

## Caspase-Glo 3/7 Activity Assay

**Cell Line:** MDA-MB-231 human breast cancer cell line was purchased from ATCC.

**Assay Format:** Caspase-Glo 3/7 luminescence assay.

**Assay Protocol:** MDA-MB-231 cells were treated with various concentrations of compound staurosporine in 384-well plate for 18 hours.

Caspase-3/7 activation was measured with Caspase-Glo 3/7 luminescence 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 4 program based on a sigmoidal dose-response equation.

