

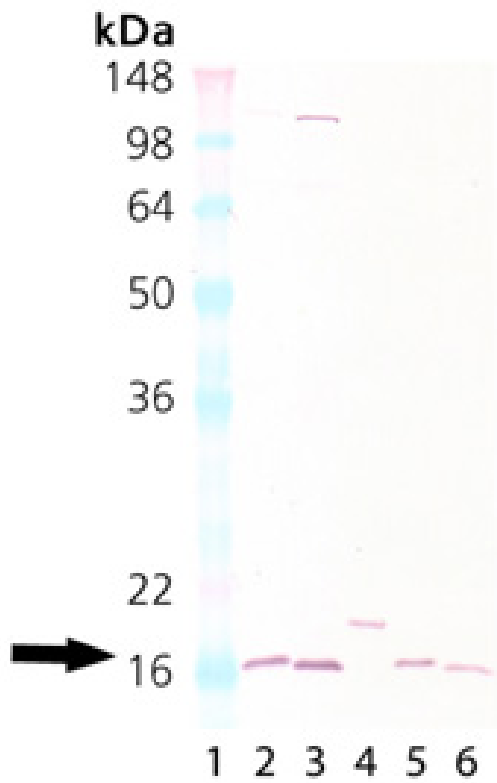
Cu/Zn SOD polyclonal antibody (biotin conjugate)

Superoxide dismutase (SOD) is responsible for the elimination of cytotoxic active oxygen by catalyzing the dismutation of the superoxide radical to oxygen and hydrogen peroxide. There are three SOD isoenzymes in mammalian cells: extracellular SOD (EC SOD, SOD-3); copper and zinc-containing SOD (Cu/Zn SOD, SOD-1); and manganese-containing SOD (Mn SOD, SOD-2). The Cu/Zn form binds Cu and Zn ions and exists as a soluble 32kDa homodimer in the cytosol. Mutations in the SOD-1 gene accounts for approximately 20% of amyelotrophic lateral sclerosis (ALS) or 'Lou Gehrig disease,' a neurodegenerative disorder characterized by the death of motor neurons in the brain, brainstem, and spinal cord that results in fatal paralysis.

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

Manuals, SDS & CofA

[View Online »](#)



Western Blot Analysis of Cu/Zn SOD: Lane 1: MWM
Lane 2: Rat liver tissue lysate, Lane 3: Mouse liver tissue lysate, Lane 4: HeLa, (cell lysate) (Prod No. ADI-LYC-HL100), Lane 5: Brain (mouse), (tissue extract) (Prod No. ADI-LYT-MB100), Lane 6: Brain (rat), (tissue extract) (Prod No. ADI-LYT-RB100).

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Superoxide Dismutase (Cu-Zn), Superoxide dismutase 1, SOD1

Application IF, IHC (PS), IP, WB

Application Notes Detects a band of ~23kDa (human) or ~19kDa (mouse and rat) by Western blot.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

Host Rabbit

Immunogen Native human Cu/Zn SOD.

Purity Detail Protein A-affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000, colorimetric) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Clam, Human, Mouse, Rat

UniProt ID P00441

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-eu@enzolifesciences.com)

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