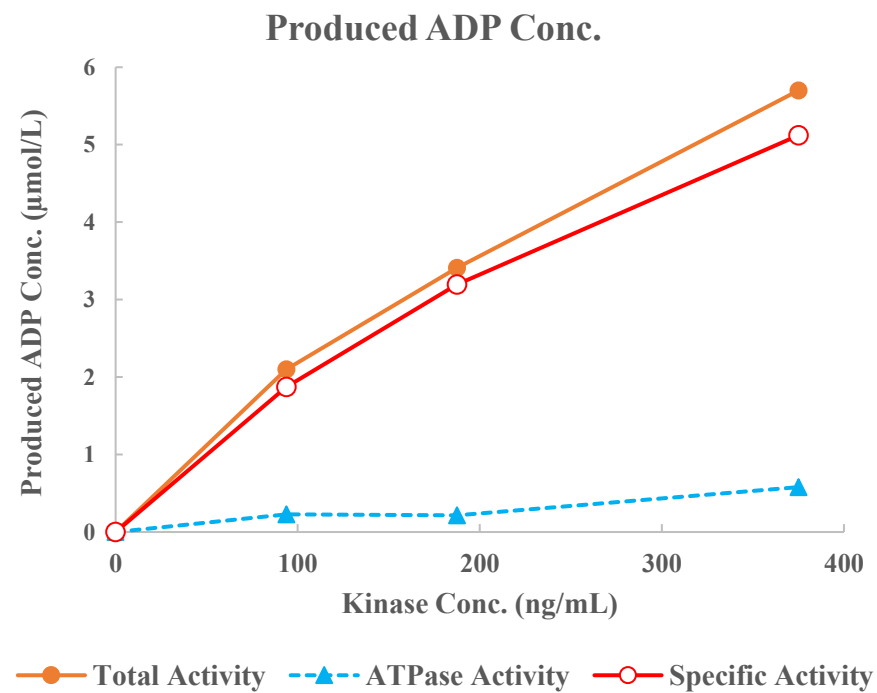
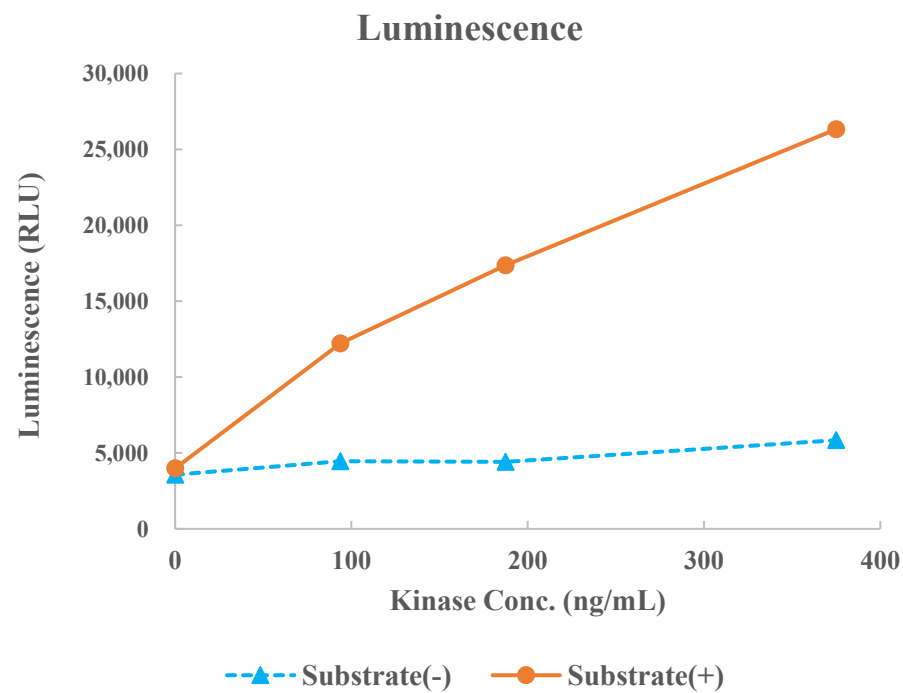


MAP4K2



Kinase	MAP4K2, 07-111
Substrate	Biotin-S6K2 peptide, BTSU14
Substrate Conc.	10 μmol/L
ATP Conc.	100 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour