

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 »](#)

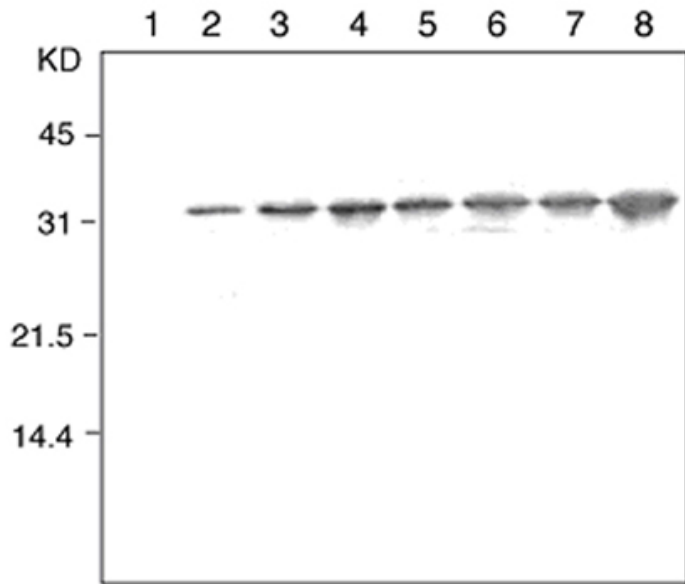
Ordering Information

[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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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