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: 75

View Online »

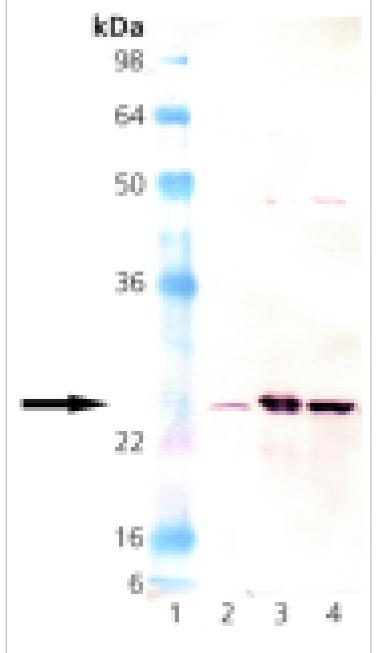
Ordering Information

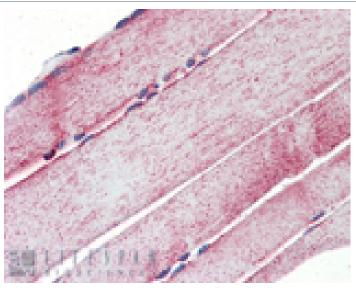
Order Online »

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

Manuals, SDS & CofA

View Online »

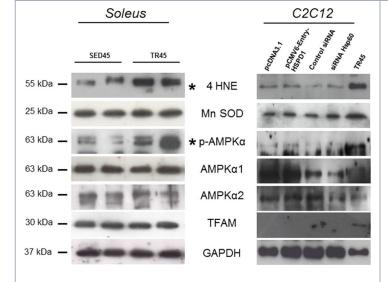




Immunohistochemistry analysis of human skeletal muscle tissue stained with Mn SOD, pAb at 10µg/ml.

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



PGC1 a1 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 α1 (113 kDa), 4 HNE (55 kDa), Mn SOD (25 kDa), p-AMPKα (63 kDa), AMPKα1 (63 kDa), AMPKα2 (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 µg of proteins were loaded in each lane; GAPDH (37 kDa) was used as the loading control. Data are presented as the means ± SD. ∂ significantly different from TR45 mice (P < 0.001). AU: Arbitrary Unit. (B) representative western blots of C2C12 cells and relative levels (bars) of PGC1 α1 (113 kDa), 4 HNE (55 kDa), Mn SOD (25 kDa), p-AMPKα (63 kDa), AMPKα1 (63 kDa), AMPKα2 (63 kDa), TFAM (30 kDa) in C2C12 myoblast transfected with pCMV6-Entry-HSPD1 vector (pcDNA3.1 was used as a negative control). 80 µg 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 ± 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 α 1 + AMPK α 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 α1 expression. *Sci Rep* (2016)

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 Manganese superoxide dismutase

Application IHC, WB

Application NotesDetects 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/ConditionsWestern 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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com