

## Mn SOD Polyclonal Antibody

(Catalog # A-6408)

### Background

Superoxide dismutase (SOD) is an antioxidant enzyme that is responsible for eliminating cytotoxic active oxygen by catalyzing the dismutation of the superoxide radical to oxygen and hydrogen peroxide. Three SOD isoenzymes in mammalian cells have been identified. They include: EC SOD (extracellular SOD), Cu/Zn SOD (copper and zinc-containing SOD) and Mn SOD (manganese-containing SOD). Mn SOD is a tumor necrosis factor (TNF)-inducible enzyme that guards cells against TNF-mediated apoptosis caused by superoxide anion detoxification and the subsequent regulation of apoptosis through cytochrome c release and the modulation of the redox state of the mitochondria. Mn SOD has been associated with tumor suppression in human breast cancer.

### Description

Rabbit polyclonal antibody to Mn SOD.

### Concentration

1mg/ml

### Formulation

Liquid. In sodium phosphate buffer, pH 7.0, containing 15mg/ml BSA, 0.09% sodium azide, and 50% glycerol.

### Immunogen

Native rat Mn SOD

### Specificity

Human, Mouse, Rat, Bovine, Chicken, Clam, Dog, Drosophila, Guinea pig, Hamster, Monkey, Pig, Rabbit, Sheep, Xenopus

### Isotype

IgG

### Purification

Immunoaffinity purified

### Storage

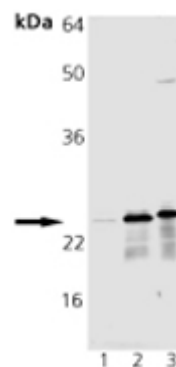
Store at -20°C. Avoid multiple freeze/thaw cycles

### Application

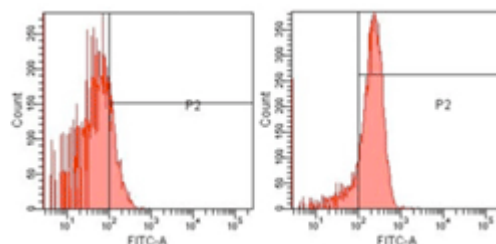
EIA, FC (1:100), IHC, IP (1:100), WB (1:1000, colorimetric)

### Ordering Information

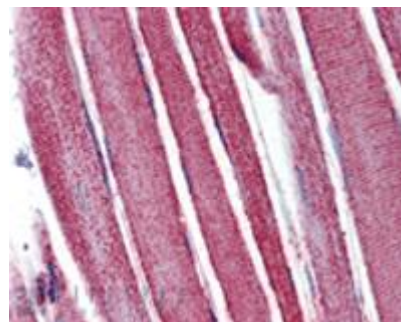
Products	Size	Cat. No.
Mn SOD Polyclonal Antibody	200 µg	A-6408-200



**Fig. 1.** Western blot analysis of Mn SOD Polyclonal Antibody: Lane 1: HeLa Cell Lysate, Lane 2: Rat Brain Tissue Extract, Lane 3: Mouse Brain Tissue Extract.



**Fig. 2.** Flow cytometry analysis of U251 cells using Mn SOD Polyclonal Antibody (right) and control rabbit IgG (left).



**Fig. 3.** Immunohistochemistry analysis of human skeletal muscle tissue stained with Mn SOD Polyclonal Antibody at 10µg/ml.

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

Epigentek Group Inc. • USA Tel: 1-877-374-4368 • Int. Tel: 1-718-484-3990 • Fax: 1-718-484-3990 • info@epigentek.com • www.epigentek.com

2015-02-02-A-6408