

## SOD3 Protein

Recombinant mature form protein expressed in Sf9 cells

### Catalog # S28-31G

Lot # L2218-6

### Product Description

Recombinant full-length human SOD3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_003102](#).

### Gene Aliases

EC-SOD

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

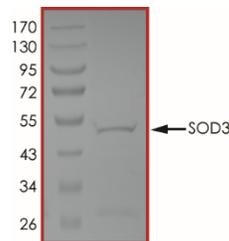
### Scientific Background

Extracellular superoxide dismutase [Cu-Zn] (SOD3) protects the extracellular space from toxic effect of reactive oxygen intermediates by converting superoxide radicals into hydrogen peroxide and oxygen. Although the function of SOD3 as a regulator of cellular growth has been well established, the role of SOD3 in tumorigenesis is still unclear.

### References

1. S. Marklund, (1973). "A novel superoxide dismutase of high molecular weight from bovine liver," Acta Chemica Scandinavica, 27,1458-1460.
2. Laukkanen MO. (2016). Extracellular Superoxide Dismutase: Growth Promoter or Tumor Suppressor? Oxid Med Cell Longev. 2016:3612589..
3. <http://www.uniprot.org/uniprot/P08294>.

### Purity



The purity of SOD3 was determined to be **>70%** by densitometry. Approx. MW **52kDa**.

## SOD3 Protein

Recombinant mature form protein expressed in Sf9 cells

Catalog #	S28-31G
Lot #	L2218-6
Purity	>70%
Concentration	0.05 µg/µl
Stability	1yr at $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**