

# EpiQuik™ Global Mono-Methyl Histone H3-K9 Quantification Kit (Colorimetric)

Base Catalog # P-3030



**PLEASE READ THIS ENTIRE USER GUIDE BEFORE USE**

The EpiQuik™ Global Mono-Methyl Histone H3-K9 Quantification Kit (Colorimetric) is suitable for specifically measuring global histone H3-K9 mono-methylation using a variety of mammalian cells (human, mouse, etc.) including fresh and frozen tissues, cultured adherent and suspension cells.

110 Bi County Blvd. Ste. 122, Farmingdale, NY 11735  
Tel: 1-877-374-4368 ■ Fax: 1-718-484-3956 ■ E-mail: [info@epigentek.com](mailto:info@epigentek.com) ■ Web: [www.epigentek.com](http://www.epigentek.com)  
© Epigentek Group Inc. All rights reserved. Products are for research use only.

## KIT CONTENTS

Components	48 assays P-3030-48	96 assays P-3030-96
C1 (10X Wash Buffer)	10 ml	20 ml
C2 (Antibody Buffer)	6 ml	12 ml
C3 (Detection Antibody, 1 mg/ml)*	5 $\mu$ l	10 $\mu$ l
C4 (Color Developer)	5 ml	10 ml
C5 (Stop Solution)	3 ml	6 ml
Standard Control (100 $\mu$ g/ml)*	10 $\mu$ l	20 $\mu$ l
8-Well Sample Strips (with Frame)	4	9
8-Well Standard Control Strips	2	3
User Guide	1	1

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

## SHIPPING & STORAGE

Upon receipt, (1) Store **C3** and **standard control** at  $-20^{\circ}\text{C}$ ; (2) Store **C5** at room temperature away from light; (3) Store **all other components** at  $4^{\circ}\text{C}$  away from light. The components of the kit are stable for up to 6 months from date of shipment, when stored properly.

**Note:** Check if buffers, **C1** and **C2**, contain salt precipitates before using. If so, warm (at room temperature or  $37^{\circ}\text{C}$ ) and shake the buffers until the salts are re-dissolved.

## MATERIALS REQUIRED BUT NOT SUPPLIED

- Orbital shaker
- Pipettes and pipette tips
- Reagent reservoir
- Microplate reader

## GENERAL PRODUCT INFORMATION

**Usage Limitation:** The *EpiQuik*<sup>™</sup> Global Mono-Methyl Histone H3-K9 Quantification Kit (Colorimetric) is for research use only and is not intended for diagnostic or therapeutic application.

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

**Safety:** Suitable lab coat, disposable gloves, and eye protection are required when working with the kit.

**Product Updates:** Epigentek reserves the right to change or modify any product to enhance its performance and design.

**Intellectual Property:** The *EpiQuik*<sup>™</sup> Global Mono-Methyl Histone H3-K9 Quantification Kit (Colorimetric) and methods of use contain proprietary technologies by Epigenetek. *EpiQuik*<sup>™</sup> is a trademark of Epigenetek Group Inc.

## A BRIEF OVERVIEW

Epigenetic activation or inactivation of genes play a critical role in many important human diseases, especially in cancer. A major mechanism for epigenetic inactivation of the genes is methylation of CpG islands in genome DNA caused by DNA methyltransferases. Histone methyltransferases (HMTs) control or regulate DNA methylation through chromatin-dependent transcription repression or activation. HMTs transfer 1-3 methyl groups from S-adenosyl-L-methionine to the lysine and arginine residues of histone proteins. ESET, G9a, SUV39-h1, SUV39-h2, SETDB1, Dim-5, and Eu-HMTase, are histone methyltransferases that catalyze methylation of histone H3 at lysine 9 (H3-K9) in mammalian cells. Mono-methyl H3-K9 is enriched in certain euchromatic domains, which have been postulated to be transcriptionally silent. The global H3-K9 mono-methylation can also be changed by inhibition or activation of HMTs. Thus, quantitative detection of global mono-methyl histone H3-K9 would provide useful information for better understanding epigenetic regulation of gene activation/silencing and for developing HMT-targeted drugs. The *EpiQuik*<sup>™</sup> Global Mono-Methyl Histone H3-K9 Quantification Kit (Colorimetric) provides a tool for measuring global mono-methylation of histone H3-K9. The kit has the following features:

- Quick and efficient procedure, which can be finished within 2.5 hours.
- Innovative colorimetric assay without the need for radioactivity, electrophoresis, or chromatography.
- Specifically captures mono-methylated H3-K9 with the detection limit as low as 2 ng/well and detection range from 20 ng-5  $\mu$ g/well of histone extracts.
- The control is conveniently included for the quantification of the amount of mono-methylated H3-K9.
- Strip microplate format makes the assay flexible: manual or high throughput.
- Simple, reliable, and consistent assay conditions.

