

Methylamp™ RNA Bisulfite Conversion Kit

Base Catalog # P-9003

PLEASE READ THIS ENTIRE USER GUIDE BEFORE USE

Uses: The converted RNA obtained with the Methylamp™ RNA Bisulfite Conversion Kit is suitable for various downstream RNA methylation analyses including methylation specific RT-PCR, MS-HRM, and bisulfite-sequencing including pyrosequencing and deep-sequencing.

Input RNA: The amount of RNA for each bisulfite reaction can be 5 ng-1 µg. For an optimal reaction, the input RNA amount should be 200-500 ng. When using the Methylamp™ RNA Bisulfite Conversion Kit for methylation-specific RT-PCR with very small amounts of input RNA (<10 ng), the number of PCR cycles should be greater than 45. The yield of RNA purified after bisulfite conversion depends on the amount of input RNA, nature of RNA, and source of the starting material.

Starting Material: RNA isolated from various tissue or cell samples can be used as starting material.

Precautions: To avoid cross-contamination, the following precautions are necessary for handling EpigenTek F-Spin Columns: Carefully pipette the sample or solution into the F-Spin Column. Use aerosol-barrier pipette tips and always change pipette tips between liquid transfers. Always cap the F-Spin Columns before placing them in a microcentrifuge. Wear gloves throughout the entire procedure. In case of contact between gloves and sample, change gloves immediately.

KIT CONTENTS

Component	50 reactions Cat. #P-9003-050	Storage Upon Receipt
Conversion Buffer	8 ml	RT
Conversion Powder	5 vials	RT
NA Binding Solution	13 ml	RT
F-Spin Column*	50	RT
F-Collection Tube	50	RT
Desulphonation Solution	300 μ l	RT
Control Primer-F (10 μM)	10 μ l	-20°C
Control Primer-R (10 μM)	10 μ l	-20°C
Elution Buffer	1.2 ml	RT

* Always cap spin columns before placing them in the microcentrifuge.

SHIPPING & STORAGE

The kit is shipped on frozen ice packs at 4°C.

Upon receipt: (1) Store the following components at -20°C immediately: **Control Primer-F** and **Control Primer-R**. Store all other components at room temperature.

The kit is stable for up to 6 months from the shipment date, when stored properly.

MATERIALS REQUIRED BUT NOT SUPPLIED

- Thermocycler with heated lid*

**Since the bisulfite reaction is not overlaid with mineral oil, only thermal cyclers with heated lids are suitable for this procedure.*

- Desktop centrifuge (up to 14,000 rpm)
- Pipette and pipette tips
- 0.2 ml PCR tubes
- 1.5 ml microcentrifuge tubes
- 100% ethanol
- RNA sample

GENERAL PRODUCT INFORMATION

Quality Control: Each lot of Methylamp™ RNA Bisulfite Conversion Kit is tested against predetermined specifications to ensure consistent product quality. EpigenTek guarantees the performance of all products in the manner described in our product instructions.

Product Warranty: If this product does not meet your expectations, simply call our technical support unit or your regional distributor. We also encourage you to contact us if you have any suggestions about product performance or new applications and techniques.

Safety: Suitable lab coat, disposable gloves, and proper eye protection are required when working with this product.

Product Updates: EpigenTek reserves the right to change or modify any product to enhance its performance and design. The information in this User Guide is subject to change at any time without notice. Be sure to use the latest User Guide for this kit which can be accessed online at www.epigentek.com/datasheet.

Usage Limitation: The Methylamp™ RNA Bisulfite Conversion Kit is for research use only and is not intended for diagnostic or therapeutic application.

Intellectual Property: The Methylamp™ RNA Bisulfite Conversion Kit and methods of use contain proprietary technologies by EpigenTek.

A BRIEF OVERVIEW

DNA cytosine methylation occurs by the covalent addition of a methyl group at the 5-carbon of the cytosine ring by DNA methyltransferases, resulting in 5-methylcytosine (5-mC). This process has been well studied and is generally associated with repression of gene expression. It was also observed that in humans, 5-mC occurs in various RNA molecules including tRNAs, rRNAs, mRNAs and non-coding RNAs (ncRNAs). At least 10,275 5-mC candidate sites were discovered in mRNAs and ncRNAs, which covered 10.6% of the total cytosine residues in the transcriptome. 5-mC seems to be enriched in some classes of ncRNA, but relatively depleted in mRNAs. However, the majority (83%) of their candidate sites were found in mRNAs. Within these transcripts 5-mC appears to be depleted within protein coding sequences but enriched in 5' and 3' UTRs. Two different methyltransferases, NSUN2 and Dnmt2 are known to catalyze the 5-mC modification in eukaryotic RNAs. Recent data strongly suggest that RNA cytosine methylation affects the regulation of various biological processes such as RNA stability and mRNA translation. Furthermore, loss of 5-mC in vault RNAs causes aberrant processing into Argonaute-associated small RNA fragments that can function as microRNAs. Thus, impaired processing of vault ncRNA may contribute to the etiology of human disorders related to NSUN2-deficiency.

Bisulfite conversion of RNA followed by RT-PCR amplification, cloning, and sequencing yields reliable information about RNA cytosine methylation states. By treating RNA with bisulfite, cytosine residues are deaminated to uracil while leaving 5-methylcytosine intact:

	<u>Unmethylated RNA</u>	<u>Methylated RNA</u>
<i>Original Sequence</i>	C-C-U-C-G-A-C-U	C-C-U- ^M C-G-A- ^M C-U
<i>After Bisulfite Conversion</i>	U-U-U-U-G-A-U-U	U-U-U- ^M C-G-A- ^M C-U

To effectively and efficiently prepare converted RNA for use in various downstream analyses, EpigenTek developed the Methylamp™ RNA Bisulfite Conversion Kit. The kit is specifically optimized and validated for bisulfite conversion of RNA and has the following advantages and features:

- Fast and convenient protocol that can be finished in 3 hours.
- Completely converts unmethylated cytosine into uracil (>99.9%) with no or negligible inappropriate/error conversion of methylcytosine to thymine (<0.1%) when the indicated range of sample RNA is used.
- Powerful protection against RNA degradation, with over 90% of RNA loss prevented.
- Included control primers are specific against bisulfite-converted RNA and can be used to test if the bisulfite conversion has been properly achieved.
- Low amount of input RNA can be used for bisulfite conversion - as low as 5 ng per reaction.
- Simple, reliable, and consistent reaction conditions.

PRINCIPLE & PROCEDURE

The Methylamp™ RNA Bisulfite Conversion Kit contains all reagents required for fast bisulfite conversion on a RNA sample. The unique conversion mix solution contains powerful RNA protection reagents to prevent chemical and thermophilic degradation, thus leading to an accelerated conversion of all cytosines to uracil with negligible methylcytosine deamination. The non-toxic RNA capture solution enables RNA to tightly bind to the column filter, so that converted RNA cleaning can be carried out on the column to effectively remove residual bisulfite and salts.



Fig 1. Schematic procedure of the Methylamp™ RNA Bisulfite Conversion Kit to obtain converted RNA.

Unconverted	G A U C C U U ^m C G A U G U C G G
Bisulfite Converted	G A U U U U U ^m C G A U G U U G G
cDNA Sequencing	G A T T T T T ^m C G A T G T T G G

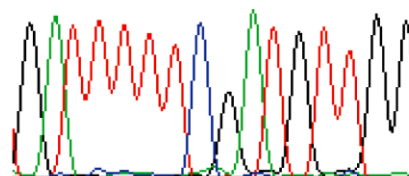


Fig 2. RNA bisulfite sequencing analysis: RNA isolated from MCF-7 cells is subject to bisulfite treatment, cDNA synthesis and sequencing after PCR amplification using 28S rRNA primers specific for bisulfite converted RNA. Unmethylated cytosines (#4, 5, 14) are converted to uracil and detected as thymine while methylated cytosine (#8) remains the same.

ASSAY PROTOCOL

For the best results, please read the protocol in its entirety prior to starting your experiment.

Starting Materials

Input RNA Amount: RNA amount can range from 5 ng to 1 µg per reaction. An optimal amount is 200–500 ng per reaction. Starting RNA may be in water or in a buffer such as TE. RNA should be high quality and relatively free of DNA. DNase I can be used to remove DNA and RNA should be eluted in RNase-free water.

RNA Storage: RNA should be stored at -20°C or -80°C until use.

1. Working Buffer and Solution Preparation

a. *Prepare Conversion Solution:*

Add 1.4 ml of **Conversion Buffer** and 40 µl of **Desulphonation Solution** to 1 vial of **Conversion Powder** to generate conversion solution. Mix by inverting and shaking the vial repeatedly for 3-4 min (trace amount of undissolved **Conversion Powder** may remain, which is normal as **Conversion Powder** is saturated in solution).

b. *Prepare 70% ethanol by adding 3 ml of distilled water to 7 ml of 100% ethanol.*

c. *Prepare 90% ethanol by adding 1ml of distilled water to 9 ml of 100% ethanol.*

d. *Prepare working desulphonation buffer:*

First dilute **Desulphonation Solution** at a 1:12 ratio by adding 5 µl of **Desulphonation Solution** to 55 µl of distilled water. Next, add 2 µl of **Diluted Desulphonation Solution** to every 1 ml of 90% ethanol, and mix.

2. RNA Bisulfite Conversion

a. Add 100 µl of the conversion solution to a PCR tube followed by adding 2-10 µl of RNA sample.

Prepared conversion solution can be stored at -20°C for up to 2 weeks without significant loss of efficiency. For the best results, the mixed solution should be used immediately.

b. Tightly close the PCR tubes and place them in a thermal cycler with heated lid. Program and run the thermal cycler :

65°C 5 min

60°C 90 min

Hold 4°C up to 16 h

Meanwhile, insert the number of **F-Spin Columns** into **F-Collection Tubes** as needed by your experiment.

3. Converted RNA Clean-Up

- a. Add 250 μ l of **NA Binding Solution** to each column. Then transfer the samples from each PCR tube (from Step 2b) to each column containing the **NA Binding Solution**. Centrifuge at 12,000 rpm for 1 min. Remove columns from collection tubes and discard the flowthrough. Place columns back into collection tubes.
- b. Add 200 μ l of 70% ethanol solution to each column. Centrifuge at 12,000 rpm for 1 min.
- c. Add 200 μ l of the working desulphonation buffer (**Desulphonation Solution** and 90% ethanol mixture according to Step 1d) to each column. Allow columns to sit for 30 min at room temperature, then centrifuge at 12,000 rpm for 1 min. Remove columns from collection tubes and discard the flowthrough. Place columns back into collection tubes.
- d. Add 200 μ l of 90% ethanol to each column. Centrifuge at 12,000 rpm for 1 min. Remove columns from collection tubes and discard the flowthrough. Place columns back into collection tubes. Add 200 μ l of 90% ethanol to each column again and centrifuge at 12,000 rpm for 1 min.
- e. Insert each column into a new 1.5 ml tube. Add 10-20 μ l of **Elution Buffer** directly to each column's filter membrane. Centrifuge at 12,000 rpm for 1 min to elute converted RNA.

Converted RNA is now ready for use, or storage at or below -20°C for up to 2 months. As the bisulfite-treated RNA is not stable, we recommend performing cDNA synthesis from bisulfite-treated RNA before next application or storage. See appendix I for cDNA synthesis.

Appendix

1. cDNA Synthesis

You can use your method of choice for cDNA synthesis. For your convenience, EpigenTek offers EpiNext™ Hi-Fi cDNA Synthesis Kit (Cat # 9004) which is optimized and validated for synthesis of cDNA from bisulfite RNA.

- a. Add the following in a 0.2-ml PCR tube on ice:

Component	Amount
Bisulfite-converted RNA (200-500 ng)	10 μ l
Random primer (50 μ M)	1 μ l
10 mM dNTP mix	1 μ l

- b. Heat in a thermocycler (no heated lid) at 65°C for 3 minutes. Place on ice immediately for at least 1 minute.

