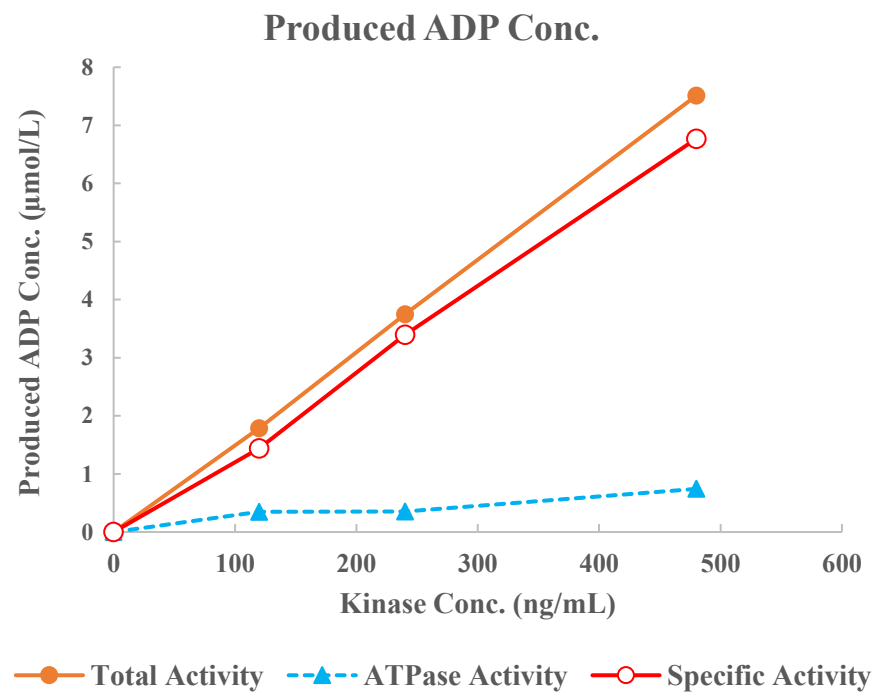
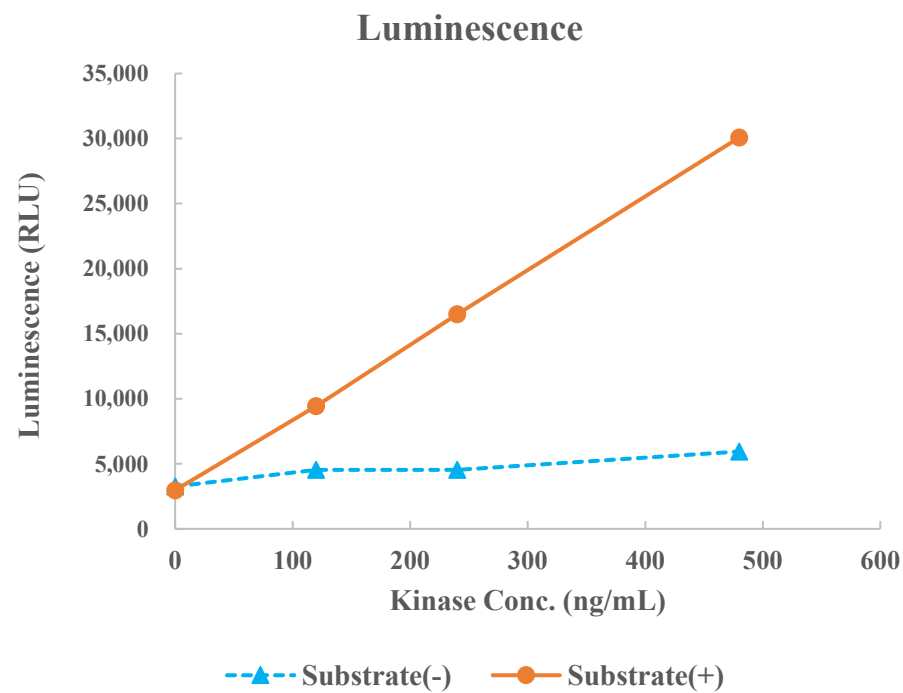


p38 α (MAPK14)



Kinase	p38 α (MAPK14), 04-152
Substrate	Biotin-Modified Erktide, BTSU03
Substrate Conc.	10 μ mol/L
ATP Conc.	150 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour