



Glutathione Peroxidase Activity Kit

For use with mammalian cells, tissues, erythrocytes,
and plasma.

Instruction Manual ADI-900-158

Sufficient Reagents for 480 tests in 5 x 96-well plates



Product Manual

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NONINFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Product Manual

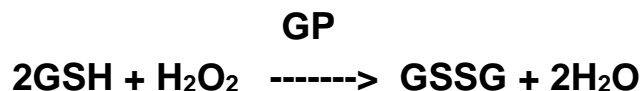
TABLE OF CONTENTS

| | |
|-----------------------------|----|
| Introduction..... | 4 |
| Materials Supplied | 5 |
| Storage..... | 6 |
| Other Materials Needed..... | 6 |
| Reagent Preparation | 7 |
| Sample Handling | 8 |
| Assay Procedure | 11 |
| Trouble Shooting | 15 |
| References | 16 |
| Contact Information | 20 |

INTRODUCTION

The Glutathione peroxidase activity kit is a complete kit for measuring the activity of glutathione-dependent peroxidases in cells, tissues, erythrocytes and plasma.

Glutathione Peroxidase is a tetramer of four identical subunits, with a molecular weight of 84,000 daltons. It requires selenium as a cofactor and contains a selenocysteine amino acid residue in the active site of each monomer that participates in the actual mechanism of the enzyme. Glutathione peroxidase (GP) is found in mammalian cells and helps to prevent lipid peroxidation of cell membranes by consuming free peroxide in the cell. The enzyme catalyzes the following reaction:



Glutathione Reductase (GR) then reduces the oxidized glutathione to complete the cycle:



Where GSH represents reduced monomeric glutathione, and GSSG represents oxidized glutathione. The oxidation of NADPH to NADP⁺ is accompanied by a decrease in absorbance at 340 nm (A₃₄₀). The rate of decrease in the A₃₄₀ is directly proportional to the Glutathione Peroxidase activity in the sample.

PRINCIPLE

1. Reaction mix, Glutathione Peroxidase, and samples are added to the appropriate wells of a 96-well plate.
2. Cumene Hydroperoxide is added to the wells to initiate the reaction.
3. The plate is transferred to a plate reader and absorbance readings are taken at 340 nm every 30 seconds or 1 minute for 10 to 15 minutes.

MATERIALS SUPPLIED

1. **Clear Microtiter Plate**

Five Plates of 96 Wells, Catalog No. 80-1639
Clear uncoated solid plates.

2. **10X Assay Buffer**

20 mL, Catalog No. 80-1662

3. **GSH + NADPH**

10 vials, Catalog No. 80-1661

4. **Glutathione Reductase**

1.1 mL, Catalog No. 80-1665

5. **Glutathione Peroxidase**

800 μ L, Catalog No. 80-1664

Glutathione Peroxidase in a 50% glycerol storage buffer

6. **Cumene Hydroperoxide**

12 mL, Catalog No. 80-1663

STORAGE

All components of this kit are stable at -20°C. All kit components are stable at their recommended storage temperatures until the kit expiration date.

OTHER MATERIALS NEEDED

1. PBS pH 7.4
2. Distilled water
3. Protease inhibitors such as phenylmethylsulfonyl fluoride (PMSF), Sigma P7626 or equivalent
4. Peroxide free Triton X-100 or Nonidet P-40 for cell extract preparation.
5. Microtubes, 0.5 and 1.5 mL
6. 15 mL conical tubes (adherent and suspension cell preparation)
7. 50 mL conical tubes (tissue preparation)
8. Precision pipettes for volumes between 1-200 μL and 100 to 1000 μL
9. Multichannel pipettor for volumes between 1 - 50 μL and 50 μL – 200 μL
10. Microplate reader or spectrophotometer capable of reading at 340 nm and taking readings every 30 seconds or 1 minute for ten to 15 minutes and exporting data to an Excel spreadsheet.
11. Microcentrifuge for processing samples
12. Sonicator or Homogenizer

REAGENT PREPARATION

1. Assay Buffer

Dilute the 10X Assay Buffer to 1X (1:10) with distilled water. The 1X Assay Buffer is stable at 4°C for up to 2 weeks. Note that 1X Assay Buffer contains 0.1 mg/mL BSA and this should be taken into account when calculating the protein concentration of your samples.

