

PRODUCT INFORMATION



Thioredoxin Reductase 1 (human, recombinant)

Item No. 39624

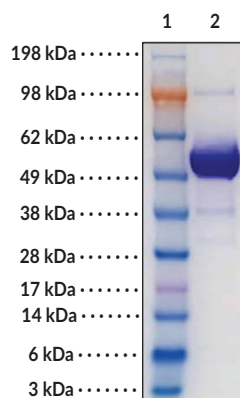
Overview and Properties

Synonyms:	NADPH-dependent Thioredoxin Reductase, TrxR1, Txnrd1
Source:	Active recombinant human TrxR1 expressed in <i>E. coli</i>
Amino Acids:	499 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: TrxR1

Coomassie stained SDS-PAGE Analysis
of 10 μg TrxR1.

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, 08/30/2023

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 1 (TrxR1) is an oxidoreductase encoded by the *TXNRD1* gene in humans and a member of the antioxidant thioredoxin system, which is involved in the maintenance of cellular thiol redox homeostasis.¹⁻³ 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 TrxR1.^{1,4} TrxR1 is ubiquitously expressed, localizes to the cytoplasm, and is regulated by the antioxidant transcription factor Nrf2.^{2,6} TrxR1 catalyzes the NADPH-dependent reduction of oxidized thioredoxin (Trx), restoring the disulfide reductase function of Trx, which regulates redox-sensitive transcription factors, such as NF- κ B and p53, and has roles in apoptosis and cell signaling.^{3,7} Genome-wide deletion of *Txnrd1* is embryonic lethal in mice.⁸ Increased serum TrxR1 activity is associated with reduced progression-free survival in patients with non-small cell lung cancer (NSCLC).⁹ Cayman's Thioredoxin Reductase 1 (human, recombinant) protein can be used for enzyme activity assay.

References

1. Mustacich, D. and Powis, G. Thioredoxin reductase. *Biochem. J.* **346**(Pt. 1), 1-8 (2000).
2. Cebula, M., Schmidt, E.E., and Arnér, E.S.J. TrxR1 as a potent regulator of the Nrf2-Keap1 response system. *Antioxid. Redox Signal.* **23**(10), 823-853 (2015).
3. 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).
4. Fritz-Wolf, K., Kehr, S., Stumpf, M., *et al.* Crystal structure of the human thioredoxin reductase-thioredoxin complex. *Nat. Commun.* **2**, 383 (2011).
6. 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).
7. 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).
8. Bondareva, A.A., Capecchi, M.R., Iverson, S.V., *et al.* Effects of thioredoxin reductase-1 deletion on embryogenesis and transcriptome. *Free Radic. Biol. Med.* **43**(6), 911-923 (2007).
9. Chen, G., Chen, Q., Zeng, F., *et al.* The serum activity of thioredoxin reductases 1 (TrxR1) is correlated with the poor prognosis in EGFR wild-type and ALK negative non-small cell lung cancer. *Oncotarget* **8**(70), 115270-115279 (2017).

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