

METTL14 antibody (pAb)

Catalog Nos: 65746, 65946, 65747

RRID: AB_3216383

Application(s): IP, WB

Reactivity: Human

Volumes: 100 µl, 50 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Isotype: IgG

Molecular Weight: 55kDa

Background: N6-methylated adenine (m6A) is prevalently present in nearly all RNA types and can be found in all organisms from bacteria to humans. It preferentially appears around stop codons and within long internal exons in mammalian messenger RNAs. m6A plays an important role in the efficiency of mRNA splicing, processing, translation efficiency, editing and mRNA stability. m6A also takes place in other RNA molecules, such as primary miRNA (pri-miRNAs).

METTL14 (methyltransferase-like 14) forms a stable N6-methyltransferase heterodimer complex with METTL3, which catalyzes the generation of m6A modification on mammalian nuclear RNAs. METTL3 is the catalytically active subunit, while **METTL14** plays a structural role critical for substrate recognition.

Immunogen: This antibody was raised against a human peptide containing methyltransferase 14.

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

Application Notes:

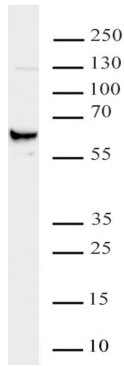
Applications Validated by Active Motif:

WB: 1:500 - 1:2000 dilution

IP: 4 µl per IP

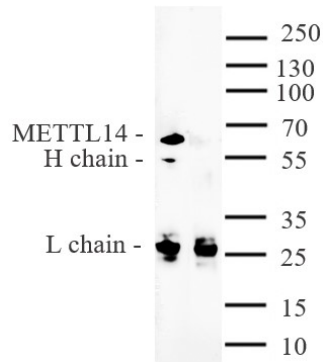
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



METTL14 antibody (pAb) tested by Western blot.

Nuclear extract of HepG2 (20 μ g / lane) was probed with METTL14 antibody (pAb) at a dilution of 1:1000.



METTL14 antibody (pAb) tested by Immunoprecipitation

500 μ l of HeLa cell nuclear extract was immunoprecipitated with 4 μ l of METTL14 antibody (Lane 1) or Rabbit IgG (Lane 2). The immunoprecipitated protein was detected by western blot using METTL14 antibody at 0.5 μ g/ml. On the gel, antibody heavy chains (50 kDa) and light chains (25 kDa) are visible.