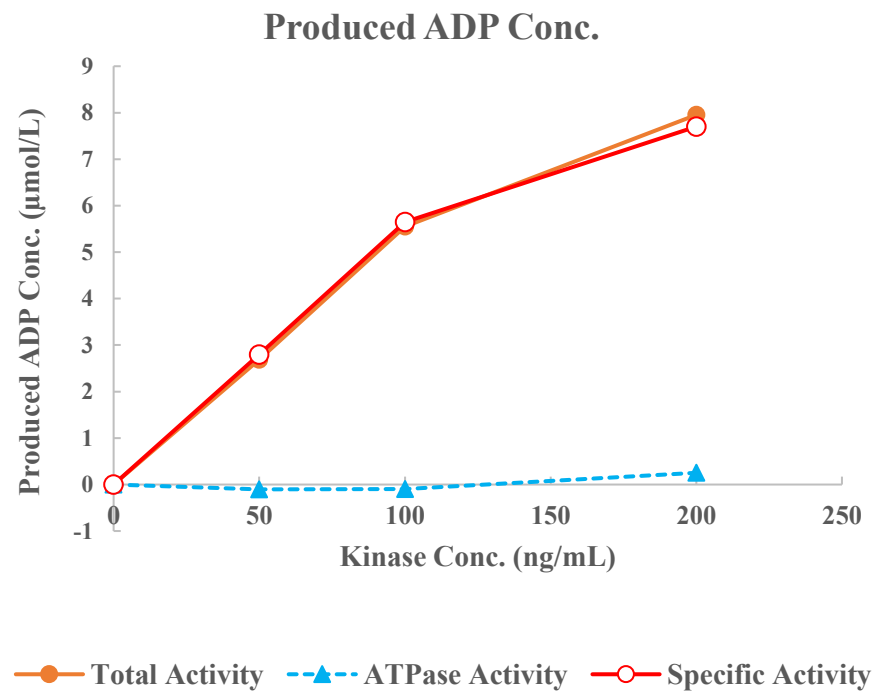
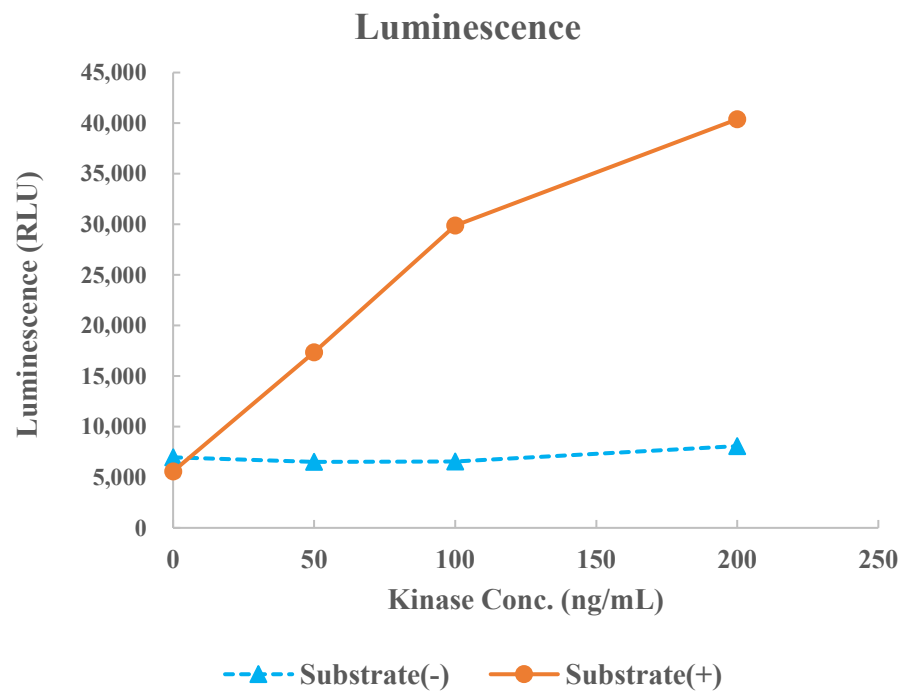


FGFR4 [V550L]



Kinase	FGFR4 [V550L], 08-526
Substrate	Biotin-CSKtide, BTSU02
Substrate Conc.	10 μmol/L
ATP Conc.	150 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour