

# EpiQuik™ Tissue Chromatin Immunoprecipitation Kit

Base Catalog # P-2003

## PLEASE READ THIS ENTIRE USER GUIDE BEFORE USE

The *EpiQuik*™ Tissue Chromatin Immunoprecipitation Kit is suitable for combining the specificity of immunoprecipitation with qualitative and quantitative PCR, MS-PCR, DNA sequencing, and southern blot as well as DNA microarray.

Like using other ChIP kits, if you use the *EpiQuik*™ ChIP kits, choice of good antibody (i.e. IP proven) is required for precipitating the fixed protein/DNA complexes.

## 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	96 reactions
	P-2003-1	P-2003-2	P-2003-3
CP1 (Wash Buffer)	28 ml	2 x 28 ml	4 x 28 ml
CP2 (Antibody Buffer)	15 ml	30 ml	2 x 30 ml
CP3 (Lysis Buffer)	4 ml	6 ml	10 ml
CP4 (ChIP Dilution Buffer)	4 ml	6 ml	10 ml
CP5 (DNA Release Buffer)	2 ml	2 x 2 ml	6 ml
CP6 (Reverse Buffer)	2 ml	2 x 2 ml	6 ml
CP7 (Binding Buffer)	5 ml	8 ml	15 ml
CP8 (Elution Buffer)	0.6 ml	1.2 ml	2 ml
Homogenizing Buffer	5 ml	10 ml	16 ml
Protease Inhibitor Cocktail (100X)*	25 $\mu$ l	50 $\mu$ l	100 $\mu$ l
Non-immune IgG (1 mg/ml)*	10 $\mu$ l	15 $\mu$ l	25 $\mu$ l
Anti-RNA Polymerase II (1 mg/ml)*	5 $\mu$ l	7 $\mu$ l	11 $\mu$ l
Proteinase K (10 mg/ml)*	25 $\mu$ l	50 $\mu$ l	100 $\mu$ l
Control Primers (GAPDH)			
Forward (20 $\mu$ M)*	10 $\mu$ l	15 $\mu$ l	20 $\mu$ l
Reverse (20 $\mu$ M)*	10 $\mu$ l	15 $\mu$ l	20 $\mu$ l
8-Well Assay Strips (with Frame)	3	6	12
8-Well Strip Caps	3	6	12
F-Spin Column	30	50	100
F-Collection Tube	30	50	100

\* Spin the solution down to the bottom prior to use.

## SHIPPING & STORAGE

The kit is shipped in three parts: the first and second parts at ambient room temperature and the third part on frozen ice packs at 4°C.

Upon receipt: (1) Store the following components at 4°C: **Protease Inhibitor Cocktail, Non-immune IgG, Anti-RNA Polymerase II, Proteinase K, Control Primers** 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.

## MATERIALS REQUIRED BUT NOT SUPPLIED

- Variable temperature waterbath
- Vortex mixer
- Desktop centrifuge (up to 14,000 rpm)
- Sonicator
- Dounce homogenizer
- Orbital shaker
- Pipettes and pipette tips
- 1.5 ml microcentrifuge tubes
- 15 ml conical tube
- Antibody of interest for chromatin immunoprecipitation
- 37% formaldehyde
- Glycine solution
- 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> ChIP kits are for research use only and are not intended for diagnostic or therapeutic application.

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

## A BRIEF OVERVIEW

Protein-DNA interaction play a critical role for cellular functions such as signal transduction, gene transcription, chromosome segregation, DNA replication and recombination, and epigenetic silencing. Identifying the genetic targets of DNA binding proteins and knowing the mechanisms of protein-DNA interaction is important for understanding cellular process.

Chromatin Immunoprecipitation (ChIP) offers an advantageous tool for studying protein-DNA interactions. Unlike other methods such as EMASA, DNA microarrays, and report gene assays, which analyze direct interactions between protein and DNA *in vitro*, ChIP can detect that a specific protein binds to the specific sequences of a gene in living cells. A typical ChIP includes the following steps: (1) formaldehyde cross-link of chromatin; (2) shearing of chromatin; (3) immunoprecipitation; (4) reversal of cross-link; and (5) DNA purification.

Requirement for using ChIP to analyze protein-DNA interaction in solid tissues and tumors is rapidly increasing. There are several methods used for chromatin immunoprecipitation, however, most of these methods available so far are very time consuming, labor intensive, low throughput, and furthermore, not specifically designed for solid tissues and tumors.

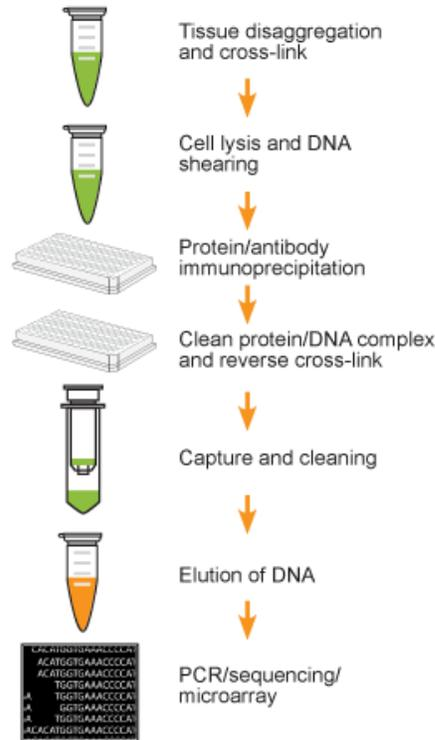
The *EpiQuik*<sup>™</sup> ChIP kits use a proprietary and unique procedure/composition to investigate protein-DNA interaction in solid tissues and tumors. The *EpiQuik*<sup>™</sup> ChIP kit series have the following features:

- The fastest procedure available, which can be finished within 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 Chromatin Immunoprecipitation Kit contains all reagents required for carrying out a successful chromatin immunoprecipitation from mammalian tissues. Particularly, this kit includes a positive control antibody (RNA polymerase II), a negative control Non-immune IgG, and GAPDH primers that can be used as a positive control to demonstrate the efficacy of the kit reagents and protocol. RNA polymerase II is considered to be enriched in the GAPDH gene promoter that is expected to be undergoing transcription in most growing mammalian cells and can be immunoprecipitated by RNA polymerase II, but not by Non-immune IgG. In this ChIP, cells are cross-linked with formaldehyde and chromatin is extracted. The chromatin is then sheared and added into the microwell immobilized with affinity antibodies. Cross-linked DNA is released from antibody-captured protein-DNA complex, reversed, 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 Chromatin 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, 1.25 M Glycine Solution.
2. Ensure that all buffers are in clear solution. Shake or vortex if these buffers precipitate.

### Antibody Binding to the Assay Strip Wells

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 strip wells once with 120  $\mu$ 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, 1  $\mu$ l of **Anti-RNA Polymerase II** as the positive control, and 2-3  $\mu$ g of your antibody of interest.
3. Cover the strip wells with Parafilm M and incubate at room temperature for 60-90 minutes. Meanwhile, prepare the cell extracts as described in the next steps.

### Tissue Disaggregation and *In Vivo* Cross-Link

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 the sample into small pieces (1-2 mm<sup>3</sup>) with a scalpel or scissors.

2. Transfer tissue pieces to a 15 ml conical tube. Prepare the **Cross-Link Solution** by adding *formaldehyde* to culture medium (final concentration is 1%. Ex: add 270  $\mu$ l of 37% *formaldehyde* to 10 ml of culture medium). Add 1 ml of **Cross-Link Solution** per every 40 mg of tissue and incubate at room temperature for 15-20 minutes on a rocking platform.
3. Add 1 ml of 1.25 M *Glycine solution* per every 9 ml of **Cross-Link Solution**, then mix and centrifuge at 800 rpm for 5 minutes. Discard the supernatant. Wash cells with 10 ml of ice-cold PBS once by centrifugation at 800 rpm for 5 minutes. Discard the supernatant.
4. 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 10-20 strokes.
5. 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.

### Cell Lysis and DNA Shearing

1. Add **CP3** containing **Protease Inhibitor Cocktail (PIC)** (Ex: 10  $\mu$ l of **PIC** to each 1 ml of **CP3**) to re-suspend the disaggregated tissue pellet (100  $\mu$ l/20 mg of tissue). Transfer cell suspension to a 1.5 ml vial (500  $\mu$ l maximum for each vial) and incubate 10 minutes on ice and vortex occasionally.
2. Shear DNA by sonication. Usually, sonicate 4 to 5 pulses of 15 to 20 seconds each at level 2 using a Branson Microtip probe, followed by 30 to 40 seconds rest on ice between each pulse. (The conditions of cross-linked DNA shearing can be optimized based on tissues and sonicator equipment. If desired, remove 5  $\mu$ l 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 at 4°C.

