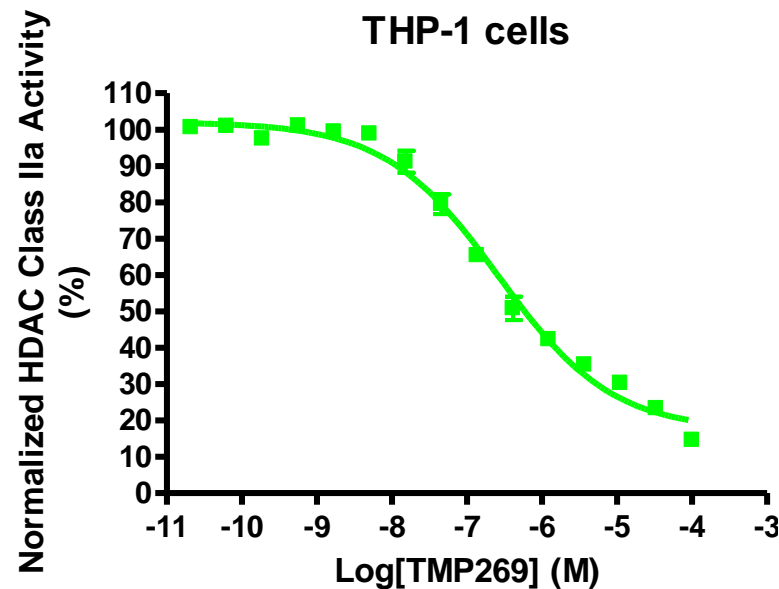


HDAC-Glo Class IIa Assay

THP-1 Cells: Human acute monocytic leukemia cell line from ATCC.

Assay format: HDAC-Glo Class IIa cell based assay.

Assay Protocol: THP-1 cells were seeded into each well of the 384-well plate. The cells were treated with various concentrations of compound TMP269 for 1 hour. HDAC-Glo Class IIa Detection Reagent was added to each well of the plate. Luminescence signal was recorded by Envision 2104 Multilabel Reader. The IC50 (EC50) value was calculated and the IC50 curve was plotted using the GraphPad Prism 4 program based on a sigmoidal dose-response equation.



	TMP269 (M)
HILLSLOPE	-0.5747
EC50	2.641e-007