

## Product datasheet

# Thioredoxin Reductase Assay Kit (Colorimetric) ab83463

[6 References](#) [5 Images](#)

### Overview

---

<b>Product name</b>	Thioredoxin Reductase Assay Kit (Colorimetric)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Urine, Cell culture extracts, Other biological fluids, Tissue Extracts
<b>Assay type</b>	Enzyme activity
<b>Assay time</b>	0h 40m
<b>Product overview</b>	Thioredoxin Reductase Assay Kit (ab83463) is a specific assay for detecting Thioredoxin Reductase (TrxR) activity.

In the thioredoxin reductase assay protocol, TrxR catalyzes the reduction of DTNB to TNB<sup>2-</sup> in the presence of NADPH, which generates a strong yellow color (OD<sub>max</sub> = 412 nm).

Other enzymes present in crude biological samples such as glutathione reductase and glutathione peroxidase can also reduce DTNB. In order to measure TrxR-only activity, a TrxR specific inhibitor is used in a separate reaction to determine TrxR specific activity. The difference between total DTNB reduction in the sample and DTNB reduction in the sample in presence of TrxR inhibitor is the value of specific TrxR activity in the sample.

Thioredoxin reductase assay protocol summary:

- add samples and standards to wells
- add reaction mix
- analyze with a microplate reader over 20 min

<b>Notes</b>	Thioredoxin reductase (TrxR, EC 1.8.1.9) is a ubiquitous mammalian enzyme that catalyzes the NADPH-dependent reduction of the redox protein thioredoxin, as well as of other endogenous and exogenous compounds such as selenite, lipid hydroperoxides and hydrogen peroxide.
--------------	---

<b>Platform</b>	Microplate reader
-----------------	-------------------

### Properties

---

<b>Storage instructions</b>	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	Identifier	100 tests
DTNB	Red	1 vial
NADPH	Blue	1 vial
TNB Standard	Brown	1 vial
TrxR Assay Buffer	WM	1 x 25ml
TrxR Inhibitor	Clear	1 vial
TrxR Positive Control	Green	1 vial

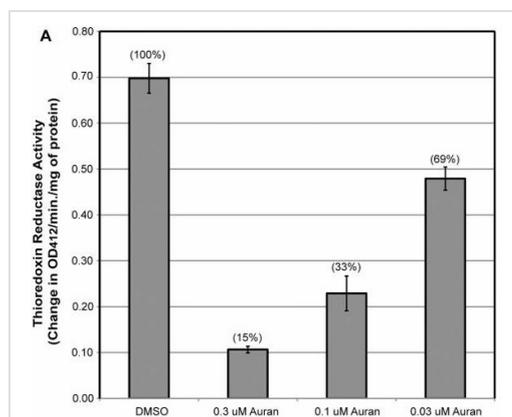
## Relevance

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.

## Cellular localization

TXNRD1: Cytoplasmic. TXNRD2: Cytoplasmic. Nuclear. Microsome. Endoplasmic reticulum. TXNRD3: Mitochondrial.

## Images



Bulman CA et al. PLoS One. 9:e0003534 (2015)

Functional Studies - beta Thioredoxin reductase

Assay Kit (ab83463)

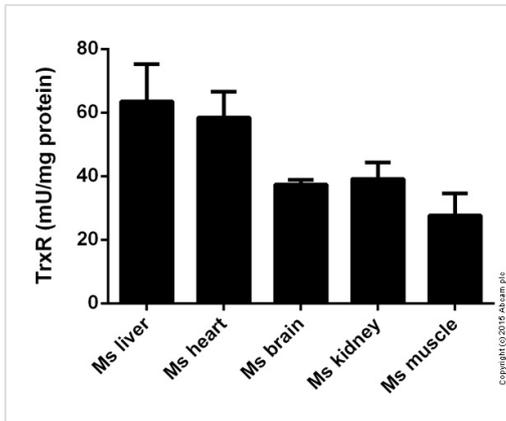
Image from Bulman CA et al., PLoS One. 2015;9(2):e0003534. Fig 4(A); doi: 10.1371/journal.pntd.0003534.

