
SOD1 Polyclonal Antibody

(Catalog #A70918)

Background

The protein encoded by this gene binds copper and zinc ions and is one of two isozymes responsible for destroying free superoxide radicals in the body. The encoded isozyme is a soluble cytoplasmic protein, acting as a homodimer to convert naturally-occurring but harmful superoxide radicals to molecular oxygen and hydrogen peroxide. The other isozyme is a mitochondrial protein. Mutations in this gene have been implicated as causes of familial amyotrophic lateral sclerosis. Rare transcript variants have been reported for this gene.

Description

SOD1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.05% proclin300, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P00441

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 2-154 of human SOD1 (NP_000445.1).

Storage

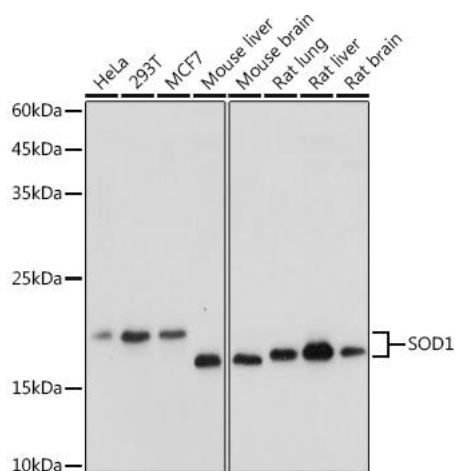
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

SOD1; ALS; ALS1; HEL-S-44; IPOA; SOD; hSod1; homodimer; superoxide dismutase 1

Application

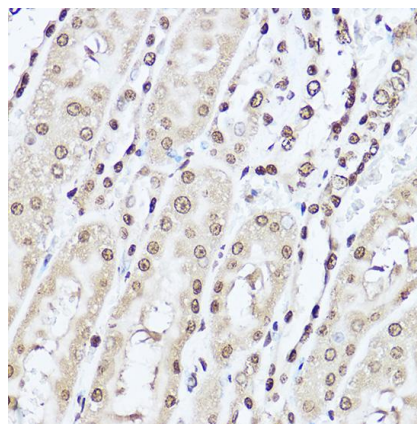
WB, IHC, IF; Recommended Dilution: WB 1:500-1:1000, IHC 1:50-1:200, IF 1:50-1:200



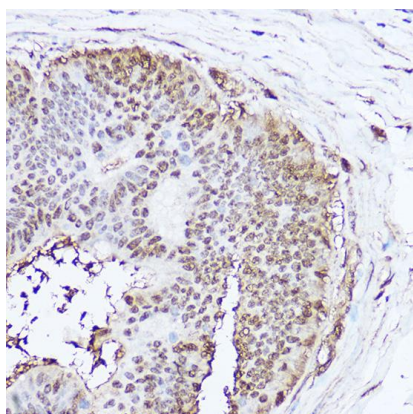
Western blot analysis of extracts of various cell lines, using at SOD1 Polyclonal Antibody 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25µg per lane.

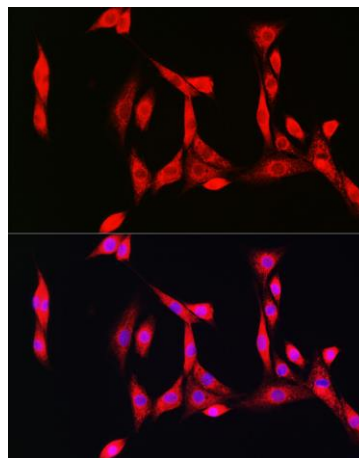
Blocking buffer: 3% nonfat dry milk in TBST.



Immunohistochemistry analysis of paraffin-embedded rat kidney using SOD1 Polyclonal Antibody at dilution of 1:100 (40x lens). Perform high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunohistochemistry analysis of paraffin-embedded human colon carcinoma using SOD1 Polyclonal Antibody at dilution of 1:100 (40x lens). Perform high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunofluorescence analysis of NIH/3T3 cells using SOD1 Polyclonal Antibody at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.