



**PRODUCT INFORMATION &  
MANUAL**

**Acid Phosphatase/ACP  
Activity Assay Kit  
(Colorimetric)  
*NBP3-25847***

For research use only.  
Not for diagnostic or therapeutic  
procedures.

[www.novusbio.com](http://www.novusbio.com) - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - [technical@novusbio.com](mailto:technical@novusbio.com)

Novus kits are guaranteed for 6 months from date of receipt

# Acid Phosphatase/ACP Activity Assay Kit (Colorimetric)

Catalog No: NBP3-25847

Method: Colorimetric method

Specification: 96T (Can detect 40 samples without duplication)

Measuring instrument: Microplate reader

Sensitivity: 0.2 U/L

Detection range: 0.2–50 U/L

Average intra-assay CV (%): 2.8

Average inter-assay CV (%): 4.5

- ▲ This kit is for research use only.
- ▲ Instructions should be followed strictly, changes of operation may result in unreliable results.
- ▲ Please kindly provide us the lot number (on the outside of the box) of the kit for more efficient service.

## General information

### ▲ Intended use

This kit can be used to measure acid phosphatase (ACP) activity in serum (plasma), tissue samples.

### ▲ Background

Acid phosphatase (ACP) is a kind of acidic hydrolytic enzyme with high content in lysosomes which catalyzes the hydrolysis of phosphate monoester to phosphoric acid under acidic conditions. ACP is ubiquitous in nature, from low organisms such as escherichia coli and yeast to higher animal and plant tissues, as well as body fluids, human liver, and prostate.

### ▲ Detection principle

Disodium p-nitrobenzene phosphate (p-NPP), a widely used phosphatase chromogenic substrate, can form p-nitrophenol under the action of acid phosphatase. Under alkaline conditions, p-nitrophenol is yellow and has a maximum absorption peak at 405 nm. The darker of the yellow product is, the higher of the ACP activity is. Therefore, the activity of ACP can be calculated by measuring the OD value at 405 nm.

### ▲ Kit components & storage

Item	Component	Specification	Storage
Reagent 1	Buffer Solution	20 mL × 1 vial	-20°C , 12 months
Reagent 2	Substrate	Powder × 3 vials	-20°C , 12 months, shading light
Reagent 3	Standard	Powder × 1 vial	-20°C , 12 months, shading light
Reagent 4	Chromogenic Agent	24 mL × 1 vial	-20°C , 12 months
	Microplate	96 wells	No requirement
	Plate Sealer	2 pieces	
<p>Note: The reagents must be stored strictly according to the preservation conditions in the above table. The reagents in different kits cannot be mixed with each other.</p>			

## ▲ Materials prepared by users

### Instruments

Micropipettor, Microplate reader (400-415 nm, optimum wavelength: 405 nm), 37°C Incubator

### Reagents

Double distilled water, PBS (0.01 M, pH 7.4)

### ▲ Safety data

Some of the reagents in the kit contain dangerous substances. It should be avoided to touch the skin and clothing. Wash immediately with plenty of water if touching it carelessly. All the samples and waste material should be treated according to the relevant rules of laboratory's biosafety.

### ▲ Precautions

Before the experiment, please read the instructions carefully, and wear gloves and work clothes.

## ▲ The key points of the assay

1. Reagent 2 working solution and standard should be stored with shading light.
2. Reagent 2 working solution should be used up within 1 day.

## Pre-assay preparation

### ▲ Reagent preparation

1. Bring all reagents to room temperature before use.

2. The preparation of reagent 2 working solution:

Dissolve reagent 2 with 1.6 mL of reagent 1. The prepared solution can be stored at  $-20^{\circ}\text{C}$  with shading light for 24 hours.

3. The preparation of 10 mmol/L standard stocking solution:

Dissolve reagent 3 with 5 mL of double distilled water and mix fully. The prepared solution can be aliquoted into smaller quantities and stored at  $-20^{\circ}\text{C}$  for 7 days with shading light.

4. The preparation of 0.5 mmol/L standard:

Mix the 10 mmol/L standard stocking solution and reagent 1 at the ratio of 1: 19. Prepare the fresh needed amount before use and store it with shading light for detection.

### ▲ Sample preparation

1. Serum (Plasma): Detect the sample directly.

2. Tissue sample: Weigh the tissue accurately. Add PBS (0.01 M, pH 7.4) in a weight (g): volume (mL) ratio of 1: 9, homogenize mechanically in ice water bath to break tissue fully. Then centrifuge at 10000 g for 10 min at  $4^{\circ}\text{C}$  and collect the supernatant for measurement. Meanwhile, determine the protein concentration of supernatant (E-BC-K318-M).

## ▲ Dilution of sample

It is recommended to take 2~3 samples with expected large difference to do pre-experiment before formal experiment and dilute the sample according to the result of the pre-experiment and the detection range (0.2–50 U/L ).

