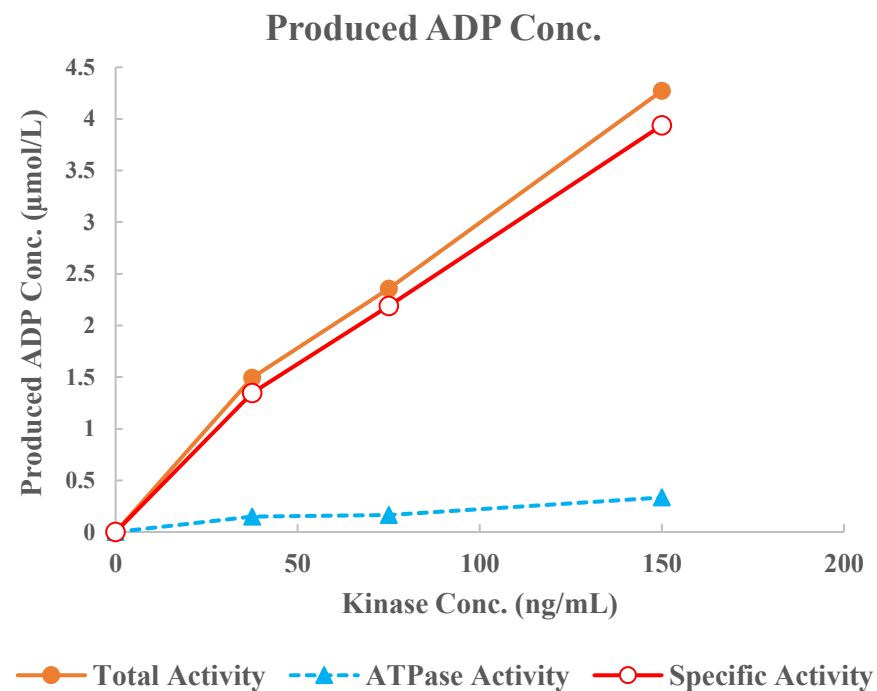
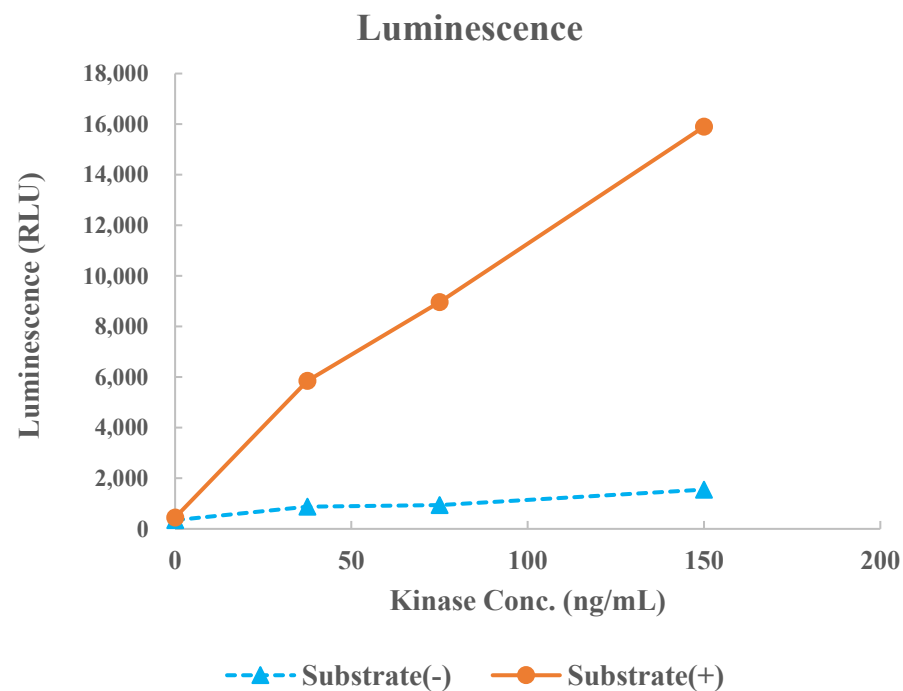


HGK(MAP4K4)



Kinase	HGK(MAP4K4), 07-137
Substrate	Biotin-Moesin-derived peptide, BTSU18
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour