

PRODUCT INFORMATION & MANUAL

Amylase (alpha/beta)
Activity Assay Kit
(Colorimetric)
NBP3-25912

For research use only.

Not for diagnostic or therapeutic procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

Amylase (αlpha/beta) Activity Assay Kit (Colorimetric)

Catalog No: NBP3-25912

Method: Colorimetric method

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

Measuring instrument: Microplate reader

Sensitivity: 0.008 U/mL

Detection range: 0.01-0.37 U/mL

Average intra-assay CV (%): 2.8

Average inter-assay CV (%): 3.6

Average recovery rate (%): 97

- ▲ 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 measure α -amylase and β -amylase activity in serum, plasma, saliva, tissue samples.

▲ Background

Amylase is a general term for hydrolase and glycogen enzymes, which generally acts on soluble starch, amylose, glycogen, etc., and is widely present in animals, plants and microorganisms. It is one of the most important enzymes in all industrial enzymes.

▲ Detection principle

The reducing sugar reacts with 3,5-dinitrosalicylic acid under heating conditions to produce a brown-red substance. Amylase activity can be calculated by measuring the OD value at 540 nm.

▲ Kit components & storage

Item	Component	Specification	Storage
Reagent 1	Substrate	10 mL×1 vial	2-8°C , 12 months
Reagent 2	Chromogenic Agent	20 mL×1 vial	2-8℃ , 12 months, shading light
Reagent 3	10 mg/mL Standard	1.5 mL×1 vial	2-8°C , 12 months
	Microplate	96 wells	No requirement
	Plate Sealer	2 pieces	

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.

▲ Materials prepared by users



✓ Instruments

Test tubes, Vortex Mixer, Centrifuge, Water bath, Microplate reader (540 nm)

▲ 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. For measuring the OD value, if there is precipitation, centrifuge at 4000 g for 5 min at room temperature and take the supernatant for determination.
- 2. When the absolute OD value is more than 0.8, sample should be diluted appropriately.

Pre-assay preparation

▲ Reagent preparation

- 1. Bring all reagents to room temperature before use. Before the experiment, preheat reagent 1 and reagent 2 at 40°C for 10 min.
- 2. If there is precipitation in reagent 1, heat it at 70-80°C in water bath until dissolved. Cool down to 40°C with fresh water before use.
- 3. If there is yellow precipitation in reagent 2, heat it at 70-80°C in water bath until dissolved. Cool down to 40°C with fresh water before use.

▲ Sample preparation

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

2. Tissue sample:

Accurately weigh 0.1 g tissue, add 0.9 mL of double distilled water and mechanical homogenate the sample in ice water bath. Collect the tissue homogenate, stand at room temperature for 15 min and oscillate per 5 min, then centrifuge at 3000 g for 10 min at room temperature, then take the supernatant and add double distilled water to a final volume of 10 mL and it is the prepared sample.

3. Saliva sample:

Use a sterile container to collect saliva samples. Remove particulates by centrifugation for 10 min at 3000 g. It is recommended to take fresh saliva sample.

▲ 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 (0.01-0.37 U/mL).

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

Sample type	Dilution factor		
Human plasma	10-15		
Human serum	10-15		
Human saliva	15-25		
Mouse serum	15-25		
Rat serum	15-25		
1% Corn grain tissue homogenate	2-5		

Note: The diluent is double distilled water.

Assay protocol

▲ Plate set up

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

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

▲ Detailed operation steps

1. The preparation of standard curve

Dilute 10 mg/mL standard with double distilled water to a serial concentration. The recommended dilution gradient is as follows: 0, 0.2, 0.4, 0.6, 0.8, 1.0, 1.2, 1.4 mg/mL. Reference is as follows:

Number	Standard concentrations (mg/mL)	10 mg/mL Standard (μL)	Double distilled water (µL)
Α	0	0	1000
В	0.2	20	980
С	0.4	40	960
D	0.6	60	940
Е	0.8	80	920
F	1.0	100	900
G	1.2	120	880
Н	1.4	140	860

2. The measurement of standard

- 1) Take 1.5 mL EP tube and number the tubes from A to H in duplication, add 75 µL of standard solution with different concentrations to the corresponding tubes.
- 2) Add 75 µL of reagent 1 to each tube.
- 3) Add 150 µL of reagent 2 to each tube.
- 4) Mix fully and incubate at 95°C for 5 min. Cool the tubes with running water and take 250 μL of supernatant to the microplate. Measure the OD value of each well with microplate reader at 540 nm.

