

# EpiQuik™ Tissue Methylated DNA Immunoprecipitation Kit

Base Catalog # P-2020

## PLEASE READ THIS ENTIRE USER GUIDE BEFORE USE

The EpiQuik™ Tissue MeDIP Kit can be used to immunoprecipitate methylated DNA from a broad range of species including human, rat, and mouse.

The EpiQuik™ Tissue MeDIP Kit is suitable for combining the specificity of methylated DNA immunoprecipitation with qualitative and quantitative PCR, and southern blot as well as DNA microarray.

## KIT CONTENTS



**Important Information:** The amount of components supplied in this kit is designed for reaction count, not sample count, such as negative IgG controls and input DNA. Thus, experiments with samples to be paired with both IgG and input may require additional columns or components to be purchased separately. Please calculate the necessary volumes based on the below kit contents and protocol prior to starting the experiment.

Components	24 reactions	48 reactions
	P-2020-24	P-2020-48
CP1 (Wash Buffer)	28 ml	2 x 28 ml
CP2 (Antibody Buffer)	15 ml	30 ml
CP3 (Lysis Buffer)	4 ml	6 ml
CP4 (ChIP Dilution Buffer)	4 ml	6 ml
CP5 (DNA Release Buffer)	2 ml	2 x 2 ml
CP6 (Reverse Buffer)	2 ml	2 x 2 ml
CP7 (Binding Buffer)	5 ml	8 ml
CP8 (Elution Buffer)	0.6 ml	1.2 ml
Homogenizing Buffer	5 ml	10 ml
Non-immune IgG (1 mg/ml)*	10 $\mu$ l	15 $\mu$ l
Anti-5-Methylcytosine (1 mg/ml)*	25 $\mu$ l	50 $\mu$ l
Proteinase K (10 mg/ml)*	25 $\mu$ l	50 $\mu$ l
8-Well Assay Strips (with Frame)	3	6
8-Well Strip Caps	3	6
F-Spin Column	30	50
F-Collection Tube	30	50

\* For maximum recovery of the products, centrifuge the original vial after thawing prior to opening the cap.

## SHIPPING & STORAGE

The kit is shipped in two parts, one part at ambient room temperature, and the second part on frozen ice packs at 4°C.

Upon receipt: (1) Store the following components at 4°C: **Non-immune IgG, Anti-5-Methylcytosine, Proteinase K** and **8-Well Assay Strips**. (2) Store **all other components** at room temperature.

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

**Note:** Avoid repeated thawing and re-freezing of temperature sensitive components. It is recommended that you aliquot accordingly ahead of time.

## MATERIALS REQUIRED BUT NOT SUPPLIED

- Variable temperature waterbath
- Vortex mixer
- Desktop centrifuge (up to 14,000 rpm)
- Dounce homogenizer
- Sonicator
- Orbital shaker
- Pipettes and pipette tips
- 1.5 ml microcentrifuge tubes
- 15 ml conical tube
- TE buffer (pH 8.0)
- Ethanol (96-100%)

## GENERAL PRODUCT INFORMATION

**Quality Control:** EpigenTek guarantees the performance of all products in the manner described in our product instructions.

**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](http://www.epigentek.com/datasheet).

**Usage Limitation:** The *EpiQuik*<sup>™</sup> Tissue MeDIP Kit is for research use only and is not intended for diagnostic or therapeutic application.

**Intellectual Property:** The *EpiQuik*<sup>™</sup> MeDIP kits and method of use contain proprietary technologies by EpigenTek. *EpiQuik*<sup>™</sup> is a trademark of EpigenTek Group Inc.

## A BRIEF OVERVIEW

The core mechanism for epigenetic alterations of genomic DNA is hypermethylation of CpG islands in specific genes and global DNA hypomethylation. Region-specific DNA methylation is mainly found in 5'-CpG-3'dinucleotides within the promoters or in the first exon of genes, which is an important pathway for the repression of gene transcription in diseased cells. Global DNA

hypomethylation is likely caused by methyl-deficiency due to a variety of environmental influences. It has been demonstrated that alterations in DNA methylation are associated with many diseases, and especially with cancer.

Highly specific isolation of methylated DNA should provide an advantage for convenient and comprehensive identification of methylation status of normal and diseased cells, such as cancer cells, which may lead to the development of new diagnostic and therapeutic methods of cancer. Several methods have been used for enriching methylated DNA, including methyl-CpG binding domain (MBD) based methylated DNA affinity column and methylated DNA immunoprecipitation. However, these methods so far are considerably time consuming, labor intensive, have low throughput, and particularly, need purified DNA as starting material.

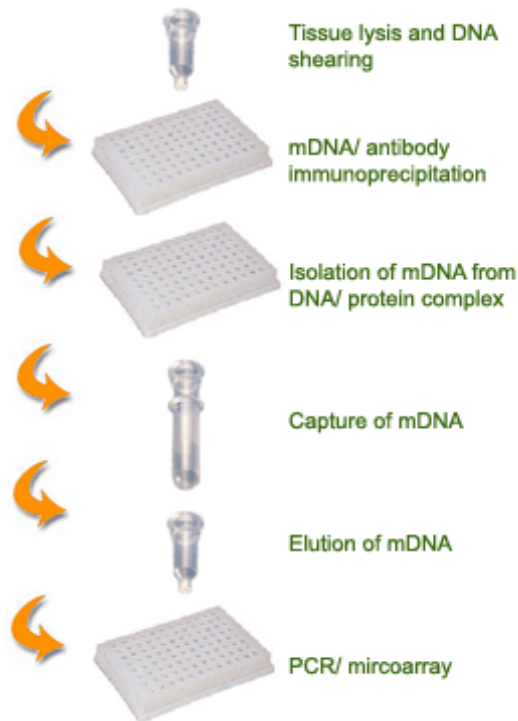
The *EpiQuik*<sup>™</sup> Tissue Methylated DNA Immunoprecipitation (MeDIP) Kit uses a proprietary and unique procedure/composition to enrich methylated DNA from various mammalian tissues. In the assay, an antibody specific to methyl cytosine is used to immunoprecipitate methylated genomic DNA. The immunoprecipitated methylated fractions can then be used for a standard DNA detection. The *EpiQuik*<sup>™</sup> Tissue Methylated DNA Immunoprecipitation Kit has the following features:

- Directly immunoprecipitate methylated fractions of DNA from tissue lysates.
- Highly efficient enrichment of methylated DNA: > 95%.
- The fastest procedure available, which can be finished within 3.5 hours.
- Strip microplate format makes the assay flexible: manual or high throughput.
- Columns for DNA purification are included: save time and reduce labor.
- Compatible with all DNA amplification-based approaches.
- Simple, reliable, and consistent assay conditions.

## PRINCIPLE & PROCEDURE

The *EpiQuik*<sup>™</sup> Tissue Methylated DNA Immunoprecipitation (MeDIP) Kit contains all reagents required for carrying out a successful methylated DNA immunoprecipitation directly from mammalian tissues. Particularly, this kit includes a ChIP-grade 5-methylcytosine antibody and a negative control Non-immune IgG. DNA in the cells is extracted, sheared, and added into the microwell immobilized with the antibody. DNA is released from the antibody-DNA complex, and purified through the specifically designed Fast-Spin Column. Eluted DNA can be used for various down-stream applications.

Schematic Procedure for Using the EpiQuik™ Tissue Methylated DNA Immunoprecipitation Kit



## PROTOCOL

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

Before starting, perform the following:

1. Prepare the following required solutions (not included): 90% Ethanol; 70% Ethanol; 1X TE Buffer (pH 8.0).
2. Ensure that all buffers are in clear solution. Shake or vortex if these buffers precipitate.

### Antibody Binding to the Assay Plate

1. Determine the number of strip wells required. Leave these strips in the plate frame (remaining unused strips can be placed back in the bag. Seal the bag tightly and store at 4°C). Wash the strip wells once with 150  $\mu$ l of **CP1**.
2. Add 100  $\mu$ l of **CP2** to each well and then add the antibodies: 1  $\mu$ l of **Non-immune IgG** as the negative control, and 0.5-1  $\mu$ l of **Anti-5-Methylcytosine** for the sample.
3. Cover the strip wells with Parafilm M and incubate at room temperature for 60 minutes. Meanwhile, prepare the cell extracts as described in the next steps.

### Cell Collection and Lysis

*Tissue Disaggregation:*

1. Place the tissue sample into a 60 or 100 mm plate. Remove unwanted tissue such as fat and necrotic material from the sample. Weigh the sample and cut it into small pieces (1-2 mm<sup>3</sup>) with a scalpel or scissors.
2. Transfer tissue pieces to a Dounce homogenizer. Add 1 ml of the **Homogenizing Buffer** per every 200 mg of tissue and disaggregate tissue pieces by 10-30 strokes.
3. Transfer homogenized mixture to a 15 ml conical tube and centrifuge at 3000 rpm for 5 minutes at 4°C. If total mixture volume is less than 2 ml, transfer mixture to a 2 ml vial and centrifuge at 5000 rpm for 5 minutes at 4°C. Remove supernatant.

### Tissue Lysis and DNA Shearing

1. Add **CP3** to re-suspend the disaggregated tissue pellet (100 µl/20 mg of tissue). Transfer the solution to a 1.5 ml vial (500 µl maximum for each vial) and incubate at room temperature for 10 minutes and vortex occasionally.
2. Shear DNA by sonication. Usually, sonicate 4-5 pulses of 15-20 seconds each at level 2 using a Branson Microtip probe, followed by 30-40 seconds rest on ice between each pulse. (The conditions of DNA shearing can be optimized based on cells and sonicator equipment. If desired, remove 5 µl of sonicated cell lysate for agarose gel analysis. The length of sheared DNA should be between 200-1000 bp.)
3. Pellet cell debris by centrifuging at 14,000 rpm for 10 minutes.

### Methylated DNA Immunoprecipitation

1. Transfer clear supernatant to a new 1.5 ml vial (supernatant can be stored at -80°C at this step). Dilute required volume of supernatant with **CP4** at a 1:1 ratio (ex: add 100 µl of **CP4** to 100 µl of supernatant).
2. Remove 5 µl of the *diluted supernatant* to a 0.5 ml vial. Label the vial as "input DNA" and place on ice.
3. Remove the incubated antibody solution and wash the strip wells three times with 150 µl of **CP2** by pipetting in and out.
4. Transfer 100 µl of the *diluted supernatant* to each strip well. Cover the strip wells with Parafilm M and incubate at room temperature (22-25°C) for 90-120 minutes on an orbital shaker (50-100 rpm).
5. Remove supernatant. Wash the wells six times with 150 µl of **CP1**. Allow 2 minutes on an orbital shaker (50-100 rpm) for each wash. Wash the wells once (for 2 minutes) with 150 µl of *1X TE Buffer*.

### Methylated DNA Isolation/Purification

1. Add 1 µl of **Proteinase K** to each 40 µl of **CP5** and mix. Add 40 µl of **CP5** containing **Proteinase K** to the samples (including the "input DNA" vial). Cover the sample wells with strip caps and incubate at 65°C in a waterbath for 15 minutes.
2. Add 40 µl of **CP6** to the samples; mix, and re-cover the wells with strip caps and incubate at 65°C in a waterbath for 30 minutes. Also add 40 µl of **CP6** to the vial containing supernatant as "input DNA." Mix and incubate at 65°C for 30 minutes.
3. Place a spin column into a 2 ml collection tube. Add 150 µl of **CP7** to the samples and transfer mixed solution to the column. Centrifuge at 12,000 rpm for 20 seconds.