Activity of endogenous *Brugia* thioredoxin reductase from soluble worm lysates following incubation with 1% DMSO or 0.3  $\mu$ M, 0.1  $\mu$ M, or 0.03  $\mu$ M of auranofin *in vitro*. Percentages indicate the percent activity of TrxR compared to DMSO controls. Thioredoxin reductase activity was significantly reduced ( $p < 0.05$ ) to 15%, 33% and 69% of endogenous activity, respectively, compared to the activity in DMSO-treated worms.

Thioredoxin reductase activity of worm lysates was assayed using female *B. malayi* treated *in vitro* with either 0.3  $\mu$ M, 0.1  $\mu$ M, or 0.03  $\mu$ M auranofin or 1% DMSO. After 5 hours of treatment, worm motility was measured using the Worminator, and then worms (24 in each group) were pooled, washed three times in PBS, and lysed by douncing in a glass homogenizer in assay buffer (ab83463) with 1 mM PMSF. The crude lysates were centrifuged at 10,000 rcf for 15 minutes at 4°C to pellet insoluble material. The total protein concentrations of soluble lysates were measured using the Bradford assay. The soluble lysates were incubated for 20 minutes in assay buffer or assay buffer with a proprietary thioredoxin reductase specific inhibitor before adding a specific substrate, DTNB (5, 5'-dithiobis (2-nitrobenzoic) acid), and measuring activity at 20 second intervals for 40 minutes using the SpectraMax Plus Microplate Reader (Molecular Devices, Sunnyvale, CA) at  $\lambda = 412$  nm. Lysates

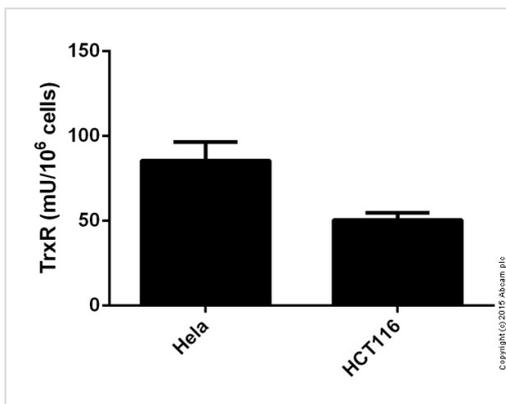
were tested in duplicate. TrxR activity was calculated based on the linear amount of TNB produced per minute per mg of total protein and adjusted for background activity from enzymes other than TrxR in the lysates.

Thioredoxin reductase measured in mouse tissue lysates showing activity (mU) per mg protein of sample tested



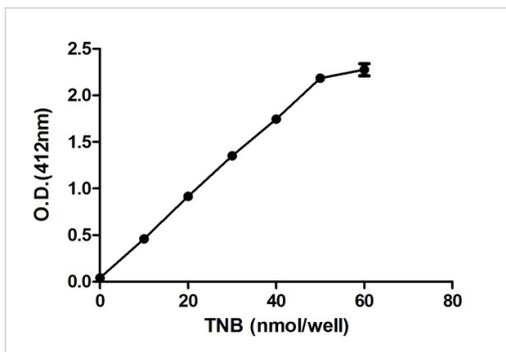
Functional Studies - beta Thioredoxin reductase Assay Kit (ab83463)

Thioredoxin reductase measured in cell lysates showing activity (mU) per 1 mln of cells tested

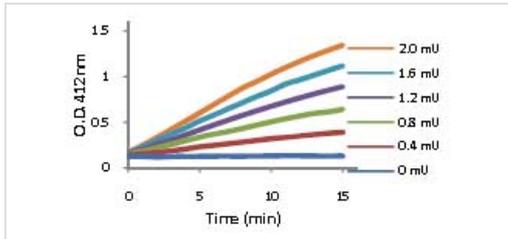


Functional Studies - beta Thioredoxin reductase Assay Kit (ab83463)

Standard curve (colourimetric) : mean of duplicates (+/-SD) with background readings subtracted



Functional Studies - beta Thioredoxin reductase Assay Kit (ab83463)



Thioredoxin reductase Kinetic Data using ab83463.

Functional Studies - Thioredoxin Reductase Assay  
Kit (ab83463)

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

### Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors