

GST Colorimetric Activity Assay Kit

(Catalog #K263-100; 100 assays; Store kit at -20°C)

I. Introduction:

Glutathione S-transferase (GST) is a family of enzymes that play an important role in detoxification of xenobiotics. GST catalyzes the formation of the thiol group of glutathione to electrophilic xenobiotics. It utilizes glutathione to scavenge potentially toxic compounds including those produced as a result of oxidative stress and is part of the defense mechanism against the mutagenic, carcinogenic and toxic effects of such compounds. The *GST Colorimetric Activity Assay Kit* is based upon the GST-catalyzed reaction between GSH and the GST substrate, CDNB (1-chloro-2,4-dinitrobenzene, which has the broadest range of isozyme detectability (e.g., alpha-, mu-, pi-, and other GST isoforms). Under certain conditions, the interaction between glutathione and CDNB is totally dependent on the presence of active GST.



The GST-catalyzed formation of GS-DNB produces a dinitrophenyl thioether which can be detected by spectrophotometer at 340 nm. One unit of GST activity is defined as the amount of enzyme producing 1 μmol of GS-DNB conjugate/min under the conditions of the assay. The kit can detect GST activity in crude cell lysate or purified protein fraction, and also quantitate GST-tagged fusion protein. Detect limit: Active GST < 4 mU.

II. Kit Contents:

Component	K263-100	Color code	Part
	100 assays	Cap color	Number
GST Assay Buffer	10 ml	NM	K263-100-1
GST Sample Buffer	25 ml	WM	K263-100-2
GST Substrate (CDNB)	1 ml	Red	K263-100-3
Glutathione (GSH, lyophilized)	2 x 17 mg	Green	K263-100-4
GST Positive Control (0.625 $\mu\text{g}/\mu\text{l}$)	20 μl	Yellow	K263-100-5

III. Reagent Preparation and Storage Conditions:

GSH: Add 550 μl of GST Sample Buffer to each vial just before use. One vial is sufficient for 50 assays. Remaining solution can be kept at -20°C for 1 week.

CDNB: This vial contains an ethanol solution of 1-chloro-2,4-dinitrobenzene (CDNB) and should be stored at -20°C.

IV. Sample Preparation Guideline:

A. Cell Sample Preparation:

1. Collect cells by centrifugation. For adherent cells, use a rubber policeman scrapping the cells to collect.
2. Homogenize or sonicate the cells in GST Sample Buffer.

3. Centrifuge at 10,000 x g for 15 minutes at 4°C.
4. Collect supernatant and use for the assay. The remaining sample should be stored at -80°C, stable for at least 1 month.

B. Tissue Sample Preparation:

1. Prior to dissection, perfuse tissue with PBS containing heparin (0.15 mg/ml) to remove red blood cells and clots.
2. Homogenize tissue in GST Sample Buffer (100 mg/0.5 ml).
3. Centrifuge at 10,000 x g for 15 minutes at 4°C.
4. Collect supernatant and use for the assay. The remaining sample should be stored at -80°C, stable for at least 1 month.

C. Plasma and Erythrocyte Sample Preparation:

1. Centrifuge an anticoagulant treated blood at 1000 x g for 10 min at 4°C.
2. Transfer the top plasma layer (without disturbing the white puffy layer) to a new tube and store on ice for assay or store at -80°C for future use. The plasma should be stable for 1 month.
3. Remove the white buffy layer and discard (leukocytes).
4. Lyse the erythrocytes (red blood cells) in 4 times its volume of ice-cold GST sample Buffer.
5. Centrifuge at 10,000 x g for 15 min at 4°C.
6. Transfer supernatant (erythrocyte lysate) to a new tube, and use it for the GST assay. The remaining samples should be stored at -80°C for future use, stable for at least one month.

D. Preparation of Bacterially Expressed GST-Fusion Protein Sample:

1. Collect bacteria by centrifugation. Freeze/thaw the pellet two times, then sonicate in GST Sample Buffer.
2. Centrifuge at 10,000 x g for 15 min at 4°C.
3. Transfer supernatant to a new tube, and use it for the GST assay. The remaining samples should be stored at -80°C for future use, stable for at least one month.

V. GST Assay Protocol:

1. Prepare samples in a total 100 μl volume with GST Sample Buffer, including a negative control with 100 μl of GST Sample buffer only and a positive control with 2 μl of GST and 98 μl of Sample Buffer.

