

## Amplite™ Fluorimetric Alkaline Phosphatase Assay Kit

*\*Near Infrared Fluorescence\**

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 11954 (500 assays)	Keep in freezer, Avoid exposure to light	Fluorescence microplate readers

### Introduction

Alkaline phosphatase is widely used in various biological assays (in particular, immunoassays) and ELISA-based diagnostics. AAT Bioquest's Amplite™ Fluorimetric Alkaline Phosphatase Assay kit uses our SunRed™-based substrate. The weakly fluorescent SunRed™ phosphate is sensitive to phosphatase-induced hydrolysis, giving the SunRed™ fluorophore that possesses intense red fluorescence. Upon phosphatase-induced hydrolysis, the SunRed™ phosphate solution has its absorption blue-shifted more than 100 nm. The maximum absorption of SunRed™ fluorophore at 633 nm makes this substrate an ideal NIR probe that can be readily detected with many fluorescence instrument systems often equipped with Cy5 settings.

Based on the near infrared fluorescence of SunRed™ fluorophore, the signal can be easily read by a fluorescence microplate reader at Ex/Em = ~620/660 nm. The kit has been used for the high throughput screening of protein phosphatase inhibitors due to its low interference from biological sample. It can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation without a separation step.

### Kit Components

Components	Amount
Component A: SunRed™ Substrate (light sensitive)	1 vial
Component B: Assay Buffer	1 bottle (25 mL)
Component C: Alkaline Phosphatase Standard	1 vial (lyophilized powder, 10 units)

### Assay Protocol for One 96-well Plate

#### Brief Summary

**Prepare assay reaction mixture (50 µL) → Add alkaline phosphatase standards and/or test samples (50 µL) → Incubate at RT or 37 °C for 30 to 120 minutes → Monitor fluorescence intensity at Ex/Em = 620/660 nm**

*Note: Thaw all the kit components at room temperature before starting the experiment.*

#### 1. Prepare 250X SunRed™ Substrate stock solution:

Add 100 µL of double sterile H<sub>2</sub>O into the vial of SunRed™ Substrate (Component A). The stock solution should be used promptly. Any remaining solution should be made in single used aliquots and frozen at -20 °C.

*Note: Avoid repeated freeze and thaw cycles.*

#### 2. Prepare assay reaction mixture:

Prepare assay reaction mixture according to the following table and keep from light.

**Table 1.** Assay reaction mixture for one 96-well plate (2X)

Components	Volume
250X SunRed™ Substrate stock solution (from Step 1)	20 µL
Assay Buffer (Component B)	5 mL
Total volume	5 mL

*Note: Prepare fresh reaction mixture for each experiment.*

#### 3. Prepare serially diluted alkaline phosphatase standards (0 to 100 mU/mL):

3.1 Add 100 µL of distilled H<sub>2</sub>O with 0.1% BSA (H<sub>2</sub>O-0.1% BSA) to alkaline phosphatase standard (Component C, 10 units) to generate a 100 units/mL alkaline phosphatase standard solution.

*Note: The alkaline phosphatase standard solution is not stable. Unused standard solution should be aliquoted and stored at -20 °C. Avoid repeated freeze and thaw cycles.*

3.2 Add 10 µL of 100 units/mL alkaline phosphatase standard solution (from Step 3.1) to 990 µL of H<sub>2</sub>O-0.1% BSA to generate a 1,000 mU/mL alkaline phosphatase standard solution.

3.3 Take 100 µL of 1,000 mU/mL alkaline phosphatase standard solution (from Step 3.2) to perform 1:3 serial dilutions to get 300, 100, 30, 10, 3, 1, 0.3, and 0 mU/mL serially diluted alkaline phosphatase standards.

3.4 Add serially diluted alkaline phosphatase standards and/or alkaline phosphatase containing test samples into a solid black 96-well microplate as described in Tables 2 and 3.

*Note: Prepare cells or tissue samples as desired. Discard unused serially diluted alkaline phosphatase standards.*

