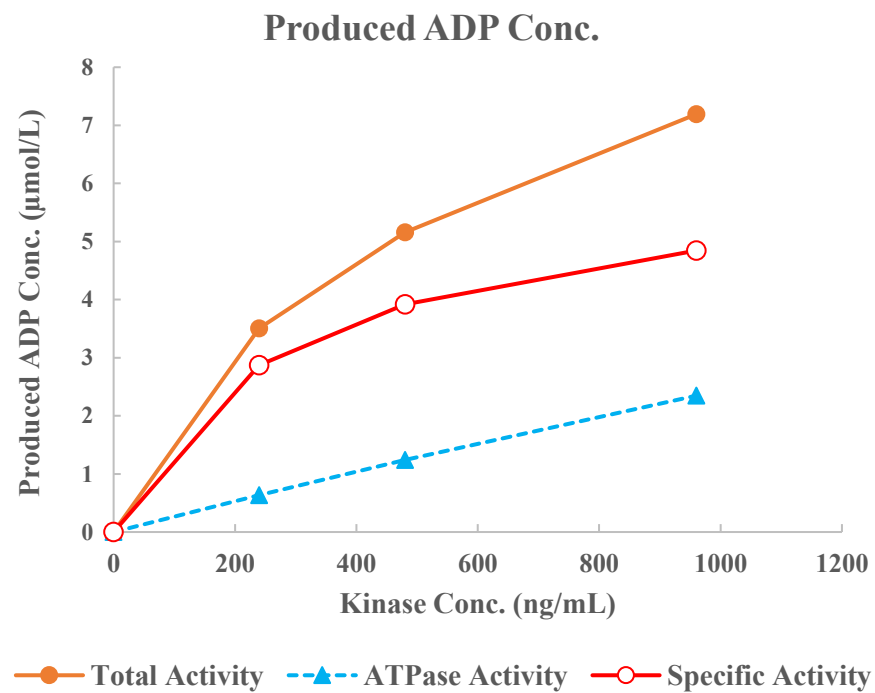
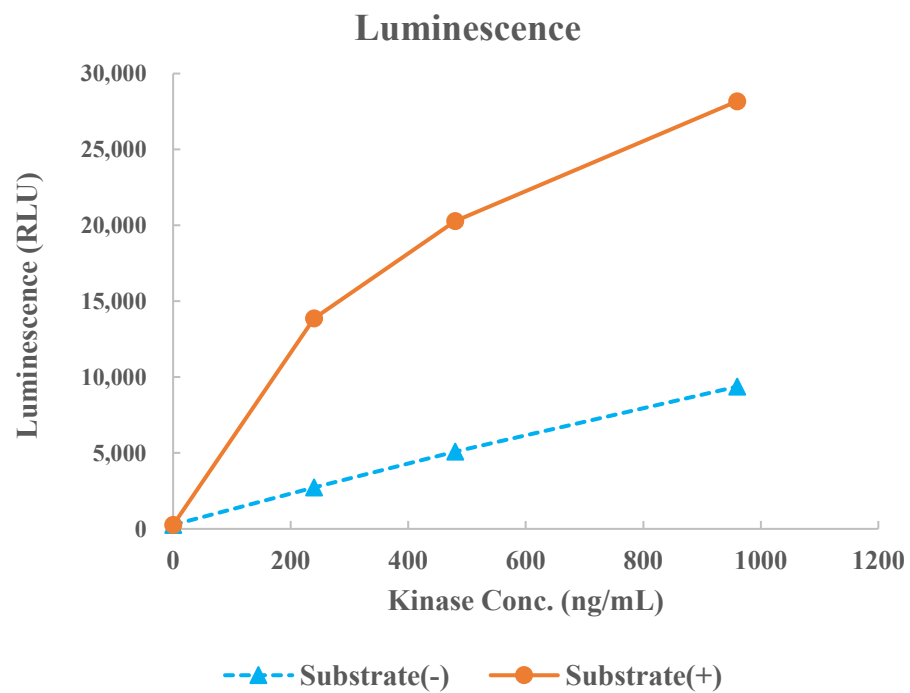


EGFR [C797S/L858R]



Kinase	EGFR [C797S/L858R], 08-563
Substrate	Biotin-Srcptide, BTSU01
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg, Mn
Metal Conc.	5, 1 mmol/L
Additives	-
Reaction Time	1 hour