



Glutathione S-Transferase Assay Kit

Item No. 703302

www.caymanchem.com

Customer Service 800.364.9897

Technical Support 888.526.5351

1180 E. Ellsworth Rd • Ann Arbor, MI • USA

TABLE OF CONTENTS

| | | |
|------------------------------|----|-----------------------------------|
| GENERAL INFORMATION | 3 | Materials Supplied |
| | 4 | Safety Data |
| | 4 | Precautions |
| | 4 | If You Have Problems |
| | 4 | Storage and Stability |
| | 4 | Materials Needed but Not Supplied |
| INTRODUCTION | 5 | Background |
| | 5 | About This Assay |
| PRE-ASSAY PREPARATION | 6 | Reagent Preparation |
| | 8 | Sample Preparation |
| ASSAY PROTOCOL | 10 | Plate Set Up |
| | 11 | Performing the Assay |
| ANALYSIS | 12 | Calculations |
| | 13 | Performance Characteristics |
| RESOURCES | 14 | Interferences and Inhibitors |
| | 16 | Troubleshooting |
| | 17 | References |
| | 18 | Plate Template |
| | 19 | Notes |
| | 19 | Warranty and Limitation of Remedy |

GENERAL INFORMATION

Materials Supplied

| Item Number | Item | Quantity |
|-------------|------------------------------------|----------|
| 703310 | GST Assay Buffer (2X) | 1 vial |
| 703312 | GST Sample Buffer (2X) | 1 vial |
| 703314 | GST Control | 1 vial |
| 703316 | GST Glutathione | 1 vial |
| 703318 | GST CDNB | 1 vial |
| 400014 | 96-Well Plate (Colorimetric Assay) | 1 plate |
| 400012 | 96-Well Cover Sheet | 1 cover |

If any of the items listed above are damaged or missing, please contact our Customer Service department at (800) 364-9897 or (734) 971-3335. We cannot accept any returns without prior authorization.



WARNING: THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

Safety Data

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent *via* email to your institution.

Precautions

Please read these instructions carefully before beginning this assay.

If You Have Problems

Technical Service Contact Information

Phone: 888-526-5351 (USA and Canada only) or 734-975-3888
Fax: 734-971-3641
Email: techserv@caymanchem.com
Hours: M-F 8:00 AM to 5:30 PM EST

In order for our staff to assist you quickly and efficiently, please be ready to supply the lot number of the kit (found on the outside of the box).

Storage and Stability

This kit will perform as specified if stored at -20°C and used before the expiration date indicated on the outside of the box.

Materials Needed But Not Supplied

1. A plate reader capable of measuring absorbance at 340 nm
2. Adjustable pipettors and a repeating pipettor
3. A source of pure water; glass distilled water or HPLC-grade water is acceptable

INTRODUCTION

Background

Glutathione S-transferases (GSTs) are ubiquitous multifunctional enzymes, which play a key role in cellular detoxification. The enzymes protect cells against toxicants by conjugating them to glutathione, thereby neutralizing their electrophilic sites, and rendering the products more water-soluble.¹ The glutathione conjugates are metabolized further to mercapturic acid and then excreted.¹ Based on their sequence homology, substrate specificity and immunological cross-reactivity, GSTs have been grouped into species-independent classes of isozymes.²⁻⁵ These classes are comprised of both cytosolic and microsomal enzymes.²⁻⁵

About This Assay

Cayman's Glutathione S-Transferase Assay Kit measures total GST activity (cytosolic and microsomal) by measuring the conjugation of 1-chloro-2,4-dinitrobenzene (CDNB) with reduced glutathione.⁶ The conjugation is accompanied by an increase in absorbance at 340 nm. The rate of increase is directly proportional to the GST activity in the sample. The Cayman GST Assay Kit can be used to measure GST activity in plasma, erythrocyte lysates, tissue homogenates, and cell lysates. Cytosolic and microsomal GST activity can be assayed separately following the procedure outlined under **Sample Preparation** (see page 8).

Reagent Preparation

1. GST Assay Buffer (2X) - (Item No. 703310)

Dilute 10 ml of GST Assay Buffer concentrate with 10 ml of HPLC-grade water. This final Assay Buffer (100 mM potassium phosphate, pH 6.5, containing 0.1% Triton X-100) should be used in the assay. When stored at 4°C, this diluted Assay Buffer is stable for at least two months. Equilibrate the diluted Assay Buffer to 25°C before using in the assay.

2. GST Sample Buffer (2X) - (Item No. 703312)

Dilute 5 ml of GST Sample Buffer concentrate with 5 ml of HPLC-grade water. This final Sample Buffer (100 mM potassium phosphate, pH 6.5, containing 0.1% Triton X-100, and 1 mM glutathione) should be used to dilute the GST samples prior to assaying. When stored at 4°C, this diluted Sample Buffer is stable for at least one month.

3. GST Control - (Item No. 703314)

This vial contains a lyophilized powder of equine liver GST. Prior to use, reconstitute the enzyme with 0.5 ml of Milli-Q water. A 20 µl aliquot of this reconstituted enzyme per well results in an increase of approximately 0.05 absorbance unit/minute under the standard assay conditions as described in **Performing the Assay** (see page 11). The reconstituted enzyme can be stored at -20°C for about two weeks. Avoid multiple freeze/thaw cycles.

4. GST Glutathione - (Item No. 703316)

This vial contains a solution of reduced glutathione and should be stored at -20°C when not being used. The reagent is ready to use as supplied.

5. GST CDNB - (Item No. 703318)

This vial contains an ethanolic solution of 1-chloro-2,4-dinitrobenzene (CDNB) and should be stored at -20°C when not being used. The reagent is ready to use as supplied. Equilibrate the reagent to 25°C before using in the assay.

Sample Preparation

The procedures listed below for tissue homogenates and cell lysates will result in assaying total GST activity (cytosolic and microsomal). To separate the two enzymes, centrifuge the 10,000 x g supernatant at 100,000 x g for 60 minutes at 4°C. The resulting 100,000 x g supernatant will contain cytosolic GST and the pellet will contain microsomal GST. Suspend the microsomal pellet in cold buffer (*i.e.*, 100 mM potassium phosphate, pH 7.0, containing 2 mM EDTA). If not assaying on the same day, freeze the samples at -80°C.

Tissue Homogenate

1. Prior to dissection, perfuse or rinse tissue with a PBS (phosphate buffered saline) solution, pH 7.4, to remove any red blood cells and clots.
2. Homogenize the tissue in 5-10 ml of cold buffer (*i.e.*, 100 mM potassium phosphate, pH 7.0, containing 2 mM EDTA) per gram tissue.
3. Centrifuge at 10,000 x g for 15 minutes at 4°C.
4. Remove the supernatant for assay and store on ice. If not assaying on the same day, freeze the sample at -80°C. The sample will be stable for at least one month.

Cell Lysate

1. Collect cells by centrifugation (*i.e.*, 1,000-2,000 x g for 10 minutes at 4°C). For adherent cells, do not harvest using proteolytic enzymes; rather use a rubber policeman.
2. Homogenize or sonicate cell pellet in 1-2 ml of cold buffer (*i.e.*, 100 mM potassium phosphate, pH 7.0, containing 2 mM EDTA).
3. Centrifuge at 10,000 x g for 15 minutes at 4°C.
4. Remove the supernatant for assay and store on ice. If not assaying on the same day, freeze the sample at -80°C. The sample will be stable for at least one month.

