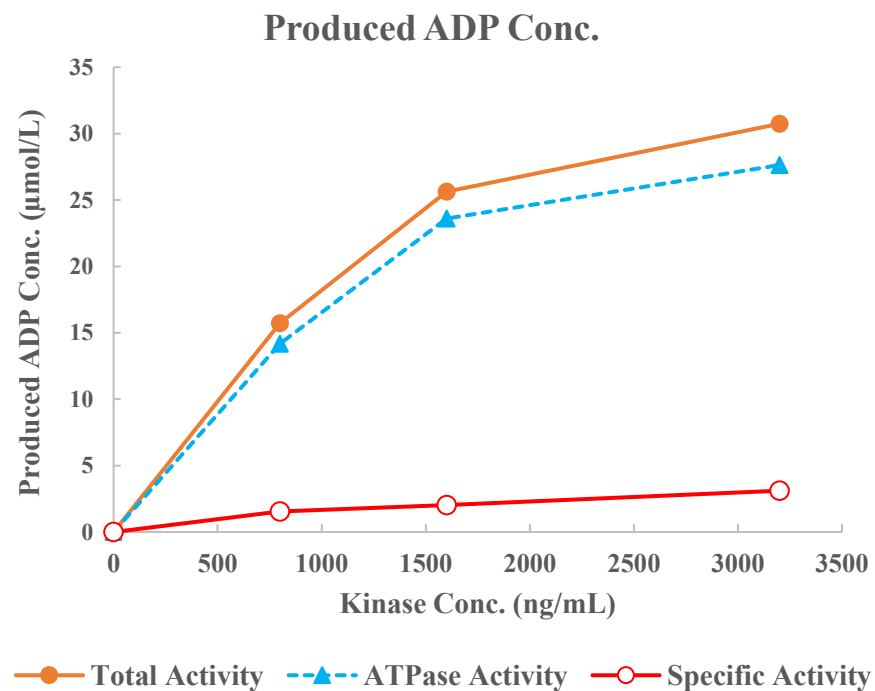
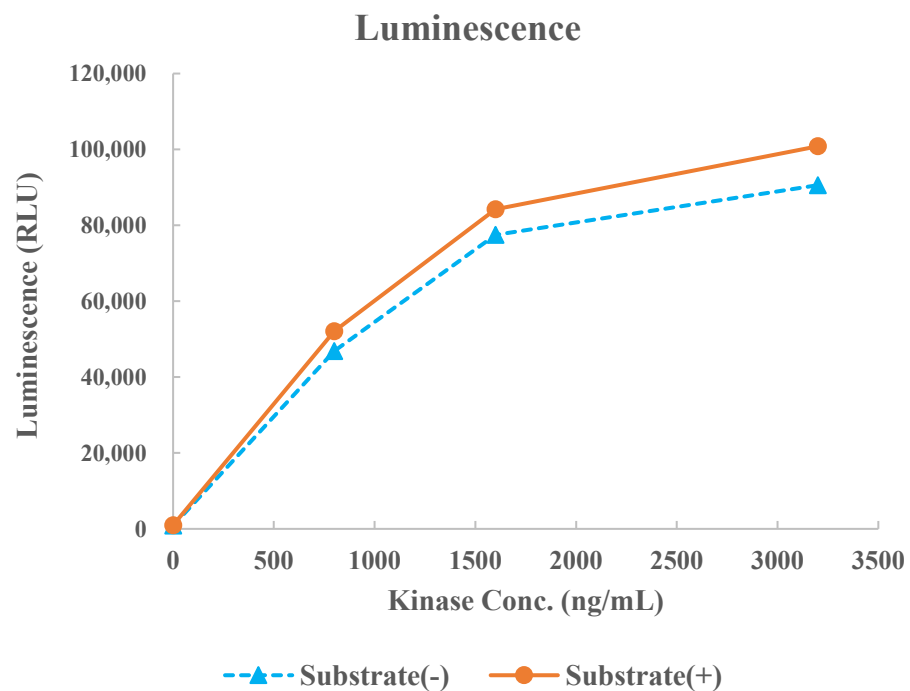


JNK3(MAPK10)



Kinase	JNK3(MAPK10), 04-150
Substrate	Biotin-Modified Erktide, BTSU03
Substrate Conc.	10 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour