

## Mycolight™ Live Bacteria Fluorescence Imaging Kit

Catalog number: 22409

Unit size: 100 Tests

Component	Storage	Amount
Component A: MycoLight™ 520	Freeze (<-15 °C), Dessicated, Avoid Light	1 vial
Component B: 10X Signal Enhancer	Freeze (<-15 °C), Avoid Light	1 vial (1 mL)
Component C: Assay Buffer	Freeze (<-15 °C), Avoid Light	1 bottle (10 mL)
Component D: DMSO	Freeze (<-15 °C), Avoid Light	1 vial (100 µL)

### OVERVIEW

The MycoLight™ Live Bacteria Fluorescence Imaging Kit provides an easy and convenient way for visualizing live bacteria through fluorescent microscope. MycoLight™ 520 is non-fluorescent esterase substrate that diffuse 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. Compare to the commonly used esterase substrate CFDA and CFDA-AM, the 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. Add MycoLight™ 520 and Signal Enhancer.
4. Incubate bacteria samples with MycoLight™ 520 and Signal Enhancer at 37°C for 5-10 minutes or room temperature for 60 minutes in dark.
5. Analyze sample by fluorescence microscope with FITC filter sets.

#### Important

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

### KEY PARAMETERS

Instrument:	Fluorescence microscope
Excitation:	488 nm
Emission:	530 nm
Recommended plate:	Black wall/clear bottom
Instrument specification(s):	FITC filter

### 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 mL of DMSO (Component D) into the 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.

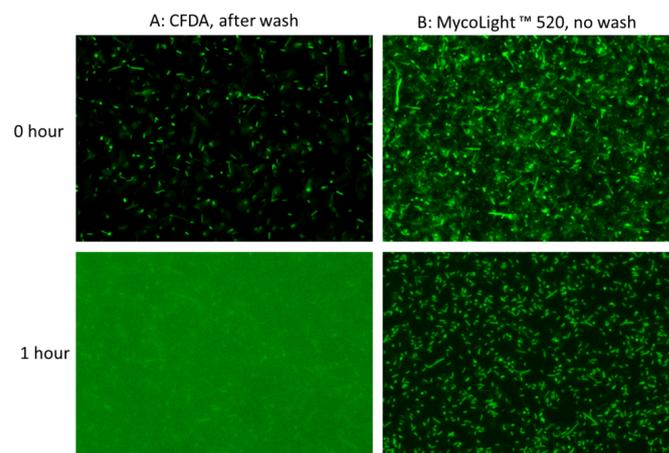
**Note** Determine the concentration of the bacterial culture before starting the treatment.

**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. Add 1 µL of the 100X MycoLight™ 520 stock solution and 10 µL of 10X Signal Enhancer (Component B) to 90 µL of the bacterial sample in Assay Buffer.
4. Mix well and incubate in dark for 5-10 min at 37°C or 60 min at RT for optimum staining results.
5. Monitor fluorescence of bacteria with a fluorescent microscope through FITC (Ex/Em = 488/530 nm) channel.

**Note** Same protocol can also be used for microplate reader assays.

### EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.** Fluorescence images of *E. coli* stained with CFDA or MycoLight™ Live Bacteria Fluorescence Imaging Kit. CFDA requires washing steps before imaging to minimize background, while no washing is needed using this kit Cat#22409. The staining efficiency of Mycolight™ 520 is much higher than CFDA as more bacteria show green fluorescence. The signal of Mycolight™ 520 remains in cells after 1 hour of staining while CFDA leaks out readily. Same amount of bacteria were presented in each sample and fluorescence images were taken under the same exposure time.

**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.