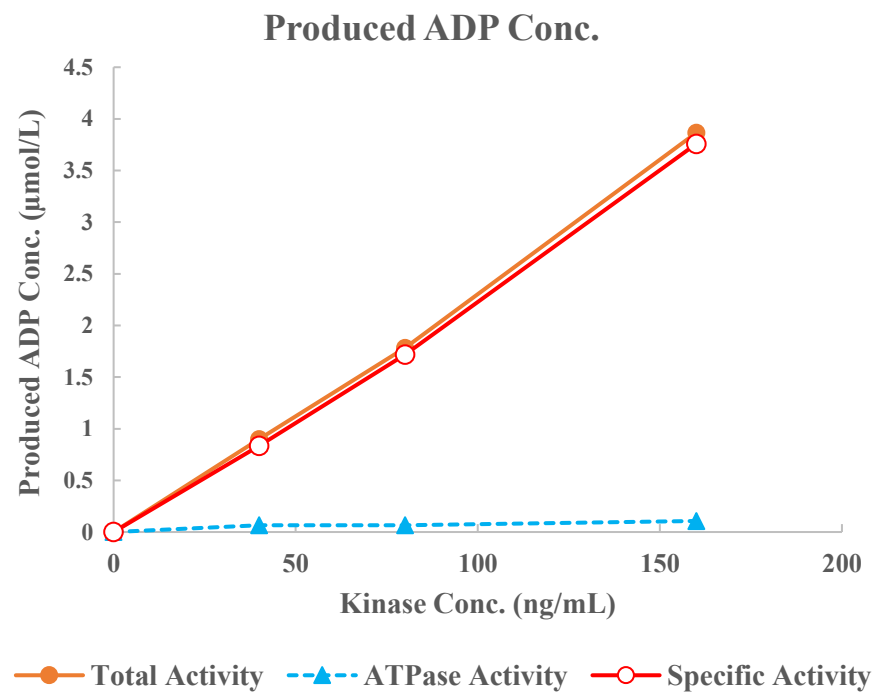
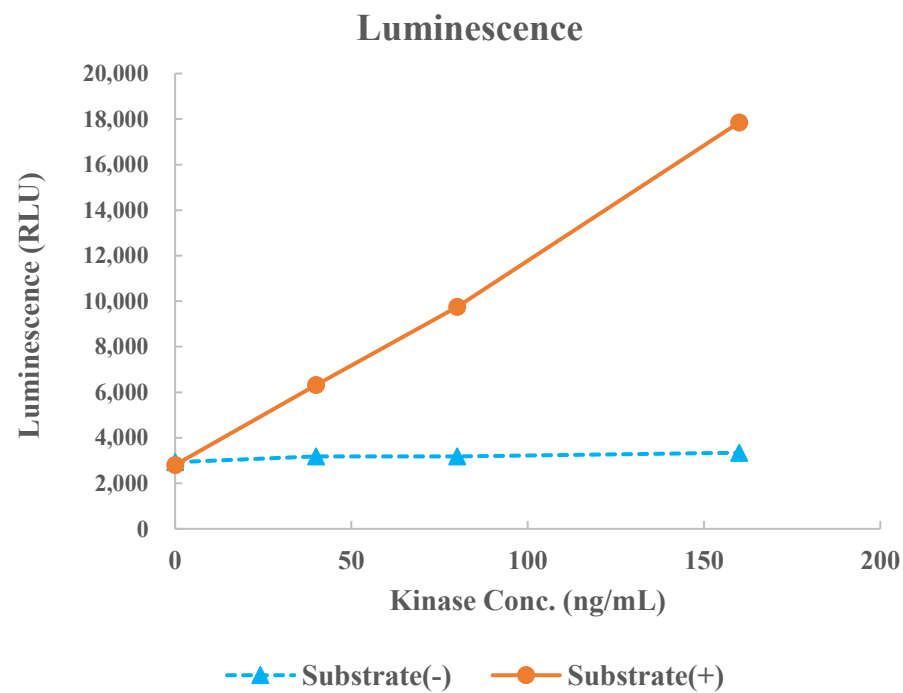


Erk2(MAPK1)



Kinase	Erk2(MAPK1), 04-143
Substrate	Biotin-Modified Erktide, BTSU03
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour