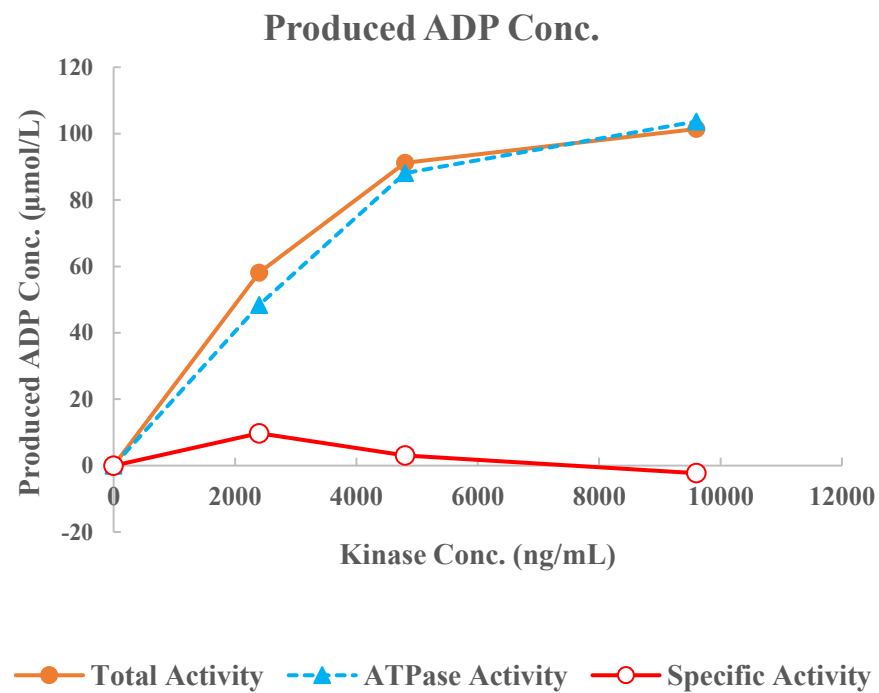
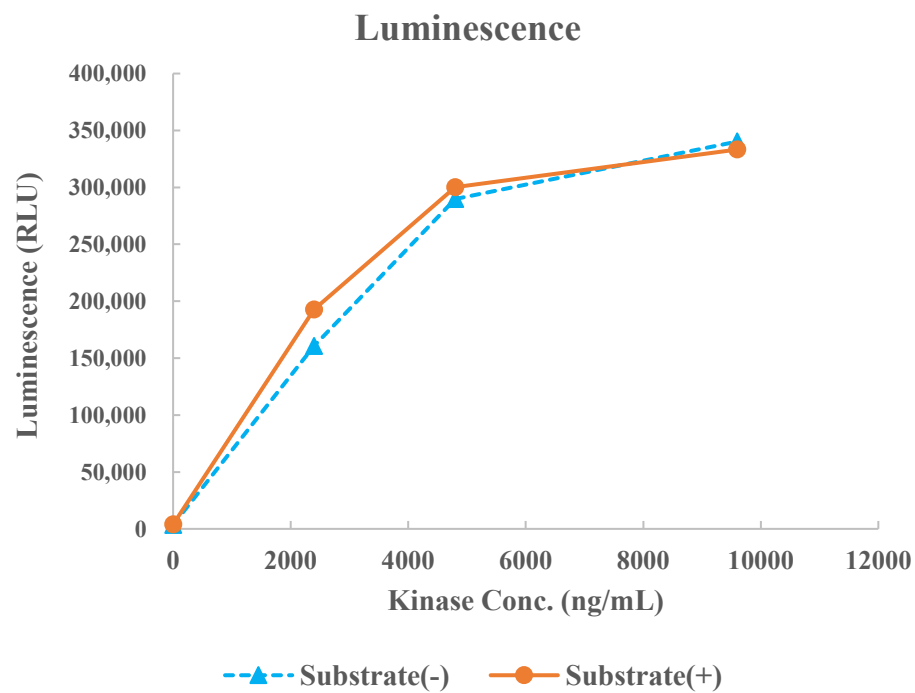


JNK1(MAPK8)



Kinase	JNK1(MAPK8), 04-163
Substrate	Biotin-Modified Erktide, BTSU03
Substrate Conc.	10 μmol/L
ATP Conc.	100 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour