

POLR2B antibody (pAb)

Catalog Nos: 61557, 61558

RRID: AB_2614965

Isotype: IgG

Application(s): IP, WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 135 kDa

Background: **POLR2B** (polymerase (RNA) II (DNA directed) polypeptide B, 140 kDa) DNA-dependent RNA polymerase catalyzes the transcription of DNA into RNA using the four ribonucleoside triphosphates as substrates. Second largest component of RNA polymerase II which synthesizes mRNA precursors and many functional non-coding RNAs. Proposed to contribute to the polymerase catalytic activity and forms the polymerase active center together with the largest subunit. Pol II is the central component of the basal RNA polymerase II transcription machinery. It is composed of mobile elements that move relative to each other. RPB2 is part of the core element with the central large cleft, the clamp element that moves to open and close the cleft and the jaws that are thought to grab the incoming DNA template (By similarity).

Immunogen: This antibody was raised against a peptide within the N-terminal region of human POLR2B.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

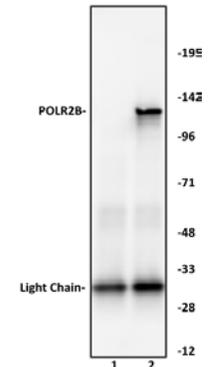
IP: 5 - 10 μ l per IP

WB*: 1:500 - 1:1,000 dilution

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

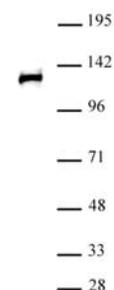
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



POLR2B antibody (pAb) tested by Immunoprecipitation.

10 μ l of POLR2B antibody was used to immunoprecipitate POLR2B from 200 μ g of HEK293 nuclear cell extract (lane 2). 10 μ l of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the POLR2B antibody at a dilution of 1:500.



POLR2B antibody (pAb) tested by Western blot.

Nuclear extract of HEK293 cells (30 μ g) probed with POLR2B antibody (pAb) at a dilution of 1:1,500.