













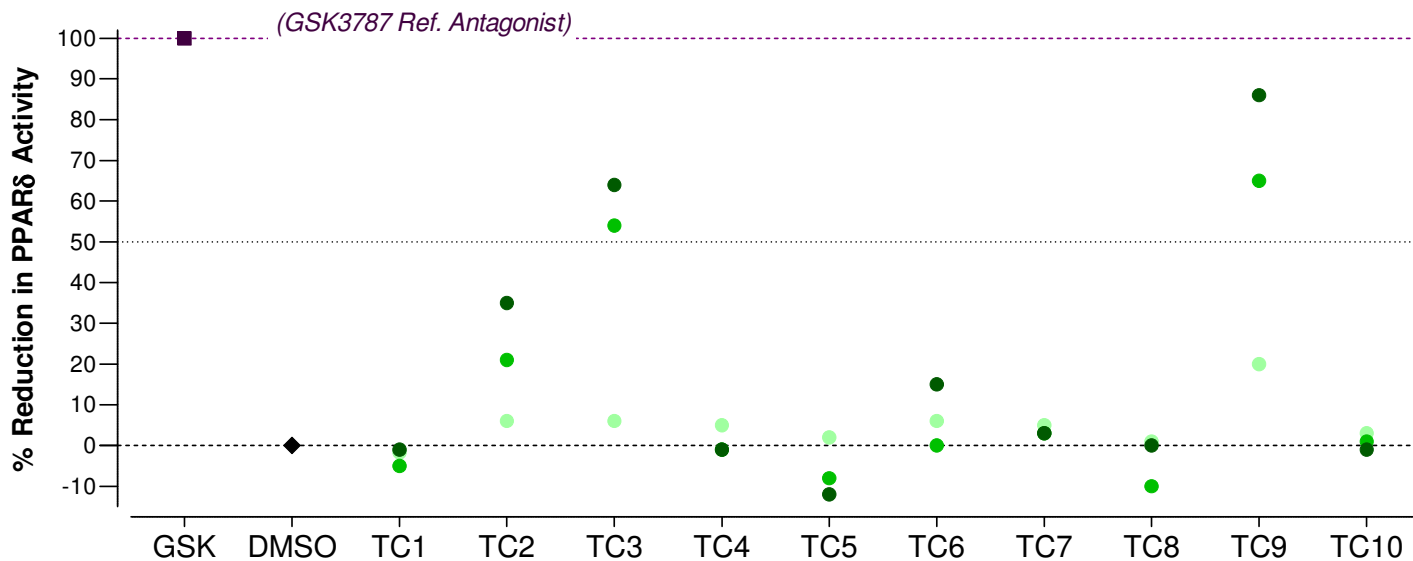








Example of Dot-plot format data presentation: analyses of 1 – 3 concentrations of test compounds.



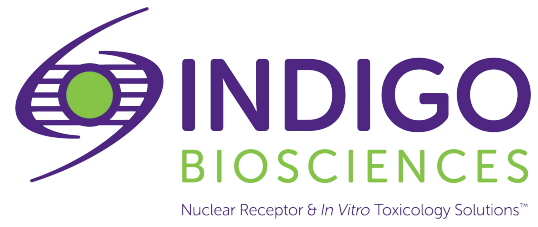
**Human PPARδ antagonist assay.**

Test compounds ● 5 μM, ● 1.67 μM, and ● 0.55 μM.

◆ 0.1% DMSO = 0% inhibition. ■ 40 μM GSK3787 (reference antagonist).

PPARδ Reporter cells are co-treated with ~EC<sub>80</sub> of GW0742 as the challenge agonist.

Percent Inhibition is:  $[100 * (1 - (RLU^{Test\ Cmpd} / RLU^{Vehicle\ control}))]$ .



*(~ End of Report ~)*