

SND1 antibody (mAb)

Catalog Nos: 61473, 61474

RRID: AB_2793651

Clone: SN4-1

Isotype: IgG1

Application(s): ICC, IF, IHC, IP, WB

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 115 kDa

Background: **SND1** (staphylococcal nuclease and tudor domain containing 1) functions as a transcriptional coactivator and a modulator of RNA metabolism. As a transcriptional coactivator, it acts as a bridging factor between STAT6 and the basal transcription factors, plays a role in PIM1 regulation of MYB activity and is a transcriptional coactivator for the Epstein-Barr virus nuclear antigen 2 (EBNA2). SND-1 has been identified as part of the RISC complex (RNA-induced silencing complex) and interacts with Ago-2. It is involved in the degradation of hyper-edited inosine containing microRNA precursors as part of the RNA interference pathway.

Immunogen: This antibody was raised against a recombinant protein corresponding to amino acids 505 - 674 of human SND1.

Buffer: Purified IgG in PBS with 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.