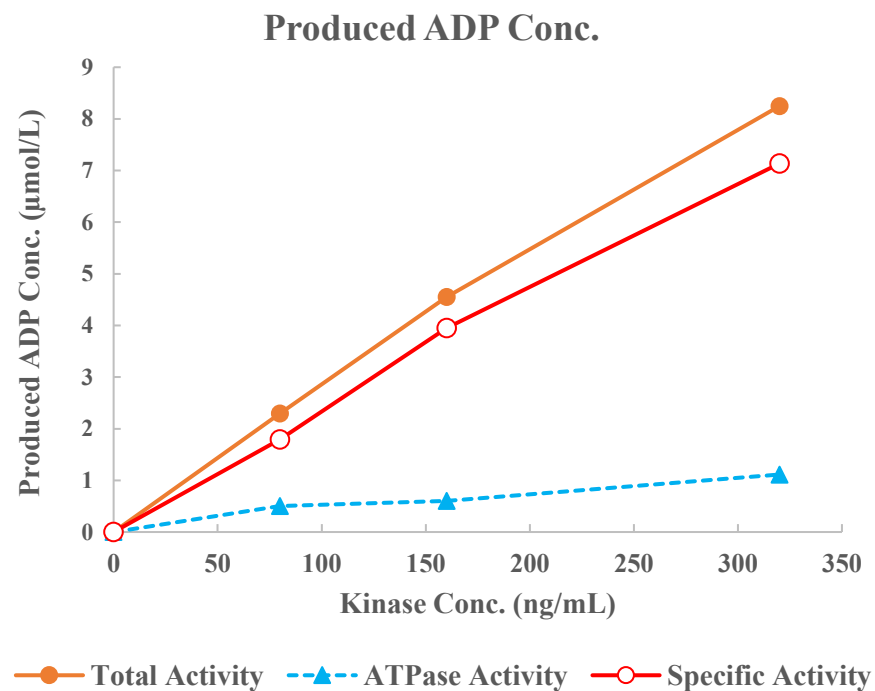
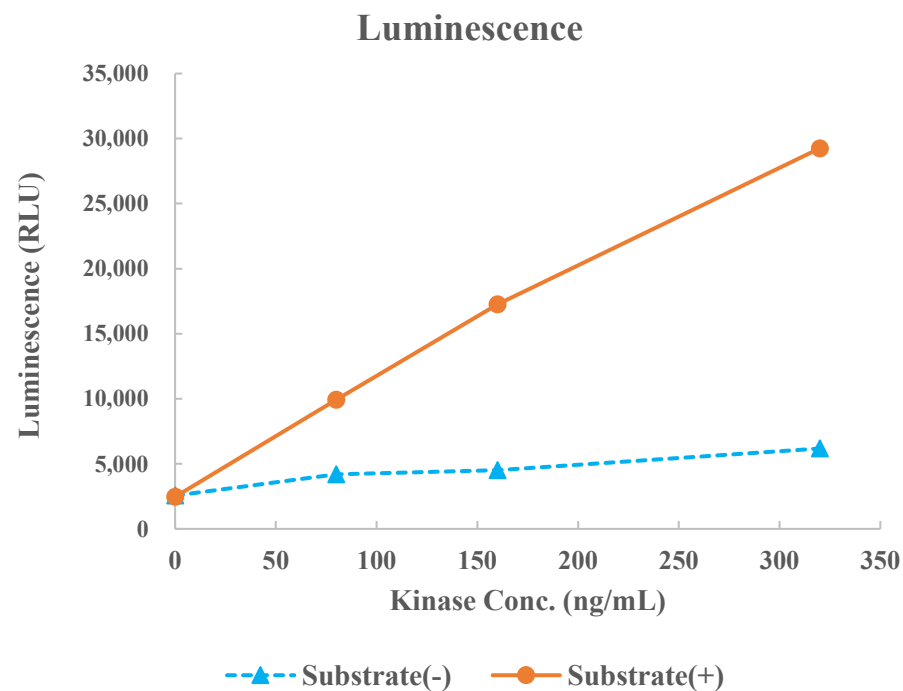


p38 β (MAPK11)



Kinase	p38 β (MAPK11), 04-153
Substrate	Biotin-Modified Erktide, BTSU03
Substrate Conc.	10 μ mol/L
ATP Conc.	75 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour