

# Thioredoxin Reductase Assay Kit

Catalog Number KA0883

100 assays

Version: 04

Intended for research use only

www.abnova.com



# **Table of Contents**

ntroduction3	5
Background	}
General Information4	ŀ
Materials Supplied	ŀ
Storage Instruction	ŀ
Assay Protocol5	)
Reagent Preparation	5
Assay Procedure	5
Data Analysis7	,
Calculation of Results7	7
Resources	\$
Troubleshooting	3



### Introduction

#### **Background**

Thioredoxin reductase (TrxR) (EC 1.8.1.9) is a ubiquitous enzyme which is involved in many cellular processes such as cell growth, p53 activity, and protection against oxidation stress, etc. The mammalian TrxR reduces thioredoxins as well as non-disulfide substrates such as selenite, lipoic acids, lipid hydroperoxides, and hydrogen peroxide. The Thioredoxin Reductase Assay Kit provides a convenient colorimetric assay for detecting TrxR activity in various samples. In the assay TrxR catalyzes the reduction of 5, 5'-dithiobis (2-nitrobenzoic) acid (DTNB) with NADPH to 5-thio-2-nitrobenzoic acid (TNB<sup>2-</sup>), which generate a strong yellow color ( $\lambda$ max = 412 nm). Since in crude biological samples other enzymes, such as glutathione reductase and glutathione peroxidase, can also reduce DTNB, therefore, TrxR specific inhibitor is utilized to determine TrxR specific activity. Two assays are performed: the first measurement is of the total DTNB reduction by the sample, and the second one is the DTNB reduction by the sample in the presence of the TrxR specific inhibitor. The difference between the two results is the DTNB reduction by TrxR.



# **General Information**

#### Materials Supplied

List of component

Component	Amount
TrxR Assay Buffer	25 mL
TNB Standard (lyophilized)	1 vial
DTNB (lyophilized)	1 vial
NADPH (lyophilized)	1 vial
TrxR Positive Control	1 vial
TrxR Inhibitor (Iyophilized)	1 vial

#### Storage Instruction

Store the kit at -20°C, protect from light. Warm Assay Buffer to room temperature before use. Briefly centrifuge vials prior to opening. Read the entire protocol before performing the assay.



# **Assay Protocol**

#### **Reagent Preparation**

- TNB Standard: Dissolve lyophilized TNB standard into 0.5 mL Assay Buffer to generate 5 mM TNB Standard. The TNB standard solution is stable for 1 week at 4°C or 2 month at -20°C.
- DTNB Solution: Dissolve DTNB into 0.9 mL Assay Buffer, sufficient for 100 assays. The DTNB solution is stable for 1 week at 4°C or 2 month at -20°C.
- NADPH: Dissolve one vial with 0.22 mL dH<sub>2</sub>O; sufficient for 100 assays. The solution is stable for 1 week at 4°C or 2 month at -20°C.
- TrxR Positive Control: Reconstitute with 90 µL Assay Buffer to generate ~0.2 mU/µL TrxR; it is stable for 1 day at 4°C or 2 month at -20°C.
- TrxR Inhibitor: Dissolve TrxR Inhibitor into 1.2 mL Assay Buffer, sufficient for 100 assays. The TrxR Inhibitor solution is stable for 2 month at -20°C.
- Ensure that the Assay Buffer is at room temperature before use. Keep samples, NADPH, TrxR inhibitor, TrxR Positive Control on ice during the assay.

#### Assay Procedure

- 1. TNB Standard Curve: Add 0, 2, 4, 6, 8, 10 μL of the TNB Standard into 96-well plate in duplicate to generate 0, 10, 20, 30, 40, 50 nmol/well standard. Bring the final volume to 100 μL with Assay Buffer.
- Sample and Positive Control Preparations: Take 20 mg Tissues or 2 x 10<sup>6</sup> Cells and homogenize in 100-200 µL cold Assay Buffer on ice (It is recommended to add Protease Inhibitor Cocktail to the buffer); Centrifuge at 10,000 x g for 15 min at 4°C; Collect the supernatant for assay and store on ice.
- 3. Serum can be tested directly. Determine the protein concentration of the supernatant using the Bradford Reagent. Keep samples at -80°C for storage.
- 4. Assay Procedure: Add 2-50 μL sample or 10 μL TrxR positive control into each well, adjust volume to 50 μL with assay buffer. 2 sets of samples should be tested as with or without TrxR Inhibitor. Add 10 μL of TrxR Inhibitor to one set of the sample for testing background enzyme activity, and add 10 μL of Assay Buffer to the other set of sample for testing total DTNB reduction, mix well.
- Reaction Mix: Mix enough reagents for the number of assays to be performed. For each well, prepare a total 40 µL Reaction Mix:
  - 30 μL Assay Buffer 8 μL DTNB Solution
  - 2 µL NADPH
- 6. Add 40 μL of the Reaction Mix to each test sample, mix well. Measure O.D.412 nm at T<sub>1</sub> to get A<sub>1t</sub> and A<sub>1l</sub>, measure O.D.412 nm again at T<sub>2</sub> after incubating the reaction at 25°C for 20 min (The incubate time can vary depend on the sample concentration) to get A<sub>2t</sub> and A<sub>2l</sub>, protect from light. The O.D. of TNB<sup>2-</sup> generated by TrxR is ΔA<sub>412 nm</sub> = (A<sub>2t</sub> A<sub>1t</sub>) (A<sub>2t</sub>-A<sub>1t</sub>).



