

Product datasheet

Thioredoxin Reductase Assay Kit (Colorimetric) ab83463

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

Overview

Product name	Thioredoxin Reductase Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Urine, Cell culture extracts, Other biological fluids, Tissue Extracts
Assay type	Enzyme activity
Assay time	0h 40m
Product overview	<p>Thioredoxin Reductase Assay Kit (ab83463) is a specific assay for detecting Thioredoxin Reductase (TrxR) activity.</p> <p>In the thioredoxin reductase assay protocol, TrxR catalyzes the reduction of DTNB to TNB^{2-} in the presence of NADPH, which generates a strong yellow color (ODmax = 412 nm).</p> <p>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.</p> <p>We have tested the GR positive control from ab83461 in the conditions of ab83463, and we do not observe activity of the GR positive control.</p> <p>Thioredoxin reductase assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add reaction mix - analyze with a microplate reader over 20 min
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K763 Thioredoxin Reductase Activity Colorimetric Assay Kit. K763-100 is the same size as the 100 test size of ab83463.</p> <p>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.</p>

Platform

Microplate reader

Properties

Storage instructions

Store at -20°C. Please refer to protocols.

Components	100 tests
DTNB	1 vial
NADPH I	1 vial
Thioredoxin Reductase Assay Buffer	1 x 25ml
Thioredoxin Reductase Inhibitor	1 vial
Thioredoxin Reductase Positive Control	1 vial
TNB Standard	1 vial

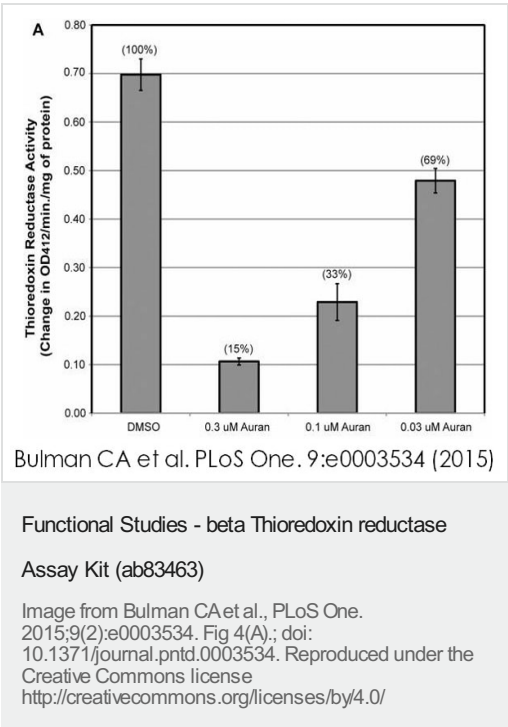
Relevance

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.

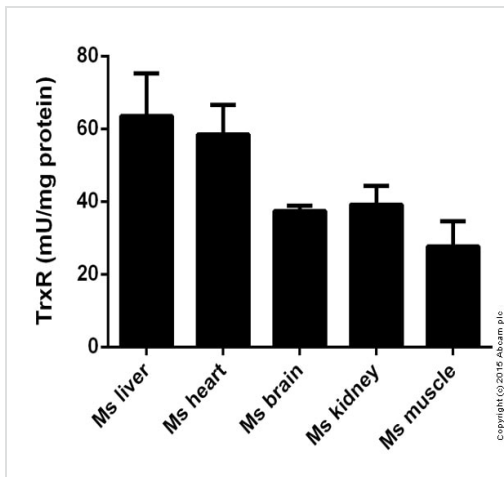
Images



Activity of endogenous *Brugia* thioredoxin reductase from soluble worm lysates following incubation with 1% DMSO or 0.3 μ M, 0.1 μ M, or 0.03 μ 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.

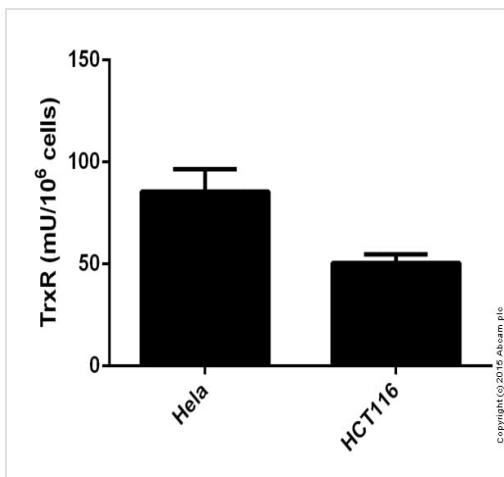
Thioredoxin reductase activity of worm lysates was assayed using female *B. malayi* treated *in vitro* with either 0.3 μ M, 0.1 μ M, or 0.03 μ M auranofin or 1% DMSO. After 5 hours of treatment, worm motility was measured using the Worminator, and then worms (24 in each group) were pooled, washed three 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.



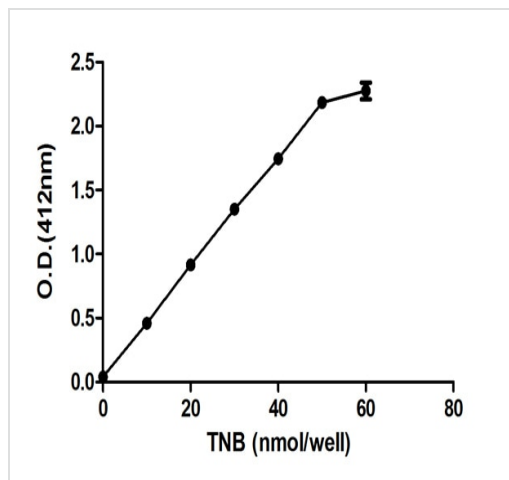
Thioredoxin reductase measured in mouse tissue lysates showing activity (mU) per mg protein of sample tested

Functional Studies - beta Thioredoxin reductase
Assay Kit (ab83463)



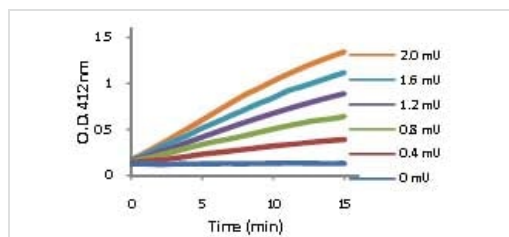
Thioredoxin reductase measured in cell lysates showing activity (mU) per 1 mln of cells tested

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. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

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