
SOD1 Protein

(Catalog No. E70000)

Description

Recombinant Human Cu/Zn Superoxide Dismutase is a single monomeric non-glycosylated polypeptide chain containing 154 amino acids.

Molecular Weight (kDa)

15.9 kDa

Formulation

20mM Tris-HCl pH-7.5 and 10% 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 95.0% as determined by (a) Analysis by RP-HPLC.

Application Notes

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

Ordering Information:

| Products | Size | Cat. No. |
|--------------|-------|-----------|
| SOD1 Protein | 25 µg | E70000 -1 |

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