



**PRODUCT INFORMATION &
MANUAL**

**Glutathione S-Transferase/
GST Activity Assay Kit
(Colorimetric)
*NBP3-24499***

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

Glutathione S-Transferase/GST Activity Assay Kit (Colorimetric)

Catalog No: NBP3-24499

Method: Colorimetric method

Specification: 100 Assays (Can detect 96 samples without duplication)

Instrument: Spectrophotometer

Sensitivity: 1 U/L

Detection range: 1-79 U/L

Average intra-assay CV (%): 1.9

Average inter-assay CV (%): 4.3

Average recovery rate (%): 105

- ▲ 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 the glutathione-S-transferase (GSH-ST) activity in serum, plasma, tissue and cell samples.

▲ Background

According to the localization of cell, glutathione-S-transferase (GSH-ST) is divided into three subfamilies: cytoplasmic GSH-ST, microsomal GSH-ST and peroxisome/mitochondrial GSH-ST. GSH-ST is a kind of detoxifying enzyme. Maintain cell integrity by catalyzing the binding of reduced glutathione to electron-friendly substance.

▲ Detection principle

GST can catalyze the binding of reduced glutathione (GSH) to dinitrobenzene (CDNB) and the product have an absorption peak at 340 nm. The activity of GSH-ST can be calculated by measuring the increasing rate of absorbance at 340 nm.

▲ Kit components & Storage

Item	Component	Specification	Storage
Reagent 1	Extracting Solution	60 mL × 2 vials	2-8°C , 12 months
Reagent 2	Buffer Solution	50 mL × 2 vials	2-8°C , 12 months
Reagent 3	Powder	2 vials	2-8°C , 12 months

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

Spectrophotometer (340 nm), Micropipettor, Incubator, Vortex mixer

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.

▲ Precautions

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

▲ The key points of the assay

1. The reagent 3 application solution and cuvette should be preheat at 37°C for 10 min .
2. The reaction time must be accurate.

Pre-assay preparation

▲ Reagent preparation

Reagent 3 application solution: Dissolve a vial of powder with 5 mL double distilled water and mix fully. Prepared solution can be stored at 2-8°C for 3 days.

▲ 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 the results of the samples are linear within 5 min.

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

Sample type	Dilution factor
Human serum (plasma)	/
10% Rat liver tissue homogenization	150-200
10% Rat lung tissue homogenization	8-12
10% Rat kidney tissue homogenization	10-15
10% Plant tissue homogenization	/

Assay protocol

▲ Detailed operating steps

1. Preheat the needed reagent 3 application solution and cuvette at 37°C for 10 min. Set the spectrophotometer to 340 nm and set the spectrophotometer to zero with double distilled water.
2. **Blank tube:** add 0.1 mL of reagent 1 to a 2 mL EP tube.
Sample tube: add 0.1 mL of sample to a 2 mL EP tube.
3. Add 0.9 mL of reagent 2 and 0.1 mL of reagent 3 application solution to each tube, mix fully and record the time immediately. Measure the absorbance at 340 nm at 20 s (A_1) and 320 s (A_2), respectively. Calculate the $\Delta A_{\text{blank or sample}} = A_2 - A_1$.

▲ Summary operation table

Preheat the Reagent 3 application solution and cuvette at 37°C for 10 min.		
	Blank tube	Sample tube
Reagent 1 (mL)	0.1	
Sample (mL)		0.1
Reagent 2 (mL)	0.9	0.9
Reagent 3 application solution (mL)	0.1	0.1
Mix fully and record the time immediately. Measure the absorbance at 340 nm at 20 s (A_1) and 320 s (A_2), respectively. Calculate the $\Delta A_{\text{blank or sample}} = A_2 - A_1$.		

