

Mn SOD (mouse), (recombinant) (His- tag)

Recombinant mouse Mn SOD (His-tag) that detoxifies superoxide radicals and protects mitochondria from oxidative stress.

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). Mn SOD is an 80 kDa tetramer that contains Mn ions and resides in the mitochondrial matrix. Mn SOD is a tumor necrosis factor (TNF)-inducible enzyme that protects cells from TNF-mediated apoptosis via superoxide anion detoxification and the subsequent regulation of apoptosis through cytochrome c release and modulation of the redox state of the mitochondria. Overexpression of this enzyme protects neurons from NMDA and nitric oxide-induced neurotoxicity.

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

kDa

750
148

98

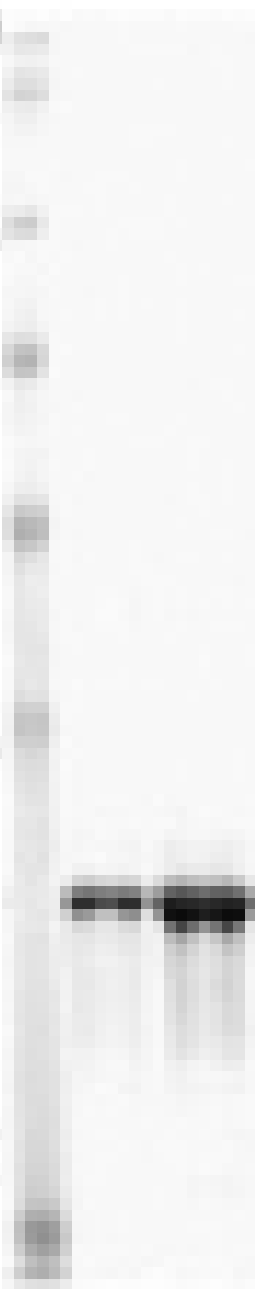
64

50

36

22

16



1 2 3

SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 2 μ g, Lane 3: 5 μ g of Mn SOD (SOD2) Recombinant Mouse Protein.

kDa

750
148

98

64

50

36

22

16



1 2 3

Western Blot analysis: Lane 1: MWM, Lane 2: Mouse Brain Tissue Extract, Lane 3: Recombinant Mn SOD Mouse Protein, probed with Mn SOD pAb (Prod. No. ADI-SOD-111).

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Superoxide Dismutase, SOD2
Application Notes	Activity Assay (superoxide dismutase). Western blot control.
Formulation	Liquid. In 50mM TRIS-HCl, pH 7.5, containing 1.5 % glycerol and 80mM sodium chloride.
MW	~25kDa
Purity	≥90% (SDS-PAGE; Western blot)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> . Mouse Mn SOD is fused at the N-terminus to a His-tag.
UniProt ID	P09671

Last modified: October 9, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com