3. The measurement of samples

- 1) Sample tube: Add 75 µL of sample to the corresponding tubes. Control tube: Add 75 µL of sample to the corresponding tubes.
- 2) Sample tube: Add 75 μL of reagent 1 to the corresponding tubes.
 Control tube: Add 75 μL of double distilled water to the corresponding tubes.
- 3) Incubate the sample tubes and control tubes at 40°C water bath for 5 min.
- 4) Add 150 µL of reagent 2 to each tube.
- 5) Mix fully and incubate at 95°C for 5 min. Cool the tubes with running water and take 250 μL of supernatant to the microplate. Measure the OD value of each well with microplate reader at 540 nm.

▲ Summary operation table

The measurement of standard

	Standard tubes
Standard solution with different concentrations (µL)	75
Reagent 1 (μL)	75
Reagent 2 (µL)	150

Mix fully and incubate at 95°C for 5 min. Cool the tubes and take 250 μ L of supernatant to the microplate. Measure the OD value of each well.

The measurement of sample

	Control tubes	Sample tubes				
Sample (µL)	75	75				
Double distilled water (µL)	75					
Reagent 1 (µL)		75				
Incubate the sample tubes and control tubes at 40°C water bath for 5 min.						
Reagent 2 (µL)	150	150				
Mix fully and incubate at 95°C for 5 min. Cool the tubes with running water and take 250 µL of supernatant to the microplate. Measure the OD value 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.

Serum (plasma) and other saliva sample:

Definition: The production of 1 mg reducing sugar catalyzed by 1 mL of sample per minute that is defined as an enzyme activity unit.

$$(\alpha+\beta)$$
 Amylase activity(U/mL) = $(\Delta A - b) \div a \times V_3 \div t \div V_2 \times f$

Tissue sample:

1) Calculate according to the protein concentration of the sample

Definition: The production of 1 mg reducing sugar catalyzed by 1 mg of tissue protein per minute that is defined as an enzyme activity unit.

$$(\alpha+\beta)$$
 Amylase activity (U/mgprot) = $(\Delta A - b) \div a \times V_3 \div t \div V_2 \div C_{pr}$

2) Calculate according to the fresh weight of sample

Definition: The production of 1 mg reducing sugar catalyzed by 1 g of tissue per minute that is defined as an enzyme activity unit.

$$(α+β)$$
 Amylase activity (U/g fresh weight) = $(ΔA - b) \div a \times V_3 \div t \div w \times \frac{V_1}{V_2} \times f$

Note:

- y: OD_{Standard} OD_{Blank}. (OD_{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.
- f: Dilution factor of sample before tested.
- $\Delta A:\ OD_{\text{Sample}} OD_{\text{Control}}.$
- V_1 : The volume of prepared tissue sample in sample preparation step (10 mL).
- V_2 : The volume of sample added to the reaction (0.075 mL).
- V_3 : The volume of enzymatic reaction (the volume of sample + the volume of reagent 1 = 0.15 mL).
- t: The time of enzymatic reaction (5 min).
- w: The weight of tissue sample (0.1 g).
- C_{pr} : Concentration of protein in sample (mgprot/mL).

Appendix I Data

▲ Example analysis

For rat serum, take 10 μ L of rat serum, dilute with double distilled water for 25 times and carry the assay according to the operation table.

The results are as follows:

standard curve: y=0.8729x-0.0112, the average OD value of the sample is 0.299, the average OD value of the control is 0.162, and the calculation result is: $(\alpha+\beta)$ Amylase activity (U/mL)

 $=(0.299-0.162+0.0112)\div0.8729\times0.15\div5\div0.075\times25=1.70 \text{ U/mL}$

Appendix II References

1. Bhatt B, Prajapati V, Patel K, Trivedi U. Kitchen waste for economical amylase production using Bacillus amyloliquefaciens KCP2. Biocatalysis and Agricultural Biotechnology. 2020: 101654.