

Mn 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: 79

[View Online »](#)

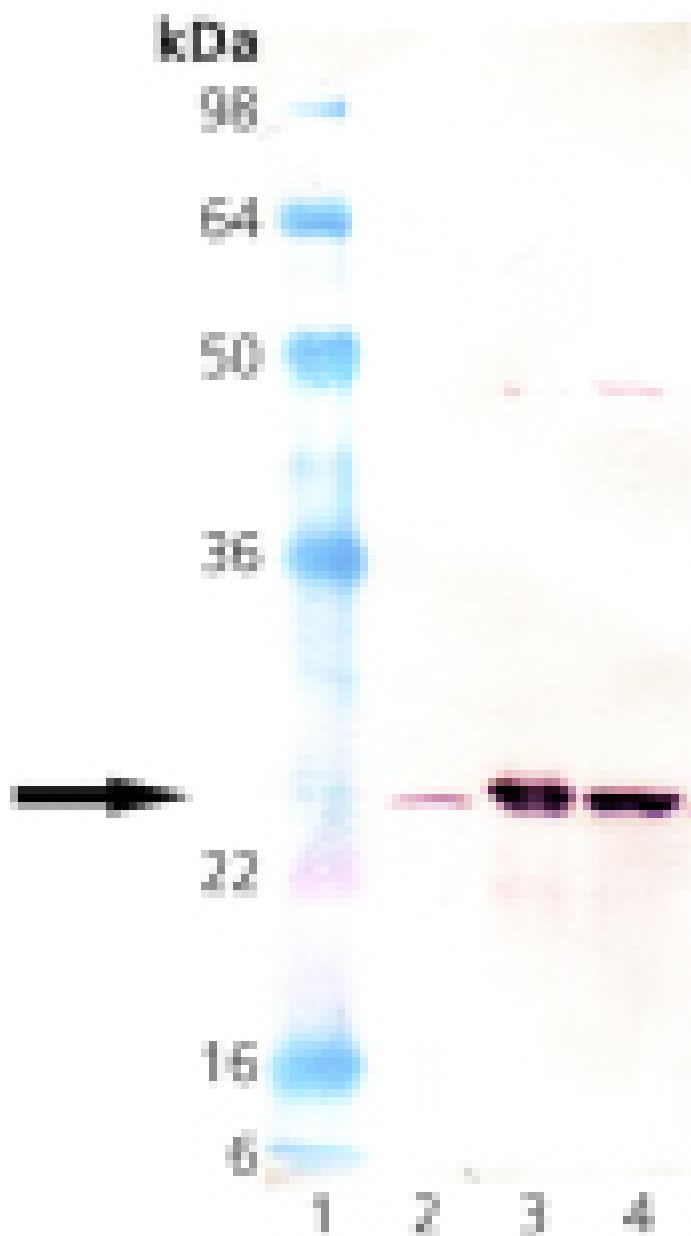
Ordering Information

[Order Online »](#)

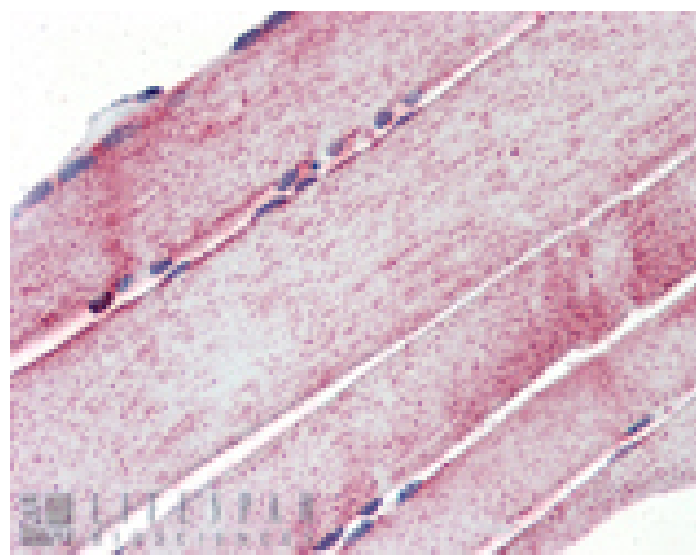
ADI-SOD-110-D	50µg
ADI-SOD-110-F	200µg

Manuals, SDS & CofA

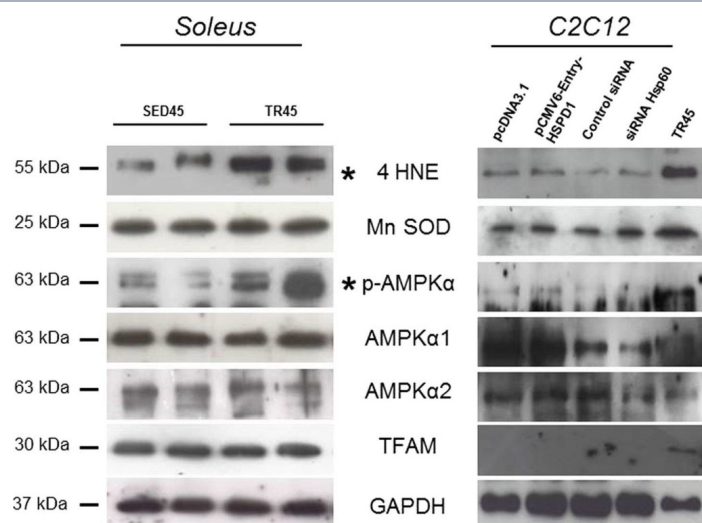
[View Online »](#)



Western blot analysis of Mn SOD: Lane 1: MW Marker, Lane 2: HeLa Cell Lysate, Lane 3: Mouse Brain Tissue Extract, Lane 4: Rat Brain Tissue Extract



Immunohistochemistry analysis of human skeletal muscle tissue stained with Mn SOD, pAb at 10µg/ml.



PGC1 $\alpha 1$ levels increase in the soleus in trained mice and in transfected C2C12 cells upon transfection with pCMV-Entry-HSPD1 vector. (A) representative western blots of soleus and relative expression levels (bars) of PGC1 $\alpha 1$ (113 kDa), 4 HNE (55 kDa), Mn SOD (25 kDa), p-AMPK α (63 kDa), AMPK $\alpha 1$ (63 kDa), AMPK $\alpha 2$ (63 kDa), TFAM (30 kDa) in soleus of sedentary (SED45, open bar, $n = 8$) and trained (TR45, shaded bar, $n = 8$) mice at 45 days. 80 μ g of proteins were loaded in each lane; GAPDH (37 kDa) was used as the loading control. Data are presented as the means \pm SD. Δ significantly different from TR45 mice ($P < 0.001$). AU: Arbitrary Unit. (B) representative western blots of C2C12 cells and relative levels (bars) of PGC1 $\alpha 1$ (113 kDa), 4 HNE (55 kDa), Mn SOD (25 kDa), p-AMPK α (63 kDa), AMPK $\alpha 1$ (63 kDa), AMPK $\alpha 2$ (63 kDa), TFAM (30 kDa) in C2C12 myoblast transfected with pCMV6-Entry-HSPD1 vector (pcDNA3.1 was used as a negative control). 80 μ g of proteins were loaded in each lane; GAPDH (37 kDa) was used as the loading control and TR45 was used as positive control. Data are presented as the means \pm SD. #significantly different from pCMV6-Entry-HSPD1 ($P < 0.0001$). AU: Arbitrary Unit. *significant results. (C) relative expression levels of 4 HNE, p-AMPK α , p-AMPK α /(AMPK $\alpha 1$ + AMPK $\alpha 2$) in the soleus. Open bars, sedentary mice; shaded bars, trained mice. AU: Arbitrary Unit.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Skeletal muscle Heat shock protein 60 increases after endurance training and induces peroxisome proliferator-activated receptor gamma coactivator 1 $\alpha 1$ expression. *Sci Rep* (2016)

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Product Details

Alternative Name	Manganese superoxide dismutase
Application	IHC, WB
Application Notes	Detects a band of ~25kDa by Western blot.
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
Gene/Protein Identifier	NP_000627 (RefSeq)
Host	Rabbit
Immunogen	Human Mn SOD protein.
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	Bovine, Dog, Guinea pig, Hamster, Human, Monkey, Mouse, Rabbit, Rat, Sheep, Xenopus
UniProt ID	P04179
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



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)