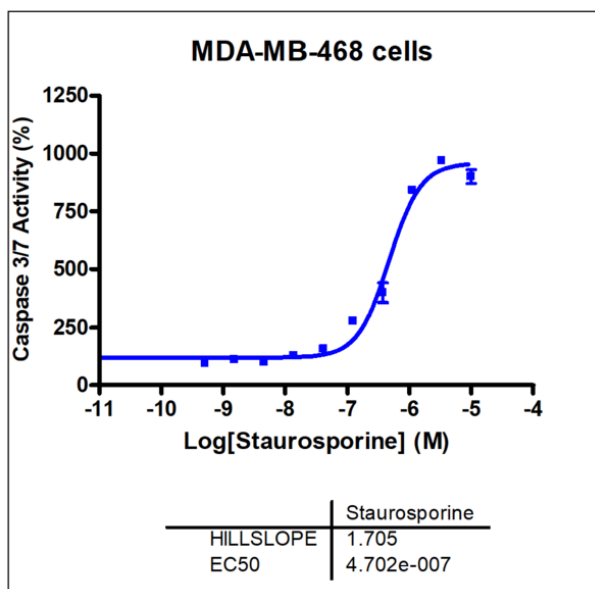


Caspase-Glo 3/7 Assay

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

Assay Properties

Cell Line	MDA-MB-468 cell line from ATCC				
Assay Format	384-well plate format using Promega's Caspase-Glo® 3/7 Assay System				
Assay Protocol	MDA-MB-468 cells were seeded into the wells of 384-well plates. The cells were treated with reference compound Staurosporine for 72 hours. Luminescence was measured 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	Luminescent signal				
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compound</th> <th>IC50 (M)</th> </tr> </thead> <tbody> <tr> <td>Staurosporine</td> <td>4.702e-007</td> </tr> </tbody> </table>	Compound	IC50 (M)	Staurosporine	4.702e-007
Compound	IC50 (M)				
Staurosporine	4.702e-007				
Screening Location	Malvern, PA, USA				



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/caspase-glo_37_assay_malvern

