

Cellular Phospho-Histone Assay

[Cell-based epigenetic assay](#) Phospho-Histone H3 ELISA Assay

Target Overview

Cell Line HT-29 human colon cell line from ATCC

Assay Properties

Assay Format ELISA

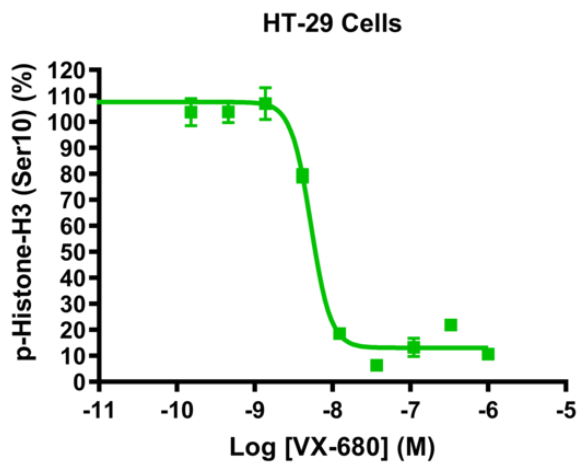
Assay Protocol HT-29 cells were seeded into the wells of 96-well plate. The cells were pre-treated with reference compound VX-680 (starting at 1 μ M, 10-dose with 3-fold dilution) for 90 minutes and then treated with phosphatase inhibitor Calyculin A for 15 minutes. Cells were lysed with cell lysis buffer and cell lysates were collected. Phospho-Histone H3 (serine 10) was measured with ELISA 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 Phospho-Histone H3

Reference Compound IC50s

Compound	IC50 (μ M)
VX-680	0.005288

Screening location Malvern, PA, USA



	VX-680
HillSlope	-3.393
EC50	5.288e-009

Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/phospho-histone_elisa_cellular_epigenetic_malvern