2. GSH + NADPH

Each vial of lyophilized GSH + NADPH contains sufficient reagent for 50 wells. Bring the number of vials required to room temperature and add 110 µL of 1X Assay Buffer to each vial. Swirl contents gently to dissolve and store on ice. The reconstituted reagent should be used within 4 hours. Discard any unused portion after use. Do not refreeze.

3. Reaction Mix

Make a 10X Reaction Mix (sufficient for 50 wells) as follows:

| | |
|---------------------------|--------|
| Glutathione Reductase | 110 µL |
| Reconstituted GSH + NADPH | 110 µL |
| 1X Assay Buffer | 880 µL |

Mix well and store on ice until added to the wells. Reaction Mix should be used within 4 hours of preparation.

4. Cumene Hydroperoxide

Thaw the cumene hydroperoxide and bring to room

temperature. Aliquot into 1.1 mL portions and freeze at -20°C. Each aliquot is sufficient for 50 wells. Note that the Cumene Hydroperoxide should be used within 1 hour at room temperature. Store on ice if in use for longer periods of time.

SAMPLE HANDLING

A. Preparation of Cell Extracts

1. **Non-adherent cells:** Centrifuge 2×10^6 to 1×10^7 non-adherent cells at 400 x g for 10 minutes at 4°C. Discard the supernatant. Suspend the cell pellet in 1 mL of ice-cold 1X PBS and transfer to a pre-chilled 1.5 mL microtube. Centrifuge at 10,000 x g for 10 seconds at 4°C. Discard the supernatant. Suspend the cell pellet in 5 pellet volumes of cold 1X Assay Buffer containing 0.4 mM PMSF, other protease inhibitors, and 1% Triton X-100 (peroxide-free).

2. **Adherent cells:** Wash the adherent cells with 1X PBS. Adherent cells may be harvested by scraping in 5 mL of ice-cold 1X PBS. Transfer to a pre chilled 15 mL tube. Centrifuge at 400 x g for 10 minutes at 4°C and discard the supernatant. Resuspend the cell pellet in 1 mL ice-cold 1X PBS and transfer to a pre chilled 1.5 mL microtube. Centrifuge at 10,000 x g for 12 seconds at 4°C. Discard the supernatant. Resuspend the cell pellet in 5 pellet volumes of ice-cold 1X Assay Buffer containing 0.4 mM PMSF, other protease inhibitors, and 1% Triton X-100 (peroxide-free).

3. Incubate the cell suspensions on ice, with periodic vortexing, for 30 minutes.

4. Microcentrifuge the disrupted cell suspension at 10,000 x g for 10-20 minutes at 4°C to remove insoluble material. Recover the supernatant to a fresh tube pre-chilled on ice.

5. Determine the protein concentration of the cleared cell lysate. Snap-freeze the cleared cell extract in small aliquots and store at -80°C . Avoid repeated freezing and thawing of the extract. The frozen cell extracts will be stable for 1 month.

B. Preparation of Plasma

1. Collect blood using an anticoagulant such as heparin, citrate, or EDTA.

2. Centrifuge at $1,500 \times g$ for 10 minutes at 4°C . Pipet off the upper yellow plasma layer without disturbing the white buffer layer (the white interface between the pelleted red blood cells and the plasma).

3. Store the plasma on ice for up to 4 hours, or freeze in small aliquots at -80°C . The frozen samples will be stable for at least 1 month.

C. Preparation of Red Blood Cell Lysates

1. Follow the directions for preparing plasma, above. Remove the buffy coat and discard.

2. Wash the red blood cells with 1X PBS at 4°C . Centrifuge at $1500 \times g$ for 10 minutes at 4°C . Discard the supernatant. Repeat this step once more.

3. Lyse the red blood cells in 4-10 volumes of 4°C deionized water by repeated gentle vortexing or mixing over a 10 minute period.

4. Centrifuge at $10,000 \times g$ for 15 minutes at 4°C . Collect the supernatant.

5. Determine the protein concentration of the erythrocyte ly-

sate. Store on ice for up to 4 hours, or freeze in small aliquots at -80°C . The frozen samples will be stable for at least one month.

