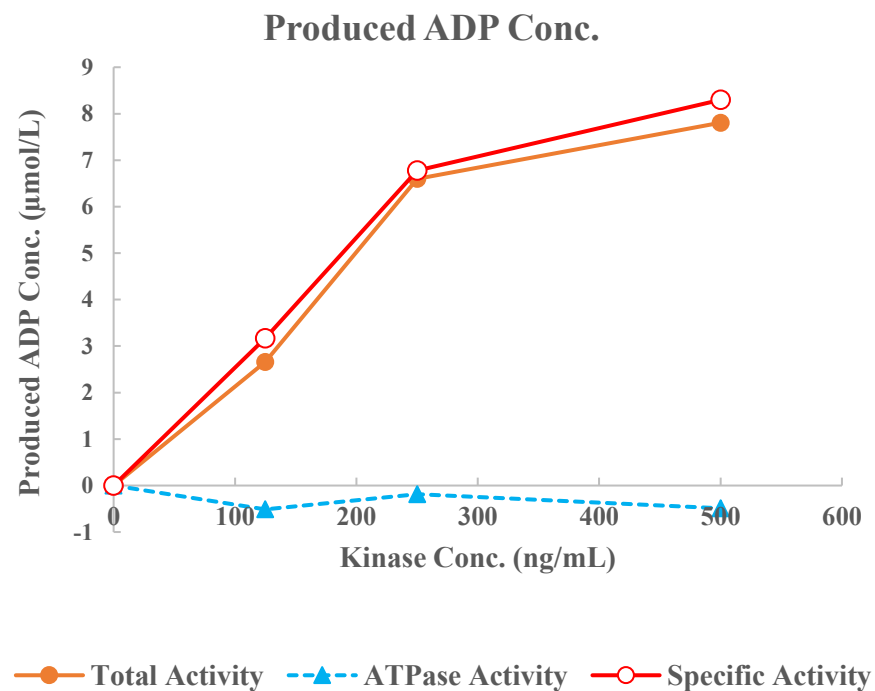
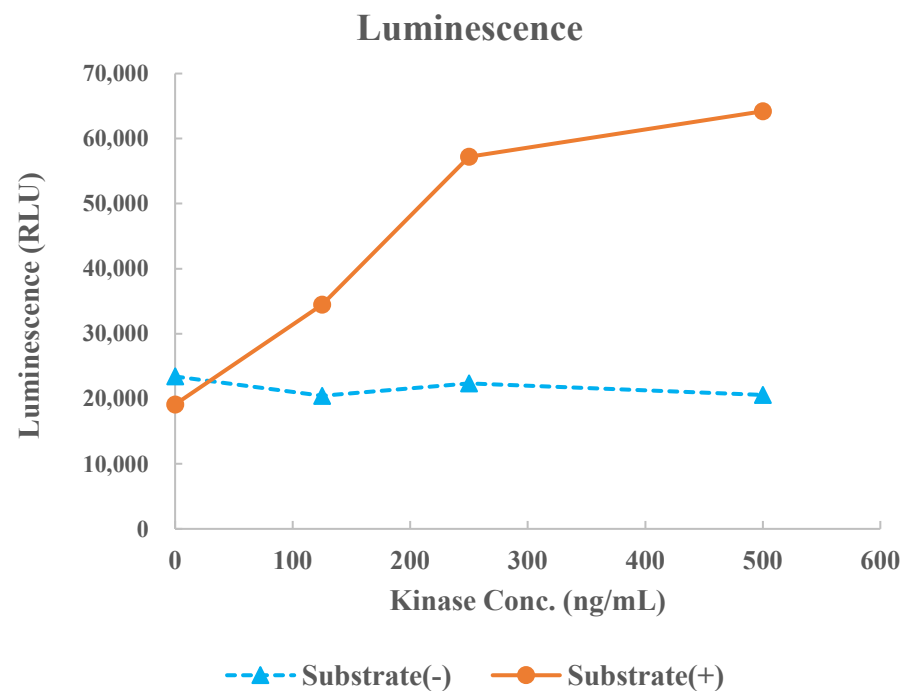


MNK1(MKNK1)



Kinase	MNK1(MKNK1) , 02-145
Substrate	Biotin-SRPKtide, BTSU70
Substrate Conc.	10 μmol/L
ATP Conc.	450 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour