

PRODUCT INFORMATION



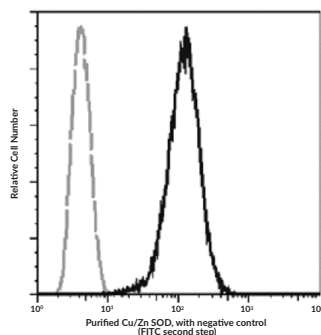
Cu/Zn SOD Monoclonal Antibody (Clone 11)

Item No. 37087

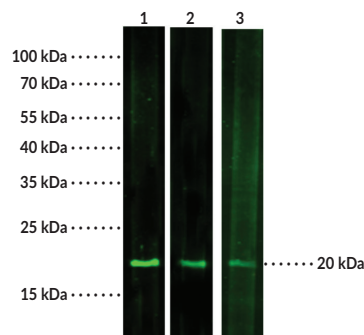
Overview and Properties

Contents: This vial contains 50, 100, or 200 μ l of protein A-affinity purified monoclonal antibody.
Synonyms: Cu/Zn Superoxide Dismutase, SOD1
Immunogen: Recombinant human Cu/Zn SOD
Cross Reactivity: (+) Cu/Zn SOD
Species Reactivity: (+) Human
Form: Liquid
Storage: -80°C (as supplied)
Stability: ≥ 1 year
Storage Buffer: 0.2 μ m filtered solution in PBS
Clone: 11
Host: Mouse
Isotype: IgG2b
Applications: Flow cytometry (FC), Immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution is 1:100-1:500 for FC, 1:500-1:1,000 for WB, and 0.2-1 μ l/mg of lysate for IP. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images

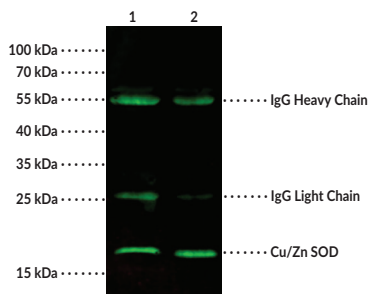


Flow cytometric analysis of human Cu/Zn SOD expression in HeLa cells. The cells were treated according to the manufacturer's manual (BD Pharmingen™), then stained with purified Cu/Zn SOD Monoclonal Antibody (Clone 11) followed by a FITC-conjugated secondary antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.



Lane 1: Jurkat whole cell lysate (30 μ g)
Lane 2: HeLa whole cell lysate (30 μ g)
Lane 3: HepG2 whole cell lysate (30 μ g)

WB of Cu/Zn SOD Monoclonal Antibody (Clone 11) at a dilution of 1:500.



Lane 1: Jurkat whole cell lysate
Lane 2: HepG2 whole cell lysate

WB of Cu/Zn SOD IP samples (500 μ g whole cell lysate and 500 μ l IgG per IP).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/13/2024

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Cu/Zn Superoxide dismutase (Cu/Zn SOD) is an antioxidant enzyme that protects cells from oxidative stress by scavenging superoxide anions.¹ It exists as a homodimer that is stabilized by a disulfide bond between the subunits. Each monomer is composed of a Greek key β -barrel and an active site channel made up of a β 4/ β 5 loop, which interacts with the dimer interface and active site zinc, and an electrostatic loop, which facilitates superoxide entry into the copper-containing active site.² Cu/Zn SOD is ubiquitously expressed and primarily localizes to the cytosol but has also been found in the nucleus, peroxisomes, lysosomes, and the intermembrane space of mitochondria.^{3,4} It catalyzes the dismutation of superoxide to hydrogen peroxide and oxygen by alternating reduction and reoxidation of copper at the enzyme active site.¹ Overexpression of *SOD1* protects against increases in reactive oxygen species (ROS) and apoptosis in an *in vitro* model of renal ischemia induced by ATP depletion.⁵ *In vivo*, transgenic mice expressing the destabilized Cu/Zn SOD^{G93A} mutation exhibit impaired spinal cord mitochondrial respiration.^{6,7} Mutations in *SOD1* are associated with amyotrophic lateral sclerosis (ALS).⁸ Cayman's Cu/Zn SOD Monoclonal Antibody (Clone 11) can be used for flow cytometry (FC), immunoprecipitation (IP), and Western blot (WB) applications.

References

1. Fukai, T. and Ushio-Fukai, M. Superoxide dismutases: Role in redox signaling, vascular function, and diseases. *Antioxid. Redox Signal.* **15(6)**, 1583-1606 (2011).
2. Perry, J.J.P., Shin, D.S., Getzoff, E.D., *et al.* The structural biochemistry of the superoxide dismutases. *Biochim. Biophys. Acta* **1804(2)**, 245-262 (2010).
3. Crapo, J.D., Oury, T., Rabouille, C., *et al.* Copper,zinc superoxide dismutase is primarily a cytosolic protein in human cells. *Proc. Natl. Acad. Sci. USA* **89(21)**, 10405-10409 (1992).
4. Okado-Matsumoto, A. and Fridovich, I. Subcellular distribution of superoxide dismutases (SOD) in rat liver. Cu, Zn-SOD in mitochondria. *J. Biol. Chem.* **276(46)**, 38388-38393 (2001).
5. Liang, H.L., Arsenault, J., Mortensen, J., *et al.* Partial attenuation of cytotoxicity and apoptosis by SOD1 in ischemic renal epithelial cells. *Apoptosis* **14(10)**, 1176-1189 (2009).
6. Mattiazz, M., D'Aurelio, M., Gajewski, C.D., *et al.* Mutated human SOD1 causes dysfunction of oxidative phosphorylation in mitochondria of transgenic mice. *J. Biol. Chem.* **277(33)**, 29626-29633 (2002).
7. Stathopoulos, P.B., Rumpfheldt, J.A.O., Karbassi, F., *et al.* Calorimetric analysis of thermodynamic stability and aggregation for apo and holo amyotrophic lateral sclerosis-associated Gly-93 mutants of superoxide dismutase. *J. Biol. Chem.* **281(10)**, 6184-6193 (2006).
8. Furukawa, Y. and O'Halloran, T.V. Posttranslational modifications in Cu,Zn-superoxide dismutase and mutations associated with amyotrophic lateral sclerosis. *Antioxid. Redox Signal.* **8(5-6)**, 847-867 (2006).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM