

---

## SOD2 Protein

(Catalog No. E70001)

---

### Description

SOD2 Human Recombinant is a single, non-glycosylated, polypeptide chain containing 219 amino acids (25-222 a.a). SOD2 protein is fused to a 20 amino acid His-Tag at N-terminus.

### Molecular Weight (kDa)

24.4 kDa

### Formulation

20mM Tris HCL pH-8, & 20% glycerol.

### Source

Expressed in E.coli expression system

### Size

25 µg

### Storage Temp & Conditions

Stored desiccated -20°C. Please prevent freeze thaw cycles.

### Purification Notes

Purity: Greater than 95% as determined by SDS-PAGE.

### Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot, Enzyme activity assay

### Ordering Information:

Products	Size	Cat. No.
SOD2 Protein	25 µg	E70001 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*