

Cellular HDAC-Glo Class IIa Assay

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

Target Overview

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

Assay Properties

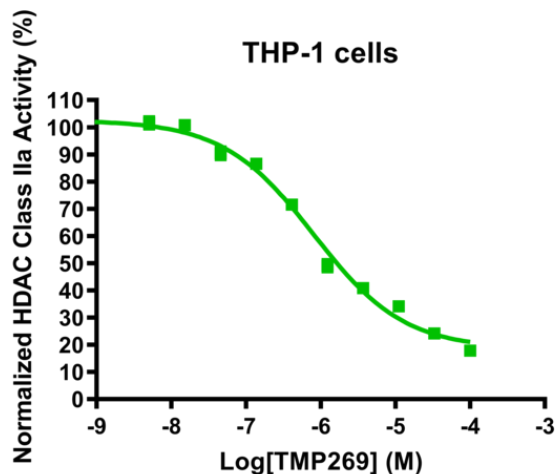
Assay Format HDAC-Glo

Assay Protocol THP-1 cells were seeded into the wells of 384-well plate. The cells were treated with reference compound TMP269 (starting at 100 μ M, 10-dose with 3-fold dilution) for 1 hour. HDAC Class IIa activity was measured with HDAC-Glo Class IIa 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 (μ M)
 TMP269 0.8095

Screening location Malvern, PA, USA



	TMP269 (M)
HillSlope	-0.7177
EC50	8.095e-007

Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/hdac-glo_classiia_cellular_epigenetic_malvern

