

EC SOD polyclonal antibody

SOD (Superoxide dismutase) 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, they are: EC SOD (extracellular SOD), Cu/Zn SOD (copper and zinc-containing SOD) and Mn SOD (manganese-containing SOD).

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

Citations: 8

[View Online »](#)

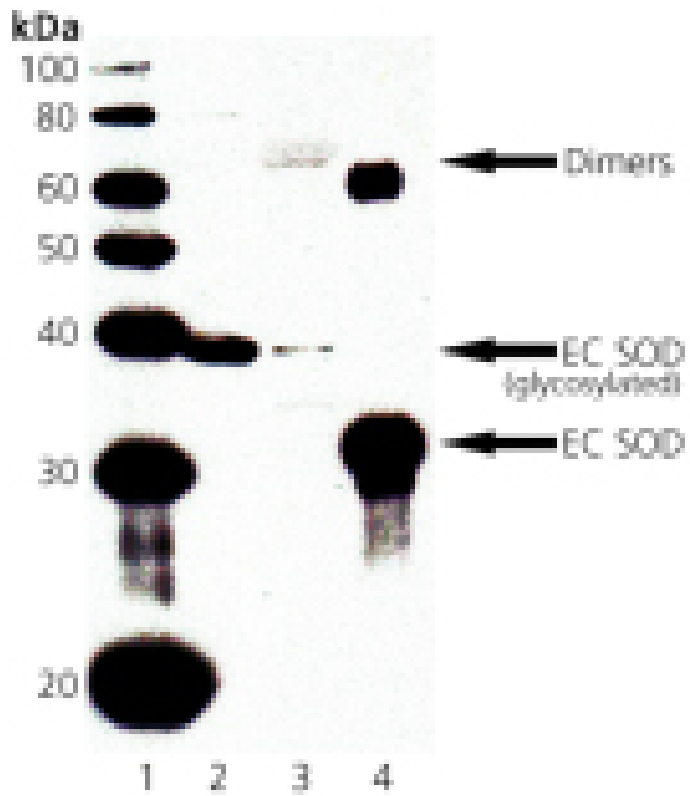
Ordering Information

[Order Online »](#)

ADI-SOD-106-F	200µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of EC SOD: Lane 1: MW Marker, Lane 2: Rat Lung Tissue Extract, Lane 3: Mouse Lung Tissue Extract, Lane 4: Human EC SOD Protein.

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, SOD3
Application	WB
Application Notes	Detects a band of ~35-40kDa (glycosylation dependent) by Western blot.
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of EC SOD. The sequence is completely conserved in human, mouse, rat and rabbit.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, ECL)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	Human, Mouse, Rat
UniProt ID	P08294
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)