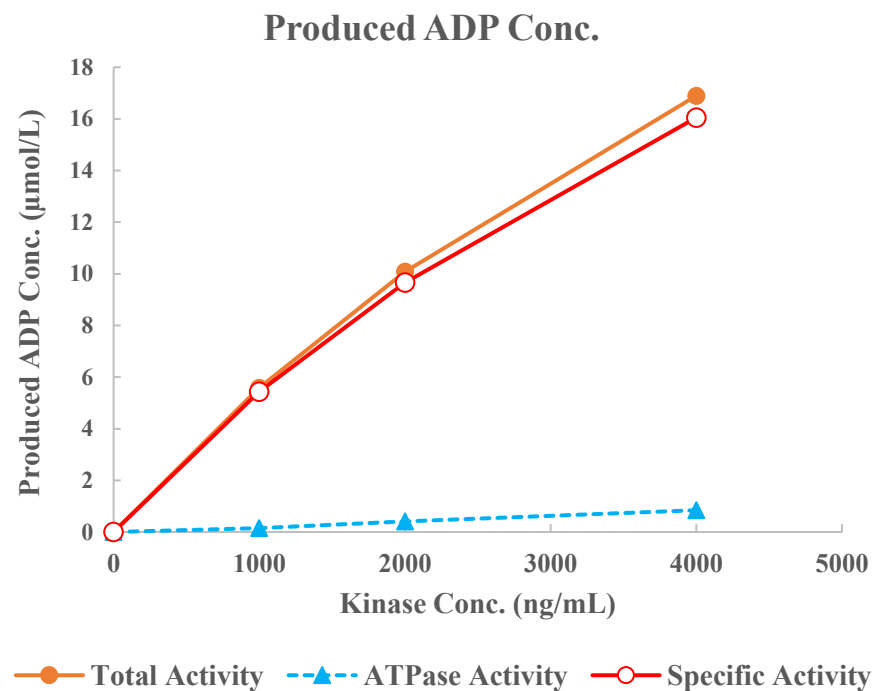
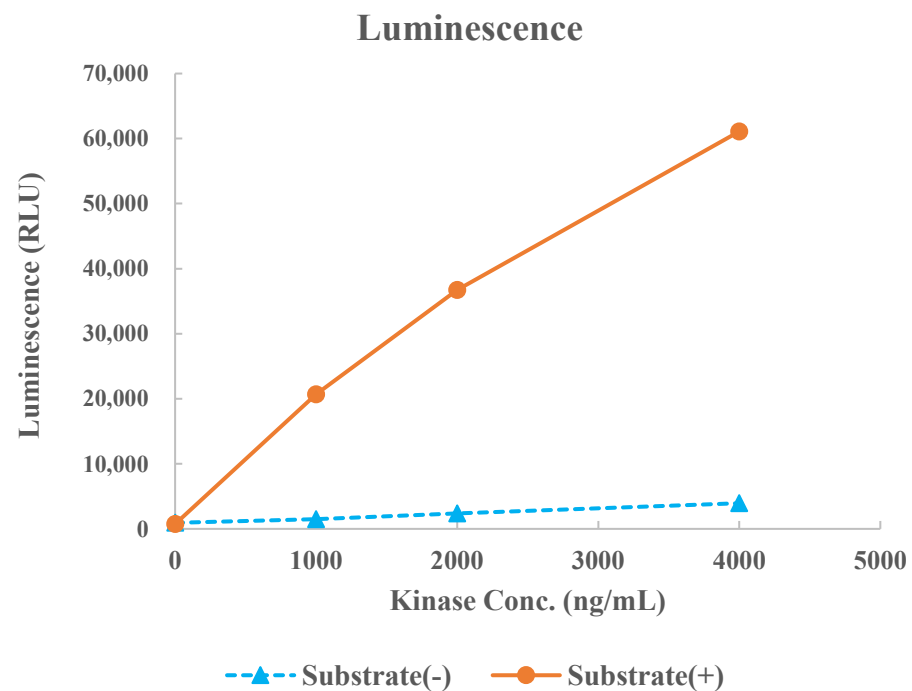


SIK3(QSK)



Kinase	SIK3(QSK), 02-130
Substrate	Biotin-AMARA peptide, BTSU28
Substrate Conc.	50 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour