

Cellular HDAC-Glo Class I/II Assay

Cell-based epigenetic assay for the detection of the relative activity of histone deacetylase (HDAC) class I/II enzymes from cells.

Target Overview

Cell Line THP-1 human acute monocytic leukemia cell line from ATCC

Assay Properties

Assay Format HDAC-Glo

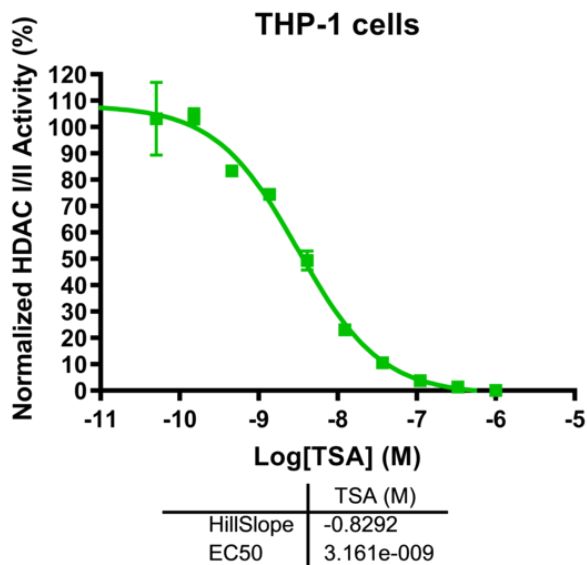
Assay Protocol THP-1 cells were seeded into the wells of 384-well plate. The cells were treated with reference compound TSA (starting at 1 μM , 10-dose with 3-fold dilution) for 1 hour. HDAC Class I/II activity was measured with HDAC-Glo Class I/II Assay on an Envision 2104 Multilabel Reader. IC50 value was calculated and IC50 curve was plotted using the GraphPad Prism program based on a sigmoidal dose-response equation.

Readout Measurement for deacetylation of the peptide substrate via luminescence

Reference Compound IC50s

Compound	IC50 (μM)
TSA	0.003161

Screening location Malvern, PA, USA



Let's Discover Together.

© Aug 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/hdac-glo_classi-ii_cellular_epigenetic_malvern

