

# PRODUCT INFORMATION



## Thioredoxin 80 (human, recombinant)

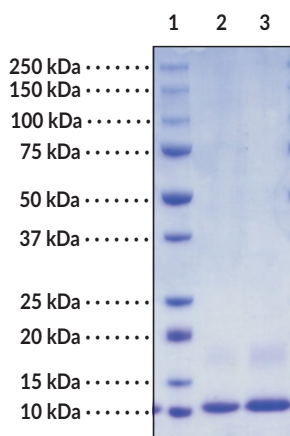
Item No. 11522

### Overview and Properties

**Synonyms:** ECEF, Eosinophil Cytotoxicity-enhancing Factor, Truncated Thioredoxin 1, Trx80, Txn80  
**Source:** Recombinant human thioredoxin 80 expressed in *E. coli*  
**Amino Acids:** 1-80  
**Uniprot No.:** P10599  
**Molecular Weight:** 9.04 kDa  
**Storage:** 4°C (as supplied)  
**Stability:** ≥6 months  
**Purity:** ≥85% estimated by SDS-PAGE  
**Supplied in:** Lyophilized in PBS, pH 7.4

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

### Image



Lane 1: MW Markers  
Lane 2: Thioredoxin 80 (2 µg)  
Lane 3: Thioredoxin 80 (4 µg)

SDS-PAGE Analysis of Thioredoxin 80.

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, 07/08/2021

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

---

Truncated thioredoxin 1 (Trx80) is a C-terminally truncated cleavage product of the oxidoreductase Trx1 that acts as a cytokine.<sup>1,2</sup> It is expressed primarily in monocytes but also in T cells, activated platelets, and cytotrophoblasts, and localizes to the cell surface.<sup>1</sup> Trx80 contains three cysteines but does not have oxidoreductase activity, and, unlike Trx, it forms homodimers. It induces proliferation of, and increases the expression of CD14, CD40, CD54, and CD86 in, monocytes, as well as enhances eosinophil-induced cytotoxicity of schistosomes and the secretion of IL-12 and IFN- $\gamma$  in peripheral blood mononuclear cells (PBMCs). It also acts as a chemoattractant for monocytes, T cells, and neutrophils. Trx80-activated monocytes inhibit growth of the bacteria *L. monocytogenes* and *B. abortus* by inducing their accumulation in lysosomes, but Trx80 increases replication of HIV in infected macrophages.<sup>3,4</sup> It inhibits amyloid- $\beta$  (A $\beta$ ) aggregation in cell-free assays and reduces the toxicity of A $\beta$  in SH-SY5Y cells.<sup>5</sup> Levels of Trx80 are reduced in postmortem brain tissue from patients with Alzheimer's disease.<sup>2</sup>

## References

---

1. Pekkari, K. and Homgren, A. Truncated thioredoxin: Physiological functions and mechanism. *Antioxid. Redox Signal.* **6(1)**, 53-61 (2004).
2. Gil-Bea, F., Akterin, S., Persson, T., *et al.* Thioredoxin-80 is a product of alpha-secretase cleavage that inhibits amyloid-beta aggregation and is decreased in Alzheimer's disease brain. *EMBO Mol. Med.* **4(10)**, 1097-1111 (2012).
3. Cortes-Bratti, X., Bassères, E., Herrera-Rodriguez, F., *et al.* Thioredoxin 80-activated-monocytes (TAMs) inhibit the replication of intracellular pathogens. *PLoS One* **6(2)**, e16960 (2011).
4. Newman, G.W., Balcewicz-Sablinska, M.K., Guarnaccia, J.R., *et al.* Opposing regulatory effects of thioredoxin and eosinophil cytotoxicity-enhancing factor on the development of human immunodeficiency virus 1. *J. Exp. Med.* **180(1)**, 359-363 (1994).
5. Mahmood, D.F.D., Abderrazak, A., El Hadri, K., *et al.* The thioredoxin system as a therapeutic target in human health and disease. *Antioxid. Redox Signal.* **19(11)**, 1266-1303 (2013).

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