

# **TREVIGEN® Instructions**

*For Research Use Only. Not For Use In Diagnostic Procedures*

---

## **Glutathione Reductase Assay Kit**

**Reagent kit for the analysis of Glutathione Reductase in  
cell extracts and tissue homogenates**

**Sufficient reagents for 100 Assays**

**Catalog# 7510-100-K**

# Glutathione Reductase Assay Kit

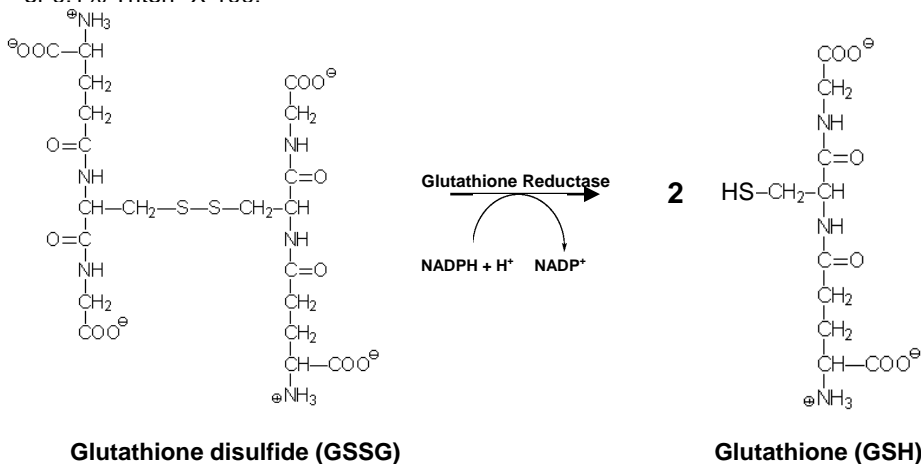
## Catalog# 7510-100-K

<b><u>Table of Contents</u></b>	<b><u>Page</u></b>
<b>I. Background</b>	<b>1</b>
<b>II. Precautions and Limitations</b>	<b>1</b>
<b>III. Materials Supplied</b>	<b>2</b>
<b>IV. Reagents/Equipment Required But Not Supplied</b>	<b>2</b>
<b>V. Reagent Preparation</b>	<b>2</b>
<b>VI. Sample Preparation</b>	<b>2</b>
<b>VII. Assay Protocol</b>	<b>3</b>
<b>VIII. Glutathione Reductase Activity Determination</b>	<b>6</b>
<b>IX. References</b>	<b>7</b>
<b>X. Related Products Available from Trevigen</b>	<b>7</b>

## I. Background

Glutathione reductase, a homodimeric flavoprotein disulfide oxidoreductase, plays an indirect but essential role in the prevention of oxidative damage within the cell by helping to maintain appropriate levels of intracellular glutathione (GSH). Glutathione reductase together with its co-factor, NADPH, catalyzes the reduction of oxidized glutathione (glutathione disulfide, GSSG) to glutathione (Figure 1). Glutathione is a water phase sulfhydryl antioxidant and antitoxin. GSH is responsible for converting hydrogen peroxide (H<sub>2</sub>O<sub>2</sub>) into water within erythrocytes (with glutathione peroxidase acting as the catalyst).

The method employed in Trevigen's Glutathione Reductase Assay Kit is a spectrophotometric assay, wherein the oxidation of NADPH to NADP<sup>+</sup> (Figure 1) is monitored as a decrease in absorbance at 340 nm. This rate of decrease in A<sub>340</sub> is directly proportional to the glutathione reductase activity in the sample because the enzyme is present at rate limiting concentrations. The unit definition for glutathione reductase activity may be expressed in terms of the oxidation of NADPH or the reduction of GSSG since their molar ratio is 1:1. One unit of glutathione reductase oxidizes 1 μmol of NADPH per minute at 25 °C, pH 7.5. Glutathione reductase activity is slightly inhibited by ammonium sulfate at concentrations greater than 100 mM, and is unaffected by 0.5 M urea, 0.1 M KCl, or 0.1% Triton<sup>®</sup> X-100.



