

HNF4A antibody (pAb)

Catalog Nos: 61189, 61190

RRID: AB_2793545

Isotype: IgG

Application(s): ICC, IF, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 60 kDa

Background: HNF4A (hepatocyte nuclear receptor 4A, NR2A1, TCF14) is an orphan nuclear receptor and activator of transcription. HNF4A is major regulator of the hepatocytic phenotype and regulates genes involved in the control of lipid homeostasis. Mutations in HNF4A can cause maturity onset diabetes of the young (MODY1). HNF4A binds DNA as a homodimer and is involved in the development of the liver, kidney and intestine.

Immunogen: This HNF4A antibody was raised against a peptide derived from the N-terminal region of human HNF4A.

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:

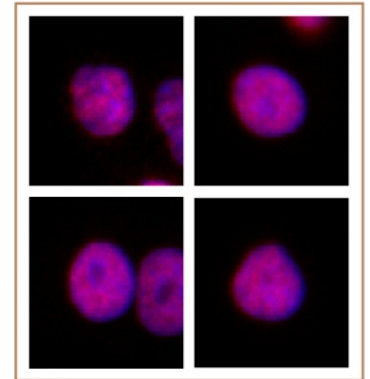
ICC/IF: 1:2,000 dilution

WB: 1:500 - 1:2,000 dilution

For optimal results, primary antibody incubations should be performed at 4°C. 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.



HNF4A antibody (pAb) tested by Immunofluorescence.

Formaldehyde fixed HeLa cells stained with HNF4A antibody at a 1:2,000 dilution.



HNF4A antibody (pAb) tested by Western blot.

Whole cell extract (20 µg) of Hep G2 cells probed with HNF4A antibody at a dilution of 1:500.