

CERTIFICATE of ANALYSIS



Thioredoxin 1 (human) Polyclonal Antibody

Item No. 11538 • Batch No. 0584926 • Expires 14SEP2023

Overview

Formulation: 100 µg antibody in 50 mM Tris-HCl, 1 mM EDTA, 15 mM sodium azide, pH 7.5
Concentration: 0.51 mg/ml
Storage/Stability: -20°C/1yr
Applications: IHC, WB
Species reactivity: (+)Human; other species not tested

Reviewed and approved by: Jeremy Bickel

Approved date: 08APR2020

A handwritten signature in black ink, appearing to read "Jeremy Bickel", written over a light blue horizontal line.

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