## PRINCIPLE & PROCEDURE

The *EpiQuik*<sup>™</sup> Global Mono-Methyl Histone H3-K9 Quantification Kit (Colorimetric) is designed for measuring global histone H3-K9 mono-methylation. In an assay with this kit, the mono-methylated histone H3 at lysine 9 is captured to the strip wells coated with an anti-monomethyl H3-K9 antibody. The captured mono-methylated histone H3-K9 can then be detected with a labeled detection antibody followed by a color development reagent. The ratio of mono-methylated H3-K9 is proportional to the intensity of absorbance. The absolute amount of mono-methylated H3-K9 can be quantitated by comparing to the standard control.



Schematic Procedure for Using the EpiQuik™ Global Mono-Methyl Histone H3-K9 Quantification Kit (Colorimetric)

## PROTOCOL

- Prepare histone extracts from cells/tissues treated or untreated by using your own successful method (acid extraction or high salt extraction).
  - For your convenience and best results, Epigentek offers the EpiQuik™ Total Histone Extraction Kit (Cat. # OP-0006) optimized for use in the EpiQuik™ modified histone quantification series.
  - Preparation of histone extracts can also be performed using the attached procedure (See Appendix). Histone extracts can be used immediately or stored at  $-80^{\circ}\text{C}$  for future use.
- 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  $4^{\circ}\text{C}$ ). Dilute **C1** with distilled water (pH 7.2-7.5) at 1:9 ratio (1 ml of **C1** + 9 ml of water).
- Add  $50\ \mu\text{l}$  of **C2** into each well. For the sample, add 50-200 ng of the histone extract into the sample wells. For standard curve, dilute **standard control** with **C2** to 1 – 100  $\text{ng}/\mu\text{l}$  at 5-7 points (e.g., 1.5, 3, 6, 12, 25, 50, and 100  $\text{ng}/\mu\text{l}$ ). Add  $1\ \mu\text{l}$  of **standard control** at the different concentrations into the standard well. For the blank, add no nuclear extracts or no standard control protein. Mix and cover the strip wells with Parafilm M and incubate at room temperature for 1-2 hours.
- Aspirate and wash the wells with  $150\ \mu\text{l}$  of **diluted C1** 3 times.
- Dilute **C3** (at 1:1000 ratio) to  $1\ \mu\text{g}/\text{ml}$  with **C2**. Add  $50\ \mu\text{l}$  of diluted **C3** to each well and incubate at room temperature for 60 min on an orbital shaker (100 rpm).
- Aspirate and wash the wells with  $150\ \mu\text{l}$  of **diluted C1** 6 times.

7. Add 100  $\mu$ l of **C4** into the wells and incubate at room temperature for 2-10 min away from light. Monitor color development in the sample and standard well (blue).
8. Add 50  $\mu$ l of **C5** to each well to stop enzyme reaction when the color in the standard wells containing the higher concentrations of standard control turns medium blue. The color should change to yellow and absorbance can be read on a microplate reader at 450 nm within 2-15 min.
9. Calculate % histone H3-K9 mono-methylation:

$$\text{Mono-methylation \%} = \frac{\text{OD (treated (tested) sample - blank)}}{\text{OD (untreated (control) sample - blank)}} \times 100\%$$

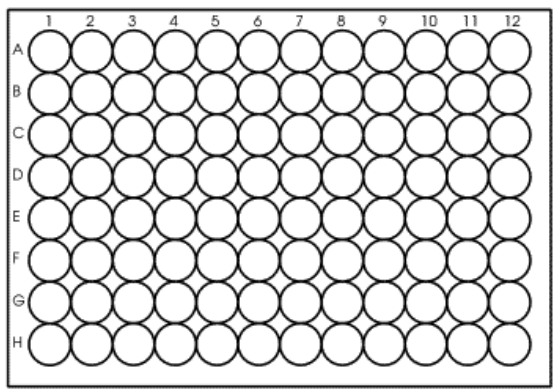
For the amount quantification, plot OD versus amount of **standard control** and determine the slope as delta OD/ng.

Calculate the amount of mono-methylated H3-K9 using the following formula:

$$\text{Amount (ng/mg protein)} = \frac{\text{OD (sample - blank)}}{\text{Protein } (\mu\text{g})^* \times \text{slope}} \times 1000$$

\* Histone extract amount added into the sample well at step 3.