Plasma and Erythrocyte Lysate

1. Collect blood using an anticoagulant such as heparin, citrate, or EDTA.
2. Centrifuge the blood at 700-1,000 x g for 10 minutes at 4°C. Pipette off the top yellow plasma layer without disturbing the white buffy layer. Store plasma on ice until assaying or freeze at -80°C. The plasma sample will be stable for at least one month.
3. Remove the white buffy layer (leukocytes) and discard.
4. Lyse the erythrocytes (red blood cells) in four times its volume of ice-cold HPLC-grade water.
5. Centrifuge at 10,000 x g for 15 minutes at 4°C.
6. Collect the supernatant (erythrocyte lysate) for assaying and store on ice. If not assaying the same day, freeze at -80°C. The sample will be stable for at least one month.

Tissue Homogenization using the Precellys 24 Homogenizer

- Prior to dissection, either perfuse or rinse tissue with phosphate buffered saline (PBS), pH 7.4, to remove any red blood cells and clots.
- Freeze organs immediately upon collection and then store at -80°C. Snap-freezing of tissues in liquid nitrogen is preferred.
- Add cold 100 mM potassium phosphate, pH 7.0, containing 2 mM EDTA.
- Homogenize tissue sample using the Precellys 24 according to appropriate settings.
- Spin the tissue homogenates at 10,000 x g for 15 minutes at 4°C.
- Collect supernatant and assay samples according to the kit booklet protocol. Samples may need to be diluted appropriately for assay and should be normalized using a protein assay.

ASSAY PROTOCOL

Plate Set Up

There is no specific pattern for using the wells on the plate. However, it is necessary to have three wells designated as non-enzymatic or background wells. The absorbance rate of these wells must be subtracted from the absorbance rate measured in the GST sample and control wells. We suggest that there be at least three wells designated as positive controls and that you record the contents of each well on the template sheet provided on page 18.

Pipetting Hints

- It is recommended that an adjustable pipette be used to deliver Assay Buffer (dilute), glutathione, and CDNB to the wells.
- Use different tips to pipette the Assay Buffer (dilute), glutathione, enzyme, and CDNB.
- Before pipetting each reagent, equilibrate the pipette tip in that reagent (*i.e.*, slowly fill the tip and gently expel the contents, repeat several times).
- Do not expose the pipette tip to the reagent(s) already in the well.

Performing the Assay

- The final volume of the assay is 200 μ l in all the wells.
 - It is not necessary to use all the wells on the plate at one time.
 - The assay temperature is 25°C.
 - Use the Assay Buffer (dilute) in the assay.
 - Monitor the increase in absorbance at 340 nm using a plate reader.
1. **Background or Non-enzymatic Wells** - add 170 μ l of Assay Buffer and 20 μ l of Glutathione to three wells.
 2. **Positive Control Wells (equine liver GST)** - add 150 μ l of Assay Buffer, 20 μ l of Glutathione, and 20 μ l of the reconstituted GST (control) to three wells.
 3. **Sample Wells** - add 150 μ l of Assay Buffer, 20 μ l of Glutathione, and 20 μ l of sample to three wells. To obtain reproducible results, the amount of GST added to the well should cause an absorbance increase between 0.012 and 0.064/min. When necessary, samples should be diluted with Sample Buffer or concentrated with an Amicon centrifuge concentrator with a molecular weight cut-off of 3,000 to bring the enzymatic activity to this level. *NOTE: The amount of sample added to the well should always be 20 μ l.*
 4. Initiate the reactions by adding 10 μ l of CDNB to all the wells you are using. Make sure to note the precise time you started and add the CDNB as quickly as possible.
 5. Carefully shake the 96-well plate for a few seconds to mix.
 6. Read the absorbance once every minute at 340 nm using a plate reader to obtain at least 5 time points.

Calculations

Determination of the Reaction Rate

1. Determine the change in absorbance (ΔA_{340}) per minute by:
 - a. Plotting the absorbance values as a function of time to obtain the slope (rate) of the linear portion of the curve (a graph is shown in Figure 1 using equine liver GST) -or-
 - b. Select two points on the linear portion of the curve and determine the change in absorbance during that time using the following equation:

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

2. Determine the rate of $\Delta A_{340}/\text{min.}$ for the background or non-enzymatic wells and subtract this rate from that of the sample wells.
3. Use the following formula to calculate the GST activity. The reaction rate at 340 nm can be determined using the CDNB extinction coefficient of $0.00503 \mu\text{M}^{-1}$. One unit of enzyme will conjugate 1.0 nmol of CDNB with reduced glutathione per minute at 25°C .

$$\text{GST Activity} = \frac{\Delta A_{340}/\text{min.}}{0.00503 \mu\text{M}^{-1}} \times \frac{0.2 \text{ ml}}{0.02 \text{ ml}} \times \text{Sample dilution} = \text{nmol/min/ml}$$

*The actual extinction coefficient for CDNB at 340 nm is $0.0096 \mu\text{M}^{-1}\text{cm}^{-1}$.⁶ This value has been adjusted for the pathlength of the solution in the well (0.524 cm).

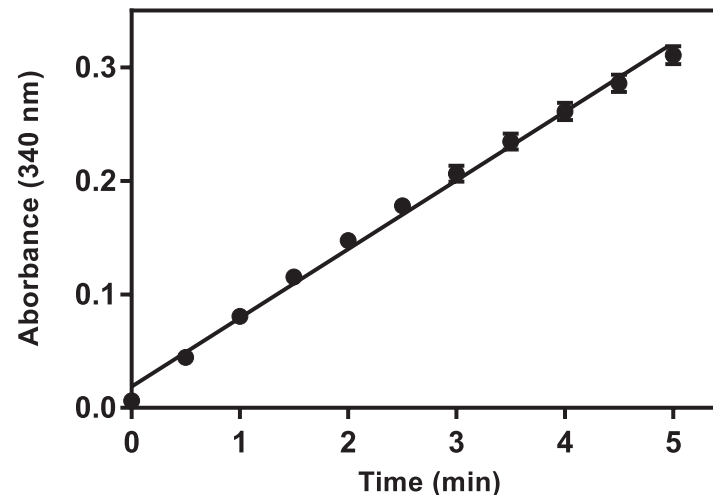


Figure 1. Activity of equine liver GST

Performance Characteristics

Precision:

When a series of eight GST samples were assayed on the same day, the intra-assay coefficient of variation was 3.6%. When a series of eight GST samples were assayed on five different days under the same experimental conditions, the inter-assay coefficient of variation was 4.8%.

Sensitivity:

The dynamic range of the assay is limited only by the accuracy of the absorbance measurement. Most plate readers are linear to an absorbance of 1.2. Samples containing GST activity between 24-128 nmol/min/ml can be assayed without further dilution or concentration. This GST activity is equivalent to an absorbance increase of 0.012 to 0.064 per minute.

Interferences and Inhibitors

- Samples that have a high intrinsic absorbance at 340 nm may exceed the absorbance maximum of the plate reader. Therefore, samples with an initial absorbance >0.7 should be diluted with Sample Buffer until the absorbance is lowered. For example, hemoglobin absorbs significantly at 340 nm, and thus erythrocyte lysates must be diluted before assaying.
- Inhibitors can be dissolved in dimethylsulfoxide or ethanol but no more than 10 µl should be added to the assay. The volume of Assay Buffer in the well should be adjusted so as to accommodate the inhibitor volume. The final volume in the wells should be 200 µl. The inhibitor should be added to the wells before the reaction is initiated with CDNB. The 100% Initial Activity well should contain GST and 10 µl of solvent (inhibitor solvent).

