

## METTL3 antibody (mAb)

**Catalog Nos:** 65687, 65688

**Clone:** 29C8

**Isotype:** IgG2a

**Application(s):** IP, WB

**Reactivity:** Human

**Volumes:** 100 µl, 10 µl

**Purification:** None

**Host:** Mouse

**Concentration:** N/A µg/µl

**Molecular Weight:** 72 kDa

**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 occurs in other RNA molecules, such as primary miRNA (pri-miRNAs).

METTL3 (methyltransferase-like 3, also known as IME4, M6A, MT-A70) forms a stable N6-methyltransferase heterodimer complex with METTL14, 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. METTL3 promotes translation by interacting with the translation initiation machinery in the cytoplasm. Its overexpression in a number of cancer cells suggests that it may participate to cancer cell proliferation by promoting mRNA translation.

**Immunogen:** This antibody was raised against a peptide within the N-terminal region of human METTL3 (N6-adenosine-methyltransferase).

**Buffer:** This antibody is provided as a hybridoma culture supernatant with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

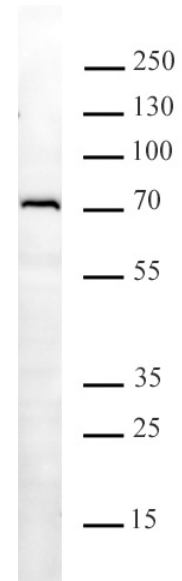
Applications Validated by Active Motif:

WB: 1:500 - 1:1000 dilution

This antibody is also available as an AbFlex® engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91333.

**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.



### METTL3 antibody (mAb) tested by Western blot

40 µg of K562 cytoplasmic extract was run on SDS-PAGE and probed with antibody at a dilution of 1:500. In addition, we recommend the addition of 0.05% Tween 20 to all blocking solutions to reduce background.