

# **Thioredoxin Reductase Assay Kit**

Item No. 10007892

## TABLE OF CONTENTS

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

## GENERAL INFORMATION

### Materials Supplied

Item Number	Item	Quantity
10009092	TrxR Assay Buffer (10X)	1 vial
10009093	Thioredoxin Reductase Control	1 vial
10009094	TrxR Inhibitor	2 vials
10009096	TrxR DTNB	1 vial
10009095	TrxR NADPH	2 vials
10009097	TrxR DMSO	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) 975-3999. We cannot accept any returns without prior authorization.



**WARNING:** This product is for laboratory research use only; not for administration to humans. Not for human or veterinary diagnostic or therapeutic use.

## Precautions

Please read these instructions carefully before beginning this assay.

For research use only. Not for human or diagnostic use.

## 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 absorbances between 405-414 nm
2. Adjustable pipettes and a repeat pipettor
3. A source of pure water; glass distilled water or HPLC-grade water is acceptable

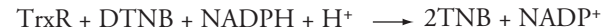
## INTRODUCTION

### Background

The thioredoxin system is one of the key cellular redox regulators. Together with the glutathione system, it controls the redox state of cysteine residues in proteins and has numerous other roles in redox regulation of cellular processes. The thioredoxin system is composed of thioredoxin reductase (TrxR) and thioredoxin. TrxRs are a family of selenium-containing pyridine nucleotide-disulfide oxidoreductases.<sup>1</sup> TrxR transfers electrons from NADPH to thioredoxin, which in turn reduces thioredoxin peroxidase, methionine sulfoxide reductase, ribonucleotide reductase, and other important redox proteins.<sup>1</sup> TrxRs are able to reduce a number of substrates other than thioredoxin, including selenite, lipid hydroperoxides, vitamin K, and hydrogen peroxide.<sup>1,2</sup> To date, two forms of mammalian TrxRs have been characterized. One is present in the cytosol (called TrxR1, TR1, or TxnRd1) and the other resides in the mitochondria (called TrxR2, TR3, or TxnRd2).<sup>3</sup> TrxRs have been implicated in playing a role in protecting against oxidative injury, cell growth and transformation, and the recycling of ascorbate.<sup>2,4</sup> Determining the involvement of TrxR in both normal and pathological cell function, as well as its drug interaction, may provide new insights into diseases and provide new treatments for cancer, AIDS, and autoimmune diseases.<sup>1,2</sup>

### About This Assay

Cayman's Thioredoxin Reductase Assay Kit provides a convenient method for detecting mammalian TrxR activity in tissue homogenates and cell lysates. It is based on the reduction of DTNB (5,5'-dithio-bis(2-dinitrobenzoic acid); Ellman's reagent) with NADPH to 5-thio-2-nitrobenzoic acid (TNB) which produces a yellow product that is measured at 405-414 nm.<sup>5,6</sup> The kit includes all reagents needed to assay mammalian TrxR activity. Measurement of TrxR activity by DTNB reduction in the absence and in the presence of aurothiomalate, a specific TrxR inhibitor included in the kit, allows for correction of non-thioredoxin reductase-independent DTNB reduction (*i.e.*, presence of glutathione).<sup>7</sup> The difference between the two results is the DTNB reduction due to TrxR activity.



## Reagent Preparation

### 1. TrxR Assay Buffer (10X) - (Item No. 10009092)

Dilute 3 ml of Assay Buffer concentrate with 27 ml of HPLC-grade water. This final Assay Buffer (50 mM potassium phosphate, pH 7.0, containing 50 mM KCl, 1 mM EDTA, and 0.2 mg/ml BSA) should be used in the assay and for reconstituting the TrxR Inhibitor and NADPH. When stored at -20°C, this diluted Assay Buffer is stable for at least six months.

### 2. Thioredoxin Reductase (control) - (Item No. 10009093)

The vial contains 100 µl of a solution of rat liver thioredoxin reductase (TrxR). The thawed enzyme should be stored on ice. The enzyme is ready to use as supplied.

