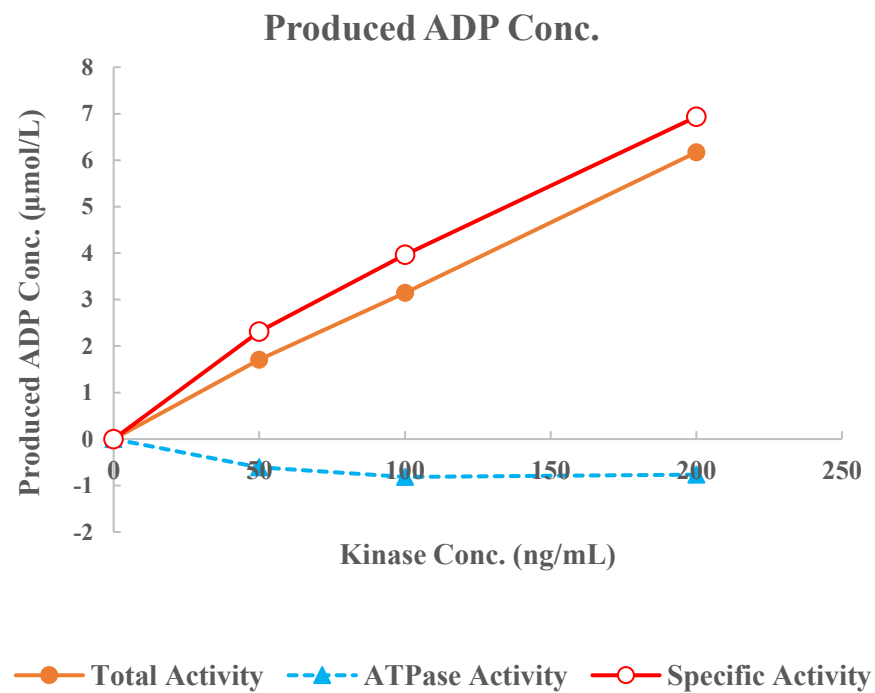
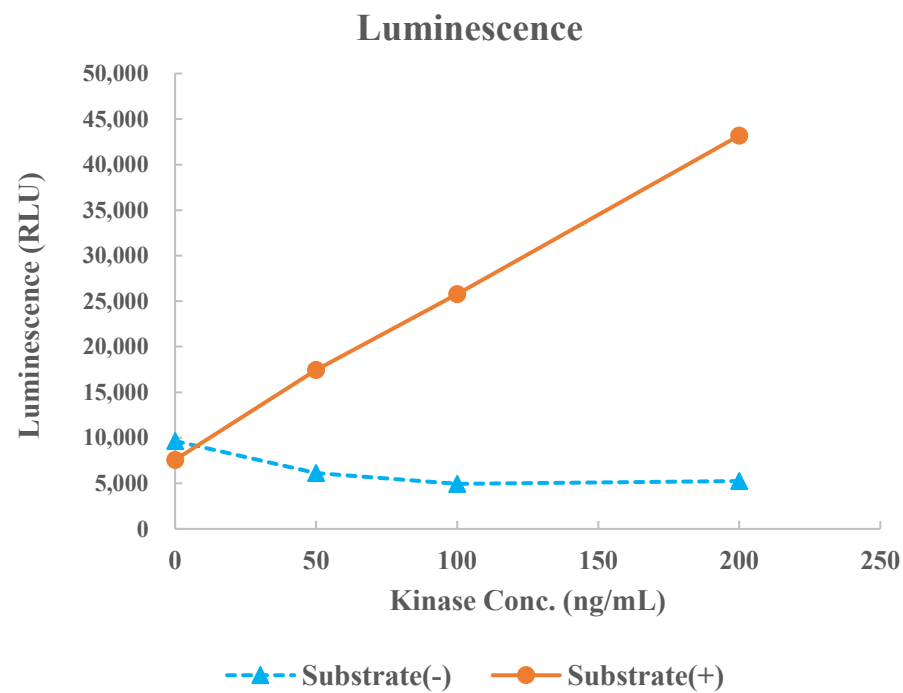


MNK2(MKNK2)



Kinase	MNK2(MKNK2), 02-146
Substrate	Biotin-SRPKtide, BTSU70
Substrate Conc.	10 μmol/L
ATP Conc.	100 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour