
HNRNPA2B1 Polyclonal Antibody

(Catalog # A73256)

This gene belongs to the A/B subfamily of ubiquitously expressed heterogeneous nuclear ribonucleoproteins (hnRNPs). The hnRNPs are RNA binding proteins and they complex with heterogeneous nuclear RNA (hnRNA). These proteins are associated with pre-mRNAs in the nucleus and appear to influence pre-mRNA processing and other aspects of mRNA metabolism and transport. While all of the hnRNPs are present in the nucleus, some seem to shuttle between the nucleus and the cytoplasm. The hnRNP proteins have distinct nucleic acid binding properties. The protein encoded by this gene has two repeats of quasi-RRM domains that bind to RNAs. This gene has been described to generate two alternatively spliced transcript variants which encode different isoforms.

Description

HNRNPA2B1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P22626

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-180 of human HNRNPA2B1 (NP_002128.1).

Storage

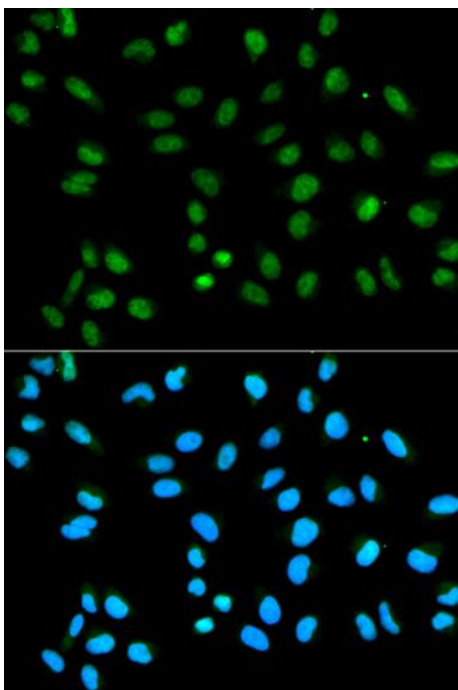
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

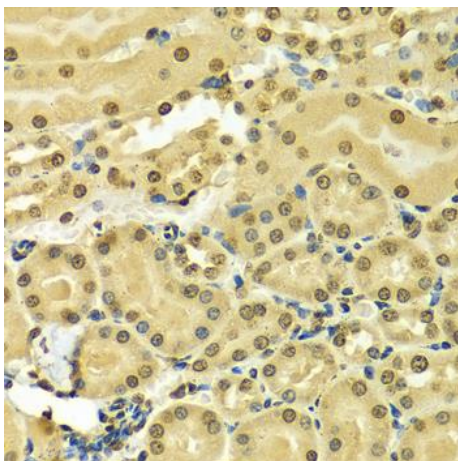
HNRNPA2; HNRNPB1; HNRPA2; HNRPA2B1; HNRPB1; IBMPFD2; RNPA2; SNRPB1; HNRNPA2B1

Application

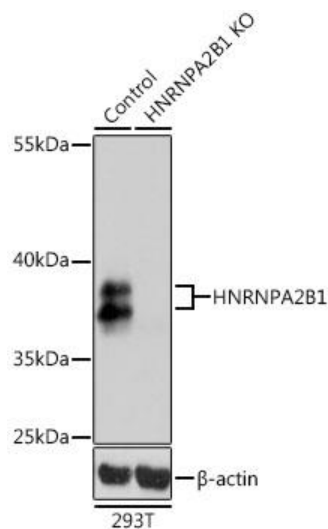
WB, IHC, IF; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:50 - 1:200



Immunofluorescence analysis of HeLa cells using HNRNPA2B1 Polyclonal Antibody. Blue: DAPI for nuclear staining.



Immunohistochemistry of paraffin-embedded mouse kidney using HNRNPA2B1 Polyclonal Antibody at dilution of 1:100 (40x lens).

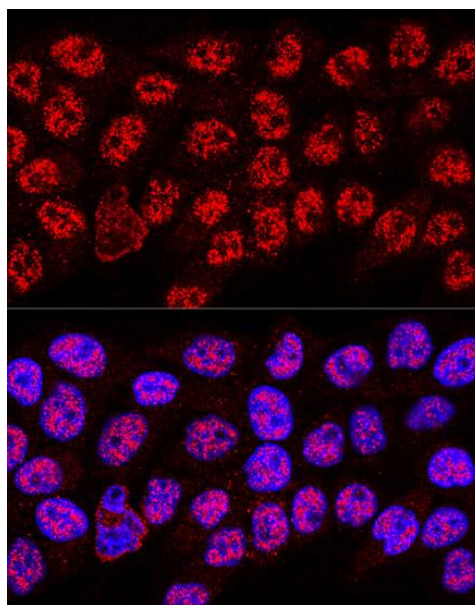


Western blot analysis of extracts from normal (control) and HNRNPA2B1 knockout (KO) 293T cells, using HNRNPA2B1 Polyclonal Antibody at 1:1000 dilution.

Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 1s.



Confocal immunofluorescence analysis of HeLa cells using HNRNPA2B1 Polyclonal Antibody at dilution of 1:400. Blue: DAPI for nuclear staining.