

# Alanine Transaminase (ALT) Activity Assay Kit (Colorimetric)

Catalog #: MA-ALT

## Introduction

Alanine transaminase (ALT), also known as alanine aminotransferase (ALAT) or serum glutamic pyruvic transaminase (sGPT), is a homodimeric enzyme dependent on cytoplasmic pyridoxal phosphate. It plays a vital role in cellular nitrogen metabolism, amino acid metabolism, and liver gluconeogenesis. ALT catalyzes the reversible transamination between alanine and  $\alpha$ -ketoglutarate to form pyruvate and glutamate. ALT is found mainly in the liver and, to a lesser extent, in kidney, heart, muscle, and pancreas tissues. Normal serum levels of ALT are low, and an elevation in serum ALT activity is a widely used marker for liver damage.

The RayBio® Alanine Transaminase Activity Assay Kit offers a user-friendly, reliable, and highly sensitive method for quantifying ALT activity in plasma, serum, cell lysates, and other biological fluid samples. In this assay, ALT catalyzes the transfer of the amino group from L-alanine to  $\alpha$ -ketoglutarate resulting in the formation of pyruvate and L-glutamate. Simultaneously, lactate dehydrogenase catalyzes the reduction of pyruvate and the oxidation of NADH to NAD. The resultant rate of absorbance decrease is directly proportional to ALT activity.

## Storage

The entire kit may be stored at 2–8 °C for up to 6 months from the date of shipment. For prepared reagent storage, see table below.

Component	Size / Description	Storage After Preparation
Microplate	A 96-well (12 strips x 8 wells) plate	RT*
Sample Buffer	10 ml	2–8 °C
Assay Buffer	10 ml	2–8 °C
Enzyme Mix Solution	2 ml	2–8 °C
Alanine Transaminase (ALT) Positive Control	1 Vial (10 $\mu$ l)	2–8 °C

RT = room temperature

\*Return unused wells to the pouch containing desiccant pack, reseal along entire edge.

## Additional Materials Required

1. Microplate reader capable of measuring absorbance at 340 nm at 37°C
2. Precision pipettes to deliver 2  $\mu$ l to 1 ml volumes
3. Multi-channel pipettes to deliver 20  $\mu$ l to 200  $\mu$ l volumes
4. Tubes to prepare sample dilutions
5. Incubator at 37°C
6. 15 ml conical tubes

## Sample Tips and General Considerations

NOTE: Optimal methods of sample preparation will need to be determined by each researcher empirically based on researched literature and knowledge of the samples.

- If not using fresh samples, freeze samples as soon as possible after collection.
- Avoid multiple freeze-thaw cycles. If possible, sub-aliquot samples prior to initial storage.
- It is strongly recommended to add a protease inhibitor cocktail to cell and tissue lysate samples.
- Avoid sonication of 1 ml or less as this can quickly heat and denature proteins.
- Most samples will not need to be concentrated. If concentration is required, a spin column concentrator with a chilled centrifuge is recommended.

### 1. Cell lysates can be prepared as follows:

For attached cells, remove supernatant from cell culture, wash cells twice with cold 1X PBS (for suspension cells, pellet the cells by spinning down the cells at 1,000 x g for 10 min) making sure to remove any remaining PBS before adding lysis buffer. Solubilize the cells at  $2 \times 10^7$  cells/ml in lysis buffer containing protease inhibitors. Pipette up and down to resuspend cells and rock the lysates gently at 2–8 °C for 30 minutes. Transfer extracts to microfuge tubes and centrifuge at 14,000 x g for 10 minutes.

It is recommended that sample protein concentrations should be determined using a total protein assay. Lysates should be used immediately or aliquot and stored at -70 °C. Thawed lysates should be kept on ice prior to use.

General tips for preparing lysate samples can be viewed on the online Resources page of the website:

<https://www.raybiotech.com/tips-on-sample-preparation/>

### 2. Plasma samples:

Collect blood with an anticoagulant such as citrate, EDTA or oxalate and mix by inversion. Centrifuge the blood at 1000 x g at 4°C for 10 minutes. Collect plasma supernatant without disturbing the white buffy layer. Sample should be tested immediately or frozen at -80°C for storage. Typically, normal human plasma has ALT activity in the range of 8-40 U/L. No dilution/Neat is recommended for a typical plasma sample.

### 3. Serum samples:

Collect blood in a tube with no anticoagulant. Allow the blood to clot at room temperature for 30 minutes. Centrifuge at 2500 x g for 20 minutes. Remove the yellow serum supernatant without disturbing the white buffy layer. Samples should be tested immediately or frozen at -80°C for storage. Typically, normal human serum has ALT activity in the range of 8-40 U/L. No dilution/Neat is recommended for a typical serum sample.

### NOTE:

If the calculated ALT activity of the sample is higher than 500 U/L, dilute the sample in Sample Buffer and repeat the assay.

# Reagent Preparation

## A. Working Solution

Mix Assay Buffer and Enzyme Mix Solution at a volumetric ratio of 5:1 to make the Working Solution. For example, mix 9.5ml of Assay Buffer and 1.9ml of Enzyme Mix Solution to prepare sufficient Working Solution for one 96-well plate. Mix well. Protect from light. The Working Solution is stable for 14 days at 4 °C.

## B. Positive Control

Alanine Transaminase (ALT) Positive Control: pipette 1  $\mu$ L of ALT Positive Control into 99  $\mu$ L Sample Buffer to create a positive control stock, mix well. Then pipette 1  $\mu$ L of the positive control stock into 99  $\mu$ L Sample Buffer to prepare the positive control sample. Mix well.

