

## Amplite™ Colorimetric Beta-Lactamase Activity Assay Kit

Ordering Information	Storage Conditions	Instrument Platform
Cat#: 12551 (200 Assays)	Keep in freezer and protect from light	Absorbance microplate readers

### Introduction

$\beta$ -Lactamases are a large family of enzymes capable of hydrolyzing  $\beta$ -lactams.  $\beta$ -Lactam ring is the common element in all beta-lactam antibiotics including penicillin derivatives, cephalosporins, monobactams, and carbapenems. Through hydrolysis,  $\beta$ -lactamase breaks the  $\beta$ -lactam ring open, thus deactivates the molecule's antibacterial properties. Bacteria from clinical and non-clinical settings are becoming increasingly resistant to  $\beta$ -lactam antibiotics by synthesizing  $\beta$ -lactamase. To overcome this resistance,  $\beta$ -lactam antibiotics are often given with  $\beta$ -lactamase inhibitors such as clavulanic acid. Therefore, detection of  $\beta$ -lactamase activity is of central importance to assess beta-lactam antibiotics as well as to prevent antibiotics resistance. AAT Bioquest's Colorimetric Beta-Lactamase Activity Assay Kit offers a sensitive colorimetric assay for measuring  $\beta$ -lactamase activity in biological samples. The  $\beta$ -lactamase activity is detected using Nitrocefin, which changes color from yellow to red upon hydrolysis by  $\beta$ -lactamase. The assay can be performed using an absorbance microplate reader by measuring the OD ratio at the wavelength of 490 nm to 380 nm.

### Kit Components

Components	Amount
Component A: Nitrocefin	1 vial (100 $\mu$ L)
Component B: Assay Buffer	1 bottle (10 mL)
Component C: $\beta$ -Lactamase Standard	1 vial (lyophilized powder)

### Assay Protocol for One 96-Well Plate

#### Brief Summary

**Prepare  $\beta$ -lactamase assay mixture (50  $\mu$ L) → Add  $\beta$ -lactamase standards or test samples (50  $\mu$ L) → Incubate at RT for 30-60 min → Monitor absorbance increase at OD ratio of 490/380 nm**

*Note: Thaw one vial of each kit component at room temperature before starting the experiment.*

#### 1. Prepare $\beta$ -lactamase standard stock solution:

Add 100  $\mu$ L of ddH<sub>2</sub>O into the vial of  $\beta$ -lactamase standard (**Component C**) to make 50 mU/mL  $\beta$ -lactamase standard stock solution.

*Note: The unused  $\beta$ -lactamase standard stock solution should be divided into single use aliquots and stored at -20 °C.*

#### 2. Prepare serial dilutions of $\beta$ -lactamase standard:

2.1 Add 10  $\mu$ L of  $\beta$ -lactamase standard stock solution (50 mU/mL, from Step 1) into 990  $\mu$ L 1×PBS buffer to generate 500  $\mu$ U/mL  $\beta$ -lactamase standard solution.

*Note: Diluted  $\beta$ -lactamase standard solution is unstable, and should be used promptly.*

2.2 Take 200  $\mu$ L of 500  $\mu$ U/mL  $\beta$ -lactamase standard solution to perform 1:2 serial dilutions in PBS to get approximately 250, 125, 62.5, 31.3, 15.6, 7.8 and 0  $\mu$ U/mL serial dilutions of  $\beta$ -lactamase standard.

2.3 Add serial dilutions of  $\beta$ -lactamase standard and  $\beta$ -lactamase containing test samples into a 96-well clear bottom microplate as described in Tables 1 and 2.

**Table 1** Layout of  $\beta$ -lactamase standards and test samples in a 96-well clear bottom microplate

BL	BL	TS	TS	....	....														
Lac 1	Lac 1	....	....	....	....														
Lac 2	Lac 2																		
Lac 3	Lac 3																		
Lac 4	Lac 4																		
Lac 5	Lac 5																		
Lac 6	Lac 6																		
Lac 7	Lac 7																		

*Note: Lac =  $\beta$ -Lactamase Standards, BL=Blank Control, TS=Test Samples.*

**Table 2** Reagent composition for each well

<b>β-Lactamase Standard</b>	<b>Blank Control</b>	<b>Test Sample</b>
Serial Dilutions*: 50 μL	1×PBS Buffer : 50 μL	50 μL

\*Note: Add the serially diluted β-lactamase standards from approximately 500 to 7.8 μU/mL into wells from Lac 1 to Lac 7 in duplicate.

### 3. Prepare β-lactamase assay mixture:

Add 50 μL nitrocefin stock solution (**Component A**) into 5mL of **Component B**, and mix well to make β-lactamase assay mixture (**Component A+B**).

Note: This β-lactamase assay mixture is enough for one 96-well plate. The β-lactamase assay mixture is not stable, prepare fresh for each use.

### 4. Run β-lactamase assay:

4.1 Add 50 μL of β-lactamase assay mixture (from Step 3) to each well of β-lactamase standard, blank control, and test samples (see Step 2.3) to make the total volume of 100 μL/well.

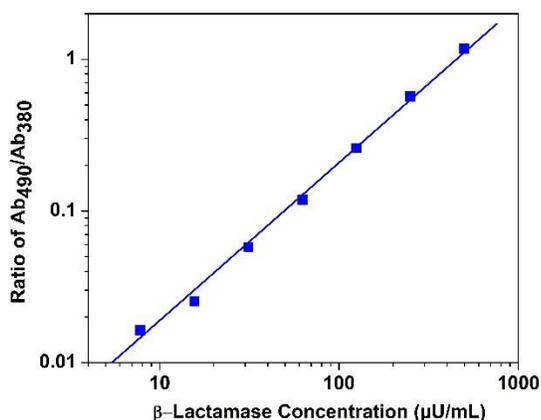
Note: For a 384-well plate, add 25 μL of sample and 25 μL of β-lactamase assay mixture into each well.

4.2 Incubate the reaction at room temperature for 30-60 minutes, protected from light.

4.3 Monitor the absorbance increase with an absorbance plate reader at OD ratio of 490/380 nm.

### Data Analysis

The absorbance reading in blank wells (with PBS and β-lactamase assay mixture only) is used as a control, and is subtracted from the values of those wells with the β-lactamase standards and test samples. A β-lactamase standard curve is shown in Figure 1. Calculate the β-lactamase concentrations of the samples according to the β-lactamase standard curve.



**Figure 1.** β-Lactamase dose response was measured with Amplite™ Colorimetric Beta-Lactamase Assay Kit (Cat#12551) on a 96-well clear bottom plate using a SpectraMax reader (Molecular Devices). As low as 15 μU/mL β-lactamase can be detected with 30 minutes incubation.

### References

1. Wilke MS, Lovering AL, Strynadka NC. (2005) β-Lactam antibiotic resistance: a current structural perspective. *Current Opinion in Microbiology* 8(5): 525-533.
2. Drawz SM, Bonomo RA. (2010). Three Decades of β-Lactamase Inhibitors. *Clinical Microbiology Reviews* 23 (1): 160–201.
3. Fisher JF, Meroueh SO, Mobashery S. (2005). Bacterial Resistance to β-Lactam Antibiotics: Compelling Opportunism, Compelling Opportunity. *Chemical Reviews* 105 (2): 395–424.