**Figure 1.** Reduction of glutathione disulfide (GSSG) by glutathione reductase and NADPH.

## II. Precautions and Limitations

1. For Research Use Only. Not for use in diagnostic procedures.
2. The physical, chemical and toxicological properties of the provided products may not yet have been fully investigated, therefore, Trevigen recommends the use of gloves, lab coats, and eye protection while using these chemical reagents. Trevigen assumes no liability for damage resulting from handling or contact with these products. MSDS are available on request.

### III. Materials Supplied

<u>Component</u>	<u>Quantity</u>	<u>Storage</u>	<u>Catalog#</u>
Glutathione Reductase*	500 µl	4 °C	7510-100-01
25X Reaction Buffer	20 ml	4 °C	7510-100-02
NADPH**	10 tubes	-20 °C	7510-100-03
GSSG Solution	6 ml	4 °C	7510-100-04
25X Sample Dilution Buffer	20 ml	4 °C	7510-100-05
25X Tissue Homogenization Buffer	40 ml	4 °C	7510-100-06

\* 0.5 U/ml in 3.6 M (NH<sub>4</sub>)<sub>2</sub>SO<sub>4</sub>, 50 mM potassium phosphate, pH 7.5

\*\* sufficient amount of reagent for 10 assays and one 5 point standard curve (each point in duplicate) per tube using 1 ml assay volumes.

### IV. Reagents/Equipment Required but not Supplied

#### Equipment

1. Spectrophotometer to read absorbance at 340 nm, preferably with a temperature-regulated cuvette holder set at 25°C.
2. Quartz cuvettes with 1 cm path length and at least a 1 ml volume
3. Pipettor
4. Pipette tips
5. Pasteur pipettes and bulb
6. Centrifuge (for cell lysis)
7. Dounce Homogenizer
8. Timer
9. Vortex

#### Reagents

1. Deionized H<sub>2</sub>O (dH<sub>2</sub>O)
2. 1 X PBS

### V. Reagent Preparation

1. NADPH: Add 525 µL dH<sub>2</sub>O to one vial of NADPH and vortex to dissolve. May be kept for up to 24 hours at 0°C - 4°C protected from light.
2. 1 X Tissue Homogenization Buffer: Prepare necessary volume from the 25 X stock solution as needed with cold deionized water.

### VI. Sample Preparation

**Note:** The final concentration of protein (from the clarified lysate/ homogenate) in each sample should be adjusted to 0.1 - 0.75 mg/ mL initially. Samples may be further diluted with 1 X Sample Dilution Buffer if necessary.

#### A. Cell Lysate Preparation

1. Detach adherent cells using a rubber policeman. Centrifuge at 250 x g for 10 minutes at 4°C. Wash the cells once with cold 1 X PBS.

2. Suspend the cell pellet in cold 1 X Tissue Homogenization Buffer. Homogenize the suspension using a Dounce homogenizer or a sonicator.
3. Centrifuge at 8,500 x g for 10 minutes at 4°C. Adjust protein concentration as needed (see note pg. 2). Transfer the supernatant to a clean tube and place on ice (if performing the assay the same day), or store at -80°C for future use.

### **B. Erythrocyte Lysate Preparation**

1. Centrifuge blood at 2,500 x g for 5 minutes at 4°C. (discard supernatant)
2. Wash the red blood cell pellet twice with 1 X PBS.
3. Lyse the erythrocyte pellet in 4 pellet volumes of cold deionized H<sub>2</sub>O. Mix well by pipetting.
4. Centrifuge the suspension at 8,500 x g for 10 minutes at 4°C. Adjust protein concentration as needed (see note pg. 2). Transfer the supernatant to a clean tube and place on ice (if performing the assay the same day) or store at -80°C for future use.

### **C. Tissue Homogenate Preparation\***

1. Wash tissue 3 times with 1X PBS and then homogenize tissue in 5-10 mL of 1X Tissue Homogenization Buffer per gram of tissue.
2. Centrifuge the suspension at 8,500 x g for 10 minutes at 4°C. Adjust protein concentration as needed (see note pg. 2). Transfer the supernatant to a clean tube and place on ice (if performing the assay the same day), or store at -80°C for future use.

## **VII. Assay Protocol**

The assay is performed at room temperature and in duplicate. Keep cell lysates, tissue homogenates, and the glutathione reductase solution on ice; all other reagents should be brought to room temperature before use.

**Note: immediately prior to use, briefly vortex each reagent. Perform each reaction separately.**

### **A. Blank**

Include a blank to account for the endogenous oxidation of NADPH over the period of the reaction.

1. Add the following components in order to a quartz cuvette:

	Blank
dH <sub>2</sub> O	930 µL
25X Reaction Buffer	40 µL
GSSG Solution	20 µL

2. Mix thoroughly by repeated pipetting with a clean Pasteur pipette.

3. Place each cuvette individually into a spectrophotometer set at 340 nm and 25°C. Adjust the absorbance reading to zero.
4. Add 10 µL of the NADPH solution prepared in Section V to the cuvette, mix thoroughly by repeated pipetting with a clean Pasteur pipette, and start a timer or a stopwatch. Record the absorbance every 60 seconds for 5 minutes.

### B. Sample

1. Add the following components in order to a quartz cuvette:

	Sample
dH <sub>2</sub> O	930 µL - X µL
25X Reaction Buffer	40 µL
Sample	X µL
GSSG Solution	20 µL

2. Mix thoroughly by repeated pipetting with a clean Pasteur pipette.
3. Place each cuvette individually into a spectrophotometer set at 340 nm and 25°C. Adjust the absorbance reading to zero.
4. Add 10 µL of the NADPH solution prepared in Section V to the cuvette, mix thoroughly by repeated pipetting with a clean Pasteur pipette, and start a timer or a stopwatch. Record the absorbance every 60 seconds for 5 minutes.

### C. Standard Curve

The kit contains sufficient Glutathione Reductase to generate ten 5-point standard curves with each point done in duplicate. The Glutathione Reductase has an activity of 0.5 U/mL (1 unit will cause the oxidation of 1 µmole of NADPH in one minute at 25°C at pH 7.5). The concentrations of Glutathione Reductase (GR) to be used are: 1, 3, 5, 7, and 10 mU/mL.

1. In the table below are the reagent volumes required for each of the five standards. Follow the procedure for preparing these solutions starting with step C-2 below:

	1	2	3	4	5
dH <sub>2</sub> O	928 µL	924 µL	920 µL	916 µL	910 µL
25X Reaction Buffer	40 µL	40 µL	40 µL	40 µL	40 µL
GSSG Solution	20 µL	20 µL	20 µL	20 µL	20 µL
GR solution	2 µL	6 µL	10 µL	14 µL	20 µL
NADPH Solution	10 µL	10 µL	10 µL	10 µL	10 µL
<i>GR concentration (mU/mL)</i>	1	3	5	7	10

2. To a quartz cuvette add the required amounts of dH<sub>2</sub>O, 25X Reaction buffer, and GSSG solution in that order.
3. Mix thoroughly by repeated pipetting with a clean Pasteur pipette.

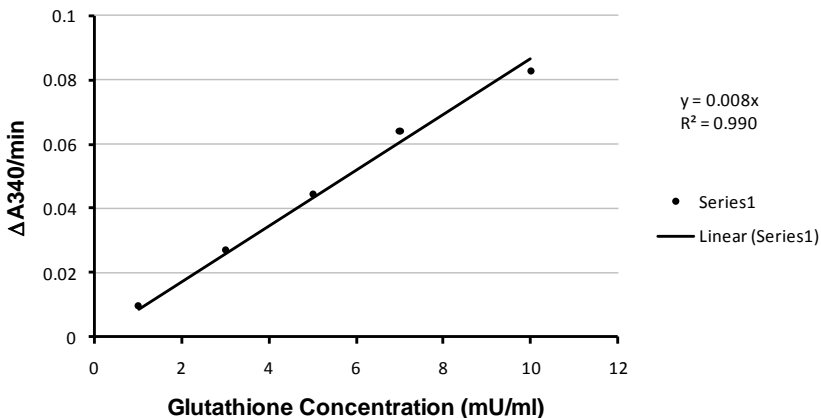
4. Add the required amount of glutathione reductase solution according to the Table in step C-1 and mix thoroughly by repeated pipetting with a clean Pasteur pipette.
5. Place the cuvette into a spectrophotometer set at 340 nm and adjust the absorbance reading to zero.
6. Add 10  $\mu$ L of the NADPH solution prepared in Section V to the cuvette, mix thoroughly by repeated pipetting with a clean Pasteur pipette, and start a timer or a stopwatch.
7. Record the absorbance every 60 seconds for 5 minutes.