▲ Calculation

1. Serum (plasma) and other liquid sample:

Definition: The amount of GSH-ST in 1 mL of sample that catalyze the combination of 1 μmol of CDNB and GSH at 37°C per minute is defined as 1 unit.

$$\text{GST activity (U/mL)} = \frac{\Delta A}{\epsilon \times d} \times 10^6 \div t \times \frac{V_1}{V_2} \times f$$

2. Tissue and cells sample:

Definition: The amount of GSH-ST in 1 mg of tissue protein that catalyze the combination of 1 μmol of CDNB and GSH at 37°C per minute is defined as 1 unit.

$$\text{GST activity (U/mgprot)} = \frac{\Delta A}{\epsilon \times d} \times 10^6 \div t \times \frac{V_1}{V_2} \times f \div C_{pr}$$

Note:

ΔA : $\Delta A_{\text{Sample}} - \Delta A_{\text{Blank}}$

ϵ : molar extinction coefficient of the product, $9.6 \times 10^3 \text{ L/mol/cm}$

d : optical path of the cuvette, 1 cm

10^6 : $1 \text{ mol} = 10^6 \mu\text{mol}$

V_1 : the total volume of the reaction system, (1.1 mL = 0.0011 L)

V_2 : the volume of sample added into the reaction system, 0.1 mL

t : reaction time, 5 min

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

f : dilution factor of sample before test

Appendix I Data

▲ Example analysis

Dilute 10% rat kidney tissue homogenate with reagent 1 at the ratio of 1:14, take 0.1 mL of diluted sample and carry the assay according to the operation table.

The results are as follows:

The average OD value of the sample (A_1) is 0.464, the average OD value of the sample (A_2) is 0.729, the average OD value of the blank (A_1) is 0.414, the average OD value of the blank (A_2) is 0.437, the concentration of protein in sample is 8.51 mgprot/mL, and the calculation result is:

$$\Delta A = (0.729 - 0.464) - (0.437 - 0.414) = 0.242$$

$$\begin{aligned} \text{GST activity (U/mgprot)} &= \frac{0.242}{1 \times 9.6 \times 10^3} \times 10^6 \div 5 \times \frac{0.0011\text{L}}{0.1\text{mL}} \times 15 \div 8.51\text{mgprot/mL} \\ &= 0.098 \text{ U/mgprot} \end{aligned}$$

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. If not detected on the same day, the serum can be stored at -80°C for a month.

▲ Plasma

Take fresh blood into the tube which has anticoagulant (Heparin is recommended), centrifuge at 700-1000 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. If not detected on the same day, the plasma can be stored at -80°C for a month.

▲ Tissue sample

Take 0.02-1g fresh tissue to wash with PBS (0.01 M, pH 7.4) at 2-8°C . Absorb the water with filter paper and weigh. Homogenize at the ratio of the volume of homogenized medium (2-8°C) (mL): the weight of the tissue (g) =9:1, then centrifuge the tissue homogenate for 10 min at 10000 g at 4°C . Take the supernatant to preserve it on ice for detection. Meanwhile, determine the protein concentration of supernatant. If not detected on the same day, the tissue sample (without homogenization) can be stored at -80°C for a month.

▲ Cells

Collect the cells and wash the cells with PBS (0.01 M, pH 7.4) for 1~2 times. Centrifuge at 1000 g for 10 min and then discard the supernatant and keep the cell sediment. Add homogenization medium at a ratio of cell number (10^6): homogenization medium (μL) =1: 300-500. Sonicate or grind with hand-operated in ice water bath. Centrifuge at 10000 g for 10 min, then take the supernatant and preserve it on ice for detection. Meanwhile, determine the protein concentration of supernatant. If not detected on the same day, the cells sample (without homogenization) can be stored at -80°C for a month.

Note:

1. Homogenized medium: Reagent 1.
2. Homogenized method:
 - (1) Hand-operated: Weigh the tissue and mince to small pieces (1 mm^3), 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.)
 - (3) Ultrasonication: Treat the cells with ultrasonic cell disruptor (200 W, 2 s/ time, interval for 3 s, the total time is 5 min).

Appendix III References

1. Dasari S, Gonuguntla S, Ganjaji M S, et al. Genetic polymorphism of glutathione S-transferases: Relevance to neurological disorders.[J]. Pathophysiology, 2018.
2. Remmerie B, Vandenbroucke K, Smet L D, et al. Expression, purification, crystallization and structure determination of two glutathione S-transferase-like proteins from *Shewanella oneidensis*[J].Acta Crystallographica, 2008, 64(6): 548-553.