

# EpiQuik™ HDAC Activity/Inhibition Assay Kit (Colorimetric)

Base Catalog # P-4002

## PLEASE READ THIS ENTIRE USER GUIDE BEFORE USE

The EpiQuik™ HDAC Activity/Inhibition Assay Kit (Colorimetric) is very suitable for measuring HDAC activity/inhibition from a broad range of species including mammalian cells/tissues, plants, and bacteria.

## KIT CONTENTS

Components	48 assays P-4002-48	96 assays P-4002-96
H1 (10X Wash Buffer)	14 ml	28 ml
H2 (HDAC Assay Buffer)	1.5 ml	3 ml
H3 (Biotinylated HDAC Substrate)*	50 $\mu$ l	100 $\mu$ l
H4 (HDAC Inhibitor, 0.5 mM)*	50 $\mu$ l	100 $\mu$ l
H5 (HDAC Assay Standard, 20 $\mu$ g/ml)*	25 $\mu$ l	50 $\mu$ l
H6 (Capture Antibody, 100 $\mu$ g/ml)*	25 $\mu$ l	50 $\mu$ l
H7 (Detection Antibody, 200 $\mu$ g/ml)*	10 $\mu$ l	20 $\mu$ l
H8 (Developing Solution)	6 ml	12 ml
H9 (Stop Solution)	3 ml	6 ml
8-Well Assay Strip (with Frame)	6	12

\* 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: the first part at ambient room temperature and the second part on frozen ice packs at 4°C.

Upon receipt: (1) Store **H3, H4, H5, and H7** at –20°C away from light; (2) Store **H1, H6, H8, and 8-Well Assay Strips** at 4°C away from light; (3) Store **all other components** at room temperature. The kit is stable for up to 6 months from the date of shipment, when stored properly.

**Note:** Check if wash buffer, **H1**, contains salt precipitates before using. If so, warm (at room temperature or 37°C) and shake the buffer until the salts are re-dissolved.

## MATERIALS REQUIRED BUT NOT SUPPLIED

- ☐ Orbital shaker
- ☐ Pipettes and pipette tips
- ☐ Microplate reader
- ☐ 1.5 ml microcentrifuge tubes

## GENERAL PRODUCT INFORMATION

**Usage Limitation:** The EpiQuik™ HDAC Activity/Inhibition Assay 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.

**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).

**Intellectual Property:** *EpiQuik*<sup>™</sup> is a trademark of EpigenTek, Inc.

## A BRIEF OVERVIEW

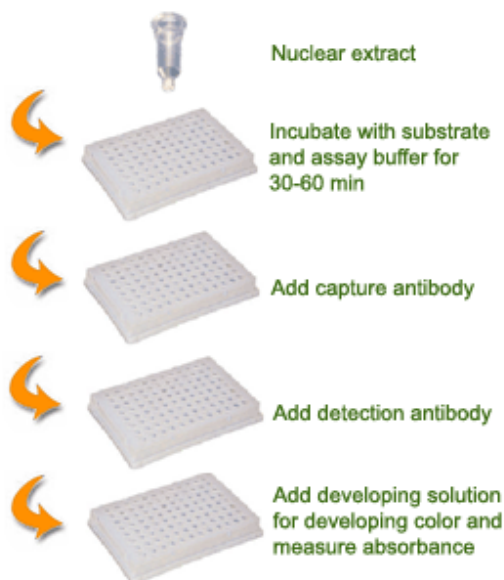
Histone deacetylases (HDACs) play a critical role in transcriptional repression of the gene expression in eukaryotic cells through catalyzing the hydrolytic removal of acetyl groups from histone lysine residues. HDACs are tightly involved in cell cycle regulation, cell proliferation, and in the development of human cancer. HDAC inhibition displays significant effects on apoptosis, cell cycle arrest, and differentiation in cancer cells. HDAC inhibitors are currently being developed as potential anticancer agents. There are several methods used for measuring HDAC activity/inhibition. However, most of these methods available so far are time consuming, laborious, produce radioactive waste, or cannot measure precise HDAC activity and inhibitory effects of inhibitors.

The *EpiQuik*<sup>™</sup> HDAC Activity/Inhibition Assay Kit (Colorimetric) uses a proprietary and unique procedure to measure HDAC activity/inhibition with the following features:

- Fast procedure, which can be finished within 3 hours.
- Innovative colorimetric assay without the use of radioactivity, extraction, or chromatography.
- Direct measurement of HDAC activity and inhibition without the use of lysyl endopeptidase, thereby avoiding the false inhibitory effect on HDACs.
- Strip microplate format makes the assay flexible: manual or high throughput analysis.
- Simple, reliable, and consistent assay conditions.

## PRINCIPLE & PROCEDURE

The *EpiQuik*<sup>™</sup> HDAC Activity/Inhibition Assay Kit (Colorimetric) is designed for measuring total HDAC activity/inhibition. In an assay with this kit, the unique acetylated histone substrate is stably captured on the strip wells. Active HDACs bind to and deacetylate histone substrate. The remaining un-deacetylated substrate can be recognized with a high affinity acetylated histone antibody. The ratio or amount of the un-deacetylated histone, which is inversely proportional to HDAC enzyme activity, can then be colorimetrically quantified through an ELISA-like reaction.



