

Product Information



Cu/Zn SOD (human) Polyclonal Antibody

Catalog No. 10011388

Contents:	This vial contains affinity-purified antibody at 1 mg/ml in PBS, pH 7.0, containing 0.1% sodium azide and 50% glycerol
Synonym:	Cu/Zn Superoxide Dimutase, SOD1
Antigen:	Human Cu/Zn SOD
Host:	Rabbit
Cross Reactivity:	(+) human, murine, bovine, monkey, coral, canine, hamster, porcine, rabbit, ovine, and rat Cu/Zn SOD. Detects a 23 kDa (human) and 19 kDa (other species) proteins corresponding to the molecular mass of Cu/Zn SOD on SDS-PAGE immunoblots.
Stability:	≥1 year at -20°C
Applications:	Western blot (WB), immunoprecipitation, EIA, and immunohistochemistry. ^{1,2} The recommended starting concentration for WB is 0.2 µg/ml, which was found to be sufficient for detection of Cu/Zn SOD in 20 µg of HeLa cell lysate. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Superoxide dismutase (SOD) is an endogenously produced intracellular enzyme present in almost every cell in the body.³ It works by catalyzing the dismutation of the superoxide radical O_2^- to O_2 and H_2O_2 , which are then metabolized to H_2O and O_2 by catalase and glutathione peroxidase.^{4,5} In general, SODs play a major role in antioxidant defense mechanisms.⁶

There are two main types of SOD in mammalian cells. One form, SOD1, contains Cu and Zn ions as a homodimer and exists in the cytoplasm. The two subunits of 16 kDa each are linked by two cystines forming an intra-subunit disulphide bridge.³ The second form, SOD2, is a manganese-containing enzyme and resides in the mitochondrial matrix. It is a homotetramer of 80 kDa. The third form, SOD3 or EC-SOD, is like SOD1 in that it contains Cu and Zn ions, however it is distinct in that it is a homotetramer, with a mass of 30 kDa and it exists only in the extra-cellular space.¹ SOD3 can also be distinguished by its heparin-binding capacity.⁷

References

1. Wispé, J.R., Clark, J.C., Burhans, M.S., *et al.* Synthesis and processing of the precursor for human manganese-superoxide dismutase. *Biochim. Biophys. Acta* **994**(1), 30-36 (1989).
2. Katsuki, H., Tomita, M., Takenaka, C., *et al.* Superoxide dismutase activity in organotypic midbrain-striatum co-cultures is associated with resistance of dopaminergic neurons to excitotoxicity. *J. Neurochem.* **76**, 1336-1345 (2001).
3. Furukawa, Y. and O'Halloran, T.V. Posttranslational modifications in Cu,Zn-superoxide dismutase and mutations associated with amyotrophic lateral sclerosis. *Antioxidants & Redox Signaling* **8**(5-6), 847-867 (2006).
4. Bannister, J.V., Bannister, W.H., and Rotilio, G. Aspects of the structure, function, and applications of superoxide dismutase. *Crit. Rev. Biochem. Mol. Biol.* **22**(2), 111-180 (1987).
5. Hassan, H.M. Biosynthesis and regulation of superoxide dismutases. *Free Radic. Biol. Med.* **5**(5-6), 377-385 (1988).
6. Gao, B., Flores, S.C., Leff, J.A., *et al.* Synthesis and anti-inflammatory activity of a chimeric recombinant superoxide dismutase: SOD2/3. *Am. J. Physiol. Lung Cell Mol. Physiol.* **284**, L917-L925 (2003).
7. Adachi, T., Ohta, H., Yamada, H., *et al.* Quantitative analysis of extracellular-superoxide dismutase in serum and urine by ELISA with monoclonal antibody. *Clin. Chim. Acta* **212**(3), 89-102 (1992).

Related Products

Superoxide Dismutase Assay Kit - Cat. No. 706002 • Hydrogen Peroxide (urinary) Assay Kit - Cat. No. 706011 • Catalase Assay Kit - Cat. No. 707002 • Antioxidant Assay Kit - Cat. No. 709001 • Cu/Zn SOD (rat) Polyclonal Antibody - Cat. No. 10011387 • Mn (rat) SOD Polyclonal Antibody - Cat. No. 10011389 • Mn SOD (human) Polyclonal Antibody - Cat. No. 10011390 • GRP90 Monoclonal Antibody (Clone 9G10) - Cat. No. 10011424 • DNA/RNA Damage Monoclonal Antibody (Clone 15A3) - Cat. No. 10011446

WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/17/2008

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com