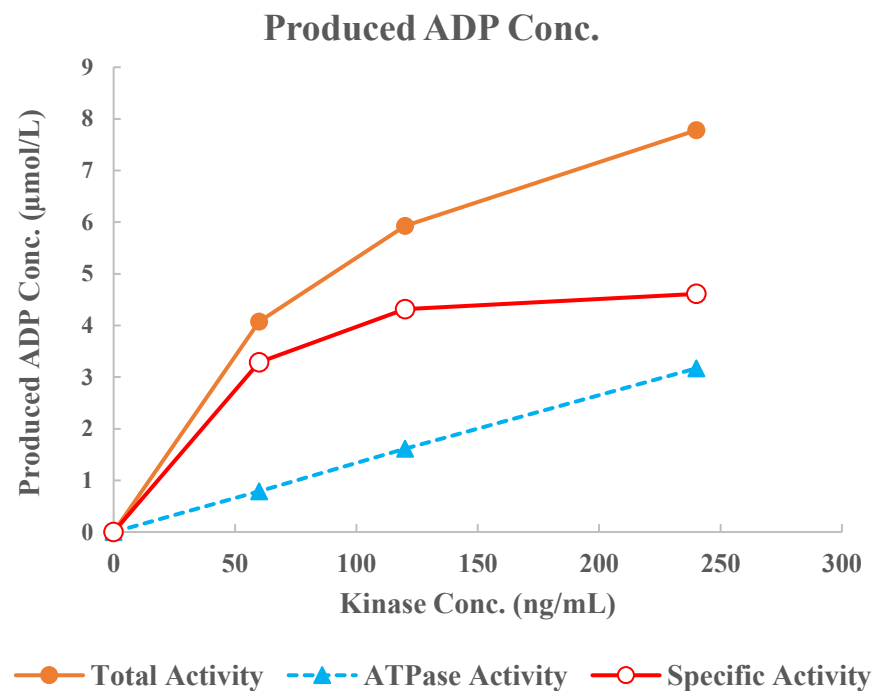
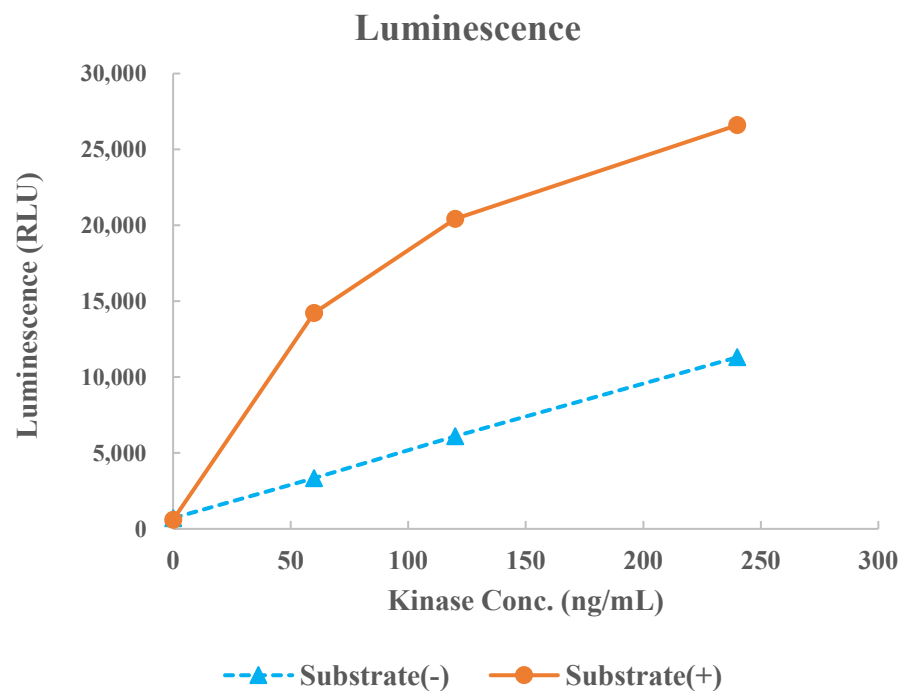


HPK1(MAP4K1)



Kinase	HPK1(MAP4K1), 07-410
Substrate	Biotin-S6K2 peptide, BTSU14
Substrate Conc.	10 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	1.25 mmol/L
Additives	-
Reaction Time	1 hour