Note: It is essential to read  $A_{1t}$ ,  $A_{1l}$ ,  $A_{2t}$  and  $A_{2l}$  in the reaction linear range. It will be more accurate if you read the reaction kinetics. Then choose  $A_{1t}$ ,  $A_{1l}$ ,  $A_{2t}$  and  $A_{2l}$  in the reaction linear range.



# **Data Analysis**

#### **Calculation of Results**

Plot the TNB standard Curve. Apply the  $\Delta A_{412nm}$  to the TNB standard curve to get B nmol of TNB (TNB amount generated between T1 and T2 in the reaction wells).

TrxR Activity =  $\frac{B}{(T2-T1) \times V}$  X Sample Dilution Factor = nmol/min/ml = mU/mL

Where: B is the TNB amount from TNB standard Curve (in nmol).

T1 is the time of the first reading  $(A_{1t} \text{ and } A_{1l})$  (in min).

T2 is the time of the second reading  $(A_{2t} \text{ and } A_{2l})$  (in min).

V is the pretreated sample volume added into the reaction well (in mL).

TrxR Unit Definition: One unit of TrxR is the amount of enzyme that generates1.0 µmol of TNB per minute at 25°C. The oxidation of 1 mole of NADPH to NADP will generate 2 mole TNB finally, therefore, 1 TNB unit equals 0.5 NADP unit.





# Resources

# Troubleshooting

#### GENERAL TROUBLESHOOTING GUIDE:

Problems	Cause	Solution
Assay not working	Use of ice-cold assay buffer	Assay buffer must be at room
	Omission of a step in the protocol	temperature
	Plate read at incorrect wavelength	• Refer and follow the data sheet
	Use of a different 96-well plate	precisely
		• Check the wavelength in the data
		sheet and the filter settings of the instrument
		• Fluorescence: Black plates (clear
		bottoms) ; Luminescence: White
		plates ; Colorimeters: Clear plates
Samples with erratic	Use of an incompatible sample type	• Refer data sheet for details about
readings		incompatible samples
	• Samples prepared in a different buffer	• Use the assay buffer provided in the kit
		or refer data sheet for instructions
	Cell/ tissue samples were not	• Use Dounce homogenizer (increase
	completely homogenized	the number of strokes); observe for
		lysis under microscope
	• Samples used after multiple free-thaw	• Aliquot and freeze samples if needed
	cycles	to use multiple times
	• Presence of interfering substance in	Troubleshoot if needed
	the sample	
	• Use of old or inappropriately stored	• Use fresh samples or store at correct
	samples	temperatures until use



Lower/ Higher	•	Improperly thawed components	•	Thaw all components completely and
readings in Samples				mix gently before use
and Standards	•	Use of expired kit or improperly stored	•	Always check the expiry date and store
		reagents		the components appropriately
	•	Allowing the reagents to sit for	•	Always thaw and prepare fresh
		extended times on ice		reaction mix before use
	•	Incorrect incubation times or	•	Refer datasheet & verify correct
		temperatures		incubation times and temperatures
	•	Incorrect volumes used	•	Use calibrated pipettes and aliquot
				correctly
Readings do not	•	Use of partially thawed components	•	Thaw and resuspend all components
follow a linear pattern				before preparing the reaction mix
for Standard curve	•	Pipetting errors in the standard	•	Avoid pipetting small volumes
	•	Pipetting errors in the reaction mix	•	Prepare a master reaction mix
				whenever possible
	•	Air bubbles formed in well	•	Pipette gently against the wall of the
				tubes
	•	Standard stock is at an incorrect	•	Always refer the dilutions in the data
		concentration		sheet
	•	Calculation errors	•	Recheck calculations after referring the
				data sheet
	•	Substituting reagents from older kits/	•	Use fresh components from the same
		lots		kit
Unanticipated results	•	Measured at incorrect wavelength	•	Check the equipment and the filter
				setting
	•	Samples contain interfering	•	Troubleshoot if it interferes with the kit
		substances		
	•	Use of incompatible sample type	•	Refer data sheet to check if sample is
				compatible with the kit or optimization
				is needed
	•	Sample readings above/below the	•	Concentrate/ Dilute sample so as to be
		linear range		in the linear range
			~	

Note: The most probable list of causes is under each problem section. Causes/ Solutions may overlap with other problems.