

# EpiQuik™ Global Mono-Methyl Histone H3-K79 Quantification Kit (Colorimetric)

Base Catalog # P-3054



## PLEASE READ THIS ENTIRE USER GUIDE BEFORE USE

The EpiQuik™ Global Mono-Methyl Histone H3-K79 Quantification Kit (Colorimetric) is suitable for specifically measuring global histone H3-K79 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.

Page 1

Printed 2017-08-17  
P-3054

## KIT CONTENTS

Components	48 assays P-3054-48	96 assays P-3054-96
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 should be stable for 6 months 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™ Global Mono-Methyl Histone H3-K79 Quantification Kit (Colorimetric) is for research use only and is not intended for diagnostic or therapeutic application

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

**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.

**Intellectual Property:** The *EpiQuik*™ Global Mono-Methyl Histone H3-K79 Quantification Kit (Colorimetric) and methods of use contain proprietary technologies by Epigentek. *EpiQuik*™ is a trademark of Epigentek Group Inc.

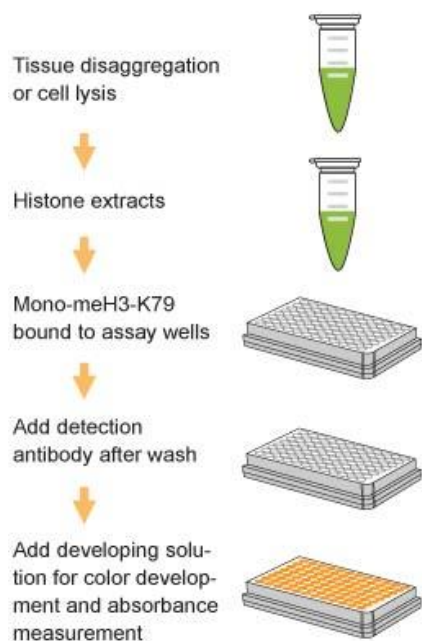
## A BRIEF OVERVIEW

Epigenetic activation or inactivation of genes plays 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. Dot1 is a histone methyltransferase that catalyzes methylation of histone H3 at lysine 79 (H3-K79) in mammalian cells. Intergenic H3-K79 mono-methylation is a mark of enhancer regions and may provide unique binding surfaces along the chromatin fiber to stabilize interactions between enhancers and their target promoters. The global H3-K79 mono-methylation can be changed by inhibition or activation of HMTs. Thus quantitative detection of global mono-methyl histone H3-K79 would provide useful information for better understanding epigenetic regulation of gene activation and for developing HMT-targeted drugs. The *EpiQuik*™ Global Mono-Methyl Histone H3-K79 Quantification Kit (Colorimetric) provides a tool for measuring mono-methylation of histone H3-K79. The kit has the following features:

- Quick and efficient procedure, which can be finished within 2.5 hours.
- Innovative colorimetric assay with no need for radioactivity, electrophoresis, and chromatography.
- Specifically capturing mono-methylated H3-K79 with the detection limit as low as 2 ng/well and detection range from 20 ng-5 µg/well of histone extracts.
- The control is conveniently included for quantification of the amount of mono-methylated H3-K79.
- Strip microplate format makes the assay flexible: manual or high throughput.
- Simple, reliable, and consistent assay conditions.

## PRINCIPLE & PROCEDURE

The *EpiQuik*™ Global Mono-Methyl Histone H3-K79 Quantification Kit (Colorimetric) is designed for measuring global histone H3-K79 mono-methylation. In an assay with this kit, the mono-methylated histone H3 at lysine 79 is captured to the strip wells coated with anti-monomethyl H3-K79 antibody. The captured mono-methylated histone H3-K79 can be then detected with a labeled detection antibody followed by color development reagent. The ratio of mono-methylated H3-K79 is proportional to the intensity of absorbance. The absolute amount of mono-methylated H3-K79 can be quantified by comparing to the standard control.



Schematic Procedure for Using the *EpiQuik™* Global Tri-Methyl Histone H3-K79 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 the 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 be also performed using the attached procedure (see the appendix). Histone extracts can be used immediately or stored at  $-80^{\circ}\text{C}$  for future use.
- Determine the number of the strip wells required. Leave these strips in the plate frame (remaining unused strips can be put 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 a 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 ng/ $\mu\text{l}$  for 5-7 points (ie: 1.5, 3, 6, 12, 25, 50, and 100 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 60-90 min.
- Aspirate and wash the wells with  $150\ \mu\text{l}$  of **diluted C1** 3 times.

5. 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).
6. Aspirate and wash the wells with 150  $\mu\text{l}$  of **diluted C1** 6 times.
7. Add 100  $\mu\text{l}$  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  $\mu\text{l}$  of **C5** to each well to stop enzyme reaction when 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-K79 mono-methylation:

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

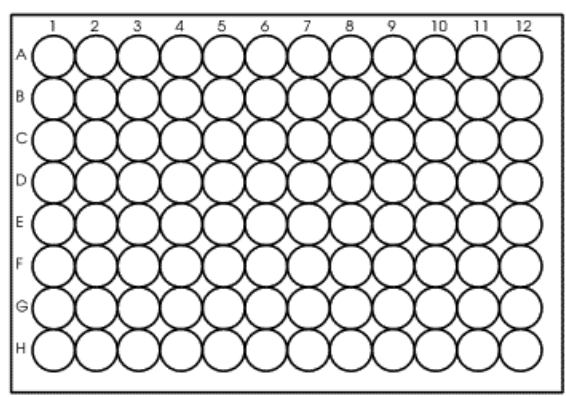
For the amount quantification, plot OD versus amount of **standard control** and determine the slope as delta OD/ng.

Calculate the amount of mono-methylated H3-K79 using the following formula:

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

\* 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 the duplicates or triplicates.

## Appendix

### Histone Extraction Protocol

1. For tissues (treated and untreated). Weigh the sample and cut the sample into small piece (1-2 mm<sup>3</sup>) with a scalpel or scissors. Transfer tissue piece 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 3,000 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 10000 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 the reagents are added in the proper order and if any steps of the procedure may have been 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 the "standard control wells," or is added insufficiently.

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

### No Signal for Only the Sample

The protein sample is not properly extracted.

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

The protein amount is added into well insufficiently.

Ensure extract contains a sufficient amount of proteins.

Protein extracts are stored incorrectly.

Ensure the protein 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.

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

Contaminated by the standard control.

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

Overdevelopment.

Decrease development time in Step 7.

## RELATED PRODUCTS

P-3055	<i>EpiQuik</i> <sup>™</sup> Global Mono-Methyl Histone H3-K79 Quantification Kit (Fluorometric)
P-3056	<i>EpiQuik</i> <sup>™</sup> Global Di-Methyl Histone H3-K79 Quantification Kit (Colorimetric)
P-3057	<i>EpiQuik</i> <sup>™</sup> Global Di-Methyl Histone H3-K79 Quantification Kit (Fluorometric)
P-3058	<i>EpiQuik</i> <sup>™</sup> Global Tri-Methyl Histone H3-K79 Quantification Kit (Colorimetric)
P-3059	<i>EpiQuik</i> <sup>™</sup> Global Tri-Methyl Histone H3-K79 Quantification Kit (Fluorometric)
P-3060	<i>EpiQuik</i> <sup>™</sup> Global Pan-Methyl Histone H3-K79 Quantification Kit (Colorimetric)
P-3061	<i>EpiQuik</i> <sup>™</sup> Global Pan-Methyl Histone H3-K79 Quantification Kit (Fluorometric)

