

# 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](#).

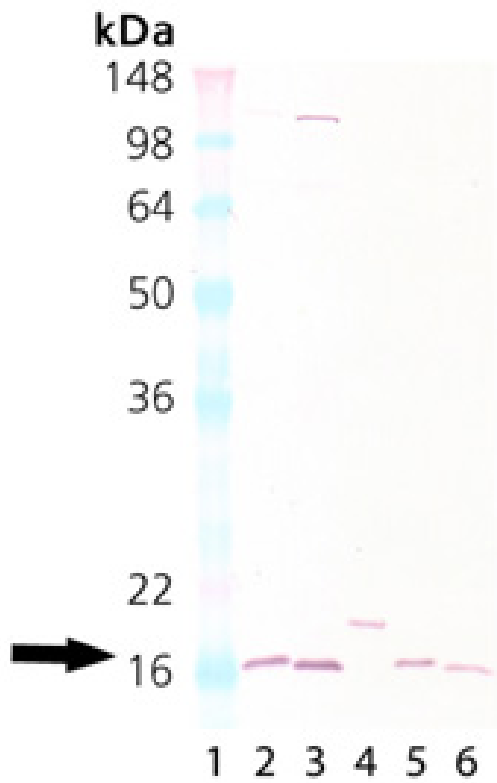
## Ordering Information

[Order Online »](#)

ADI-SOD-100B-0050	50µg
ADI-SOD-100B-0200	200µg

## 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)