

CERTIFICATE of ANALYSIS



Thioredoxin Reductase 1 (C. elegans, recombinant)

Item No. 39628 • Batch No. 0824699 • Expires 19AUG2027

Overview

Formulation:	TE buffer with 50% glycerol
Concentration:	6.6 mg/mL *Buffer
Storage/stability:	-80°C/1yr
Unit definition:	One unit is defined as the amount of enzyme required to reduce 1 μ mol DTNB per minute in 0.5 ml standard DTNB assay with 2.5 mM DTNB and 0.3 NADPH in TE buffer (50 mM Tris-HCl, 2 mM EDTA, pH 7.5)
Purity:	$\geq$ 95% estimated by SDS-PAGE

Reviewed and approved by: Justine Albertson

Approved date: 19AUG2026

A handwritten signature in black ink, appearing to read "J. Albertson", written over a faint, circular, textured background.

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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