- c. Add the following to the tube on ice:

Component	Amount
5X RT Reaction Buffer	4 μ l
0.1M DTT	2 μ l
RNase Inhibitor	1 μ l
RT Enzyme Mix	1 μ l
Total Volume	20 μl

Vortex the sample briefly to mix and collect by centrifugation. Incubate as follows: 42°C for 45 min followed by 80°C for 5 min (no heated lid).

Store the cDNA synthesis reaction at -20°C, or proceed directly to next application such as methylation specific PCR (see Appendix 2 “Working with Methylation Specific qPCR”) or bisulfite-sequencing.

2. Working with Methylation Specific qPCR

When working with methylation specific-PCR, we recommend using the Methylamp™ MS-qPCR Fast Kit (Cat # P-1028) which contains a hot start polymerase system and has been optimized to decrease the overall methylation specific-qPCR amplification time. The master mix is provided at a 2X concentration for easier preparation of PCR reactions requiring only the addition of primers and templates. With this kit, the MS-qPCR can be finished in as short as 70 min.

Prepare the PCR Reactions

Component	Size (μ l)	Final Concentration
Methylamp Master Mix (2X)	10 μ l	1X
Forward Primer	1 μ l	0.4-0.5 μ M
Reverse Primer	1 μ l	0.4-0.5 μ M
cDNA Template	1-2 μ l	50 pg-0.1 μ g
RNase-free H ₂ O	6-7 μ l	
Total Volume	20 μl	

For the negative control, use RNase-free water instead of cDNA template.

Program the PCR Reactions

Cycle Step	Temp	Time	Cycle
<i>Activation</i>	95°C	7 min	1
<i>Cycling</i>	95°C	10 sec	40-45
	55°C	10 sec	
	72°C	8 sec	
<i>Final Extension</i>	72°C	1 min	1

TROUBLESHOOTING

Problem	Possible Causes	Suggestions
RNA is poorly modified	Poor RNA quality (RNA is severely degraded).	Check if the sample RNA 260/280 ratio is between 1.9 - 2.0.
	Too little RNA or too much RNA (i.e., < 1ng or >1 µg).	Increase or decrease input RNA to within the correct range, or to the optimal amount of 200-500 ng.
	Temperature or thermal cycling condition is incorrect.	Check for appropriate temperature or thermal cycling conditions.
	Insufficient RNA clean-up.	Ensure that 2 µl of diluted Desulphonation Solution is added into every 1 ml of 90% ethanol in Step 1d.
	Kit is not stored or handled properly.	Store all components of the kit at room temperature.
Eluate contains little or no RNA	Poor input RNA quality (degraded).	Check if RNA is degraded.
	NA Binding Solution is not added into the sample.	Ensure that NA Binding Solution is added in Step 3a.
	Concentration of ethanol solution used for RNA clean-up is not correct.	Use 90% ethanol for RNA clean-up.
Poor results in downstream methylation-specific qRT-PCR	Little or no PCR product even in positive control.	Ensure that cDNA synthesis is properly performed.
		Ensure that all PCR components were added and that suitable PCR program is used (PCR cycle should be >40).
		PCR primers and probes were not appropriate or were incorrectly designed. Ensure the primer and probes are suitable for MS-PCR.
	Significant non-specific PCR products.	Failed bisulfite conversion. Ensure that all steps of the bisulfite treatment and cleanup protocol were followed and that input RNA amount is within the recommended range.
		Primers and probes are not specific for converted RNA and target genes. Check the primer and probe design.

RELATED PRODUCTS

DNA Bisulfite Modification

P-1001	Methylamp™ DNA Modification Kit
P-1010	Methylamp™ One-Step DNA Modification Kit
P-1016	Methylamp™ Whole Cell Bisulfite Modification Kit
P-1026	BisulFlash™ DNA Modification Kit

cDNA Synthesis

P-9004	EpiNext™ Hi-Fi cDNA Synthesis Kit
--------	-----------------------------------

Methylation Analysis

P-1028	Methylamp™ MS-qPCR Fast Kit
--------	-----------------------------