

# MycoLight™ Flow Cytometric Live Bacteria Assay Kit

Catalog number: 22407  
Unit size: 100 Tests

Component	Storage	Amount
Component A: MycoStain It™ 520	Freeze (<-15 °C), Minimize light exposure	2 vials
Component B: Assay Buffer	Freeze (<-15 °C)	1 bottle (50 mL)
Component C: DMSO	Freeze (<-15 °C)	1 vial (200 µL)

## OVERVIEW

The MycoLight Flow Cytometric Live Bacteria Assay Kit provides an easy and convenient method for evaluating bacterial vitality as a function of the intracellular esterase activity. MycoLight™ 520 is non-fluorescent esterase substrate that diffuses into both Gram positive and Gram-negative bacteria. Upon hydrolysis by bacterial intracellular non-specific esterase, a green fluorescent product is produced and accumulated within bacteria. Compared to the commonly used esterase substrate CFDA and CFDA-AM, this kit provides brighter and more stable signal with lower background and easier staining protocol.

## AT A GLANCE

### Protocol summary

1. Prepare 100X dye stock solution.
2. Prepare bacteria samples.
3. Incubate bacteria samples with MycoLight™ 520 at 37°C for 5-10 minutes or room temperature for 60 minutes in dark.
4. Analyze sample by flow cytometry with FITC channel.

### Important

Thaw one of each kit component at room temperature before starting the experiment.

## KEY PARAMETERS

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

## 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.

1. MycoLight™ 520 stock solution (100X):  
Add 100 µL of DMSO (Component C) into one vial of MycoLight™ 520 (Component A) to make 100X stock solution.

## SAMPLE EXPERIMENTAL PROTOCOL

1. Prepare bacteria sample with concentration in range of 10<sup>6</sup> to 10<sup>8</sup> cells/ml. Grow bacteria into late log phase in appropriate medium. Remove medium by centrifugation at 10,000 x g for 10 minutes and re-suspend the pellet in Assay Buffer (Component B).

**Note** Measure the optical density of the bacterial culture at wavelength = 600 nm (OD<sub>600</sub>) to determine the cell number. For *E. coli* culture, OD<sub>600</sub> = 1.0 equals 8 x 10<sup>8</sup> cells/ml.

2. Treat cells with test compounds as desired. Remove treatments by centrifugation at 10,000 x g for 10 minutes and re-suspend the pellet in

appropriate amount of Assay Buffer (Component B) so the concentration of bacteria in the treated sample is the same as the live.

**Note** Determine the concentration of the bacterial culture before starting the treatment.

**Note** Dead bacteria can serve as negative control, it is recommended to kill bacteria with 70% ethanol for 30 min followed by 60 min of boiling.

3. Example of Live/Dead bacterial mixture preparation: Mix seven different proportions of the bacterial suspensions as in Table 1 for a total volume of 200 µL for each sample.
4. Add 2 µL of the 100X MycoLight™ 520 stock solution to 200 µL of the bacterial sample in Assay Buffer.
5. Mix well and incubate in dark for 5-10 min at 37°C or 60 min at RT for optimum staining results.
6. Add 300 µL of Assay Buffer (Component C) to increase volume before analyzing the cells with a flow cytometer.
7. Monitor fluorescence of bacteria with a flow cytometer through FITC channel (488/530 nm).

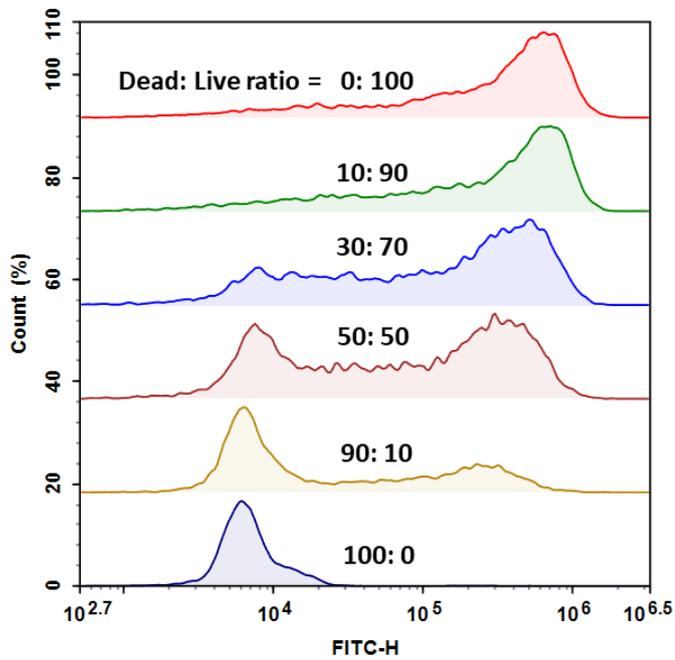
**Note** To exclude debris, it is recommended to set the threshold of the flow cytometer as the following: FSC >10,000, SSC >5,000.

**Note** The efficiency of MycoLight™ 520 is highly strain dependent and the staining conditions should be optimized accordingly.

**Table 1.** Example Volumes of Live and Dead samples to mix to achieve various Live/Dead ratios.

Ratio of Dead:Live Cells	µL of Dead Sample	µL of Live Sample
0:100	0	200
10:90	20	180
30:70	60	140
50:50	100	100
90:10	180	20
100:0	200	0

## EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.** Relative viability of *E.coli* suspension was analyzed using ACEA NovoCyte flow cytometer in FITC channel. The readings (Count(%)) were obtained from various Live/Dead *E.coli* mixtures (Table 1). The live and dead population in each mixture can be easily distinguished by the two distinct peaks.

**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.