**D. Standard Curve Calculations**

1. Determine the rate of decrease in absorbance @ 340nm per minute ( $\Delta A_{340}$  /min) for each of the standards and the blank:

$$\frac{(A_{340} \text{ @ 1 min.}) - (A_{340} \text{ @ 5 min.})}{4 \text{ min}} = \Delta A_{340}/\text{min}$$

2. Determine the average of the duplicate  $\Delta A_{340}$  /min values for each standard and blank.
3. Calculate the net rate for each standard by subtracting the rate obtained for the blank.
4. Plot GR concentration versus these new  $\Delta A_{340}$ / min. values.



**Figure 2.** Plot of glutathione reductase concentration vs. the rate of decrease in absorbance at 340 nm due to the oxidation of NADPH by GR.

### VIII. Glutathione Reductase Activity Determination

The GR activity in the sample(s) may be calculated by one of two methods.

**Method 1:** Enter average A<sub>340</sub> for standards and samples into calculation worksheet (MS Excel required) provided on Trevigen website [http://www.trevigen.com/docs/1427816753.7510-100-k\\_calculation\\_worksheet.xls](http://www.trevigen.com/docs/1427816753.7510-100-k_calculation_worksheet.xls)

**Method 2:**

1. Determine the rate of decrease in absorbance per minute ( $\Delta A_{340}/\text{min}$ ) for both samples and blanks.
2. Calculate the net rate by subtracting the blank rate from the sample rate.
3. Convert the net rate ( $\Delta A_{340}/\text{min}$ ) to concentration of NADPH consumed, which is equal to the activity of Glutathione Reductase in mU/mL. The molar extinction coefficient ( $E^M$ ) for NADPH is  $6220 \text{ M}^{-1}\text{cm}^{-1}$  and  $E^M = 6.22 \times 10^{-3} \text{ nmol/mL}$  if the pathlength is 1 cm. One unit of glutathione reductase is defined as the amount of enzyme required to catalyze the reduction of one micromole of GSSG per minute at pH 7.5 and 25°C. One molecule of NADPH is consumed per molecule of GSSG reduced. Therefore, the reduction of NADPH (measured by loss of A<sub>340</sub>) directly correlates with GSSG reduction.

1 U of Glutathione Reductase = 1  $\mu\text{mol}$  GSSG/min = 1  $\mu\text{mol}$  NADPH/min  
 or 1 mU of Glutathione Reductase = 1 nmole NADPH consumed/min

If  $E^M$  is the molar extinction coefficient for NADPH at 340 nm,

$$E^M = 6220 \text{ M}^{-1}\text{cm}^{-1} = 6.22 \times 10^{-3} \mu\text{M}^{-1}\text{cm}^{-1} = 6.22 \times 10^{-3} \text{ L}/\mu\text{mol}/\text{cm}$$

$$= 6.22 \times 10^{-3} \text{ mL}/\text{nmol} \text{ if the pathlength is 1 cm}$$

$$\frac{\Delta A_{340} / \text{min}}{(6.22 \times 10^{-3} \mu\text{M}^{-1}\text{cm}^{-1})} = \frac{\Delta A_{340} / \text{min}}{6.22 \times 10^{-3} \text{ mL}/\text{nmol}} = x \text{ nmol NADPH}/\text{mL}$$

$$= x \text{ mU}/\text{mL Glutathione Reductase}$$

4. Correct for the sample dilution in the assay and for the sample dilution performed prior to the assay.

For Example:

If the sample volume was 200  $\mu\text{L}$  and was diluted 1/50 prior to the assay,

Mean Sample Rate, $\Delta A_{340}/\text{min}$	= 0.0325 $\Delta A_{340}/\text{min}$
Mean Blank Rate, $\Delta A_{340}/\text{min}$	= 0.0005 $\Delta A_{340}/\text{min}$
Net Rate, $\Delta A_{340}/\text{min}$	= 0.0320 $\Delta A_{340}/\text{min}$ (0.0325 - 0.0005)
Glutathione Reductase Activity	= 0.0320/6.22 $\times 10^{-3}$ = 5.14 mU/mL
Assay Dilution Correction	= 1000 $\mu\text{L}/200 \mu\text{L} \times 5.14$ = 25.7 mU/ml
Sample Dilution Correction	= 50 $\times 25.7$ = 1285 mU/ml



