

N6-Methyladenosine (m6A) antibody (mAb)

Catalog Nos: 61755, 61756

RRID: AB_2793759

Clone: 17-3-4-1

Isotype: IgG1, k

Application(s): DB, IP

Reactivity: Budding Yeast, Human, Not Species Specific

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: Not Applicable

Background: N6-Methyladenosine (m6A) is an RNA modification on the N-6 position of adenosine. This modification has been found to be abundant in the 3' UTR and stop codons of mammalian mRNA. m6A is associated with miRNA binding sites suggesting a potential role in epigenetic gene regulation. FTO and ALKBH are demethylases for 6-methyladenosine while a multiprotein complex that includes METTL3 functions as the methyltransferase. Recent findings revealed that m6A is also present on metazoan DNA, suggesting a genuine epigenetic role for this modification in the context of DNA as well.

Immunogen: This antibody was raised against 6-Methyladenosine conjugated to BSA.

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

Application Notes:

For N6-Methyladenosine (m6A), we also offer an AbFlex[®] N6-Methyladenosine (m6A) Recombinant Antibody (rAb). For details, see Catalog No. 91261.

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.