HNRNPC antibody (pAb)



Catalog Nos: 65742, 65942, 65743 RRID: AB_3216376 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: HNRNPC (Heterogeneous Nuclear Ribonucleoprotein C) is a member of the subfamily nuclear ribonucleoproteins (hnRNPs) involved in pre-mRNA processing, mRNA metabolism, and transport. 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 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 can act as a tetramer and is involved in the assembly of 40S hnRNP particles. Multiple transcript variants encoding at least two different isoforms have been described for this gene (C1 and C2).

HNRNPC is a critical alternative splicing regulator that complexes with heterogeneous nuclear RNA (hnRNA) and plays an essential role in RNA maturation, stabilization, translocation, and posttranscriptional modification. Cellular homeostasis, among other cellular processes, is regulated by RNA alternative splicing. Alternative splicing disfunction contributes to many disorders, including cancer. The up regulation of HNRNPC expression has been tied to cancer progression and is thought to function by regulating RNA splicing, RNA expression, RNA stability, mRNA translation, m6A RNA methylation, and dsRNA induced interferon response.

Immunogen: Peptide containing a sequence fragment corresponding to human heterogeneous nuclear ribonucleoprotein C.

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: WB: 0.5 - 2 µg/ml

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.05% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

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.

Application Key: ChIP = Chromatin Immunoprecipitation; FACS = Flow Cytometry; IF = Immunofluorescence; IHC = Immunohistochemistry; IP = Immunoprecipitation; WB = Western Blot





HNRNPC antibody (pAb) tested by Western blot.

Detection of HNRNPC by Western blot. The analysis was performed using 20 μ g of HeLa nuclear extract and 0.5 μ g of HNRNPC antibody (pAb. A protein band of approximate molecular weight of 37 kDa was detected.