## PLATE CONFIGURATION



- **Strip 1-3 (for 96 assays) or strip 1-2 (for 48 assays):** standard wells (**green trimmed**); the standard curve can be generated with 5-8 concentration points (includes blank).
- Example amount of standard control/well: **A1:** 100 ng; **B1:** 50 ng; **C1:** 25 ng; **D1:** 12 ng; **E1:** 6 ng; **F1:** 3 ng; **G1:** 1.5 ng; **H1:** 0 ng.
- **Strip 4-12 (for 96 assays) or strip 3-6 (for 48 assays):** sample wells (**No label**).
- Each sample or standard point can be assayed in duplicates or triplicates



## Appendix

### Histone Extraction Protocol

1. For tissues (treated and untreated), weigh the sample and cut the sample into small pieces (1-2 mm<sup>3</sup>) with a scalpel or scissors. Transfer tissue pieces to a Dounce homogenizer, add TEB buffer (PBS containing 0.5% Triton X 100, 2 mM PMSF and 0.02% NaN<sub>3</sub>) at 200 mg/ml, and disaggregate tissue pieces by 50-60 strokes. Transfer homogenized mixture to a 15 ml conical tube and centrifuge at 3000 rpm for 5 min at 4°C. If total mixture volume is less than 2 ml, transfer mixture to a 2 ml vial and centrifuge at 10,000 rpm for 1 min at 4°C. Remove supernatant.

For cells (treated and untreated), harvest cells and pellet the cells by centrifugation at 1000 rpm for 5 min at 4°C. Resuspend cells in TEB buffer at 10<sup>7</sup> cells/ml and lyse cells on ice for 10 min with gentle stirring. Centrifuge at 3000 rpm for 5 min at 4°C. If total volume is less than 2 ml, transfer cell lysates to a 2 ml vial and centrifuge at 10,000 rpm for 1 min at 4°C. Remove supernatant.

2. Resuspend cell/tissue pellet in 3 volumes (approx. 200 µl/10<sup>7</sup> cells or 200 mg tissues) of extraction buffer (0.5N HCl + 10% glycerol) and incubate on ice for 30 min.
3. Centrifuge at 12,000 rpm for 5 min at 4°C and remove the supernatant fraction to new vial.
4. Add 8 volumes (approx. 0.6 ml/10<sup>7</sup> cells or 200 mg tissues) of acetone and leave at -20°C overnight.
5. Centrifuge at 12,000 rpm for 5 min and air-dry the pellet. Dissolve the pellet in distilled water (30-50 µl/10<sup>7</sup> cells or 200 mg tissues).
6. Quantify the protein concentration. Aliquot the extract and store the extract at -20°C or -80°C.

### TROUBLESHOOTING

#### No Signal for Both the Standard Control and the Samples

Reagents are added incorrectly.

Check if reagents are added in order and if some steps of the procedure were omitted by mistake.

Incubation time and temperature are incorrect.

Ensure the incubation time and temperature described in the protocol are followed correctly.

#### No Signal or Very Weak Signal for Only the Standard Control

The amount of standard control is not added into "standard control wells" or is added insufficiently.

Ensure sufficient amount of control is properly added to the standard control well.

### No Signal for Only the Sample

The protein sample is not properly extracted.

The protein amount is added into well insufficiently.

Protein extracts are incorrectly stored.

Ensure the procedure and reagents are correct for the nuclear protein extraction

Ensure extract contains sufficient amount of protein.

Ensure the nuclear extracts are stored at  $-20^{\circ}\text{C}$  or  $-80^{\circ}\text{C}$ .

### High Background Present for the Blank

The well is not washed sufficiently.

Contaminated by the standard control.

Overdevelopment.

Check if wash at each step is performed according to the protocol.

Ensure the well is not contaminated by adding the control protein or by using control protein contaminated tips.

Decrease development time in Step 7

## RELATED PRODUCTS

P-3003	<i>EpiQuik</i> <sup>™</sup> Histone Methyltransferase Activity/Inhibition Assay Kit (H3-K9)
P-3016	<i>EpiQuik</i> <sup>™</sup> <i>In Situ</i> Histone H3-K9 Methylation Assay Kit
P-3031	<i>EpiQuik</i> <sup>™</sup> Global Mono-Methyl Histone H3-K9 Quantification Kit (Fluorometric)
P-3032	<i>EpiQuik</i> <sup>™</sup> Global Di-Methyl Histone H3-K9 Quantification Kit (Colorimetric)
P-3033	<i>EpiQuik</i> <sup>™</sup> Global Di-Methyl Histone H3-K9 Quantification Kit (Fluorometric)
P-3034	<i>EpiQuik</i> <sup>™</sup> Global Tri-Methyl Histone H3-K9 Quantification Kit (Colorimetric)
P-3035	<i>EpiQuik</i> <sup>™</sup> Global Tri-Methyl Histone H3-K9 Quantification Kit (Fluorometric)
P-3036	<i>EpiQuik</i> <sup>™</sup> Global Pan-Methyl Histone H3-K9 Quantification Kit (Colorimetric)
P-3037	<i>EpiQuik</i> <sup>™</sup> Global Pan-Methyl Histone H3-K9 Quantification Kit (Fluorometric)

