

Caspase-Glo 3/7 Assay

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

Target Overview

Cell Line MDA-MB-468 cell line from ATCC

Assay Properties

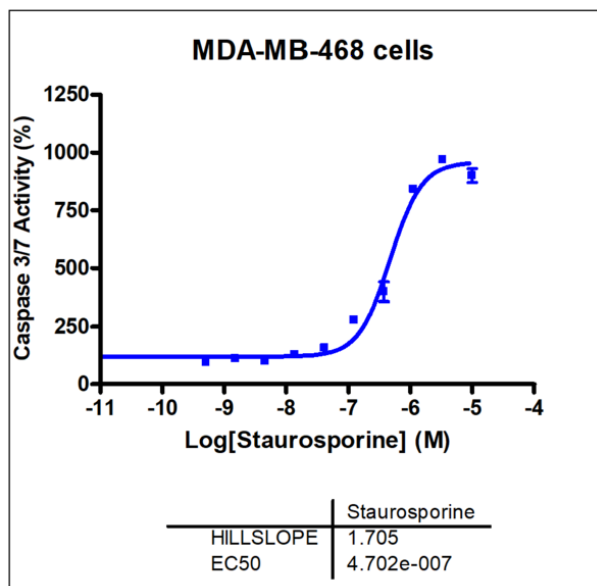
Assay Format 384-well plate format using Promega's Caspase-Glo® 3/7 Assay System

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	Compound	IC50 (M)
	Staurosporine	4.702e-007

Screening Location Malvern, PA, USA



Let's Discover Together.

