

hnRNPA2B1 antibody (pAb)

Catalog Nos: 65726, 65926, 65727**RRID:** AB_3216298**Application(s):** WB**Reactivity:** Human**Quantities:** 100 µg, 50 µg, 10 µg**Purification:** Protein A Chromatography**Host:** Rabbit**Isotype:** IgG**Molecular Weight:** 37 kDa

Background: hnRNPA2B1 (Heterogeneous Nuclear Ribonucleoprotein A2/B1, also known as RNPA2; HNRPA2; HNRPB1; SNRPB1; HNRNPA2; HNRNPB1; IBMPFD2; HNRPA2B1) 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. hnRNPA2B1 includes two different isoforms A2 and B1. It is reported that hnRNPA2B1 is a nuclear “reader” of the m6A mark, acting as an adaptor that recruits the Microprocessor complex to a subset of precursor miRNAs, facilitating their processing into mature miRNAs.

Immunogen: This hnRNPA2B1 antibody was raised against a fusion protein corresponding to human hnRNPA2B1.

Buffer: PBS containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

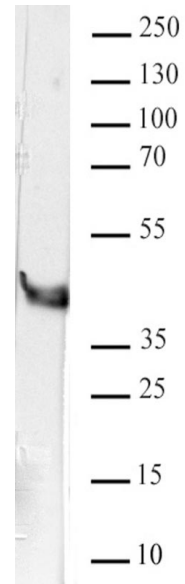
Application Notes:

Applications Validated by Active Motif:

WB: 0.5 µg/mL-2 µg/mL

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.

**hnRNPA2B1 antibody (pAb) tested by Western blot.**

REH nuclear extract (20 µg per lane) was probed with hnRNPA2B1 antibody (0.5 µg/mL).