# Assay Procedure

Positive control, samples, and Sample Buffer (used as a blank) should be assayed in duplicate or triplicate. A freshly prepared positive control should be used each time the assay is performed.

1. Set up a microplate reader or a microplate incubator at 37°C.
2. Prepare Working Solution (See Reagent Preparation, section A), and incubate it at 37°C for at least five minutes.
3. Prepare Positive Control (See Reagent Preparation, section B).
4. Pipette 10  $\mu$ L of sample, Sample Buffer (as the blank) and Positive Control into each well of the 96-well plate.
5. Transfer 100  $\mu$ L of pre-warmed Working Solution into each well (it is recommended to use a multi-channel pipette), mix and incubate at 37°C for one minute.
6. After one minute, read and record absorbance at 340nm (A1 Reading). Repeat readings every minute for the next four minutes at 37°C. (A2, A3, A4, A5 Reading).

## Calculation of Results

Calculate the mean absorbance for each set of duplicate/triplicate samples, Positive Control, and Sample Buffer blank.

Calculate the average absorbance difference per minute  $\Delta OD_{340nm}/min$  for each sample, the Positive Control, and Sample Buffer blank.

$$\Delta OD_{340nm}/min = \frac{(A1-A2) + (A2-A3) + (A3-A4) + (A4-A5)}{4}$$

Subtract the absorbance difference per minute of the Blank wells from absorbance difference per minute of the sample and positive control, this is the corrected absorbance difference per minute.

$$\Delta OD'_{340nm} /min = \Delta OD_{340nm}/min - \Delta OD_{340nm}^{Blank} /min$$

One international Unit (U/L) is defined as the amount of enzyme that catalyzes the transformation of one micromole of substrate per minute.

$$ALT \text{ activity (U/L)} = \frac{\Delta OD'_{340nm} \times 110 \mu l \times 1000}{1.836 \text{ mM}^{-1} \times 10 \mu l} \times \text{sample dilution}$$

Where:  $\Delta OD'_{340nm} /min$  = Corrected absorbance change per minute

1000 = Conversion of U/ml to U/L

$1.836 \text{ mM}^{-1}$  is the adjusted extinction coefficient for NADH at 340nm with the current path length of the solution in a 96-well plate.

$$ALT \text{ activity International System of Units (SI Units) (nkat/L)} = ALT \text{ activity (U/L)} * 16.67$$

*Note:* If the calculated ALT activity of the sample is higher than 500 U/L, dilute the sample in Sample Buffer and repeat the assay.

## A. Typical Data

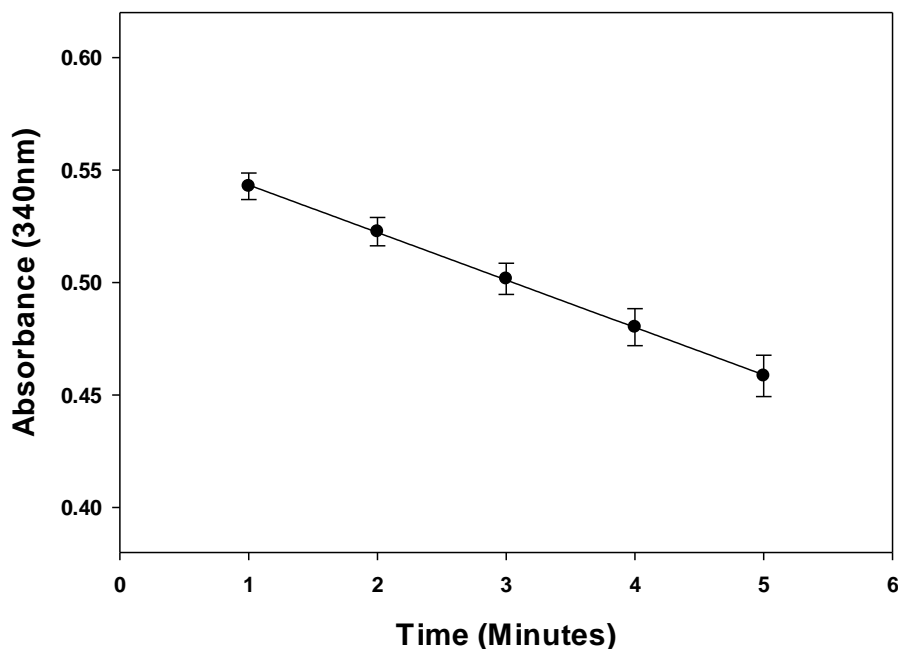


Figure 1. Example of OD value of Alanine Transaminase (ALT) Positive Control verse Time. This example is for demonstration purposes only.

	$\Delta OD_{340nm}/min$	ALT activity (U/L)
<b>Human Serum</b>	0.002	11.9841
<b>Human Plasma (EDTA)</b>	0.002	11.9841
<b>Human Plasma (Citrate)</b>	0.00175	10.4860
<b>Human Plasma (Heparin)</b>	0.0015	8.9880
<b>HepG2 cell lysate (protein concentration 4.28mg/ml)</b>	0.03083	184.7545
<b>ALT Positive Control</b>	0.02858	171.2729

Table 1. Typical data of different sample types.

## **B. Linear Range**

Up to 500 U/L ALT activity

## **C. Reproducibility**

Intra-assay Precision (Precision within an assay):

To assess intra-assay precision, 16 wells per sample (total of 4 samples) were tested on a single plate. The intra-assay coefficient of variation was found to be 8.5%.

Inter-assay Precision (Precision between assays):

To assess inter-assay precision, 4 samples were tested in separate assays (n=4). The inter-assay coefficient of variation was found to be 9.3%.

This product is for research use only.