## IX. References

1. Tietze, F., Enzymic method for quantitative determination of nanogram amounts of total and oxidized glutathione: applications to mammalian blood and other tissues. *Anal. Biochem.*, **27**, 502-522 (1969).
2. Carlberg, I., Mannervik, B., Glutathione reductase. *Meth. Enzymol.*, **113**, 485-490 (1985).
3. Smith IK, Vierheller TL, Thorne CA, Assay of glutathione reductase in crude tissue homogenates using 5,5'-dithiobis(2-nitrobenzoic acid). *Anal. Biochem.*, **175**, 408-413 (1988).
4. Dolphin, D., Poulson, R., and Avramovic, O. (Eds.), *Glutathione: Chemical, Biochemical, and Metabolic Aspects*, Volumes A and B, Wiley and Sons (1989).
5. Dringen R, Gutterer JM. Glutathione reductase from bovine brain. *Methods Enzymol.* **348**, 281-288 (2002).

## X. Related Products Available From Trevigen

Contact Trevigen for details of our unique product line for studying DNA damage and repair. All of Trevigen's kits include highly qualified enzymes, substrates, buffers, full instructions for use, and a synopsis specific for your kit.

### PARP Assay Kits:

Catalog #	Description	Size
4690-096-K	HT F Homogeneous PARP Inhibition Assay kit	96 tests
4677-096-K	HT Universal Colorimetric PARP Assay w/ Histone Coated Strip Wells	96 samples
4676-096-K	Universal Chemiluminescent PARP Assay w/Histone Coated Strip Wells	96 samples
4684-096-K	HT Universal Colorimetric PARP/Apoptosis Assay	96 samples
4685-096-K	HT Universal Chemiluminescent PARP/Apoptosis	96 samples

### DNA Damage Antibodies:

Catalog #	Description	Size
4335-MC-100	Anti-PAR polymer mAb (10HA)	100 µl
4336-BPC-100	Anti- PAR polymer polyclonal	100 µl
4338-MC-50	Anti-human/murine-PARP mAb (clone C2-10)	50 µg
4354-MC-050	Anti-8-oxo-dG (clone 2E2)	50 µl
4360-MC-100	Anti-BPDE (clone 8E11)	100 µg
4418-APC-100	Anti-Phosphorylated Histone H2AX polyclonal	100 µl

### Glutathione Kits

Catalog #	Description	Size
7511-100-K	HT Glutathione Assay Kit	384 tests
7512-100-K	HT Glutathione Peroxidase Assay Kit	480 tests
7513-500-K	HT Glutathione Reductase Assay Kit	500 tests

### CometAssay® Kits:

Catalog #	Description	Size
4250-050-K	CometAssay® Kit	50 samples
4251-050-K	CometAssay® Silver Staining Kit	50 samples
4252-040-K	CometAssay® Higher Throughput Kit	40 samples
4253-096-K	CometAssay® Kit 96 Wells	96 samples

### FLARE™ Assay Kits:

Catalog #	Description	Damage Recognized	Size
4040-100-FK	Fpg Kit	8-oxoguanine, DNA containing formamidopyrimidine moieties	75 samples
4040-100-FM			100 samples
4045-01K-FK	Endonuclease III Kit	Thymine Glycol, 5,6-dihydrothymine, urea, 5-hydroxy-6-hydrothymine, 5,6-dihydro-uracil, alloxan, 5-hydroxy-6-hydrouracil, uracil glycol, 5-hydroxy-5-methylhydantoin, 5-hydroxycytosine,5-hydroxy-uracil, methyl-taronylurea, thymine ring saturated or fragmentation product	75 samples
4045-01K-FM			100 samples
4130-100-FK	hOGG1 Kit	8-oxoguanine, DNA containing formamidopyrimidine moieties	75 samples
4130-100-FM			100 samples

**The product accompanying this document is intended for research use only and is not intended for diagnostic purposes or for use in humans.**



Please  
Recycle

**Trevigen, Inc.**

8405 Helgerman Ct. Gaithersburg, MD 20877

Tel: 1-800-873-8443 • 301-216-2800

Fax: 301-560-4973

e-mail: [info@trevigen.com](mailto:info@trevigen.com)

[www.trevigen.com](http://www.trevigen.com)