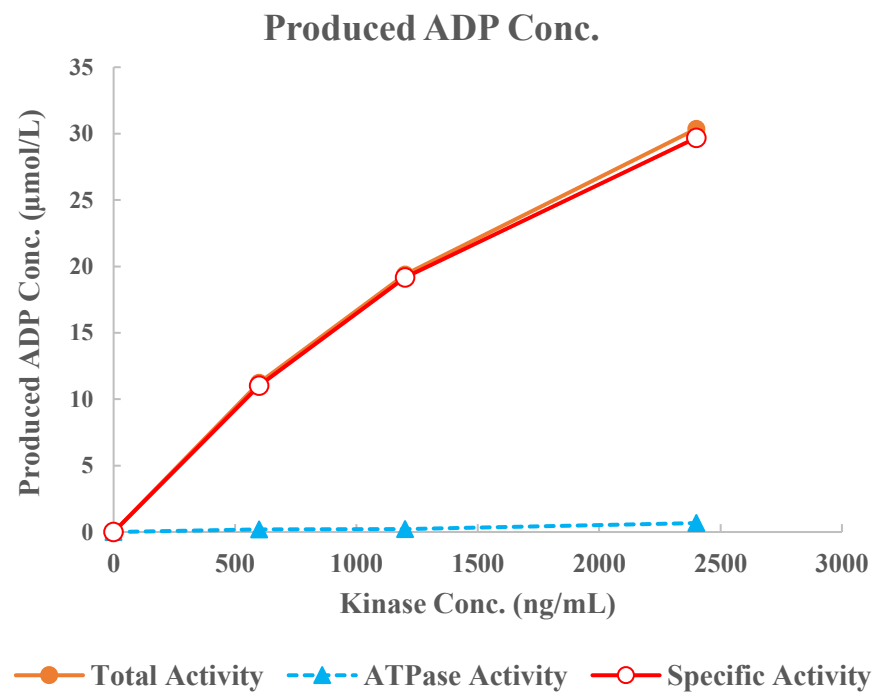
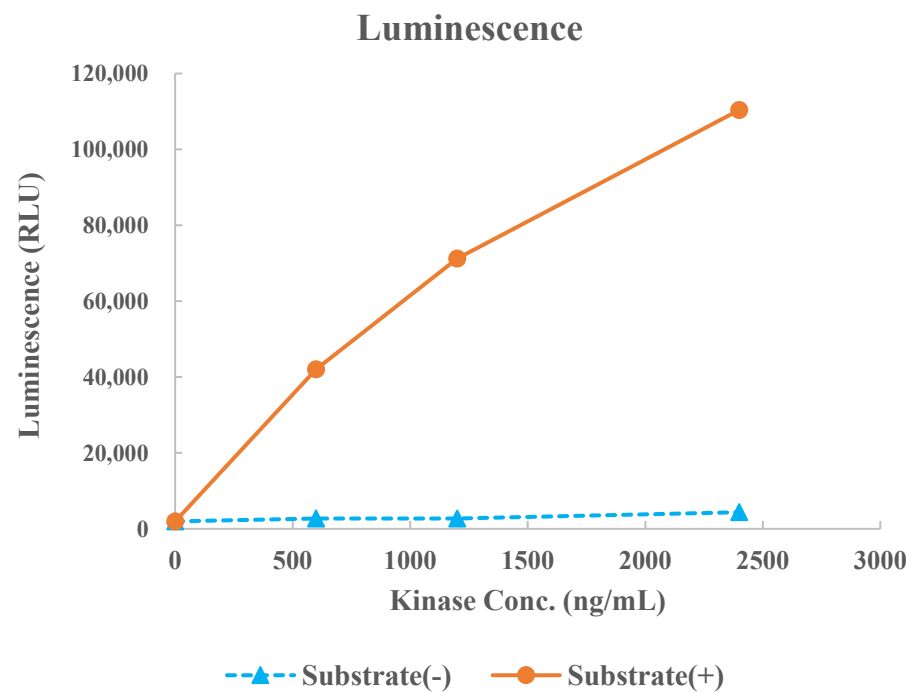


QIK(SNF1LK2)



Kinase	QIK(SNF1LK2), 02-129
Substrate	Biotin-AMARA peptide, BTSU28
Substrate Conc.	50 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour