

Amplite™ Fluorimetric Alkaline Phosphatase Assay Kit

Green Fluorescence

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 11953 (500 assays)	Keep in freezer Avoid exposure to light	Fluorescence microplate readers

Introduction

Alkaline phosphatase (ALP) (EC 3.1.3.1) is a hydrolase enzyme responsible for removing phosphate groups from many types of molecules, including nucleotides, proteins, and alkaloids. An important use of alkaline phosphatase is as a label for enzyme immunoassays. Alkaline phosphatase is a highly sensitive enzyme for ELISA, immuno-histochemical as well as Northern, Southern and Western blot applications. It is widely used in various biological assays (in particular, immunoassays) and ELISA-based diagnostics.

Our Amplite™ Fluorimetric Alkaline Phosphatase Assay Kit uses our FDP, a fluorogenic phosphatase substrate, to quantify the alkaline phosphatase activity in solutions, in cell extracts, and on solid surfaces (such as PVDF membranes). It can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation without a separation step. Its signal can be easily read by a fluorescence microplate reader at Ex/Em = ~490/525 nm. The kit provides an optimized “mix and read” assay protocol which is compatible with HTS liquid handling instruments.

Kit Components

Components	Amount
Component A: FDP (light sensitive)	1 vial
Component B: Assay Buffer	1 bottle (25 mL)
Component C: Alkaline Phosphatase Standard	1 vial (lyophilized powder, 10 units)
Component D: DMSO	1 vial (500 µL)
Component E: Stop Solution	1 bottle (25 mL)

Assay Protocol for One 96-well Plate

Brief Summary

Prepare assay reaction mixture (50 µL) → Add alkaline phosphatase standards or test samples (50 µL) → Incubate at RT or 37 °C for 10-30 minutes → Monitor fluorescence intensity at Ex/Em = 490/ 525 nm

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

1. Prepare 250X FDP stock solution:

Add 100 µL of DMSO (Component D) into the vial of FDP (Component A). The FDP stock solution should be used promptly. Any remaining solution should be aliquoted and refrozen 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

Components	Volume
250X FDP (from Step 1)	20 µL
Assay Buffer (Component B)	5 mL
Total volume	5 mL

5.6 Monitor the fluorescence increase with a fluorescence plate reader at Ex/Em = 490 ± 10/525 ± 10 nm.

Data Analysis

The fluorescence in blank wells (with equal volume of assay reaction mixture and H₂O-0.1% BSA only) is used as a control, and is subtracted from the values for those wells with alkaline phosphatase reactions. An alkaline phosphatase standard curve is shown in Figure 1.

Note: The fluorescence background increases with time due to spontaneous hydrolysis, thus it is important to subtract the fluorescence intensity value of the blank wells for each data point.

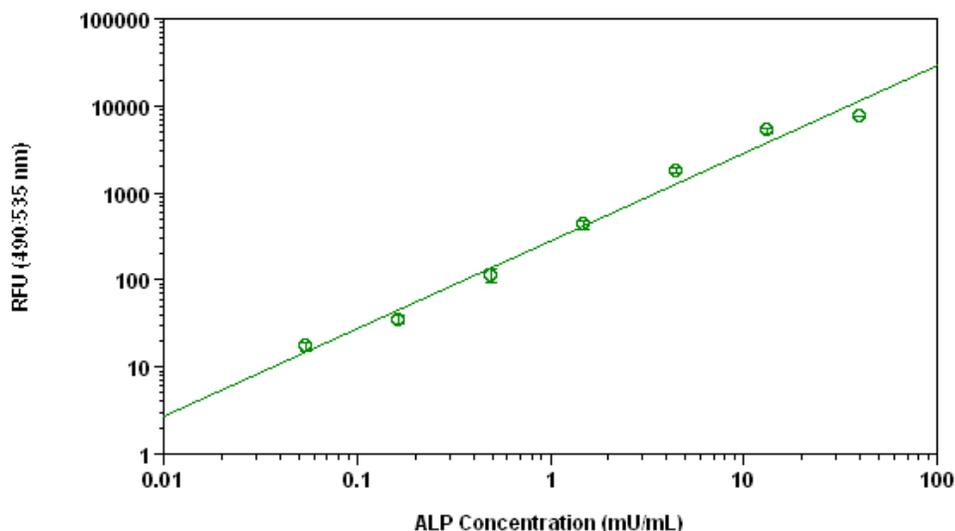


Figure 1. Alkaline phosphatase dose response was measured with the Amplite™ Fluorimetric Alkaline Phosphatase Assay Kit in a solid black 96-well plate using a Gemini microplate reader (Molecular Devices). As low as 0.01 mU/well of alkaline phosphatase can be detected with 30 minutes incubation (n=3).

References

1. Zhu X, Jiang C. (2006) 8-Quinoyl phosphate as a substrate for the fluorimetric determination of alkaline phosphatase. Clin Chim Acta.
2. Ali AT, Penny CB, Paiker JE, Psaras G, Ikram F, Crowther NJ. (2006) The effect of alkaline phosphatase inhibitors on intracellular lipid accumulation in preadipocytes isolated from human mammary tissue. Ann Clin Biochem, 43, 207.
3. Lee DH, Lim BS, Lee YK, Yang HC. (2006) Effects of hydrogen peroxide (H₂O₂) on alkaline phosphatase activity and matrix mineralization of odontoblast and osteoblast cell lines. Cell Biol Toxicol, 22, 39.
4. Ali AT, Penny CB, Paiker JE, van Niekerk C, Smit A, Ferris WF, Crowther NJ. (2005) Alkaline phosphatase is involved in the control of adipogenesis in the murine preadipocyte cell line, 3T3-L1. Clin Chim Acta, 354, 101.
5. Rieu JP, Ronzon F, Place C, Dekkiche F, Cross B, Roux B. (2004) Insertion of GPIanchored alkaline phosphatase into supported membranes: a combined AFM and fluorescence microscopy study. Acta Biochim Pol, 51, 189.
6. Palermo C, Manduca P, Gazzero E, Foppiani L, Segat D, Barreca A. (2004) Potentiating role of IGFBP-2 on IGF-II-stimulated alkaline phosphatase activity in differentiating osteoblasts. Am J Physiol Endocrinol Metab, 286, E648.

Warning: This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at info@aatbio.com if you have any questions.