4. Add 200  $\mu$ l of 70% ethanol to the column, centrifuge at 12,000 rpm for 20 seconds. Remove the column from the collection tube and discard the flowthrough.
5. Replace column to the collection tube. Add 200  $\mu$ l of 90% ethanol to the column and centrifuge at 12,000 rpm for 20 seconds.
6. Remove the column and discard the flowthrough. Replace column to the collection tube and wash the column again with 200  $\mu$ l of 90% ethanol at 12,000 rpm for 35 seconds.
7. Place the column in a new 1.5 ml vial. Add 10-20  $\mu$ l of **CP8** directly to the filter in the column and centrifuge at 12,000 rpm for 20 seconds to elute purified DNA.

Methylated DNA is now ready for use or storage at  $-20^{\circ}\text{C}$ .

**Note:** For PCR positive control (methylation) and negative control (unmethylation), the primers for highly methylated sequences of *H19ICR*, *LAP* or *XIST* and the primer for unmethylated  $\beta$ -actin or *GAPDH* sequence could be used, respectively. For conventional PCR, the number of PCR cycles may need to be optimized for the better PCR results.

**References:** Weber M et al: *Nature Genetics*, 37: 853-862, 2005.

## TROUBLESHOOTING

### Little or No PCR Product

- |  |  |
|--|--|
| 1. Insufficient tissues.                                   | Increase tissue amount (ex: > 20 mg of tissue/per reaction).   |
| 2. Insufficient tissue lysis.                              | Follow the guidelines in the protocol. Check the tissue lysis by observing a 5 $\mu$ l portion of the tissue lysate under the microscope.  |
| 3. Insufficient/too much sonication.                       | Follow the protocol instruction for obtaining the appropriate sized DNA. Keep the sample on ice during the sonication.                     |
| 4. Incorrect temperature/insufficient time for DNA release | Follow the guidelines in the protocol for appropriate temperature and time.  |
| 5. Incorrect PCR conditions.                               | Check if all PCR components are added.<br>Increase amount of DNA added to PCR reaction.<br>Increase the number of cycles for PCR reaction. |
| 6. Incorrect or bad primers                                | Ensure the designed primers are specific to the target sequence.   |
| 7. The column is not washed with 90% ethanol.              | Ensure that wash solution is 90% ethanol.  |

8. DNA is not completely passed through the filter.

Increase centrifuge time to 1 minute at steps 3 to 7 of "Methylated DNA Isolation/Purification."

### Little or No Amplification Difference Between the Sample and the Negative Control

1. Insufficient wash at each wash step.

Follow the protocol for appropriate wash.

2. Antibody is added into the well for the negative control by mistake.

Ensure antibody is added into the correct well.

3. Too many PCR cycles.

If using conventional PCR, decrease the cycles to appropriate cycle number. Differences between quantities of starting DNA can be measured generally within the linear PCR amplification phase.

## RELATED PRODUCTS

P-2002	EpiQuik™ Chromatin Immunoprecipitation (ChIP) Kit
P-2003	EpiQuik™ Tissue Chromatin Immunoprecipitation (ChIP) Kit
P-2006	EpiQuik™ Methyl-Histone H3-K9 ChIP Kit
P-2007	EpiQuik™ Methyl-Histone H3-K4 ChIP Kit
P-2008	EpiQuik™ Tissue Methyl-Histone H3-K9 ChIP Kit
P-2009	EpiQuik™ Tissue Methyl-Histone H3-K4 ChIP Kit
P-2010	EpiQuik™ Acetyl-Histone H3 ChIP Kit
P-2011	EpiQuik™ Acetyl-Histone H4 ChIP Kit
P-2012	EpiQuik™ Tissue Acetyl-Histone H3 ChIP Kit
P-2013	EpiQuik™ Tissue Acetyl-Histone H4 ChIP Kit
P-2015	EpiQuik™ Methyl-Histone H3-K27 ChIP Kit
P-2016	EpiQuik™ Tissue Methyl-Histone H3-K27 ChIP Kit
P-2019	EpiQuik™ Methylated DNA Immunoprecipitation Kit