D. Preparation of Tissue Homogenates

1. Prior to dissection, perfuse the tissue with 1X PBS plus 0.16 mg/mL heparin to remove blood components and clots.

2. Homogenize the tissue in 5-10 mL per gram of tissue of cold 1X Assay Buffer containing 0.4 mM PMSF, other protease inhibitors, and 1% Triton X-100 (peroxide-free).

3. Centrifuge at 10,000 x g for 10-20 minutes at 4°C to remove insoluble material. Recover the supernatant to a fresh tube prechilled on ice.

4. Determine the protein concentration of the cleared tissue lysate.

5. Alternative protocol: Mince the tissue in cold 1X PBS and make a single cell suspension by forcing the tissue through a stainless steel wire mesh screen using a pestle. Centrifuge the single cell suspension at 1,000 x g for 10 minutes at 4°C . Discard the supernatant. Resuspend the pellet in 5 pellet volumes of ice-cold 1X Assay Buffer containing 0.4 mM PMSF, other protease inhibitors, and 1% Triton X-100 (peroxide-free).

6. Incubate the cell suspensions on ice, with periodic vortexing, for 30 minutes.

7. Microcentrifuge the disrupted tissue single cell suspension at 10,000 x g for 10-20 minutes at 4°C to remove insoluble material. Recover the supernatant to a fresh tube prechilled on ice (If supernatant is cloudy, repeat centrifugation).

8. Determine the protein concentration of the cleared tissue lysate.

9. Store the clarified tissue extracts on ice and assay immediately, or snap- freeze the cleared tissue extract in small aliquots and store at -80°C . Avoid repeated freezing and thawing of the extract. The frozen cell extracts will be stable for 1 month.

ASSAY PROCEDURE

1. Set up your plate reader to measure absorbance at 340 nm every 30 seconds or 1 minute. Include a 10 second orbital shake prior to the initial read. Accuracy and consistency of results is dependent on maintaining a constant temperature. Set the plate chamber of your instrument to 25°C if possible. Blank your readings with respect to 200 μL of 1X Assay Buffer in one well.
2. Plan your experiment to measure each sample in triplicate and at different dilutions or amounts of protein. Some samples, particularly erythrocyte lysates, may have to be diluted significantly with 1X Assay Buffer to achieve a reasonable rate of decrease in absorbance at 340 nm. Note that the final volume of the reaction is 200 μL .
3. Set up the following reactions in a 96 well plate (per well):

| | |
|--|-------------------|
| 1X Assay Buffer: | 140 μL |
| 10X Reaction Mix: | 20 μL |
| Glutathione Peroxidase, sample, or control | 20 μL |

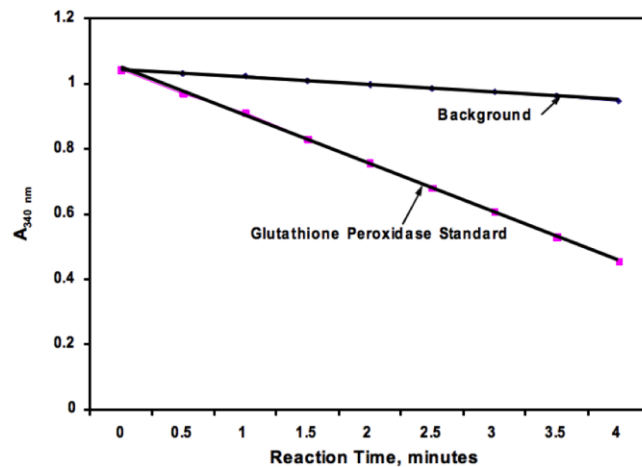
Note: The Glutathione Peroxide is provided to serve as a positive control to ensure the assay is working. It should not be used to construct a calibration curve to measure Glutathione Peroxidase concentration in samples.

4. Initiate the reactions by quickly adding 20 μ L of Cumene Hydroperoxide to each well using a multichannel pipettor.
5. Immediately begin measuring absorbance at 340 nm every 30 seconds or 1 minute over a 10-15 minute period.
6. **Controls:** Include a background set of wells where 20 μ L of 1X Assay Buffer is added instead of sample or Glutathione Peroxidase. The rate of decrease of absorbance at 340 nm in the background is subtracted from that of the samples or standard to obtain the net rate of decrease of absorbance at 340 nm for the calculation of Glutathione Peroxidase activity in your samples.