### 3. TrxR Inhibitor - (Item No. 10009094)

The vial contains a lyophilized powder of sodium aurothiomalate (ATM). Reconstitute the vial with 1 ml of diluted Assay Buffer before use. The reconstituted Inhibitor is stable for four hours. ATM is a specific thioredoxin inhibitor. Since several enzymes present in biological samples can reduce DTNB, the ATM is used to determine the reduction of DTNB specific to thioredoxin reductase. The concentration of Inhibitor used in the assay, 20 µM, will effectively remove all thioredoxin reductase activity.<sup>8</sup>

### 4. TrxR DTNB - (Item No. 10009096)

Weigh 4 mg of DTNB into another vial, add 2 ml of dimethylsulfoxide (DMSO) and vortex until dissolved. Store the reagent at room temperature in the dark and use within four hours.

### 5. TrxR NADPH - (Item No. 10009095)

The vial contains a lyophilized powder of NADPH. Reconstitute the vial with 2 ml of diluted Assay Buffer and store on ice. The reconstituted NADPH is stable for six hours.

### 6. TrxR DMSO - (Item No. 10009097)

The vial contains 10 ml of DMSO. It is ready to use as supplied. Once thawed, DMSO can be stored at room temperature for six months.

## Sample Preparation

### Tissue Homogenate

The amount of TrxR activity in animal tissues varies from organ to organ. Values range from 0.05-0.6 units per mg of protein for crude extracts.

1. Prior to dissection, either perfuse tissue or rinse tissue with a phosphate buffered saline (PBS) 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.*, 50 mM potassium phosphate, pH 7.4, containing 1 mM EDTA) per gram of 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

Typically cell culture extracts have a range of 0.4-4 units per 10<sup>8</sup> cells (0.04-0.25 units per mg of protein)

1. Collect cells (1 x 10<sup>8</sup>) 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 cell pellet in 5-10 ml of cold buffer (*i.e.*, 50 mM potassium phosphate, pH 7.4, containing 1 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.

## ASSAY PROTOCOL

### Plate Set Up

There is no specific pattern for using the wells on the plate. However, it is necessary to have two wells designated as background wells, two wells designated as positive control wells, and two background wells treated with inhibitor. We recommend that each sample also be treated with inhibitor.

A typical layout of samples to be measured in duplicate is given below (see Figure 1). We suggest that you record the contents of each well on the template sheet provided on page 19.

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

B - Background  
B+I - Background + ATM  
C - Positive Control  
S - Sample  
S+I - Sample + ATM  
★ - Additional Samples

Figure 1. Sample plate format

### Pipetting Hints

- It is recommended that an adjustable pipette be used to deliver reagents to the wells.
- 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.

### General Information

- The final volume of the assay is 200  $\mu$ l in all the wells.
- It is not necessary to use all the wells on the plate at one time.
- The assay temperature is 22°C.
- We recommend assaying samples in the presence and absence of ATM. Since several enzymes present in biological samples can reduce DTNB, ATM is used to determine the reduction of DTNB due only to TrxR activity.
- We recommend assaying samples in triplicate, but it is the user's discretion.
- Thirty samples can be assayed in triplicate or forty-five in duplicate.
- Monitor the absorbance at 405-414 nm.

## Performing the Assay

1. **Background Wells** - add 160  $\mu\text{l}$  of diluted Assay Buffer to two wells.
2. **Background + ATM Wells** - add 140  $\mu\text{l}$  of diluted Assay Buffer and 20  $\mu\text{l}$  of ATM to two wells. *NOTE: ATM does slightly react with DTNB so it is important to subtract this activity from the sample + ATM wells.*
3. **Positive Control Wells (rat liver TrxR)** - add 140  $\mu\text{l}$  of diluted Assay Buffer and 20  $\mu\text{l}$  of rat liver TrxR (control) to two wells.
4. **Sample Wells** - add 140  $\mu\text{l}$  of diluted Assay Buffer and 20  $\mu\text{l}$  of sample to two wells.
5. **Sample + ATM Wells** - add 120  $\mu\text{l}$  of diluted Assay Buffer, 20  $\mu\text{l}$  of sample, and 20  $\mu\text{l}$  of ATM to two wells.
6. Initiate the reactions by adding 20  $\mu\text{l}$  of NADPH and 20  $\mu\text{l}$  of DTNB to all the wells being used. Carefully shake the microtiter plate for 10 seconds to mix.
7. Read the absorbance once every minute at 405-414 nm using a plate reader to obtain at least five time points.

## ANALYSIS

### Calculations

#### Determination of the Reaction Rate

1. Determine the change in absorbance ( $\Delta A_{405}$ ) per minute by either:
  - a. Plotting the average absorbance values as a function of time to obtain the slope (rate) of the linear portion of the curve (a graph is shown using rat liver TrxR, see Figure 2, page 12) - 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:

$$\Delta A_{405}/\text{min.} = \frac{A_{405}(\text{Time 2}) - A_{405}(\text{Time 1})}{\text{Time 2 (min.)} - \text{Time 1 (min.)}}$$

2. Determine the rate  $\Delta A_{405}/\text{min.}$  for the background and subtract this rate from all the wells, including the background + ATM, sample, sample + ATM, and positive control wells.
3. Use the following formulas to calculate the TrxR activity. The reaction rate at 405 or 414 nm can be determined using either the DTNB extinction coefficient of  $6.35 \text{ mM}^{-1}$  (405 nm) or  $6.75 \text{ mM}^{-1}$  (414 nm). The actual extinction coefficients for DTNB at 405 nm and 414 nm are  $12.8 \text{ mM}^{-1}\text{cm}^{-1}$  and  $13.6 \text{ mM}^{-1}\text{cm}^{-1}$ , respectively. The values have been adjusted for the pathlength of the solution in the well (0.496 cm). One unit is defined as the NADPH-dependent production of 2  $\mu\text{mol}$  of 2-nitro-5-thiobenzoate per minute at 22°C.

Corrected  $\Delta A/\text{min.}$  (sample) =

$$\Delta A/\text{min (sample)} - [\Delta A/\text{min (sample + ATM)} - \Delta A/\text{min (Bkg + ATM)}]$$

TrxR Activity ( $\mu\text{mol}/\text{min}/\text{ml}$ ) =

$$\frac{\text{Corrected } \Delta A/\text{min. (sample)}}{6.35 \text{ mM}^{-1}} \times \frac{0.2 \text{ ml}}{0.02 \text{ ml}} \times \text{Sample Dilution}$$

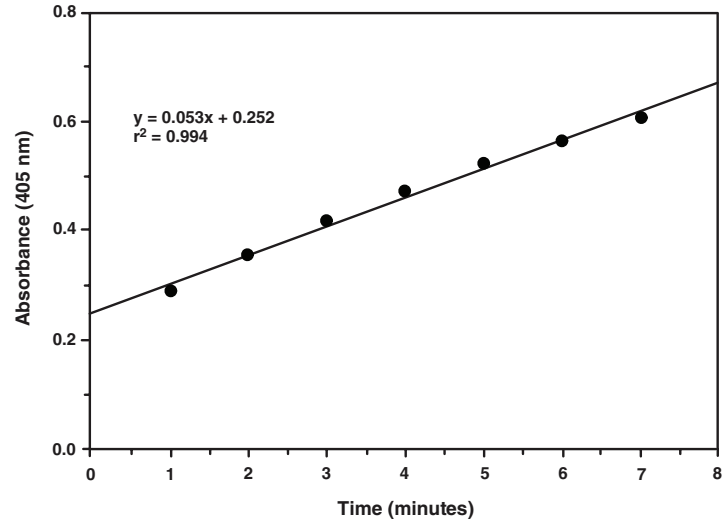


Figure 2. Activity of the rat liver TrxR

## Performance Characteristics

### Sensitivity:

Under the standardized conditions of the assay described in this booklet, the detection range of the assay is from 0.015-0.08  $\mu\text{mol}/\text{min}/\text{ml}$  of TrxR activity, which is equivalent to an absorbance increase of 0.01-0.05 per minute.

### Precision:

When a series of twenty-one thioredoxin reductase measurements were performed on the same day, the intra-assay coefficient of variation was 2.3%. When a series of twenty-one thioredoxin reductase measurements were performed on five different days under the same experimental conditions, the inter-assay coefficient of variation was 6.1%.

### Interferences

The following reagents were tested for interference in the assay.

	Reagent	Will Interfere (Yes or No)
Buffers:	Tris (pH 8)	No
	HEPES (pH 7)	No
	Phosphate (pH 7)	No
	Borate (pH 9)	Yes
Detergents:	Triton X-100 ( $\leq 1\%$ )	No
	Polysorbate 20 ( $\leq 1\%$ )	No
	CHAPS ( $\leq 1\%$ )	No
Protease Inhibitors/ Chelators:	Antipain ( $\leq 0.1$ mg/ml)	No
	PMSF ( $\leq 200$ $\mu$ M)	No
	Leupeptin ( $\leq 10$ $\mu$ g/ml)	No
	Trypsin ( $\leq 10$ $\mu$ g/ml)	No
	Chymostatin ( $\leq 10$ $\mu$ g/ml)	No
	EDTA ( $\leq 10$ mM)	No
	EGTA ( $\leq 1$ mM)	No
Solvents:	Ethanol (10 $\mu$ l)	No
	Methanol (10 $\mu$ l)	Yes
	Dimethylsulfoxide (10 $\mu$ l)	No
Others:	Glutathione	Yes
	BSA ( $\leq 1\%$ )	No
	Glycerol ( $\leq 10\%$ )	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
No activity was detected in the sample	TrxR activity was too low	Concentrate the sample using an Amicon concentrator with a molecular weight cut-off of 30,000 and re-assay
The sample rate did not change in the presence of ATM	TrxR activity was too low	Concentrate the sample using an Amicon concentrator with a molecular weight cut-off of 30,000 and re-assay
The sample starting absorbance is $> 1.0$	There maybe interference due to thiol-containing compounds	Dilute the sample and re-assay

## References

1. Mustacich, D. and Powis, G. Thioredoxin reductase. *Biochem J.* **346**, 1-8 (2000).
2. Becker, K., Gromer, S., Schirmer, R.H., *et al.* Thioredoxin reductase as a pathophysiological factor and drug target. *Eur. J. Biochem.* **267**, 6118-6125 (2000).
3. Turanov, A.A., Su, D., and Gladyshev, V.N. Mouse mitochondrial thioredoxin reductase: Characterization of alternative cytosolic forms and cellular targets. *J. Biol. Chem.* **281(32)**, 22953-22963 (2006).
4. Sun, Q.-A., Wu, Y., Zappacosta, F., *et al.* Redox regulation of cell signaling by selenocysteine in mammalian thioredoxin reductases. *J. Biol. Chem.* **274(35)**, 24522-24530 (1999).
5. Smith, A.D., Morris, V.C., and Levander, O.A. Rapid determination of glutathione peroxidase and thioredoxin reductase activities using a 96-well microplate format: Comparison to standard cuvette-based assays. *Int. J. Vitam. Nutr. Res.* **71(1)**, 87-92 (2001).
6. Hill, K.E., McCollum, G.W., and Burk, R.F. Determination of thioredoxin reductase activity in rat liver supernatant. *Anal. Biochem.* **253**, 123-125 (1997).
7. Smith, A.D., Guidry, C.A., Morris, V.C., *et al.* Aurothioglucose inhibits murine thioredoxin reductase activity *in vivo*. *J. Nutr.* **129**, 194-198 (1999).
8. Gromer, S., Arscott, L.D., Williams, C.H., Jr., *et al.* Human placenta thioredoxin reductase. Isolation of the selenoenzyme, steady kinetics, and inhibition by therapeutic gold compounds. *J. Biol. Chem.* **273(32)**, 20096-20101 (1998).

## Related Products

Antioxidant Assay Kit - Item No. 709001  
Ascorbate Assay Kit - Item No. 700420  
Catalase Assay Kit - Item No. 707002  
Glutathione Assay Kit - Item No. 703002  
Glutathione Peroxidase Assay Kit - Item No. 703102  
Glutathione Reductase Assay Kit - Item No. 703202  
Glutathione S-Transferase Assay Kit - Item No. 703302  
Hydrogen Peroxide (Urinary) Assay Kit - Item No. 706011  
8-Isoprostane EIA Kit - Item No. 516351  
Lipid Hydroperoxide Assay Kit - Item No. 705002  
iPF<sub>2α</sub>-VI EIA Kit - Item No. 516301  
Protein Carbonyl Assay Kit - Item No. 10005020  
Protein Carbonyl Fluorometric Assay Kit - Item No. 700490  
Superoxide Dismutase Assay Kit - Item No. 706002  
TBARS Assay Kit - Item No. 10009055  
Xanthine Oxidase Assay Kit - Item No. 10010895

## Warranty and Limitation of Remedy

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery. Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence. This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

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

## NOTES

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.

©04/25/2011, Cayman Chemical Company, Ann Arbor, MI, All rights reserved. Printed in U.S.A.