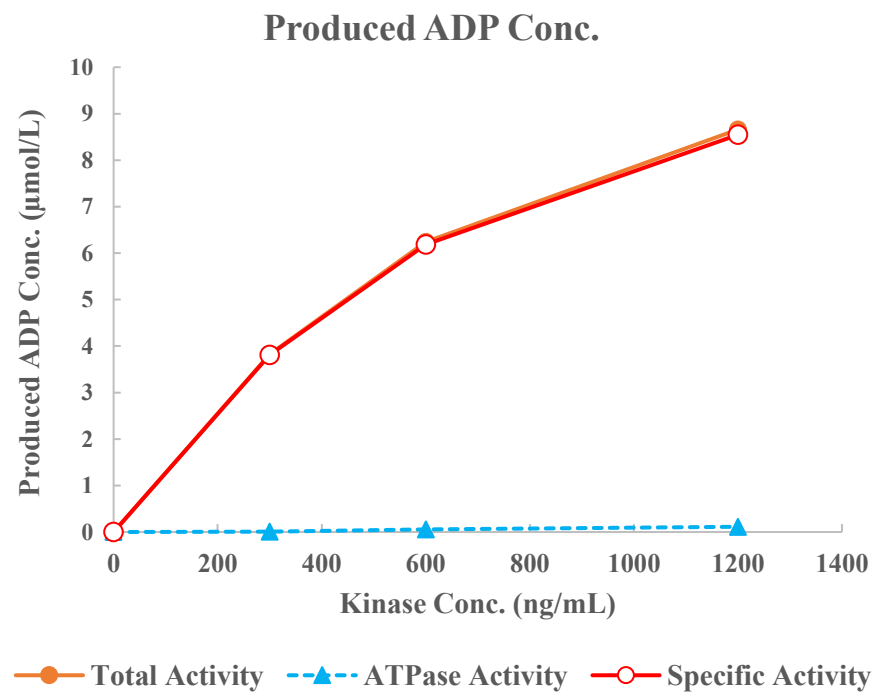
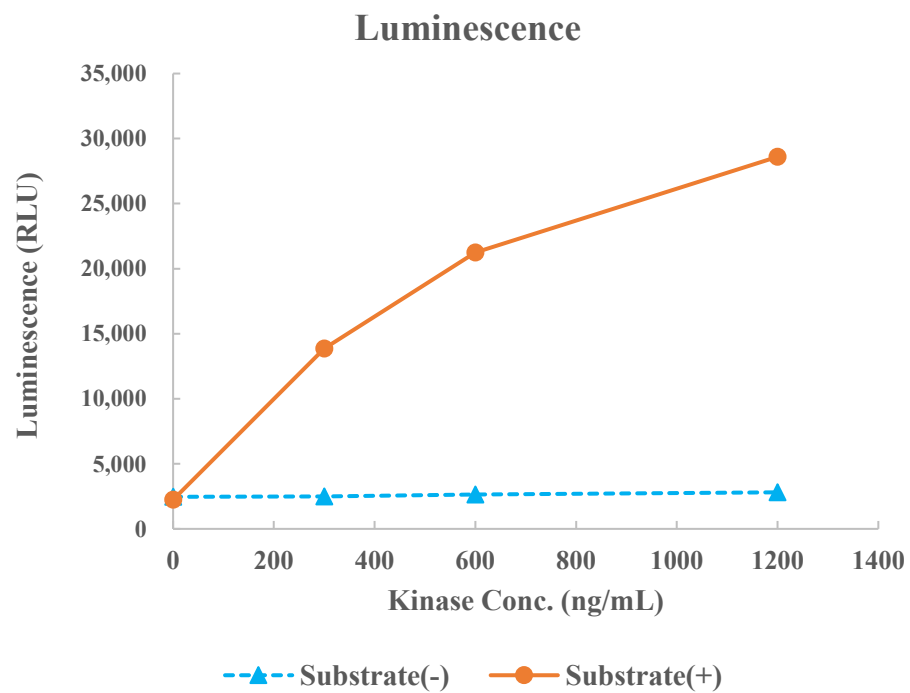


ALK [G1202R]



Kinase	ALK [G1202R], 08-544
Substrate	Biotin-Src tide, BTSU01
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour