

# Thioredoxin monoclonal antibody (2C9)

Thioredoxins act as antioxidants by reducing the proteins by cysteine thiol-disulfide exchange. Several plasmid expression vectors have been constructed that direct the synthesis of foreign polypeptides in *E. coli* as fusions with C-terminal Thioredoxin (TrxA) to offer soluble expression of normally insoluble or difficult to express proteins such as cytokines and growth factors.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

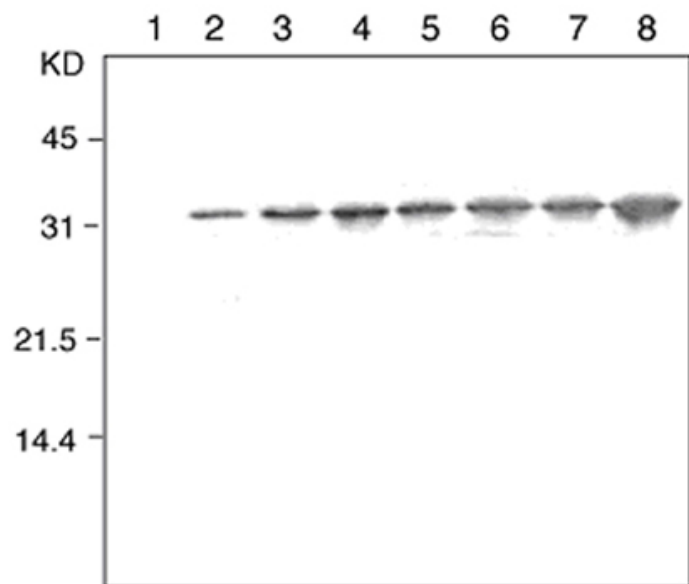
## Ordering Information

[Order Online »](#)

ADI-MSA-150-E	100µg
---------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Thioredoxin fusion protein expressed cell lysate after no induction (1), or induction for 30 min (2), 60 min (3), 90 min (4), 120 min (5), 150 min (6), 180 min (7), 210 min (8)

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name      TXN

Application      WB

Clone      2C9

Formulation      Liquid. In PBS containing 50% glycerol.

Host      Mouse

Immunogen      E. coli thioredoxin (trxA).

Isotype      IgG1

Purity Detail      Protein A affinity purified.

Recommendation Dilutions/Conditions      Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity      E. coli

Specificity      Recognizes trx and trx-tag-f thioredoxin.

Worry-free Guarantee      This antibody is covered by our [Worry-Free Guarantee](#)

.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)