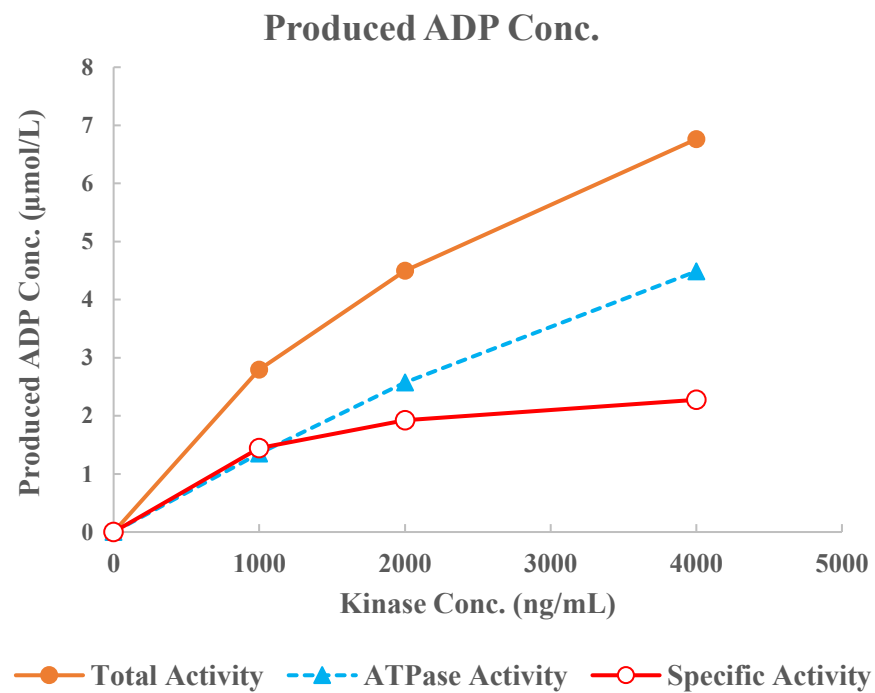
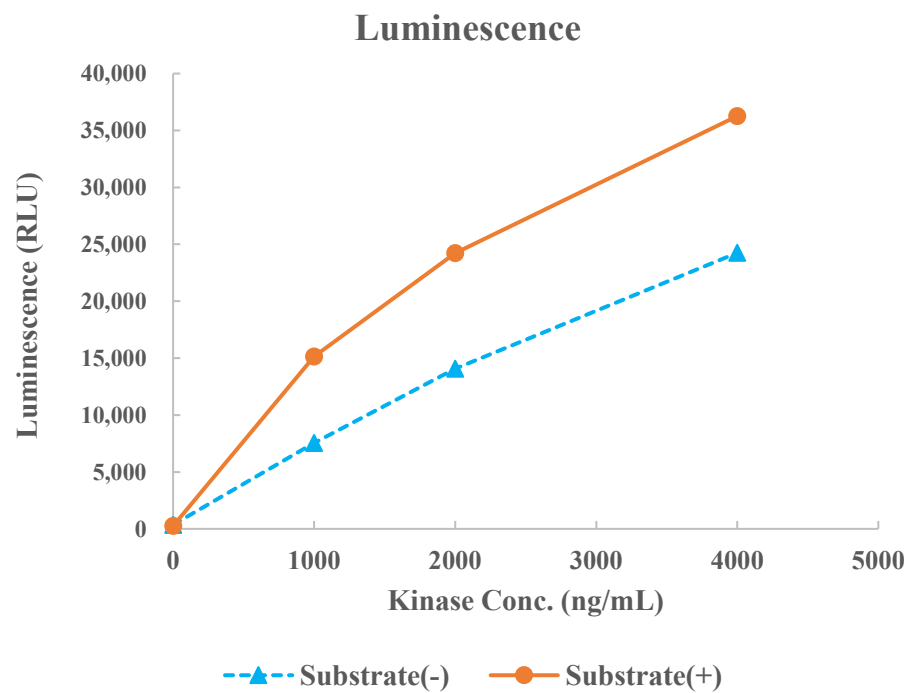


PDK1(PDPK1)



Kinase	PDK1(PDPK1), 01-132
Substrate	Biotin-T308tide, BTSU38
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	2 μmol/L PIftide, 20 μmol/L Cantharidin
Reaction Time	1 hour