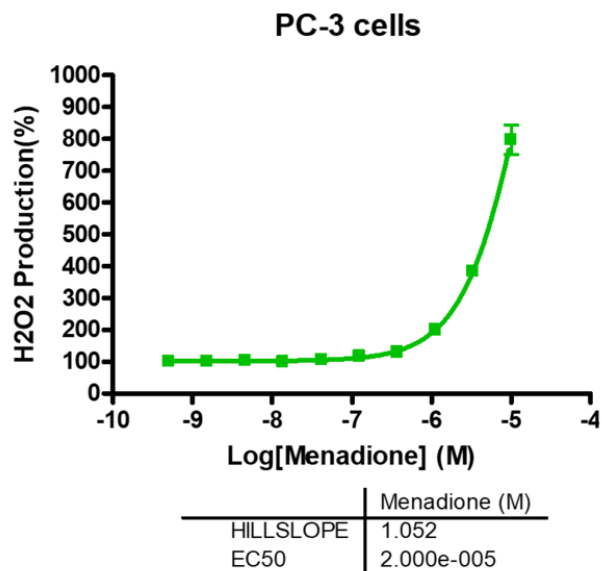


ROS-GLO H2O2 Assay

[Apoptosis assays](#) are available for single and multi-time point analysis as well as real-time live-cell imaging.

Target Overview	
Cell Line	PC-3 cell line from ATCC
Assay Properties	
Assay Format	384-well plate format using ROS-GLO H2O2 Assay System
Assay Protocol	PC-3 cells were treated with various concentrations of compound menadione in 384-well plate for 6 hours. ROS H2O2 production was measured with ROS-Glo H2O2 assay. 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 program based on a sigmoidal dose-response equation.
Readout	Luminescent signal
Reference	Compound IC50 (M)
Compound IC50s	Menadione 2.000e-005
Screening Location	Malvern, PA, USA



Let's Discover Together.

© Dec 2021 Reaction Biology. All rights reserved.

https://www.reactionbiology.com/datasheet/ros-glo_h2o2_assay_malvern