The following reagents were tested for interference in the assay.

| | Reagent | Will Interfere (Yes or No) |
|---------------------------------|-------------------------------|----------------------------|
| Buffers: | Tris | No |
| | Phosphate | No |
| Detergents: | Triton X-100 ($\leq 0.1\%$) | No |
| | Polysorbate 20 (0.1%) | Yes |
| | CHAPS (0.1%) | Yes |
| Protease Inhibitors/ Chelators: | Antipain (0.1 mg/ml) | Yes |
| | Leupeptin (10 µg/ml) | Yes |
| | Trypsin (10 µg/ml) | Yes |
| | Chymostatin (10 µg/ml) | Yes |
| | PMSF (≤ 200 µM) | No |
| | EGTA (≤ 5 mM) | No |
| | EDTA (≤ 2 mM) | No |
| Solvents: | Ethanol (10 µl) | No |
| | Methanol (10 µl) | Yes |
| | Dimethyl Sulfoxide (10 µl) | No |
| Others: | Glycerol ($\leq 10\%$) | No |
| | BSA ($\leq 1\%$) | No |

Troubleshooting

| Problem | Possible Causes | Recommended Solutions |
|--|---|---|
| Erratic values; dispersion of duplicates/triplicates | A. Poor pipetting/technique B. Bubble in the well(s) | A. Be careful not to splash the contents of the wells B. Carefully tap the side of the plate with your finger to remove bubbles |
| The initial absorbance in the wells is less than 0.1 | CDNB was not added to the wells | Make sure to add all components to the wells and re-assay |
| No increase in absorbance was observed in the sample wells | Enzyme activity was too low or glutathione was not added to the wells | Concentrate your sample by using an Amicon centrifuge concentrator with a 3,000 MW cut-off and re-assay; make sure to add components to the wells |
| Reaction rate was too fast | Too much enzyme added to well(s) | Dilute your samples with Sample Buffer and re-assay |
| The initial absorbance in the sample wells is above 0.7 | | Dilute your samples with Sample Buffer and re-assay |

References

1. Boyland, E. and Chasseaud, L.F. The role of glutathione and glutathione S-transferases in mercapturic acid biosynthesis. *Adv. Enzymol.* **32**, 173-219 (1969).
2. Mannervik, B. The isoenzymes of glutathione transferase. *Adv. Enzymol. Relat. Areas Mol. Biol.* **57**, 357-417 (1985).
3. Mannervik, B., Ålin, P., Guthenberg, C., *et al.* Identification of three classes of cytosolic glutathione transferase common to several mammalian species: correlation between structural data and enzymatic properties. *Proc. Natl. Acad. Sci. USA* **82**, 7202-7206 (1985).
4. Morgenstern, R., DePierre, J.W., and Ernster, L. Activation of microsomal glutathione S-transferase activity by sulfhydryl reagents. *Biochem. Biophys. Res. Commun.* **87**, 657-663 (1979).
5. Jakoby, W.B. The glutathione S-transferases: A group of multifunctional detoxification proteins. *Adv. Enzymol. Relat. Areas Mol. Biol.* **46**, 383-414 (1978).
6. Habig, W.H., Pabst, M.J., and Jakoby, W.B. Glutathione S-transferases. The first enzymatic step in mercapturic acid formation. *J. Biol. Chem.* **249**, 7130-7139 (1974).

| | | | | | | | | |
|----|---|---|---|---|---|---|---|---|
| 12 | | | | | | | | |
| 11 | | | | | | | | |
| 10 | | | | | | | | |
| 9 | | | | | | | | |
| 8 | | | | | | | | |
| 7 | | | | | | | | |
| 6 | | | | | | | | |
| 5 | | | | | | | | |
| 4 | | | | | | | | |
| 3 | | | | | | | | |
| 2 | | | | | | | | |
| 1 | | | | | | | | |
| | A | B | C | D | E | F | G | H |

Warranty and Limitation of Remedy

Buyer agrees to purchase the material subject to Cayman’s Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

This document is copyrighted. All rights are reserved. This document may not, in whole or part, be copied, photocopied, reproduced, translated, or reduced to any electronic medium or machine-readable form without prior consent, in writing, from Cayman Chemical Company.

©12/07/2017, Cayman Chemical Company, Ann Arbor, MI, All rights reserved. Printed in U.S.A.

