

Amplite™ Colorimetric Alanine Aminotransferase Assay Kit

Blue Color

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 13803 (200 assays)	Keep in frozen and avoid light	Absorbance microplate readers

Introduction

Alanine aminotransferase (ALT), also called serum glutamate pyruvic transaminase (GPT), is a member of transferase family. It catalyzes the reversible transfer of an α -amino group between alanine and glutamate, and is an important enzyme in amino acid metabolism. ALT is found mainly in liver and small amount in heart, muscle, and kidneys. In healthy subjects, serum ALT levels are low. However, when cells are damaged, such as acute and chronic hepatitis, obstructive jaundice, carcinoma of liver, myocardial infarction, ALT may leak into the blood stream and the ALT levels are significantly elevated. Therefore, determination of serum ALT level has great clinical and diagnostic significance.

Amplite™ Colorimetric Alanine Aminotransferase Assay Kit provides a quick and sensitive method for the measurement of ALT in various biological samples. ALT catalyzes the reaction of alanine and α -ketoglutarate to pyruvate and glutamate:



The product glutamate is measured by the generation of a blue color product through an enzyme coupled reaction cycle. The signal can be read by an absorbance microplate reader at the absorbance ratio of $A_{570\text{nm}}$ to $A_{610\text{nm}}$. With the Amplite™ Colorimetric Alanine Aminotransferase Assay Kit, as little as 10 mU/mL ALT was detected in a 100 μ L reaction volume. The assay is robust, and can be readily adapted for a wide variety of applications.

Kit Components

Components	Amount
Component A: ALT Enzyme Mixture	1 bottle (lyophilized powder)
Component B: ALT Assay Buffer	1 bottle (10 mL)
Component C: NAD	1 vial
Component D: ALT Positive Control	1 vial (10 U)

Assay Protocol for One 96-Well Plate

Brief Summary

Prepare ALT reaction mixture (50 μ L) → Add ALT standards or test samples (50 μ L) → Incubate at 37°C for 60 min to 120 min → Monitor absorbance increase at the absorbance ratio of $A_{570\text{nm}}/A_{610\text{nm}}$

Note: Thaw one bottle Component A and B at room temperature before starting the experiment.

1. Prepare serial dilutions of ALT standard (1 to 1000 mU/mL):

- 1.1 Add 100 μ L DPBS into the vial of ALT Positive Control (Component D) to make 100U/mL ALT stock solution.

Note: The unused ALT stock solution should be divided into single use aliquots and stored at -20°C.

- 1.2 Add 10 μ L of 100 U/mL ALT standard solution (from Step 1.1) into 990 μ L DPBS buffer with 0.1% BSA to generate 1 U/mL ALT standard solution.

Data Analysis

The absorbance in blank wells (with the DPBS buffer with 0.1% BSA only) is used as a control, and is subtracted from the values for those wells with the ALT reactions. An ALT standard curve is shown in Figure 1. *Note: The fluorescence background increases with time, thus it is important to subtract the intensity value of the blank wells from that of each data point.*

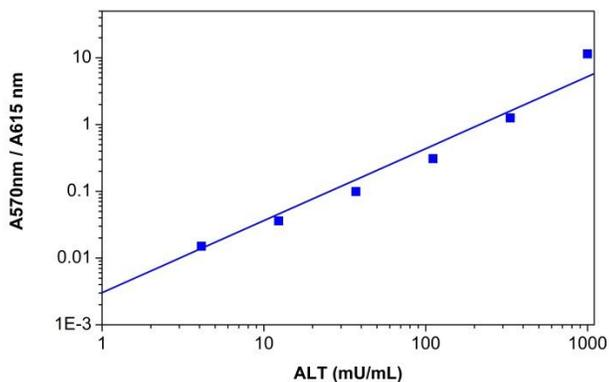


Figure 1. ALT dose response was measured with Amplite™ Colorimetric Alanine Aminotransferase Assay Kit in a 96-well black/clear bottom plate using a SpectraMax microplate reader (Molecular Devices). As low as 15.6 mU/mL ALT can be detected with 90 min incubation (n=3) at 37°C.

References

1. Hayashi H, Mizuguchi H, Miyahara I, Nakajima Y, Hirotsu K, Kagamiyama H (2003). "Conformational change in aspartate aminotransferase on substrate binding induces strain in the catalytic group and enhances catalysis". *J Biol Chem* 278 (11): 9481–9488.
2. Gaze DC (2007). "The role of existing and novel cardiac biomarkers for cardioprotection". *Curr. Opin. Invest. Drugs* 8 (9): 711–7.
3. Berg, JM; Tymoczko, JL; Stryer, L (2006). *Biochemistry*. W.H. Freeman. pp. 656–660.
4. Nalpas B, Vassault A, Charpin S, Lacour B, Berthelot P (1986). "Serum mitochondrial aspartate aminotransferase as a marker of chronic alcoholism: diagnostic value and interpretation in a liver unit". *Hepatology* 6 (4): 608–614.
5. Almo SC, Smith DL, Danishefsky AT, Ringe D (March 1994). "The structural basis for the altered substrate specificity of the R292D active site mutant of aspartate aminotransferase from *E. coli*". *Protein Eng.* 7 (3): 405–412.

Warning: This kit is only sold to our authorized distributors or end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of the 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.