Calculation of Results

A. Determination of Glutathione Peroxidase Activity.

1. Calculate the mean absorbance at each time point of the triplicate values for the samples, Glutathione Peroxidase Standard, and background.
2. Plot the mean absorbance versus time. A representative standard curve is shown in Figure 1:



3. Determine the slope from a linear portion of the curve for the sample, Glutathione Peroxidase standard, and background. Express the results as the change in absorbance per minute ($\Delta A_{340}/\text{min}$).
4. Subtract the $\Delta A_{340}/\text{min}$ for the background from that of the samples.
5. One Unit of Glutathione Peroxidase is defined as the amount of enzyme that will cause the oxidation of 1 nmole of NADPH to NADP^+ per minute at 25°C . The reaction rate can be calculated knowing the extinction coefficient of NADPH which is $0.00622 \mu\text{M}^{-1}\text{cm}^{-1}$. Since the path length



Product Manual

of the samples in the wells is 0.61 cm, the extinction coefficient is modified to $0.00379 \mu\text{M}^{-1}$. Calculate the Glutathione Peroxidase activity in your samples from the following equations

(Where Y is the volume of your samples):

Glutathione Peroxidase Activity =

$$\frac{\Delta A_{340}/\text{min}}{0.00379 \mu\text{M}^{-1}} \times \frac{0.2 \text{ mL}}{Y \text{ mL}} \times \text{Sample Dilution}$$

$$= \text{nmole}/\text{min}/\text{mL} = \text{Units}/\text{mL}$$

TROUBLE SHOOTING

| Problem | Cause | Solution |
|--|--|---|
| Erratic values. Poor reproducibility of triplicates | Poor pipetting technique | Depress plunger of your pipettor all the way. Fill with solution. Depress to first stop. Fill with solution. Depress to first stop. Repeat as often as necessary. |
| | Bubbles in well | Avoid making bubbles |
| No decrease in absorbance observed in sample wells or Glutathione Peroxidase standard. | Enzyme activity in the samples is too low. | Increase volume and/or concentration of your samples. |
| | Failure to add cumene hydroperoxide | Add cumene hydroperoxide to wells. |
| Initial absorbance is below 0.5 or the absorbance decreases very rapidly | Glutathione Peroxidase activity is too high in your samples. | Dilute your samples with 1X Assay Buffer and re-assay. |
| Initial absorbance is less than 0.1 and there is no decrease in absorbance | Failure to add GSH + NADPH to 10X Reaction Mix | Add GSH + NADPH to 10X Reaction Mix |

| Problem | Cause | Solution |
|--|----------------------------|---|
| Initial absorbance in sample wells is greater than 1.5 | Sample is too concentrated | Dilute your sample with 1X Assay Buffer and reassay |

REFERENCES

1. Ozdemir G, Ozden M, Maral H, Kuskay S, Cetinalp P, Tarkun I. Malondialdehyde, glutathione, glutathione peroxidase and homocysteine levels in type 2 diabetic patients with and without microalbuminuria. *Ann Clin Biochem.* 2005, 42:99-104.
2. Prasad NR, Menon VP, Vasudev V, Pugalendi KV. Radioprotective effect of sesamol on gamma-radiation induced DNA damage, lipid peroxidation and antioxidants levels in cultured human lymphocytes. *Toxicology.* 2005, 209:225-35.
3. Sindhu RK, Ehdai A, Farmand F, Dhaliwal KK, Nguyen T, Zhan CD, Roberts CK, Vaziri ND. Expression of catalase and glutathione peroxidase in renal insufficiency. *Biochim Biophys Acta.* 2005, 1743:86-92.
4. Szelestei T, Bahring S, Wagner Z, Aydin A, Molnar GA, Kocsis B, Nagy J, Wittmann I. Serum levels of L-arginine analogues and glutathione peroxidase and catalase gene variants in Type 2 diabetes mellitus patients.
5. Chu FF, Esworthy JH, Doroshov JH. Role of Se-dependent glutathione peroxidases in gastrointestinal inflammation and cancer. *Free Rad Biol Med.* 2004. 36:1481-95.



Product Manual

NOTES:



Product Manual

NOTES:



Product Manual

NOTES:



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com