Note: We recommend preparing several dilutions of your sample and perform duplicate wells for each measurement.

2. Add 10 μl of Glutathione to each well containing the sample or control above.
3. Prepare Substrate Mix by adding 10 μl CDNB Solution + 90 μl of GST Assay Buffer for each sample including the standard.
4. Mix well and transfer 100 μl of the Mix into each sample (including the standard) well.
5. Carefully shake the plate to start the reaction.
6. Read the absorbance once every minute at 340 nm using a plate reader to obtain at least 5 time points.

1. Calculation of GST Assay Results:

A. Determine the change in absorbance (ΔA_{340}) per minute by:

- Plotting the absorbance values as a function of time to obtain the slope (rate) of the linear portion of the curve.
- Select two points on the linear portion of the curve and determine the change in absorbance during that time, using the following equation:

$$\Delta A_{340}/\text{min} = \frac{A_{340} (\text{Time 2}) - A_{340} (\text{Time 1})}{\text{Time 2 (min)} - \text{time 1 (min)}}$$

B. Determine the rate of $\Delta A_{340}/\text{min}$ for the background wells and subtract the rate from that of the sample wells.

C. Use the following formula to calculate the GST activity. The reaction rate at 340 nm can be determined using the GS-DNB extinction coefficient at 340 nm $0.0096 \mu\text{M}^{-1}\text{cm}^{-1}$. The value has been adjusted for the path length of the solution in the well 0.524 cm).

$$\begin{aligned} \text{GST Activity} &= \frac{\Delta A_{340}/\text{min} \times 0.2 \text{ ml Reaction Volume}}{0.0096 \mu\text{mol}^{-1} \text{cm}^{-1} \times 1000 \text{ ml} \times 0.524 \text{ cm} \times A} \times \text{Sample Dilution Factor} \\ &= 0.0396 \times \Delta A_{340}/\text{min} \times \text{Sample Dilution/A} \quad (\mu\text{mol}/\text{min}) \end{aligned}$$

Where:

$0.0096 \mu\text{mol}^{-1}\text{cm}^{-1} \times 1000 \text{ ml}$ is GS-DNB extinction coefficient.

A is sample volume added in the reaction well in milliliter (ml).

0.524 cm is light path of the 0.2 ml Reaction Volume in 96 well plate (cm).

FOR RESEARCH USE ONLY! Not to be used in human.

FEATURED PRODUCTS:

Apoptosis Detection Kits & Reagents

- Annexin V Kits & Bulk Reagents
- Caspase Assay Kits & Reagents
- Mitochondrial Apoptosis Kits & Reagents
- Nuclear Apoptosis Kits & Reagents

Cell Fractionation System

- Mitochondria/Cytosol Fractionation Kit
- Nuclear/Cytosol Fractionation Kit
- Membrane Protein Extraction Kit
- Cytosol/Particulate Rapid Separation Kit
- Mammalian Cell Extraction Kit
- FractionPREP Fractionation System

Cell Proliferation & Senescence

- Quick Cell Proliferation Assay Kit
- Senescence Detection Kit
- High Throughput Apoptosis/Cell Viability Assay Kits
- LDH-Cytotoxicity Assay Kit
- Bioluminescence Cytotoxicity Assay Kit
- Live/Dead Cell Staining Kit

Cell Damage & Repair

- HDAC Fluorometric & Colorimetric Assays & Drug Discovery Kits
- HAT Colorimetric Assay Kit & Reagents
- DNA Damage Quantification Kit
- Glutathione & Nitric Oxide Fluorometric & Colorimetric Assay Kits

Signal Transduction

- cAMP & cGMP Assay Kits
- Akt & JNK Activity Assay Kits
- Beta-Secretase Activity Assay Kit

Adipocyte & Lipid Transfer

- Recombinant Adiponectin, Survivin, & Leptin
- CETP Activity Assay & Drug Discovery Kits
- Total Cholesterol Quantification Kit

Molecular Biology & Reporter Assays

- siRNA Vectors
- Cloning Insert Quick Screening Kit
- Mitochondrial & Genomic DNA Isolation Kits
- 5 Minutes DNA Ligation Kit
- 20 Minutes Gel Staining/Destaining Kit

Growth Factors and Cytokines

Monoclonal and Polyclonal Antibodies and Cytokines