

HDAC Inhibitor Drug Screening Kit

(Catalog #K340-100; 100 assays; Store kit at -20°C)

I. Introduction:

Inhibition of histone deacetylase (HDAC) has been implicated to modulate transcription and to induce apoptosis or differentiation in cancer cells. However, screening of compounds for HDAC inhibition has been difficult due to the lack of convenient tools for analyzing HDAC activity. The new HDAC Inhibitor Drug Screening Kit provides a fast, fluorescence-based method that eliminates radioactivity, extractions, or chromatography, as used in traditional assays. The new procedure requires only two easy steps, both performed on the same microtiter plate. First, your inhibitor candidates are mixed with HeLa Nuclear Extract and HDAC fluorometric substrate, which comprises an acetylated lysine side chain. Deacetylation of the substrate sensitizes the substrate, so that, in the second step, treatment with the Lysine Developer produces a fluorophore. The fluorophore can be easily analyzed using a fluorescence plate reader or a fluorometer. The assay is well suited for high throughput screening applications.

II. Kit Contents:

Component	K340-100	Color Code
	100 assays	Cap Color
HDAC Substrate [Boc-Lys(Ac)-AMC, 4 mM]	500 µl	Amber
10X HDAC Assay Buffer	1.0 ml	Green
Lysine Developer	1.0 ml	Orange
HDAC Inhibitor (Trichostatin A, 1 mM)	10 µl	Blue
HeLa Nuclear Extract (5 mg/ml)	200 µl	Red
Deacetylated Standard [Boc-Lys-AMC, 4 mM]	20 µl	Yellow

III. Histone Deacetylase Assay Protocol:

A. General Consideration:

- Read the entire protocol before beginning the procedure.
- The HeLa nuclear extract should be refreeze immediately at -70°C after each use to avoid loss of activity.
- The kit provides sufficient reagents for 100 positive control assays with the HeLa Nuclear Extract and 5 Negative Control assays with the HDAC Inhibitor, Trichostatin A.

B. HDAC Assay Protocol:

1. Dilute test samples into 85 µl (final volume) with ddH₂O in each well. For positive reading, add 85 µl ddH₂O only. For negative reading, add 83 µl of ddH₂O and 2 µl of Trichostatin.
2. Add 10 µl of the 10X HDAC Assay Buffer and 2 µl of HeLa Nuclear Extract to each well. Mix thoroughly.
3. Add 5 µl of the HDAC Fluorometric Substrate to each well. Mix thoroughly. Incubate plates at 37°C for 30 minutes (or longer if desired).

4. Stop the reaction by adding 10 µl of Lysine Developer and mix well. Incubate the plate at 37°C for 30 min.
5. Read sample in a fluorescence plate reader with Ex. = 350-380 nm and Em. = 440-460 nm. Signal should be stable for several hours at room temperature. Inhibitor activity can be expressed as inhibition of Relative Fluorescence Units.

C. Standard Curve (optional):

1. If desired, a standard curve can be prepared using the known amount of the Deacetylated Standard included in the kit. The exact concentration range of the Deacetylase Standard will vary depending on the fluorometer model, the gate setting, and the exact wavelength used. We recommend starting with a dilution range of 1-20 µM in Assay Buffer.
2. Add 90 µl each of the dilutions and also 10 µl of the 10X Assay Buffer (as zero) into a set of wells on the microtiter plate.
3. Add 10 µl of Lysine Developer to each well and incubate at 37°C for 30 min (Note: Incubation time should be kept the same for both standard and test samples.)
4. Read samples in a fluorescence plate reader or a fluorometer with Ex. = 350-380 nm and Em. = 440-460 nm.
5. Plot fluorescence signal (y-axis) versus concentration of the Deacetylated Standard (x-axis). Determine the slope as AFU/µM.
6. Based on the slope, you can determine the absolute amount of deacetylated lysine generated in your sample.

RELATED PRODUCTS:

Apoptosis Detection Kits & Reagents

- Annexin V Kits & Bulk Reagents
- Caspase Assay Kits & Reagents
- Mitochondrial Apoptosis Kits & Reagents
- Nuclear Apoptosis Kits & Reagents
- Apoptosis Inducers and Set

Cell Fractionation System

- Mitochondria/Cytosol Fractionation Kit
- Nuclear/Cytosol Fractionation Kit
- Membrane Protein Extraction Kit
- Cytosol/Particulate Rapid Separation Kit
- Mammalian Cell Extraction Kit
- FractionPREP Fractionation System

Cell Proliferation & Senescence

- Quick Cell Proliferation Assay Kit
- Senescence Detection Kit
- LDH-Cytotoxicity Assay Kit

Cell Damage & Repair

- HDAC Fluorometric & Colorimetric Assays Kits
- HAT Colorimetric Assay Kit & Reagents
- DNA Damage Quantification Kit
- Glutathione & Nitric Oxide Fluorometric & Colorimetric Assay Kits
- cAMP & cGMP Assay Kits