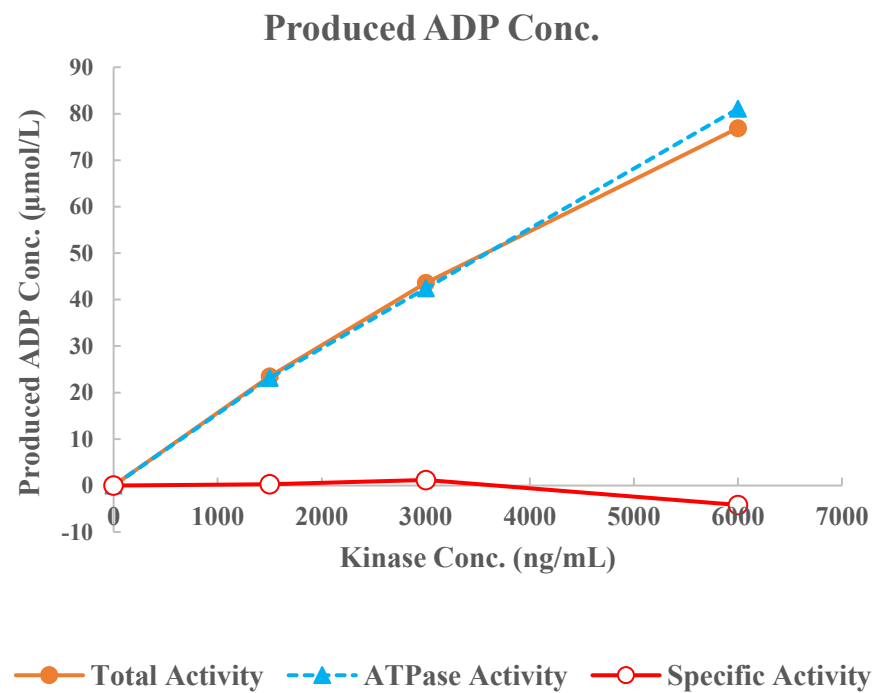
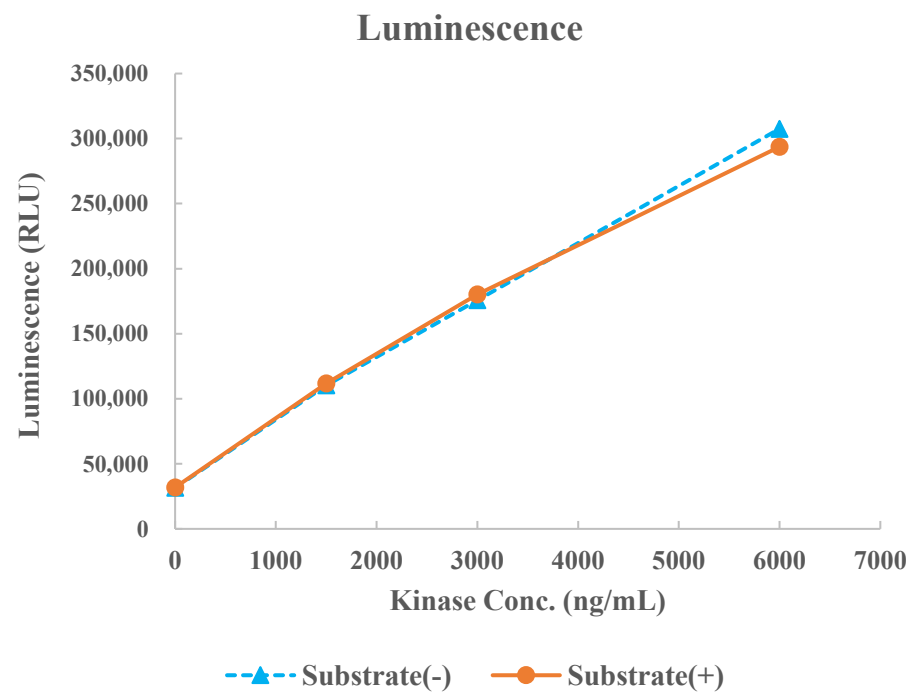


Erk5(MAPK7)



Kinase	Erk5(MAPK7), 04-146
Substrate	Biotin-EGFR-derived peptide, BTSU31
Substrate Conc.	10 μmol/L
ATP Conc.	1000 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour