

PRODUCT INFORMATION

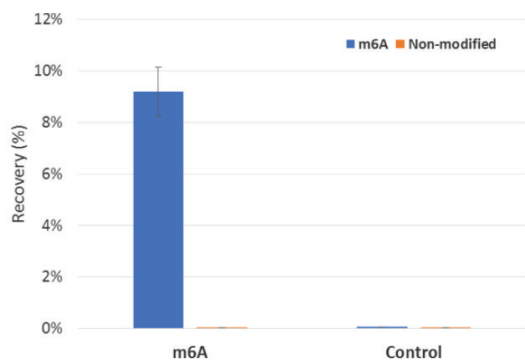
IgG4 Fc (human) Rabbit Monoclonal Antibody - Biotinylated (Clone RM362)

Item No. 32164

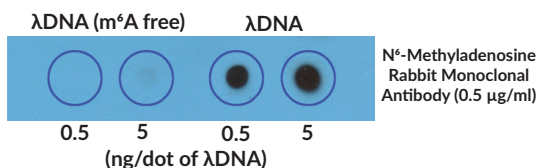
Overview and Properties

Contents:	This vial contains 50 µg of protein A-affinity purified monoclonal antibody.
Synonym:	Immunoglobulin G4
Immunogen:	Peptide from the Fc region of human IgG4
Cross Reactivity:	(+) IgG4; (-) Human IgG1, IgG2, IgG3, IgM, IgA, IgD, IgE; (-) Goat, mouse, rat IgG
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM217
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA; the recommended starting concentration is .05-0.2 µg/ml for ELISA. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



MeRIP was performed using m⁶A and control RNA from NEB (E1610). Antibodies (N⁶-Methyladenosine Rabbit Monoclonal Antibody (Clone RM362) and Polyethylene Glycol) were mixed for a short time, with Protein A beads before the addition of m⁶A labelled and non-labelled RNA in equal amounts. Captured RNA was then washed, eluted and run through RT-PCR. Resulting cDNA was then analyzed using real time PCR with specific probes to m⁶A labelled and non-labelled sequences.



DB of λDNA with or without m⁶A using N⁶-Methyladenosine Rabbit Monoclonal Antibody (Clone RM362). The membrane was pre-spotted with 5 and 0.5 ng/dot of λDNA. m⁶A-free DNA was isolated from bacteriophage λ grown in an adenine methylase-deficient *E. coli* host.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/15/2024

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

N⁶-Methyladenosine (m⁶A) is an abundant post-transcriptional RNA modification that is found in various classes of eukaryotic or viral RNA, including mRNA, rRNA, tRNA, snRNA, microRNA (miRNA), and long non-coding RNA (lncRNA).^{1,2} m⁶A modifications are enriched near stop codons of mRNA transcripts, increasing along the length of the coding region and decreasing after 3'-UTRs.^{2,3} m⁶A formation is a reversible process that occurs primarily in the nucleus.⁴ Its formation is catalyzed by a methyltransferase complex that is composed of the writers METTL3 and METTL14, which preferentially methylate RNA substrates with a GGACU domain, and the accessory protein Wilms tumor 1-associating protein (WTAP), which recruits target RNAs.² m⁶A demethylation is mediated by the erasers fat mass and obesity-associated (FTO) protein and ALKBH5. Upon export into the cytosol, m⁶A modifications are recognized by a variety of reader proteins, including YTH-domain family (YTHDF) members, which regulate RNA translation. m⁶A modifications regulate multiple steps of RNA processing, including pre-mRNA splicing, nuclear export, translation, and decay, influencing a variety of critical cellular functions, such as differentiation, neurodevelopment, immunity, and oncogenesis.^{1,2,5} Global m⁶A mRNA levels are decreased in tumor tissue isolated from patients with endometrial cancer.⁶ Cayman's N⁶-Methyladenosine Rabbit Monoclonal Antibody (Clone RM362) can be used for dot blot (DB), ELISA, and methylated RNA immunoprecipitation (MeRIP) applications.

References

1. Yang, C., Hu, Y., Zhou, B., *et al.* The role of m⁶A modification in physiology and disease. *Cell Death Dis.* **11(11)**, 960 (2020).
2. Zhang, C., Fu, J., and Zhou, Y. A review in research progress concerning m⁶A methylation and immunoregulation. *Front Immunol.* **10**, 922 (2019).
3. Meyer, K.D., Saletore, Y., Zumbo, P., *et al.* Comprehensive analysis of mRNA methylation reveals enrichment in 3' UTRs and near stop codons. *Cell* **149(7)**, 1635-1646 (2012).
4. Zaccara, S., Ries, R.J., and Jaffrey, S.R. Reading, writing and erasing mRNA methylation. *Nat. Rev. Mol. Cell Biol.* **20(10)**, 608-624 (2019).
5. Sun, T., Wu, R., and Ming, L. The role of m⁶A RNA methylation in cancer. *Biomed. Pharmacother.* **112**, 108613 (2019).
6. Liu, J., Eckert, M.A., Harada, B.T., *et al.* m⁶A mRNA methylation regulates AKT activity to promote the proliferation and tumorigenicity of endometrial cancer. *Nat. Cell Biol.* **20(9)**, 1074-1083 (2018).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM