

PRODUCT INFORMATION & MANUAL

Nitric Oxide Assay Kit (Colorimetric) NBP3-24525

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

Nitric Oxide Assay Kit (Colorimetric)

Catalog No: NBP3-24525

Method: Colorimetric method

Specification: 96T (Can detect 80 samples without duplication)

Instrument: Microplate reader

Sensitivity: 0.16 µmol/L

Detection range: 0.16-100 µmol/L

Average intra-assay CV (%): 2.4

Average inter-assay CV (%): 3.7

Average recovery rate (%): 102

- ▲ 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 for detection of nitric oxide (NO) in serum, plasma, saliva, animal and plant tissue samples.

▲ Background

The half-life of NO is extremely short, it exists in form of nitrate or nitrite produced by vascular endothelial cell, vascular smooth muscle cell, platelet, and macrophage and so on. The concentration of NO can be indirectly measured by detecting that of nitrate or nitrite.

NO react with oxygen and water to generate nitrate or nitrite which can form a kind of pale red azo compound when meet with nitrate chromogenic reagent, the absorbance of the compound can be measured to calculate the concentration of NO indirectly.

▲ Detection principle

NO is easily oxidized to form NO₂ in vivo or in aqueous solution, and a reddish azo compound is formed with the color developing agent, and the concentration of the azo compound is linearly related to the concentration of NO. The concentration of NO can be calculated indirectly by measuring the OD value at 550 nm.

$$NO_2$$
 + R_1 NH_2 H^+ $N^{\geq N^+}$ R_1 R_2 R_2 R_2

▲ Kit components & storage

Item	Component	Specification	Storage	
Reagent 1	Sulphate Solution	24 mL × 1 vial	2-8°C , 12 months	
Reagent 2	Alkali Reagent	12 mL × 1 vial	2-8°C , 12 months	
Reagent 3	Chromogenic Agent A	1.9 mL × 2 vials	2-8°C , 12 months, shading light	
Reagent 4	Chromogenic Agent B	Powder × 1 vial	2-8℃, 12 months, shading light	
Reagent 5	Acid Solution	1.3 mL × 2 vials	2-8°C , 12 months	
Reagent 6	Sodium Nitrite Standard	Powder × 2 vials	-20°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



1 Instruments

Microplate reader(540-550 nm), Micropipettor, Vortex mixer, Centrifuge



Reagents:

Double distilled water, Normal saline (0.9% NaCl), 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. Use disposable EP tubes or clean glass tubes with stopper for centrifugation.
- 2. The supernatant for assay should not contain sediment, otherwise it will affect the results.
- 3. All reagents should be prepared the day before the assay, let it fully dissolved. Please add reagents to the bottom of well vertically and slowly, avoid to add on the wall of well and generate bubble.
- 4. Serum samples can be stored for 3 days at $4^{\circ}\mathbb{C}$ and for a month at -20 $^{\circ}\mathbb{C}$.

Pre-assay preparation

▲ Reagent preparation

- 1. If there is any crystal precipitation in reagent 3, please dissolve it fully with water bath at above 60 ℃ before use.
- 2. Preparation of reagent 4 working solution
 - Dissolve a vial of reagent 4 with 3.8 mL of double distilled water fully. The prepared solution can be stored at 4°C for 2 months with shading light.
- 3. Preparation of chromogenic reagent
 - Mix the reagent 3, reagent 4 working solution and reagent 5 at a ratio of 3:3:2 fully. Prepare the fresh solution before use and it can't be used when its color gets darker.
- Preparation of 2 mmol/L Sodium Nitrite Standard
 Dissolve reagent 6 with 2 mL of double distilled water. Prepare the needed amount before use.

▲ Sample preparation

The samples should be prepared as conventional methods. Also please refer to appendix II.

▲ 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 (0.16-100 µmol/L).

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

Sample type	Dilution factor
Human serum	1
Human plasma	1
10% Mouse liver tissue homogenization	1
Rat serum	1
Rat plasma	1
10% Epipremnum aureum tissue homogenization	1

Note: The diluent is double distilled water, normal saline (0.9% NaCl) or PBS (0.01 M, pH 7.4).

Assay protocol

▲ Plate set up

	1	2	3	4	5	6	7	8	9	10	11	12
Α	Α	Α	S1	S9	S17	S25	S33	S41	S49	S57	S65	S73
В	В	В	S2	S10	S18	S26	S34	S42	S50	S58	S66	S74
С	С	С	S3	S11	S19	S27	S35	S43	S51	S59	S67	S75
D	D	D	S4	S12	S20	S28	S36	S44	S52	S60	S68	S76
Е	Е	Е	S5	S13	S21	S29	S37	S45	S53	S61	S69	S77
F	F	F	S6	S14	S22	S30	S38	S46	S54	S62	S70	S78
G	G	G	S7	S15	S23	S31	S39	S47	S55	S63	S71	S79
Н	Н	Н	S8	S16	S24	S32	S40	S48	S56	S64	S72	S80

Note: A, blank well; B-H, standard wells; S1-S80, sample wells.

▲ Detailed operation steps

1. The preparation of standard curve

Dilute 2 mmol/L sodium nitrite standard with double distilled water to a serial concentration. The recommended dilution gradient is as follows: 0, 10, 20, 30, 40, 60, 80, 100 μ mol/L. Reference is as follows:

Number	Standard concentrations (µmol/L)	2 mmol/L sodium nitrite standard (μL)	Double distilled water (µL)
Α	0	0	2000
В	10	10	1990
С	20	20	1980
D	30	30	1970
Е	40	40	1960
F	60	60	1940
G	80	80	1920
Н	100	100	1900

2. The measurement of samples

(1) Standard tubes: Take a* µL of sodium nitrite standard solution with different concentrations to 1.5 mL EP tubes.

Sample tubes: Take a* µL of sample to 1.5 mL EP tubes.

Note: a* = Sample volume= Standard volume.

For serum or plasma samples, a* is 200-300 µL.

For tissue, a* is $100-300 \mu L$.

- (2) Add 200 µL of reagent 1 and mix fully with a vortex mixer.
- (3) Add 100 μL of reagent 2 and mix fully with a vortex mixer.
- (4) Stand for 15 min at room temperature, centrifuge at 3100 g for 10 min. (If there is precipitate in the supernatant, please transfer the supernatant to a new EP tube and centrifuge again.)
- (5) Take 160 μ L of supernatant to the corresponding wells of microplate for chromogenic reaction.
- (6) Add 80 μL of chromogenic reagent to each well, oscillate for 2 min and stand at room temperature for 15 min.
- (7) Measure the OD value at 550 nm with microplate reader.

▲ Summary operation table

1. Pre-treatment

Standard	Sample
a*	
	a*
200	200
100	100
	a* 200

Mix fully and stand for 15 min at room temperature, centrifuge at 3100 g for 10 min, take 160 μL of the supernatant for chromogenic reaction.

2. Chromogenic reaction

	Standard	Sample
Supernatant (µL)	160	160
Chromogenic reagent (µL)	80	80

Mix thoroughly for 2 min, stand for 15 min at room temperature, measure the \mbox{OD} 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. If the sample have been diluted, the concentration calculated from the standard curve must be multiplied by the dilution factor. The actual concentration is the calculated concentration multiplied by dilution factor. The standard curve is: y= ax + b.

1. Serum (plasma):

NO content (
$$\mu$$
mol/L) = (ΔA_{550} -b) ÷ a × f

2. Tissue:

NO content (
$$\mu$$
mol/gprot) = (ΔA_{550} -b) ÷ a × f ÷ C_{pr}

Note:

- y: The absolute OD value of standard;
- x: The concentration of standard;
- a: The slope of standard curve;
- b: The intercept of standard curve.
- ΔA_{550} : Absolute OD (OD_{Sample} OD_{Blank})
- f: Dilution factor of sample before test.
- C_{pr}: Concentration of protein in sample (gprot/L)

Appendix I Data

▲ Example analysis

Dilute human serum for 2 times, carry the assay according to the operation table.

The results are as follows:

standard curve: y = 0.00215 x + 0.00514, the average OD value of the sample is 0.056, the average OD value of the blank is 0.035, the concentration of protein in sample is 9.23 gprot/L, and the calculation result is:

NO content(μ mol/L)= (0.056-0.035-0.00514)÷0.00215=3.20 (μ mol/L)

Appendix II Sample preparation

The following sample pretreatment methods are for reference only.

▲ Serum

Collect fresh blood and stand at 25° C for 30 min to clot the blood. Then centrifuge at 2000 g for 15 min at 4° C. Take the serum (which is the upper light yellow clarified liquid layer) to preserve it on ice for detection.

▲ Plasma

Take fresh blood into the tube which has anticoagulant, centrifuge at 1000-2000 g for 10 min at 4° C. Take the plasma (which is the upper light yellow clarified liquid layer, don't take white blood cells and platelets in the middle layer) to preserve it on ice for detection.

▲ Tissue sample

Take 0.02-1g fresh tissue to wash with PBS (0.01 M, pH 7.4) at 2-8 $^{\circ}$ C . Absorb the water with filter paper and weigh. Homogenize at the ratio of the volume of homogenized medium (2-8 $^{\circ}$ C) (mL): the weight of the tissue (g) =9:1, then centrifuge the tissue homogenate for 10 min at 10000 g at 4 $^{\circ}$ C . Take the supernatant to preserve it on ice for detection. Meanwhile, determine the protein concentration of supernatant.

▲ Saliva

Gargle with clear water, collect the saliva 30 min later, centrifuge at 10000 g for 10 min at 4° C . Take the supernatant and preserve it on ice for detection.

Blake singered

Note:

- 1. Homogenized medium: PBS (0.01 M, pH 7.4).
- 2. Homogenized method:
- (1) Hand-operated: Weigh the tissue and mince to small pieces (1 mm³), then put the tissues pieces to glass homogenized tube. Add homogenized medium into homogenized tube, place the tube into the ice bath with left hand, and insert the glass tamping rod vertically into the homogenized tube with the right hand to grind up and down for 6-8 min.
 - Or put the tissue into the mortar, and add liquid nitrogen to grind fully. Then add the homogenized medium to homogenize.
- (2) Mechanical homogenate: Weigh the tissue to EP tube, add the homogenized medium to homogenize the tissue with homogenizer instrument (60 Hz, 90s) in the ice bath. (For samples of skin, muscle and plant tissue, the time of homogenization can be properly prolonged.)

Appendix III References

- 1. Davis K L, Martin E, Turko I V, Murad F. Novel effects of nitric oxide. Annu Rev Pharmacol Toxicol, 2001, 41(1): 203-236.
- 2. Sun J, Zhang X, Broderick M, Fein H. Measurement of Nitric Oxide Production in Biological Systems by Using Griess Reaction Assay. Sensor, 2003, 3(8): 276-284