

m⁶A RNA Methylation Assay Kit (Fluorometric) ab233491

[2 References](#) [1 Image](#)

Overview

Product name	m ⁶ A RNA Methylation Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Other biological fluids, Tissue, Adherent cells, Suspension cells
Assay type	Quantitative
Sensitivity	5 pg/well
Assay time	3h 45m
Product overview	<p>m⁶A RNA Methylation Assay Kit (Fluorometric) (ab233491) is a complete set of optimized buffers and reagents to fluorometrically quantify methylated N⁶-methyladenosine (m⁶A) in RNA. It is suitable for a direct detection of m⁶A RNA methylation status using total RNA isolated from any species such as mammals, plants, fungi, bacteria and viruses.</p>

This kit contains a unique binding solution allowing RNA >70 nts to be tightly bound to the wells, which enables quantification of m⁶A from both mRNA and nc-RNA such as tRNA, rRNA and snRNA. The optimized antibody and enhancer solutions allow high specificity to m⁶A, with no cross-reactivity to unmethylated adenosine within the indicated concentration range of the sample RNA. Also included are universal positive and negative controls which are suitable for quantifying m⁶A from any species.

Notes

N⁶-methyladenosine (m⁶A) is the most common and abundant modification in RNA molecules present in eukaryotes. The m⁶A modification is catalyzed by a methyltransferase complex METTL3 and removed by the recently discovered m⁶A RNA demethylases FTO and ALKBH5, which catalyze m⁶A demethylation in an α -ketoglutarate (α -KG)- and Fe²⁺-dependent manner. It was shown that METTL3, FTO, and ALKBH5 play important roles in many biological processes, ranging from development and metabolism to fertility. m⁶A accounts for more than 80% of all RNA base methylations and exists in various species. m⁶A is mainly distributed in mRNA and also occurs in non-coding RNA such as tRNA, rRNA, and snRNA. The relative abundance of m⁶A in mRNA transcripts has been shown to affect RNA metabolism processes such as splicing, nuclear export, translation ability and stability, and RNA transcription. Abnormal m⁶A methylation levels induced by defects in m⁶A RNA methylase and demethylase could lead to dysfunction of RNA and diseases. For example, abnormally low levels of m⁶A in target mRNAs due to increased FTO activity in patients with FTO mutations, through an as-yet undefined pathway, contributes to the onset of obesity and related diseases. The dynamic and reversible chemical m⁶A modification in

RNA may also serve as a novel epigenetic marker of profound biological significance. Therefore, more useful information for a better understanding of m⁶A RNA methylation levels and distribution on RNA transcripts could benefit diagnostics and therapeutics of disease.

Platform Microplate reader

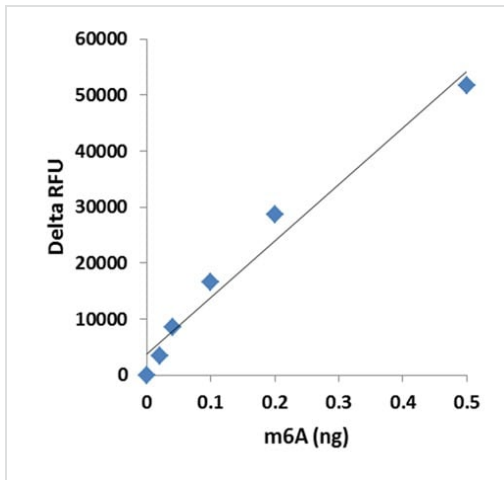
Properties

Storage instructions Please refer to protocols.

Components	1 x 48 tests	1 x 96 tests
10X Wash Buffer	1 x 14ml	1 x 28ml
8-Well Assay Strips (With Frame)	6 units	12 units
Binding Solution	1 x 5ml	1 x 10ml
Capture Antibody, 1000 X	1 x 5µl	1 x 10µl
Detector Antibody, 1000 X	1 x 6µl	1 x 12µl
Dilution Buffer	1 x 4ml	1 x 8ml
Enhancer Solution	1 x 5µl	1 x 10µl
Fluoro Developer	1 x 8µl	1 x 16µl
Fluoro Enhancer	1 x 8µl	1 x 16µl
Negative Control, 100 µg/mL	1 x 10µl	1 x 20µl
Positive Control, m6A 2 µg/mL	1 x 10µl	1 x 20µl

Relevance N⁶-Methyladenosine (m⁶A) is an abundant modification found in mRNA, tRNA, snRNA, as well as long non-coding RNA, in all species. RNA adenosine methylation is catalyzed by a multicomponent complex composed of METTL3/MT-A70, METTL14, and WTAP in mammals. METTL3 & METTL14 are responsible for the methyltransferase activity of the complex, and WTAP mediates substrate recruitment.

Images



Example of a standard curve generated with m6A RNA Methylation Assay Kit (ab233491).

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors