

PRODUCT INFORMATION



Thioredoxin Reductase 2 (human, recombinant)

Item No. 39625

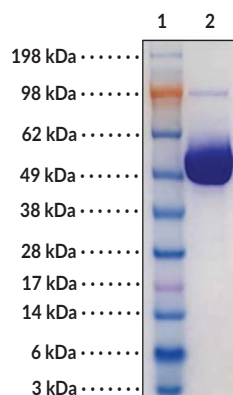
Overview and Properties

Synonyms:	NADPH-dependent Thioredoxin Reductase 2, TrxR2, Txnrd2
Source:	Active recombinant human TrxR2 expressed in <i>E. coli</i>
Amino Acids:	488 residues
Storage:	-20°C (as supplied)
Stability:	≥1 year
Purity:	≥95% estimated by SDS-PAGE
Supplied in:	TE Buffer with 50% glycerol
Protein	
Concentration:	1 mg/ml
Unit Definition:	One unit is defined as the amount of enzyme required to reduce 1 μmol DTNB per minute in 0.5 ml standard DTNB assay with 2.5 mM DTNB and 0.3 NADPH in TE buffer (50 nM Tris-HCl, 2 mM EDTA, pH 7.5)

Special Conditions: Centrifuge tube briefly before opening

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
Lane 2: TrxR2

Coomassie stained SDS-PAGE Analysis
of 10 μg TrxR2.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/31/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Thioredoxin reductase 2 (TrxR2) is an oxidoreductase encoded by the *TXNRD2* gene in humans and a member of the antioxidant thioredoxin system.¹ It exists as a homodimer and contains a dimer interface domain, FAD- and NADPH-binding domains, an N-terminal redox catalytic site, and a C-terminal selenocysteine residue, which is essential for the catalytic activity of TrxR2.^{1,2} TrxR2 is ubiquitously expressed and localizes primarily to the mitochondria.³ It catalyzes the NADPH-dependent reduction of oxidized thioredoxin 2 (Trx2), restoring the disulfide reductase function of Trx2, which reduces peroxiredoxin-3 (Prx3) dimers.^{4,5} TrxR2 has roles in the maintenance of mitochondrial integrity and redox homeostasis in the aging heart and in the regulation of chondrocyte viability, proliferation, and differentiation.² Genome-wide deletion of *Txnrd2* is embryonic lethal in mice.⁶ *TXNRD2* expression and TrxR2 protein levels are increased in bone biopsies from patients with osteosarcoma that later progress to metastatic disease.⁷ SNPs in *TXNRD2* are associated with serum markers of selenium status in patients with advanced- or high-stage, but not localized low-grade, prostate cancer.⁸ Cayman's Thioredoxin Reductase 2 (human, recombinant) protein can be used for enzyme activity assays.

References

1. Mustacich, D. and Powis, G. Thioredoxin reductase. *Biochem. J.* **346**(Pt. 1), 1-8 (2000).
2. Scalcon, V., Bindoli, A., and Rigobello, M.P. Significance of the mitochondrial thioredoxin reductase in cancer cells: An update on role, targets and inhibitors. *Free Radic. Biol. Med.* **127**, 62-79 (2018).
3. Li, W., Bandyopadhyay, J., Hwaang, H.S., et al. Two thioredoxin reductases, *trxr-1* and *trxr-2*, have differential physiological roles in *Caenorhabditis elegans*. *Mol. Cells* **34**(2), 209-218 (2012).
4. Holmgren, A. and Lu, J. Thioredoxin and thioredoxin reductase: Current research with special reference to human disease. *Biochem. Biophys. Res. Commun.* **396**(1), 120-124 (2010).
5. Watanabe, R., Nakamura, H., Masutani, H., et al. Anti-oxidative, anti-cancer and anti-inflammatory actions by thioredoxin 1 and thioredoxin-binding protein-2. *Pharmacol. Ther.* **127**(3), 261-270 (2010).
6. Conrad, M., Jakupoglu, C., Moreno, S.G., et al. Essential role for mitochondrial thioredoxin reductase in hematopoiesis, heart development, and heart function. *Mol. Cell. Biol.* **24**(21), 9414-9423 (2004).
7. Topkas, E., Cai, N., Cumming, A., et al. Auranofin is a potent suppressor of osteosarcoma metastasis. *Oncotarget* **7**(1), 831-844 (2016).
8. Méplan, C., Rohrmann, S., Steinbrecher, A., et al. Polymorphisms in thioredoxin reductase and selenoprotein K genes and selenium status modulate risk of prostate cancer. *PLoS One* **7**(11), e48709 (2012).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM