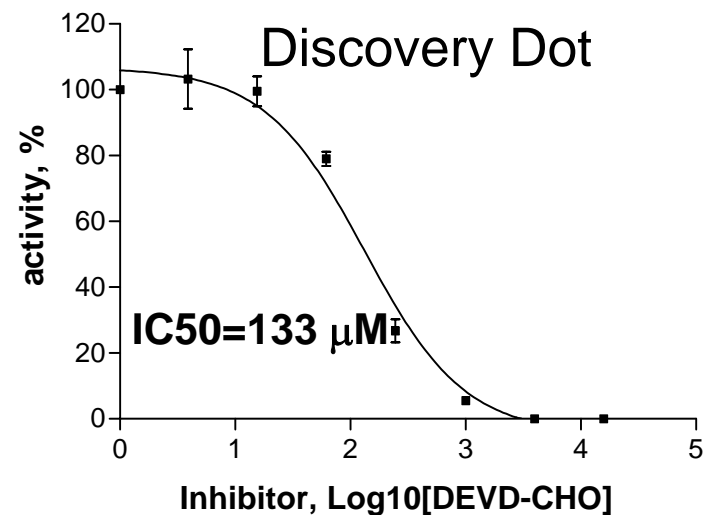
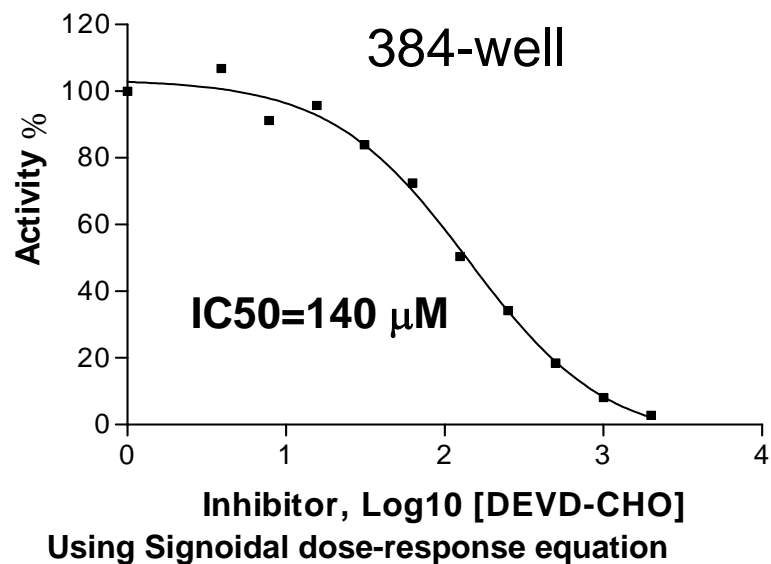


## Caspase 6 IC50: 384 Well – Discovery Dot Comparison



Caspase 6, Substrate: Z-DEVD-R110 (50  $\mu$ M ) Inhibitor: Z-DEVD-CHO