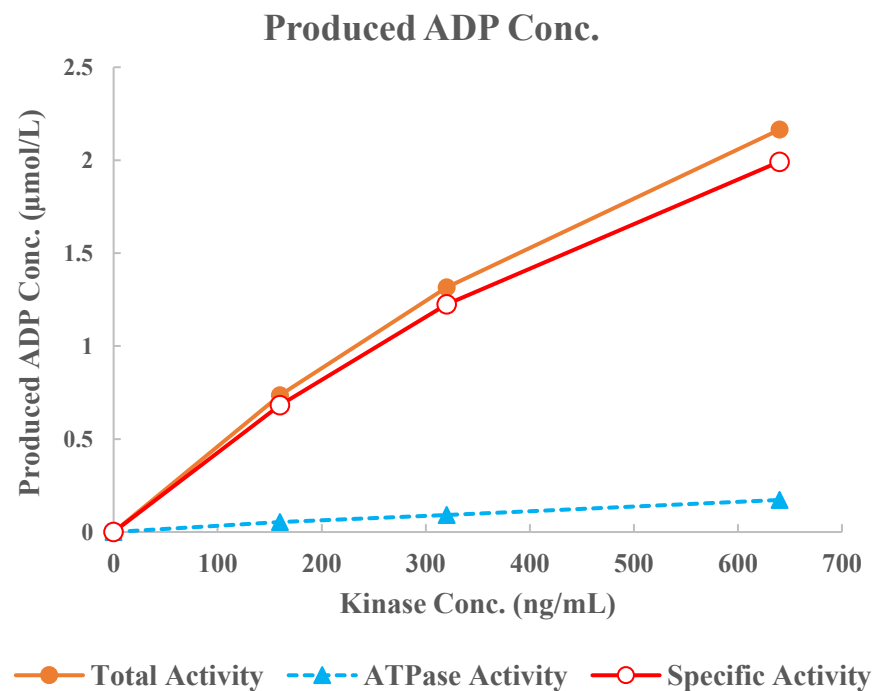
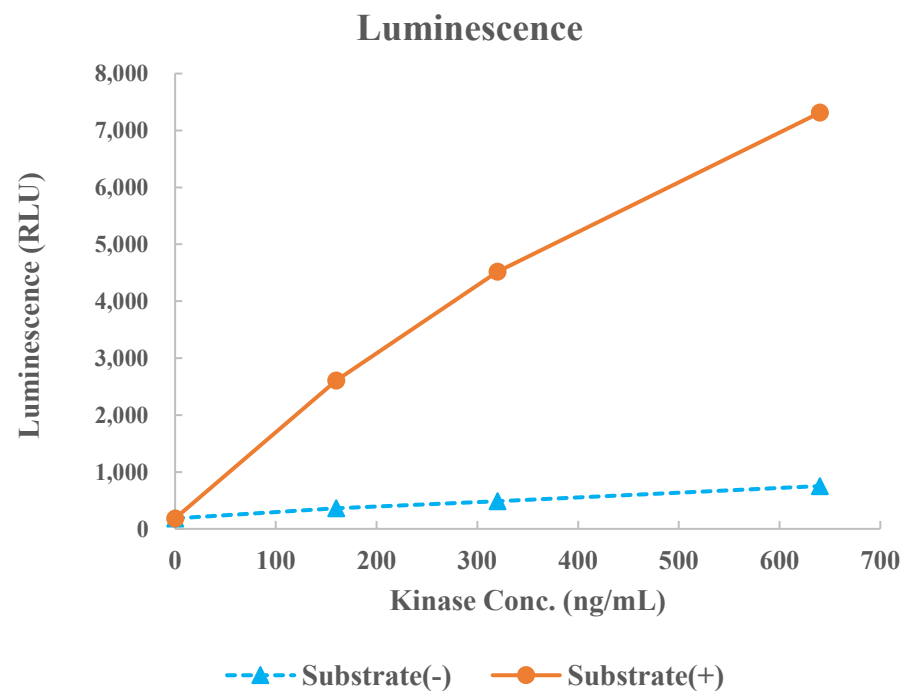


EGFR [d746-750/T790M]



Kinase	EGFR [d746-750/T790M], 08-528
Substrate	Biotin-Srcide, BTSU01
Substrate Conc.	10 μmol/L
ATP Conc.	5 μmol/L
Metal	Mg, Mn
Metal Conc.	5, 1 mmol/L
Additives	-
Reaction Time	1 hour