# EC SOD polyclonal antibody

SOD (Superoxide dismutase) is responsible for the elimination of cytotoxic active oxygen by catalyzing the dismutation of the superoxide radical to oxygen and hydrogen peroxide. There are three SOD isoenzymes in mammalian cells, they are: EC SOD (extracellular SOD), Cu/Zn SOD (copper and zinc-containing SOD) and Mn SOD (manganese-containing SOD).

This antibody is covered by our Worry-Free Guarantee.

Citations: 7

View Online »

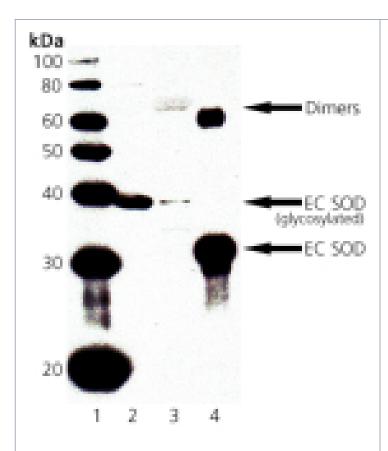
## **Ordering Information**

Order Online »

ADI-SOD-106-D	50µg
ADI-SOD-106-F	200μg

Manuals, SDS & CofA

**View Online »** 



Western blot analysis of EC SOD: Lane 1: MW Marker, Lane 2: Rat Lung Tissue Extract, Lane 3: Mouse Lung Tissue Extract, Lane 4: Human EC SOD Protein.

#### **Handling & Storage**

**Handling** Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

### Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name Superoxide Dismutase, SOD3

**Application** WB

**Application Notes**Detects a band of ~35-40kDa (glycosylation dependent) by

Western blot.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium

azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to a portion of EC SOD.

The sequence is completely conserved in human, mouse,

rat and rabbit.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000, ECL)Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

**Source** Purified from rabbit serum.

Species Reactivity Human, Mouse, Rat

UniProt ID P08294

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com