

m1A (N1-methyladenosine) RNA Enrichment and Quantification Kit

Catalog #: QMERIP-M1A

User Manual

Last Revised: October 14th, 2024

Introduction

N1-methyladenosine (m1A), the methylation of adenosine in the N1 position, is a prevalent and reversible post-transcriptional RNA modification that can occur on messenger RNA (mRNA), transfer RNA (tRNA) and ribosomal RNA (rRNA). Recent studies have revealed the crucial role of m1A RNA modification in multiple cellular processes, including RNA structural stability, folding, interactions with proteins, cell viability, impaired self-renewal ability, cell proliferation, and cell death. Methylated RNA immunoprecipitation (MeRIP) is a popular method to observe and map the location of RNA modifications in specific genes as well as transcriptome-wide.

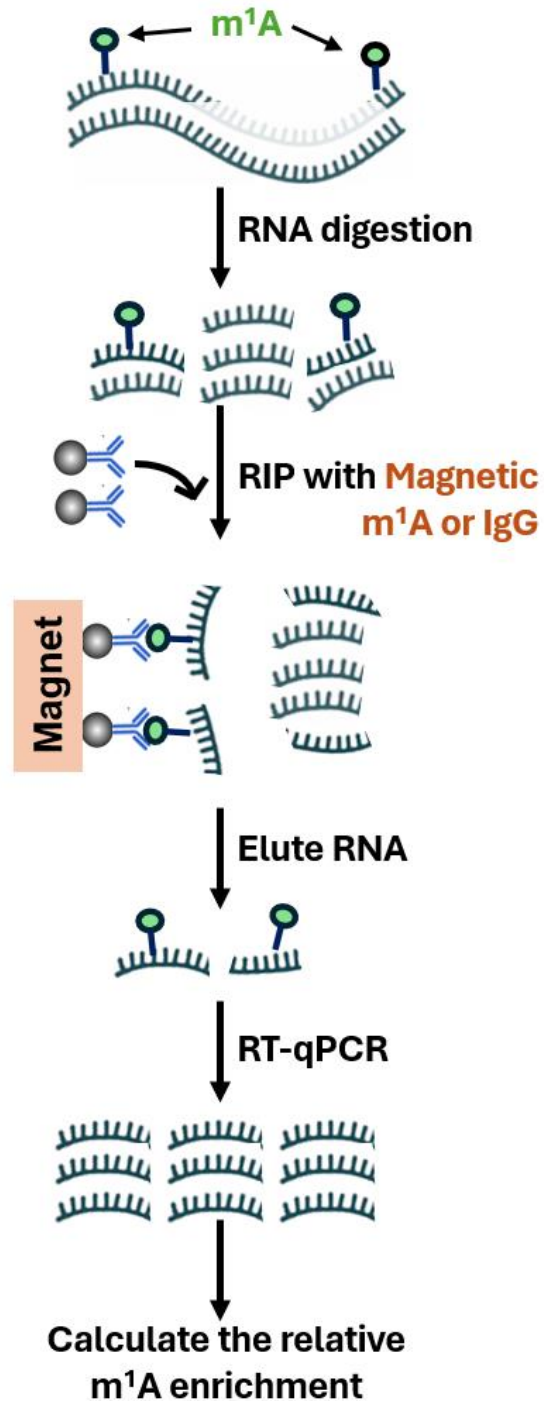
The RayBio® m1A (N1-methyladenosine) RNA Enrichment and Quantification Kit (QMERIP-M1A) identifies the abundance and enrichment location of m1A by MeRIP and quantitative real-time PCR. The enriched m1A RNA could also be used for transcriptome-wide profiling of m1A. In this assay, RNA sample is digested into fragments consisting of 100~500 nucleotides followed by incubation with an antibody against m1A conjugated to a magnetic bead. Then, the enriched RNA is released, purified, and eluted for RT-qPCR analysis. Included in the kit is a non-immune IgG control conjugated to a magnetic bead and two pairs of primer controls for HeLa cell RNA. QMERIP-M1A uses a fast and streamlined procedure so the assay can be finished in 5 hours with a hands-on time of less than 1 hour. This kit has been validated for human, mouse, and rat m¹A enrichment.

Product Use

QMERIP-M1A is for research use only. It is not approved for application in clinical or in vitro diagnostic procedures.

How It Works

Figure 1. QMERIP-M1A workflow



Storage / Stability

The entire kit can be stored at -20°C or -80°C for a period of 12 months prior to opening. After opening, the reagents are valid for at least 6 months if stored as described in the table below. Avoid repeated freeze-thaw cycles.

Kit Components

| Name | Catalog # | Size / Qty | Storage After Opening |
|---------------------------------------|------------|-----------------|-----------------------|
| RNA Digestion Buffer (10X) | RIP-DB | 1 mL / 1 tube | -20°C |
| Magnetic bead-conjugated m1A antibody | RIP-M1A | 0.3 mL / 1 tube | -20°C |
| Magnetic bead-conjugated IgG antibody | RIP-IGG | 20 µL / 1 tube | -20°C |
| RNase Inhibitor | RIP-RI | 0.6 mL / 1 tube | -20°C |
| ACTB (Negative) Primers* | RIP-NP-M1A | 20 µL / 1 tube | -20°C |
| APT5A (Positive) Primers* | RIP-PP-M1A | 20 µL / 1 tube | -20°C |
| One-Step RT Master Mix (2X) | RIP-RTM | 1 mL / 1 tube | -20°C |
| IP Buffer (5X) | RIP-IB | 30 mL / 1 tube | 4°C |
| Elution Buffer | RIP-EB | 4 mL / 1 tube | 4°C |
| 0.5 M EGTA | RIP-EGTA | 0.2 mL / 1 tube | 4°C |
| NaOH | RIP-NA | 0.6 mL / 1 tube | 4°C |
| Nuclease-Free H ₂ O | RIP-H2O | 1 mL / 1 tube | 4°C |

*Primers included in the kit have only been validated for Hela cell RNA. If running a different RNA sample type, you will need to supply your own primers for the target gene of interest

Additional Materials Required

- Fluorescence PCR instrument and compatible PCR plate or tubes
- Sterile nuclease-free pipette tips (barrier tips recommended) and microfuge tubes
- Vortex Mixer
- Magnetic Separator
- Total RNA, mRNA, tRNA or rRNA
- Primers for target gene of interest, with or without probe

Important Notes Before Starting

Please carefully read through the entire protocol before starting this assay. This protocol requires multiple steps but can be finished in one day. The hands-on time is less than 1 hour. The approximate time required for each step is provided in **Table 2** below.

Table 2. Procedure Overview & Estimated Timetable

| Protocol Step | Time Required | Stopping Points and Protocol Notes |
|-----------------|---------------|---|
| RNA Digestion | ~50 minutes | Can pause after this step and store fragmented RNA at -80°C |
| RIP and Elution | ~1.5 hours | Can pause after this step and store eluted RNA at -80°C |
| RT-qPCR | ~2.0 hours | Continue to data analysis |
| Data Analysis | ~0.5 hours | $\Delta\Delta Cq$ Calculations |

Reagent Preparation

Place all reagents on ice for 30 minutes to thaw before use. Then, mix gently with a vortex and briefly centrifuge at 1,500x g for 30 seconds to collect contents at the bottom.

RNA Sample Preparation: Extract or purify total RNA, tRNA, rRNA or mRNA from virous cells or tissues. Using high-quality and intact RNA is critical as degraded RNA can affect the quality of the assay. We recommend checking the RNA integrity by Bioanalyzer or agarose gel before starting the assay. The total RNA, tRNA, rRNA or mRNA input amount can range from 5 μ g to 50 μ g for each sample (this includes the 3 reactions per sample:

Input, m1A-IP and IgG-IP). For optimal quantification, we recommend 20 µg per sample for total RNA, and 10 µg for the other RNA types (tRNA, rRNA and mRNA).

Assay Procedure

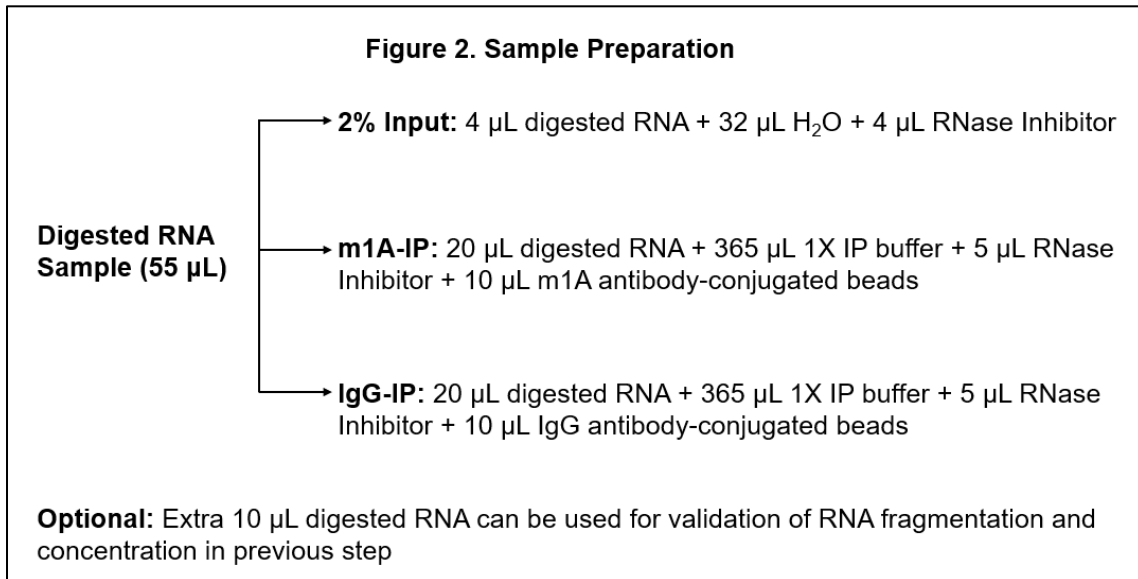
1. RNA Digestion

- 1.1. Adjust the RNA sample to a total volume of 45 µL in a 200-µL PCR tube with nuclease-free water. Add 5 µL of 10X RNA Digestion Buffer (RIP-DB). Mix well by pipetting up and down and spin down the tubes to collect contents to the bottom.
- 1.2. Incubate the RNA samples at 37°C in the thermal cycler block for 30 minutes. Immediately remove the tubes from the block and add 5 µL of 0.5 M EGTA (RIP-EGTA) to each tube. Vortex and spin down the tube and place it on ice.
- 1.3. **(Optional)** Validate digested RNA size distribution by running 0.5 µg of the digested RNA sample on a 2% (wt/vol) agarose gel or Agilent 2100 Bioanalyzer with an Agilent RNA 6000 Pico kit. The distribution of RNA fragment sizes should be centered around 100~500 bp. Validate digested RNA concentration by NanoDrop.

2. m1A RNA Immunoprecipitation (MeRIP)

- 2.1. **Sample Preparation (see Figure 2 for reference):** For each RNA sample, place three new 1.5 mL Eppendorf tubes on ice and label them “2% Input” (see below for calculation), “m1A-IP”, and “IgG-IP”. Into the “2% Input” tube, pipette 4 µL of digested RNA sample, then add 32 µL of nuclease-free H₂O and 4 µL of RNase inhibitor (RIP-RI). Put the “2% Input” tube on ice or store at -20°C until Step 3.1 of RT-qPCR. For the “m1A-IP” and “IgG-IP” tubes, pipette 20 µL of digested RNA sample into each tube.

2% Input Calculation: [(4 µL of digested RNA / 20 µL digested RNA for m1A-IP reaction / 10 (Dilution factor) x 100% = 2%)]



- 2.2. Dilute 5X IP buffer (RIP-IB) with nuclease-free H₂O to a 1X solution. Add 365 µL 1X IP buffer to the “m1A-IP” and “IgG-IP” tubes. Then, add 5 µL RNase inhibitor (RIP-RI) to each tube.
- 2.3. Thoroughly resuspend the m1A antibody-conjugated beads (RIP-M1A) and IgG antibody-conjugated beads (RIP-IGG) by end-over-end rotation or by pipetting up and down. No clumps of beads should be visible. Add 10 µL m1A antibody-conjugated beads and 10 µL IgG antibody-conjugated beads into the “m1A-IP” and “IgG-IP” tubes, respectively.
- 2.4. Incubate the “m1A-IP” and “IgG-IP” tubes with rotation at low speed for 60 minutes at room temperature.
- 2.5. Centrifuge the tubes briefly and place them on the magnetic separator for 1 minute, then carefully remove the supernatant.

Caution: Be careful not to disturb or discard the beads that contain the RNA

- 2.6. Remove the tubes from the magnet. Add 0.5 mL 1X IP Buffer to each tube and mix gently by pipetting up and down several times to completely resuspend the antibody-

conjugated beads. Place the tubes on the magnetic separator for 1 minute, then carefully remove the supernatant.

- 2.7. Repeat the above step for two additional washes. Be sure to remove the final IP Buffer leaving only the antibody-conjugated beads.
- 2.8. Remove the tubes from the magnet and place them on ice. Add 30 μ L of Elution Buffer and 5 μ L of RNase inhibitor (RIP-RI) to each tube and mix by gently pipetting up and down several times to completely resuspend the antibody-conjugated beads.
- 2.9. Centrifuge the tubes briefly to collect contents to the bottom of the tube. Place the tubes on the magnetic separator for 1 minute.
- 2.10. Transfer the supernatant containing eluted RNA fragments to new 1.5 mL microcentrifuge tubes. Take special care to not aspirate the beads, as it will increase background noise.
- 2.11. Add 5 μ L NaOH (RIP-NA) into all IP tubes. Continue to RT-qPCR or store all RNA samples at -80°C until further use.

3. RT-qPCR Assay for m1A-RIP

The m1A enriched RNA can now be detected by quantitative RT-qPCR assay. This kit includes APT5A (encodes a subunit of mitochondrial ATP synthase) as a positive control (RIP-PP-M1A) and ACTB (Actin Beta) as a negative control (RIP-NP-M1A) for m1A modification in HeLa cells. Calculation of each m1A enriched RNA sample can be performed using the comparative C_q ($\Delta\Delta\text{C}_q$) method with three PCR amplicons: 2% Input, m1A-IP and IgG-IP. In addition, we recommend verifying successful RNA retrieval by running a gel after RT-qPCR.

Note: The included APT5A and ACTB primers have only been validated for HeLa cell RNA. We do not recommend using the included primers with other RNA types.

Optional: If you are running a different RNA sample type and you want to ensure the assay is functioning properly, you can include a separate HeLa cell RNA sample with the included primers for positive and negative controls.

3.1. Prepare the RT-qPCR reaction mix as stated in **Table 3**. Use enough reagents for two extra reactions to account for loss of volume. We recommend running all RNA samples (each digested RNA sample includes 2% Input, m1A-IP and IgG-IP) in triplicate. So, each RNA sample will have 9 reactions (3 for each condition).

Calculate the number of reactions needed = [1 (2% Input) + 1 (m1A-IP) + 1 (IgG-IP)] x n (number of digested RNA samples) x 3 (triplicates) + 2 (To account for loss of volume).

Example: For 3 samples in triplicate, total reactions = (1+1+1) x 3 x 3 + 2 = 29.

Table 3: RT-qPCR Reaction Components

| Components | Volume per Reaction | Master Mix for 29 Reactions |
|-----------------------------|----------------------------|------------------------------------|
| One-Step RT Master Mix (2X) | 10 µL | 290 µL |
| Forward primer (10 µM)* | 1 µL | 29 µL |
| Reverse primer (10 µM)* | 1 µL | 29 µL |
| Nuclease-free Water | 6 µL | 174 µL |
| RNA Sample | 2 µL | Add to each tube individually |
| Total Volume | 20 µL | 522 µL |

*Primers included in the kit have only been validated for HeLa cell RNA. If running a different RNA sample type, you will need to supply your own primers for the target gene of interest.

3.2. Prepare a master mix of all components except RNA according to the table. Mix thoroughly by gently pipetting up and down. Collect liquid to the bottom of the tube by brief centrifugation.

- 3.3. Aliquot 18 μ L master mix into qPCR tubes or plate. For the best results, ensure accurate and consistent pipetting volumes and minimize bubbles.
- 3.4. Add 2 μ L RNA sample to the appropriate qPCR tubes or plate. Seal tubes with flat, optically transparent caps; seal plate with optically transparent film. Take care to properly seal plate edges and corners to prevent artifacts caused by evaporation.
- 3.5. Spin tubes or plate briefly to remove bubbles and collect liquid (1 minute at 2,500 – 3,000 rpm).
- 3.6. Program qPCR instrument with indicated thermocycling protocol (see **Table 4** below).

Use the SYBR (or Taqman™) mode setting on the qPCR instrument. Refer to **Table 4** for SYBR PCR program setup (Use the Applied Biosystems QuantStudio 5 Real-Time PCR System as an example).

Table 4. RT-qPCR program (For SYBR model)

| | Steps | Temperature (°C) | Time | Cycles |
|--------|---|-------------------------|-------------|---------------|
| Step 1 | Reverse Transcription | 55 | 10 minutes | 1 |
| Step 2 | Initial Denaturation | 95 | 3 minutes | 1 |
| Step 3 | Denature | 95 | 10 seconds | 40-45 |
| | Anneal, extend, and detect fluorescence | 60 | 60 seconds | |
| Step 4 | Melt Curve (SYBR model) | 60-95 | Various | 1 |

- 3.7. Results and Export: In the analysis settings, select automatic baseline, or select according to your lab validation. Then, export the results as an excel file.