Schematic Procedure for Using the EpiQuik™ HDAC Activity/Inhibition Assay Kit (Colorimetric)

## PROTOCOL

1. Prepare nuclear extracts by using your own successful method. For your convenience and the best results, EpigenTek offers a nuclear extraction kit (Cat. No. OP-0002-1) optimized for use in the EpiQuik™ series. Nuclear extracts can be used immediately or stored at  $-80^{\circ}\text{C}$  for future use.
2. 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 **H1** (10X Wash Buffer) with distilled water (pH 7.2-7.5) at a 1:10 ratio (ex: 1 ml of **H1** + 9 ml of distilled water).
3. Dilute **H3** at a 1:50 ratio with **diluted H1**, and add  $50\ \mu\text{l}$  of the **diluted H3** into each sample well and each control well (control well is **H3** only). Do not add diluted **H3** to the wells for the blank and standard curve. For preparation of the standard curve, add  $50\ \mu\text{l}$  of **diluted H1** into the wells (without **H3** added), followed by adding  $1\ \mu\text{l}$  of **H5** at different amounts (0.1 – 10 ng). Cover the wells with Parafilm M and incubate at room temperature for 30-45 minutes.
4. Aspirate and wash each well with  $150\ \mu\text{l}$  of **diluted H1** two times.
5. Add  $28\ \mu\text{l}$  of **H2** and  $2\ \mu\text{l}$  of nuclear extracts (4-20  $\mu\text{g}$ ) or HDAC enzymes to each strip well, except the wells for the control, blank, and standard curve. Mix, cover the strip wells, and incubate at  $37^{\circ}\text{C}$  for 45-60 minutes. For the control and standard curve, add  $2\ \mu\text{l}$  of **H2** instead of nuclear extract. For HDAC inhibition, add  $2\ \mu\text{l}$  of different amounts of **H4** or tested inhibitors, and reduce **H2** volume to  $26\ \mu\text{l}$ . For the blank, add  $30\ \mu\text{l}$  of **H2** into the blank wells.
6. Aspirate and wash each well with  $150\ \mu\text{l}$  of **diluted H1** three times.

7. Dilute **H6** (at a 1:100 ratio) to 1 µg/ml with **diluted H1**. Add 50 µl of the **diluted H6** to each strip well and incubate at room temperature for 60 minutes on an orbital shaker (50-100 rpm).
8. Aspirate and wash each well with 150 µl of **diluted H1** four times.
9. Dilute **H7** (at a 1:1000 ratio) to 0.2 µg/ml with **diluted H1**. Add 50 µl of the **diluted H7** to each strip well and incubate at room temperature for 25-30 minutes.
10. Aspirate and wash each well with 150 µl of **diluted H1** four to five times.
11. Add 100 µl of **H8** to each well and incubate at room temperature for 2-10 minutes away from light. Monitor the color development in the sample and standard wells (blue).
12. Add 50 µl of **H9** 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 minutes.
13. Calculate HDAC activity or inhibition. For simple calculation:

$$\text{HDAC activity (OD/h/ml)} = \frac{[\text{OD (control - blank)} - \text{OD (sample - blank)}]}{\text{reaction time (0.5-1 hour)}} \times \text{sample dilution}^*$$

$$\text{Inhibition \%} = \left(1 - \frac{[\text{OD (control - blank)} - \text{OD (inhibitor sample - blank)}]}{[\text{OD (control - blank)} - \text{OD (no inhibitor sample - blank)}]}\right) \times 100\%$$

For an accurate calculation, plot OD value versus amount of **H5** and determine the slope as delta OD/ng.

Calculate HDAC activity using the following formula:

$$\text{Activity (ng/h/ml)} = \frac{[\text{OD (control-blank)} - \text{OD (sample - blank)}]}{\text{slope} \times \text{reaction time (0.5-1 hour)}} \times \text{sample dilution}^*$$

*\* If there is no dilution before adding protein extracts (2 µl) into the well, the sample dilution factor should be 500 (1000:2).*

## TROUBLESHOOTING

### No Signal for the Sample

The protein sample is not properly extracted.

Ensure the protein extraction protocol is suitable for nuclear protein extraction.

The protein amount is added into well insufficiently.

Ensure extract contains a sufficient amount of protein.

The sample is not prepared from fresh cells or tissues.

The nuclear extracts from frozen cells or tissues significantly lose enzyme activity. A fresh sample should be used.

Nuclear extracts are stored incorrectly.

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

Reagents are added incorrectly.

Check if 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.

Absence of HDAC activity in the sample due to treatment.

N/A.

#### High Background Present for the Blank

The well is not washed sufficiently.

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

Overdevelopment.

Decrease development time in step 11.

## RELATED PRODUCTS

P-4005	EpiQuik™ HDAC1 Assay Kit
P-4006	EpiQuik™ HDAC2 Assay Kit
P-4007	EpiQuik™ HDAC8 Assay Kit