

PRODUCT INFORMATION & MANUAL

Polyphenol Oxidase/PPO Activity Assay Kit (Colorimetric) NBP3-25779

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

Polyphenol Oxidase (PPO) Activity Assay Kit

Catalog No: NBP3-25779

Method: Colorimetric method

Specification: 100Assays (Can detect 50 samples without duplication)

Measuring instrument: Spectrophotometer

Average intra-assay CV (%): 4.6

Average inter-assay CV (%): 9.8

- ▲ 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 detect polyphenol oxidase (PPO) activity in plant tissue samples.

▲ Background

Polyphenol oxidase (PPO) is one of the most widely distributed metalloproteinases in the nature. It is ubiquitous in plants, fungi, and insects. The activity of polyphenol oxidase can be detected even on the decaying plant residues of the soil. PPO catalyzes the formation of lignin and quinone compounds, which can prevent cells from being harmed by pathogens, and can also play a direct role in disease resistance by forming quinone substances. Therefore, through the research on the activity of PPO, it can more directly reflect the disease resistance of plant organism in the process of growth.

▲ Detection principle

Polyphenol oxidase (PPO) can catalyze phenolic compounds into quinone substances. The latter has specific absorption at 410 nm. The activity of PPO can be calculated indirectly by measuring the OD value at 410 nm.

▲ Kit components & Storage

	Component	Specification	Storage
Reagent 1	Extracting Solution	60 mL × 2 vials	2-8°C , 12 months
Reagent 2	Buffer Solution	40 mL × 2 vials	2-8°C , 12 months
Reagent 3	Substrate	20 mL × 1vial	2-8°C , 12 months, shading light

Note: The reagents must be stored strictly according to the preservation conditions in the above table.

▲ Materials prepared by users



1 Instruments

Spectrophotometer (410 nm), Tubes, Micropipette, Vortex mixer, 100°C Water bath



Reagents:

Double distilled water

▲ 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.

A Precautions

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

▲ The key points of the assay

- 1. The temperature and time of incubation at 37°C must be accurately.
- 2. The explosion-proof EP tubes are recommended to use for the 100°C water bath.
- 3. It is a normal phenomenon that suspended substance appeared in some tubes, you can centrifuge at 11000 g for 10 min at room temperature, then take the supernatant for measuring the OD value.

Pre-assay preparation

▲ Reagent preparation

- 1. Preheat the reagent 1 at 37°C for 20 min before use, and then use it after completely clarified.
- 2. Bring the reagent 2 and reagent 3 to room temperature before use.

▲ Sample preparation

1. Extraction of crude enzyme solution A

Accurately weigh the plant tissue sample, add reagent 1 according to the ratio of Weight (g): Volume (mL) =1:9. Mechanical homogenate the sample in ice water bath. Centrifuge at 11000 g for 15 min, take the supernatant for detection. Meanwhile, determine the protein concentration of supernatant (E-BC-K168-M,E-BC-K168-S).

2. Extraction of crude enzyme solution B (For control tubes)

After the crude enzyme solution A was extracted, 50% of the supernatant was taken to a new 1.5mL EP tube and boiled at 100° C for 5 min. Cool the tubes with running water and crude enzyme solution B was prepared.

▲ 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.

Sample type	Dilution factor
10% Pepper tissue homogenization	1
10% Corn tissue homogenization	1
10% Potato tissue homogenization	1
10% Ginger tissue homogenization	1
10% Apple tissue homogenization	1
10% Pear tissue homogenization	1
10% Chinese yam tissue homogenization	1

Assay protocol

▲ Detailed operating steps

- Control tube: Add 600 μL of reagent 2 into 1.5 mL EP tubes.
 Sample tube: Add 600 μL of reagent 2 into 1.5 mL EP tubes.
- 2. Add 150 µL of reagent 3 into each tubes.
- 3. Sample tube: Add 150 μ L of crude enzyme solution A into sample tubes. Control tube: Add 150 μ L of crude enzyme solution B into control tubes.
- 4. Mix fully with the vortex mixer, incubate accurately at 37°C for 3 min, incubate at 100°C water bath for 5 min immediately. Then cool the tubes to room temperature with running water.
- 5. Set spectrophotometer to zero with double distilled water and measure the OD value of each tube with 1 mL quartz cuvette at 410 nm. (The OD value of the sample tube is record as A_1 , the OD value of the control tube is record as A_2 , \triangle $A = A_1 A_2$).

▲ Summary operation table

	Sample tube	Control tube
Reagent 2 (µL)	600	600
Reagent 3 (µL)	150	150
Crude enzyme solution A (µL)	150	
Crude enzyme solution B (µL)		150

Mx fully, incubate accurately at 37°C for 3 min, incubate at 100°C water bath for 5 min immediately. Then cool the tubes to room temperature with running water. Set spectrophotometer to zero with double distilled water and measure the OD value of each tube with 1 mL quartz cuvette at 410 nm.

▲ Calculation

Definition: 0.01 OD value changed at 410 nm by 1 mg of tissue protein sample per minute in the reaction system at 37°C that is defined as an enzyme activity unit.

PPO activity(U/mgprot)=
$$\Delta A \div 0.01 \div V \div C_{pr} \div T \times f = 222.2 \times \Delta A \div C_{pr} \times f$$

Note:

 ΔA : $\Delta A = A_1 - A_2$

V: The volume of sample added to the reaction, 0.15 mL.

T: Reaction time, 3 min;

 C_{pr} : The concentration of protein in sample, mgprot/mL.

f: The dilution factor of sample before tested.

Appendix I Data

▲ Example analysis

For chinese yam tissue, take 0.1 g of chinese yam tissue, add 0.9 mL of reagent 1, then homogenize the sample in ice water bath, centrifuge at 10000 g for 10 min at 4°C , then take 0.15 mL of chinese yam tissue supernatant and carry the assay according to the operation table. The results are as follows:

the average OD value of the sample (A_1) is 0.324, the average OD value of the control (A_2) is 0.129, $\Delta A = A_1 - A_2 = 0.324 - 0.129 = 0.195, the concentration of protein in sample is 1.97 mgprot/mL, and the calculation result is:$

PPO activity (U/mgprot) =
$$\frac{222.2 \times 0.195}{1.97} = 21.99 \text{ U/mgprot}$$