer of your choice to get  $1 - 5 \times 10^5$  cells per tube (Optional).
5. Monitor the fluorescence increase using a fluorescence microscope equipped with a DAPI filter set.

**EXAMPLE DATA ANALYSIS AND FIGURES**

**Figure 1.** Fluorescence images of intracellular lipid droplets in control (Left) and Oleic Acid treated HeLa cells (Right) using Cell Navigator® Lipid Droplets Fluorescence Assay Kit. HeLa cells were incubated with 300  $\mu$ M of Oleic Acid for 24 hours to induce intracellular lipid droplets formation. After washing with PBS, the cells were labeled with 1X Droplite™ Blue. Images were acquired with fluorescence microscope using DAPI filter set.

**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.