

SREBP-1 antibody (mAb)

Catalog Nos: 39939, 39940

RRID: AB_2616606

Clone: IgG-2A4

Isotype: IgG

Application(s): ChIP, EMSA, IP, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 68 and 125 kDa

Background: SREBP-1 and SREBP-2 (Sterol regulatory element-binding protein 1 and 2) are members of the basic-helix-loop-helix-leucine zipper (bHLH-ZIP) family of transcription factors and are involved in the regulation of cholesterol homeostasis. SREBP-1 and SREBP-2 are normally membrane bound at the ER and nuclear envelope, but in sterol-depleted cells, the membrane-bound precursors are cleaved to generate a soluble isoforms that translocate to the nucleus. These active forms, which include the bHLH-ZIP domain, activate transcription of genes involved in the uptake and synthesis of cholesterol.

Immunogen: This SREBP-1 antibody was raised against a recombinant protein corresponding to amino acids 301-407 of human SREBP-1.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.