

PRODUCT INFORMATION



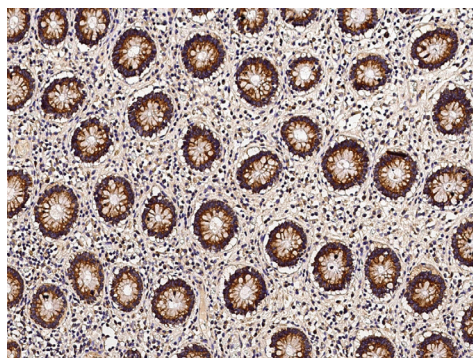
Thioredoxin 2 Rabbit Monoclonal Antibody (Clone 054)

Item No. 37075

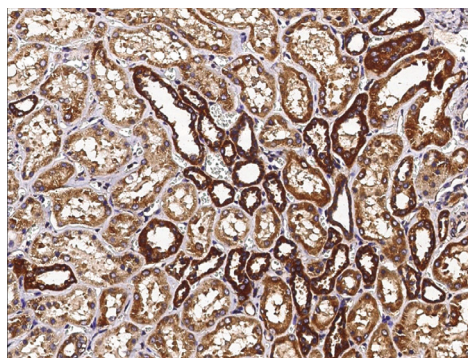
Overview and Properties

Contents:	This vial contains 50 or 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	Mitochondrial Thioredoxin, mtTrx, Trx2, Txn2
Immunogen:	Recombinant human Trx2
Cross Reactivity:	(+) Trx2
Species Reactivity:	(+) Human
Uniprot No.:	Q99757
Form:	Liquid
Storage:	-80°C (as supplied)
Stability:	≥1 year
Storage Buffer:	0.2 µm filtered solution in PBS
Clone:	054
Host:	Rabbit
Isotype:	IgG
Application:	Immunohistochemistry (IHC); the recommended starting dilution is 1:100-1:500. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Immunohistochemical staining of formalin-fixed and paraffin embedded human Trx2 in human small intestine with Thioredoxin 2 Rabbit Monoclonal Antibody (Clone 054) at 1:200 dilution.



Immunohistochemical staining of formalin-fixed and paraffin embedded human Trx2 in human kidney with Thioredoxin 2 Rabbit Monoclonal Antibody (Clone 054) at 1:200 dilution.

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, 07/19/2022

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 2 (Trx2) is a thiol-disulfide oxidoreductase and part of the antioxidant thioredoxin system that is involved in the maintenance of cellular redox homeostasis.¹ It is ubiquitously expressed, with the highest levels of expression in metabolically active tissues such as the heart, skeletal muscle, testis, kidney, adrenal gland, and brain, and is localized to the mitochondria.^{2,3} Trx2 contains two active site cysteine residues, but lacks the additional structural cysteines present in Trx1 that participate in dimerization.⁴ During the catalytic cycle, the active site cysteines are oxidized to a disulfide upon reduction of oxidized protein disulfide substrates and are subsequently restored to their reduced state by thioredoxin reductase (TrxR) and NADPH.¹ Trx2 is upregulated by TNF- α and reactive oxygen species (ROS), inhibits TNF- α -induced NF- κ B activation, and inhibits apoptosis through redox-sensitive binding and regulation of apoptosis signal-regulating kinase 1 (ASK1).^{5,6} Homozygous disruption of the gene encoding Trx2, *Txn2*, causes exencephaly and is embryonic lethal in mice.⁷ Trx2 deficiency is associated with premature neurodegeneration.⁸ Cayman's Thioredoxin 2 Rabbit Monoclonal Antibody (Clone 054) can be used for immunohistochemistry (IHC).

References

1. Miranda-Vizuete, A., Damdimopoulos, A.E., and Spyrou, G. The mitochondrial thioredoxin system. *Antioxid. Redox Signal.* **2(4)**, 801-810 (2000).
2. Spyrou, G., Enmark, E., Miranda-Vizuete, A., *et al.* Cloning and expression of a novel mammalian thioredoxin. *J. Biol. Chem.* **272(5)**, 2936-2941 (1997).
3. Rybnikova, E., Damdimopoulos, A.E., Gustafsson, J.A., *et al.* Expression of novel antioxidant thioredoxin-2 in the rat brain. *Eur. J. Neurosci.* **12(5)**, 1669-1678 (2000).
4. Smeets, A., Evrard, C., Landtmeters, M., *et al.* Crystal structures of oxidized and reduced forms of human mitochondrial thioredoxin 2. *Protein Sci.* **14(10)**, 2610-2621 (2005).
5. Hansen, J.M., Zhang, H., and Jones, D.P. Mitochondrial thioredoxin-2 has a key role in determining tumor necrosis factor- α -induced reactive oxygen species generation, NF- κ B activation, and apoptosis. *Toxicol. Sci.* **91(2)**, 643-650 (2006).
6. Zhang, R., Al-Lamki, R., Bai, L., *et al.* Thioredoxin-2 inhibits mitochondria-located ASK1-mediated apoptosis in a JNK-independent manner. *Circ. Res.* **94(11)**, 1483-1491 (2004).
7. Nonn, L., Williams, R.R., Erickson, R.P., *et al.* The absence of mitochondrial thioredoxin 2 causes massive apoptosis, exencephaly, and early embryonic lethality in homozygous mice. *Mol. Cell. Biol.* **23(3)**, 916-922 (2003).
8. Holzerova, E., Danhauser, K., Haack, T.B., *et al.* Human thioredoxin 2 deficiency impairs mitochondrial redox homeostasis and causes early-onset neurodegeneration. *Brain* **139(Pt. 2)**, 346-354 (2016).

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