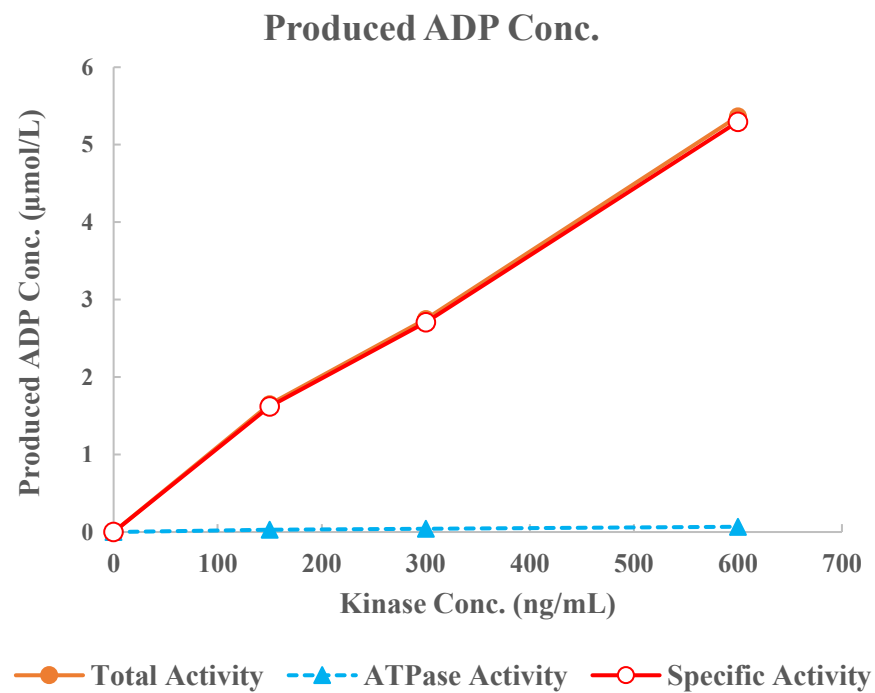
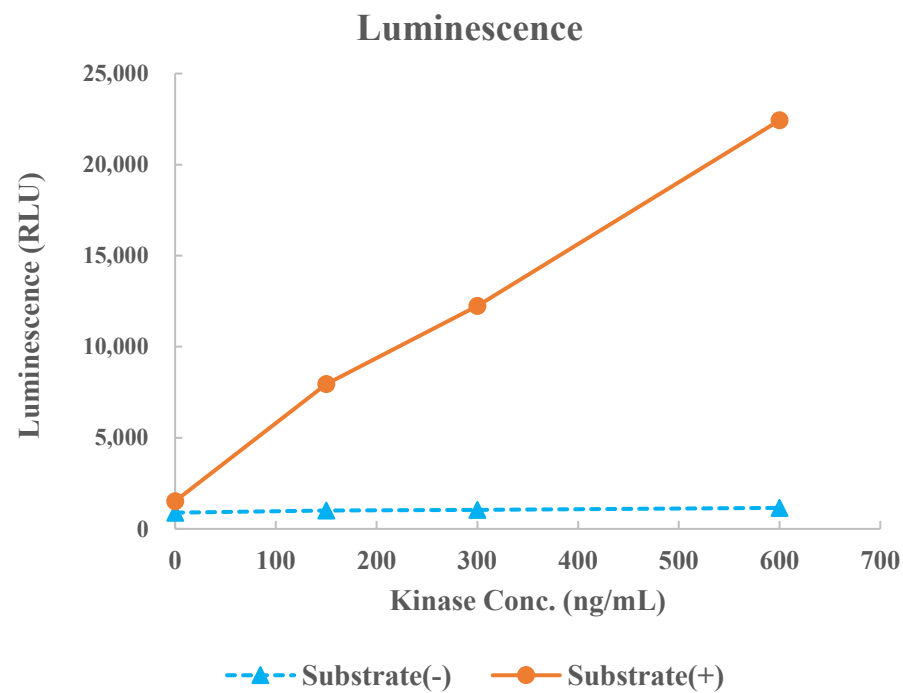


PRKX



Kinase	PRKX, 01-130
Substrate	Biotin-Kemptide, BTSU09
Substrate Conc.	50 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour