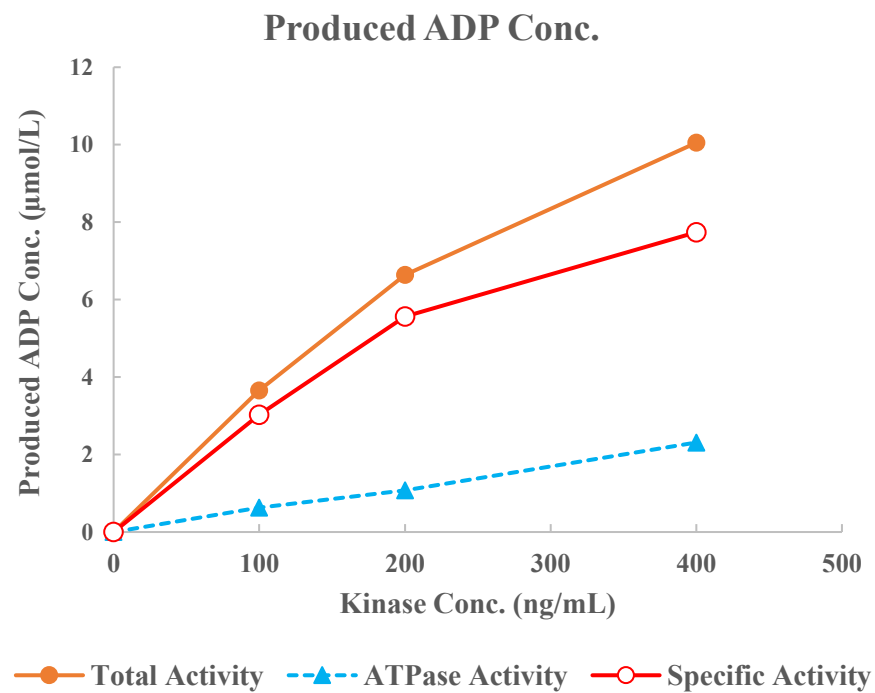
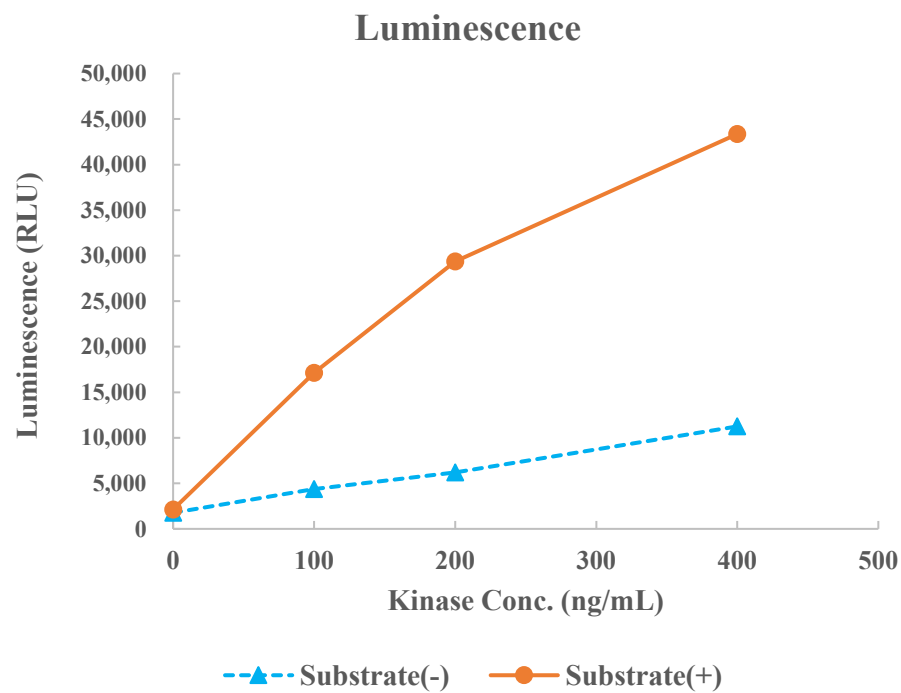


TAK1-TAB1(MAP3K7)



Kinase	TAK1-TAB1(MAP3K7), 09-019
Substrate	Biotin-LRRKtide, BTSU48
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	1.25 mmol/L
Additives	-
Reaction Time	1 hour