

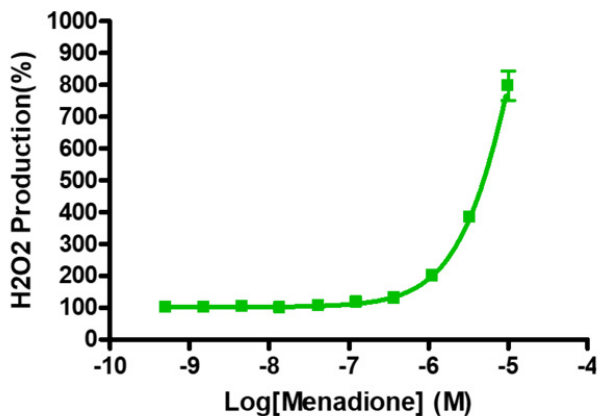
# ROS-GLO H<sub>2</sub>O<sub>2</sub> Assay

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

## Assay Properties

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

### PC-3 cells



	Menadione (M)
HILLSLOPE	1.052
EC50	2.000e-005