4. Calculate the Relative m1A Enrichment

4.1. Calculate the mean Cq value (Quantification Cycle Value) for each set of triplicates: 2% Input, m1A-IP and IgG-IP samples. Then, use the comparative $\Delta\Delta Cq$ method to analyze the relative m1A enrichment.

4.2. Normalize Cq values of the 2% Input, m1A-IP and IgG-IP samples to the average ΔCq of 2% input samples: $\text{Normalized } \Delta Cq = Cq_{IP} - Cq_{\text{Average 2\% Input}}$

4.3. Calculate the percent of 100% Input for each sample: $2^{-\Delta Cq (\text{normalized IP})} \times 50 \times 100\%$

4.4. The relative m1A enrichment measured by the m1A-IP over the IgG-IP:

$$= 2^{-\Delta\Delta Cq} = 2^{-[\Delta Cq (\text{normalized m1A-IP}) - \Delta Cq (\text{normalized IgG-IP})]}$$

or

$$= 2^{-\Delta Cq (\text{normalized m1A-IP})} \times 100\% / 2^{-\Delta Cq (\text{normalized IgG-IP})} \times 100\%$$

Normally, relative m1A enrichment ≥ 8 means significant enrichment with P value <0.05 by appropriate statistical analysis. See example calculations in Typical Data section below.

Assay Procedure Summary

1. Prepare All Reagents, Samples and Standards as Instructed
2. RNA Digestion
3. m1A RNA Immunoprecipitation (RIP)
4. RT-qPCR Assay for m1A-RIP
5. Calculate the Relative m1A Enrichment

Typical Data: Example of QMERIP-M1A Analysis:

The individual and mean Cq value for each set of triplicates samples shown in table below.

| Samples | Replicates | ACTB | | | APT5A | | |
|-----------|----------------|---------------|---------------|---------------|---------------|---------------|---------------|
| | | 2% Input | IgG-IP | m1A-IP | 2% Input | IgG-IP | m1A-IP |
| Sample #1 | Rep-1 | 16.500 | 32.731 | 32.344 | 18.137 | 30.062 | 26.075 |
| | Rep-2 | 16.763 | 34.958 | 32.559 | 18.266 | 30.239 | 26.251 |
| | Rep-3 | 16.295 | 34.411 | 32.153 | 18.668 | 29.869 | 26.827 |
| | Average | 16.520 | 34.033 | 32.352 | 18.357 | 30.057 | 26.384 |
| Sample #2 | Rep-1 | 16.915 | 34.532 | 32.301 | 19.542 | 29.730 | 25.936 |
| | Rep-2 | 17.590 | 34.776 | 32.275 | 18.833 | 29.617 | 25.946 |
| | Rep-3 | 16.751 | 34.433 | 32.988 | 19.122 | 29.571 | 25.858 |
| | Average | 17.085 | 34.580 | 32.521 | 19.166 | 29.639 | 25.913 |
| Sample #3 | Rep-1 | 17.808 | 32.433 | 32.858 | 20.084 | 31.295 | 26.700 |
| | Rep-2 | 17.082 | 32.777 | 31.860 | 19.573 | 31.424 | 26.342 |
| | Rep-3 | 17.681 | 30.081 | 32.108 | 19.839 | 30.826 | 26.280 |
| | Average | 17.524 | 31.764 | 32.276 | 19.832 | 31.182 | 26.441 |

The average of ACTB Cq _{2% Input} for 3 samples is: $(16.520 + 17.085 + 17.524) / 3 = 17.043$

The average of APT5A Cq _{2% Input} for 3 samples is: $(18.357 + 19.166 + 19.832) / 3 = 19.118$

Normalized $\Delta Cq = Cq_{IP} - Cq_{Average\ 2\% Input}$ shown in the table below.

| Samples | ACTB | | | APT5A | | |
|-----------|----------|--------|--------|----------|--------|--------|
| | 2% Input | IgG-IP | m1A-IP | 2% Input | IgG-IP | m1A-IP |
| Sample #1 | -0.523 | 16.990 | 15.309 | -0.761 | 10.939 | 7.266 |
| Sample #2 | 0.043 | 17.537 | 15.478 | 0.048 | 10.521 | 6.795 |
| Sample #3 | 0.481 | 14.721 | 15.233 | 0.714 | 12.064 | 7.323 |

The percentage of RIP by Input: $2^{-\text{normalized } \Delta Cq} \times 50 \times 100\%$ shown in the table below:

| Samples | ACTB | | | APT5A | | |
|-----------|----------|--------|--------|----------|--------|--------|
| | 2% Input | IgG-IP | m1A-IP | 2% Input | IgG-IP | m1A-IP |
| Sample #1 | 143.724 | 0.038 | 0.123 | 169.510 | 2.547 | 32.481 |
| Sample #2 | 97.093 | 0.026 | 0.110 | 96.760 | 3.402 | 45.019 |
| Sample #3 | 71.660 | 0.185 | 0.130 | 60.969 | 1.168 | 31.233 |

Result: The relative m1A enrichment = The ratio of percentage of m1A-IP to IgG-IP

The relative m1A enrichment of ACTB in Sample #1 = $0.123/0.038 = 3.237$

The relative m1A enrichment of APT5 in Sample #1 = $32.481/2.547 = 12.753$

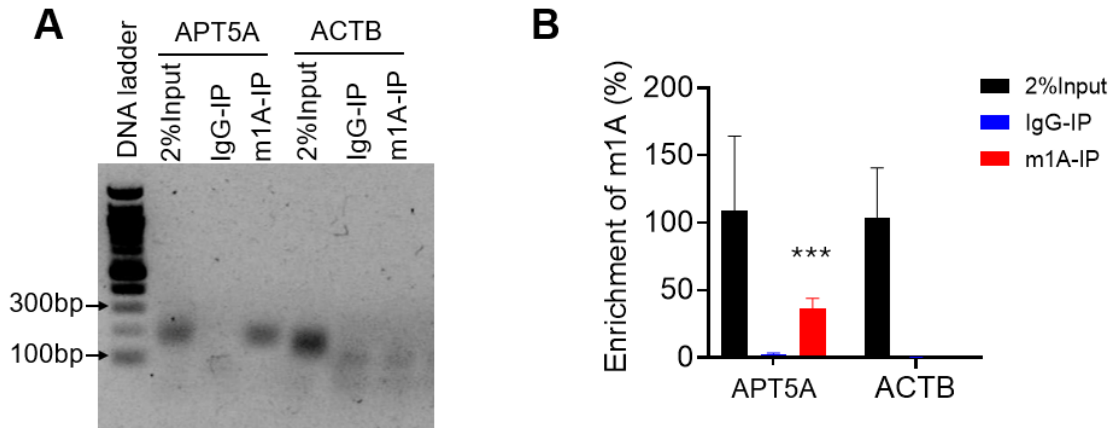


Figure 3. QMERIP-M1A enrichment test by PCR and RT-qPCR. (A) Representative images of agarose gel running after PCR amplification for the enrichment of m1A targeted APT5A and ACTB mRNA. (B) Bar graph showing the percentage enrichment of m1A targeted APT5A and ACTB mRNA using RT-qPCR assays. 2% Input RNA sample before IP was used as the positive control and IgG-IP was used as the negative control. Almost no PCR products were amplified from the m1A-IP ACTB RNA samples. Error bars, SD ($n = 3$ independent experiments). *** P -value <0.001 compared to IgG-IP was calculated using the student's t -test.

Precision

Intra-assay CV: $<5\%$

Inter-assay CV: $<20\%$

TROUBLESHOOTING

| Problem | Cause | Solution |
|---|---|--|
| Little or no enriched RNA | Insufficient amount of qualified RNA or m1A-containing RNA | Use a higher amount of RNA |
| | Poor enrichment with antibody | Increase incubation time of the antibody with RNA to overnight at 4°C |
| | Inappropriate RNA fragmenting condition | RNA digestion enzyme contained in buffer may be degraded due to improper storage. Ensure the proper storage conditions are used for this component |
| | Improper primer or probe for PCR | Redesign the primer with/without probe, check m1A occupancy in your targeting region or gene |
| | RNA degraded | Use RNase inhibitor as recommended in this protocol. Make sure that all working conditions are RNase-free and RNases are not being introduced. Use RNase-inactivating reagents to ensure the work area and materials are RNase-free |
| | Incorrect temperature and/or insufficient time during RNA release | Ensure that proper incubation times and temperatures described in the protocol are followed correctly |
| No difference in enriched RNA between m1A-IP and IgG-IP | Improper washing during RIP steps | Check if washing is performed according to the protocol. Increase wash time at each wash step or add an additional wash step |
| | Not enough fragmentation | Optimize fragmentation conditions to obtain appropriately sized fragments |