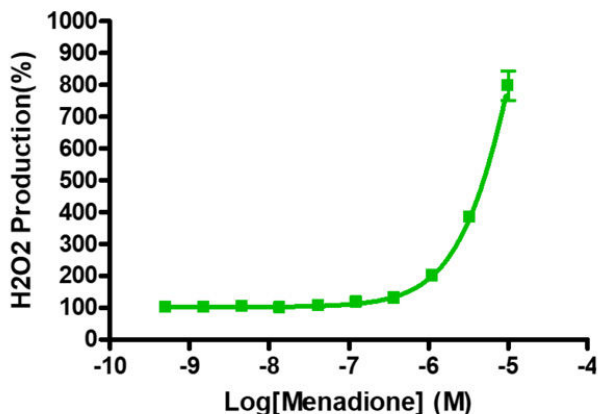


ROS-GLO H2O2 Assay

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

Assay Properties	
Cell Line	PC-3 cell line from ATCC
Assay Format	384-well plate format using ROS-GLO H ₂ O ₂ Assay System
Assay Protocol	PC-3 cells were treated with various concentrations of compound menadione in 384-well plate for 6 hours. ROS H ₂ O ₂ production was measured with ROS-Glo H ₂ O ₂ assay. Luminescence signal was recorded by Envision 2104 Multilabel Reader. The IC ₅₀ (EC ₅₀) value was calculated and the IC ₅₀ curve was plotted using the GraphPad Prism program based on a sigmoidal dose-response equation.
Readout	Luminescent signal
Reference Compound IC ₅₀ s	Compound IC₅₀ (M) Menadione 2.000e-005
Screening Location	Malvern, PA, USA
Further information	More information can be found on our website Apoptosis Assay Services for Drug Discovery

PC-3 cells



	Menadione (M)
HILLSLOPE	1.052
EC50	2.000e-005

Let's Discover Together.

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

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