The recommended dilution factor for different samples is as follows (for reference only):

Sample type	Dilution factor
10% <i>Epipremnum aureum</i> tissue homogenate	5-10
Mouse plasma	5-10
Rat plasma	5-10
Human urine	1
Human plasma	5-10
10% Rat spleen tissue homogenate	20-30
10% Rat liver tissue homogenate	20-30
10% Rat kidney tissue homogenate	20-30

**Note:** The diluent is PBS (0.01 M, pH 7.4).

## Assay protocol

### ▲ Plate set up

	1	2	3	4	5	6	7	8	9	10	11	12
A	A	A	S1	S9	S17	S25	S33	S1'	S9'	S17'	S25'	S33'
B	B	B	S2	S10	S18	S26	S34	S2'	S10'	S18'	S26'	S34'
C	C	C	S3	S11	S19	S27	S35	S3'	S11'	S19'	S27'	S35'
D	D	D	S4	S12	S20	S28	S36	S4'	S12'	S20'	S28'	S36'
E	E	E	S5	S13	S21	S29	S37	S5'	S13'	S21'	S29'	S37'
F	F	F	S6	S14	S22	S30	S38	S6'	S14'	S22'	S30'	S38'
G	G	G	S7	S15	S23	S31	S39	S7'	S15'	S23'	S31'	S39'
H	H	H	S8	S16	S24	S32	S40	S8'	S16'	S24'	S32'	S40'

Note: A-H, standard wells; S1-S40, sample wells; S1'-S40', control wells.

## ▲ Detailed operation steps

### The preparation of standard curve

Dilute 0.5 mmol/L standard solution with reagent 1 to a serial concentration. The recommended dilution gradient is as follows: 0, 0.05, 0.1, 0.2, 0.25, 0.3, 0.4, 0.5 mmol/L. Reference is as follows:

Number	Standard concentrations (mmol/L)	0.5 mmol/L standard solution ( $\mu\text{L}$ )	Reagent 1 ( $\mu\text{L}$ )
A	0	0	200
B	0.05	20	180
C	0.1	40	160
D	0.2	80	120
E	0.25	100	100
F	0.3	120	80
G	0.4	160	40
H	0.5	200	0

### The measurement of samples

1. **Standard well:** Take 40  $\mu\text{L}$  of standards with different concentrations into the standard wells  
**Sample well:** Add 40  $\mu\text{L}$  of sample into the sample wells.  
**Control well:** Add 40  $\mu\text{L}$  of sample into the control wells.
2. Add 40  $\mu\text{L}$  of reagent 1 into the standard wells and control wells.  
 Add 40  $\mu\text{L}$  of reagent 2 working solution into the sample wells
3. Mix fully with microplate reader for 3 s and incubate at 37°C for 10 min
4. Add 160  $\mu\text{L}$  of reagent 4 into each well.
5. Mix fully with microplate reader for 3 s. Measure the OD values of each well at 405 nm with microplate reader.

### ▲ Summary operation table

	Standard well	Sample well	Control well
Standards with different concentrations ( $\mu\text{L}$ )	40		
Sample ( $\mu\text{L}$ )		40	40
Reagent 1 ( $\mu\text{L}$ )	40		40
Reagent 2 Working solution ( $\mu\text{L}$ )		40	
Mix fully and incubate at 37°C for 10 min.			
Reagent 4 ( $\mu\text{L}$ )	160	160	160
Mix fully. Measure the OD values of each well.			

## ▲ Calculation

Plot the standard curve by using OD value of standard and correspondent concentration as y-axis and x-axis respectively. Create the standard curve with graph software (or EXCEL). The concentration of the sample can be calculated according to the formula based on the OD value of sample.

The standard curve is:  $y = ax + b$ .

### 1. Serum (plasma) sample

Definition: The amount of 1  $\mu\text{mol}$  p-nitrophenol produced by 1 L serum (plasma) per minute hydrolysis PNPP at 37°C is defined as 1 activity unit.

$$\text{ACP activity (U/L)} = (\Delta A - b) \div a \div T \times f \times 1000^*$$

### 2. Tissue sample

Definition: The amount of 1  $\mu\text{mol}$  p-nitrophenol produced by 1 g tissue protein per minute hydrolysis PNPP at 37°C is defined as 1 activity unit.

$$\text{ACP activity (U/L)} = (\Delta A - b) \div a \div T \times f \times 1000^* \div C_{pr}$$

#### Note:

y:  $OD_{\text{Standard}} - OD_{\text{Blank}}$  ( $OD_{\text{Blank}}$  is the OD value when the standard concentration is 0);

x: The concentration of Standard;

a: The slope of standard curve;

b: The intercept of standard curve;

$\Delta A$ : ( $OD_{\text{Sample}} - OD_{\text{Control}}$ );

f: Dilution factor of sample before test;

T: Reaction time, 10 min;

$C_{pr}$ : Concentration of protein in tissue sample, gprot/L;

1000\*: 1 mmol=1000  $\mu\text{mol}$

## Appendix I Data

### ▲ Example analysis

For rat kidney tissue, take 10% rat kidney tissue homogenate diluted for 20 times and carry the assay according to the operation table. The results are as follows:

Standard curve:  $y = 2.1298x + 0.0025$ , the average OD value of the sample is 0.470, the average OD value of the control is 0.06, the concentration of protein in sample is 9.47 gprot/L, and the calculation result is:

$$\text{ACP activity (U/gprot)} = (0.470 - 0.06 - 0.0025) \div 2.1298 \div 10 \div 9.47 \times 20 \times 1000 \\ = 40.4 \text{ U/gprot}$$

## Appendix II References

1. Bull H, Murray P G, Thomas D, et al. Acid phosphatases [J]. *Molecular Pathology*, 2002, 55(2):65.
2. Mclachlan K D. Acid phosphatase activity of intact roots and phosphorus nutrition in plants. 2. Variations among wheat roots [J]. *Crop and Pasture Science*, 1980, 31(3): 441-448.