

Information for 100 cell line panel service

Test date:

Testing will start on: February 01, May 01, July 01, September 01, and November 01

Compound Treatment Time

72 hours

Readout

CellTiter-Glo® Luminescent

Testing Format

IC50 (10-dose with 3-fold dilution in duplicate) in 384-well plate format

Compound Submission

Dissolved in DMSO or as pre-weighed powder form

Tolerable Solvent

Max 0.2 % DMSO

Reference Compound

included (staurosporine)

Report

IC50 curves and IC50 values including raw data for own analysis