### Protein/DNA Immunoprecipitation

1. Transfer supernatant to a 1.5 ml vial (supernatant can be stored at -80°C at this step). Dilute the required volume of supernatant with **CP4** at a 1:1 ratio (ex: add 100  $\mu$ l of **CP4** to 100  $\mu$ l of cell supernatant).
2. Remove 5  $\mu$ l 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  $\mu$ l of **CP2** by pipetting in and out.
4. Transfer 100  $\mu$ l of the *diluted supernatant* to each strip well. Cover the strip wells with Parafilm M and incubate at room temperature (22-25°C) for 60-90 minutes on an orbital shaker (50-100 rpm).
5. Remove supernatant. Wash the wells with 150  $\mu$ l of **CP1** six times. Allow 2 minutes on an orbital shaker (100 rpm) for each wash. Wash the wells once (for 2 minutes) with 150  $\mu$ l of 1X *TE Buffer*.

### Cross-Linked DNA Reversal/DNA Purification

1. Add 1  $\mu$ l of **Proteinase K** to each 40  $\mu$ l of **CP5** and mix. Add 40  $\mu$ l 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  $\mu$ l of **CP6** to the samples; mix, re-cover the wells with strip caps and incubate at 65°C in a waterbath for 90 minutes. Also add 40  $\mu$ l of **CP6** to the vial containing *supernatant*, labeled as "input DNA". Mix and incubate at 65°C for 90 minutes.
3. Place a spin column into a 2 ml collection tube. Add 150  $\mu$ l 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 15 seconds. Remove the column from the collection tube and discard the flowthrough.
5. Replace the 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.

DNA is now ready for use or storage at -20°C.

**Note:** If a conventional PCR or a SYBR green real time PCR is performed, control primers (110 bp, for human tissues) included in the kit can be used as the positive control. For mouse or rat tissues, Control primers may be needed to be designed by the user. For conventional PCR, the number of PCR cycles may need to be optimized for better PCR results.

## TROUBLESHOOTING

### Little or No PCR Products

- |  |   |
|--|---|
| 1. Insufficient amount of tissue.  | Increase tissue amount (ex: >10 mg of tissue/per reaction).   |
| 2. Insufficient or too much cross-linking.                                   | Check if the appropriate cross-link step is carried out according to the protocol.  |
| 3. 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. |
| 4. Insufficient/too much sonication.   | Follow the protocol instructions for obtaining the appropriate sized DNA. Keep the sample on ice during the sonication.                   |
| 5. Antibody does not bind to protein.  | Check if the subclass or isotype of the antibody is correct. Choose a ChIP or IP grade antibody.  |
| 6. Incorrect temperature/time for DNA release and reversal of cross-linking. | Follow the guidelines in the protocol for appropriate temperature and time.   |

- |  |  |
|--|--|
| 7. 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. |
| 8. Wrong or bad primers.                             | Ensure the designed primers are specific to the target sequence.   |
| 9. The column is not washed with 90% ethanol.        | Ensure that wash solution is 90% ethanol.  |
| 10. DNA is not completely passed through the filter. | Increase centrifuge time to 1 minute at steps 3 to 7 of "Cross-Linked DNA Reversal/DNA Purification."                                      |

#### **Little or No Amplification Difference Between the Positive Control and the Negative Control**

- |  |   |
|--|---|
| 1. Insufficient wash at each wash step.  | Follow the protocol for appropriate wash.   |
| 2. Positive control 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 the appropriate cycle number. Differences in starting DNA quantities may be measured within the linear PCR amplification phase. |

## **RELATED PRODUCTS**

P-2002	<i>EpiQuik</i> <sup>™</sup> Chromatin Immunoprecipitation (ChIP) Kit
P-2006	<i>EpiQuik</i> <sup>™</sup> Methyl-Histone H3-K9 ChIP Kit
P-2007	<i>EpiQuik</i> <sup>™</sup> Methyl-Histone H3-K4 ChIP Kit
P-2008	<i>EpiQuik</i> <sup>™</sup> Tissue Methyl-Histone H3-K9 ChIP Kit
P-2009	<i>EpiQuik</i> <sup>™</sup> Tissue Methyl-Histone H3-K4 ChIP Kit
P-2010	<i>EpiQuik</i> <sup>™</sup> Acetyl-Histone H3 ChIP Kit
P-2011	<i>EpiQuik</i> <sup>™</sup> Acetyl-Histone H4 ChIP Kit
P-2015	<i>EpiQuik</i> <sup>™</sup> Methyl-Histone H3-K27 ChIP Kit
P-2016	<i>EpiQuik</i> <sup>™</sup> Tissue Methyl-Histone H3-K27 ChIP Kit
P-2017	<i>EpiQuik</i> <sup>™</sup> Methyl-CpG Binding Domain Protein 2 ChIP Kit
P-2018	<i>EpiQuik</i> <sup>™</sup> Tissue Methyl-CpG Binding Domain Protein 2 ChIP Kit