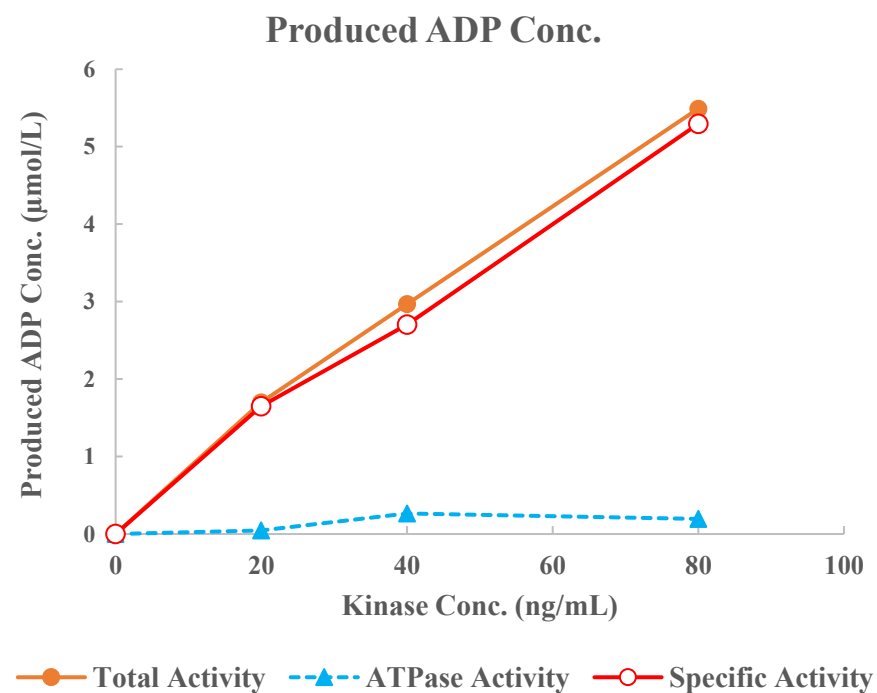
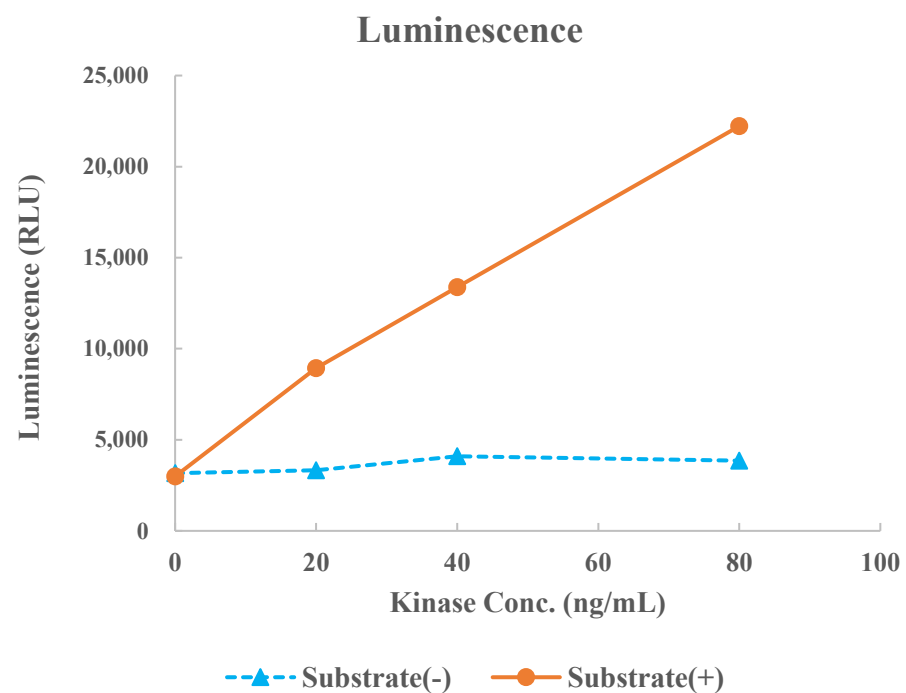


AKT2



Kinase	AKT2, 01-102
Substrate	Biotin-Crosstide, BTSU19
Substrate Conc.	10 μmol/L
ATP Conc.	100 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour