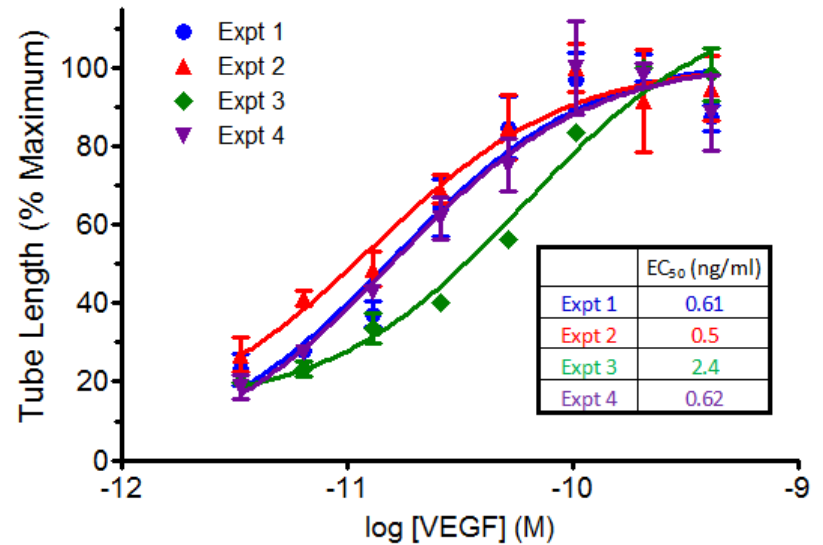


VEGF concentration response using 24 wells of a 96-well plate (8 concentrations n=3).



Resulting EC50 taken at endpoint for 4 successive experiments