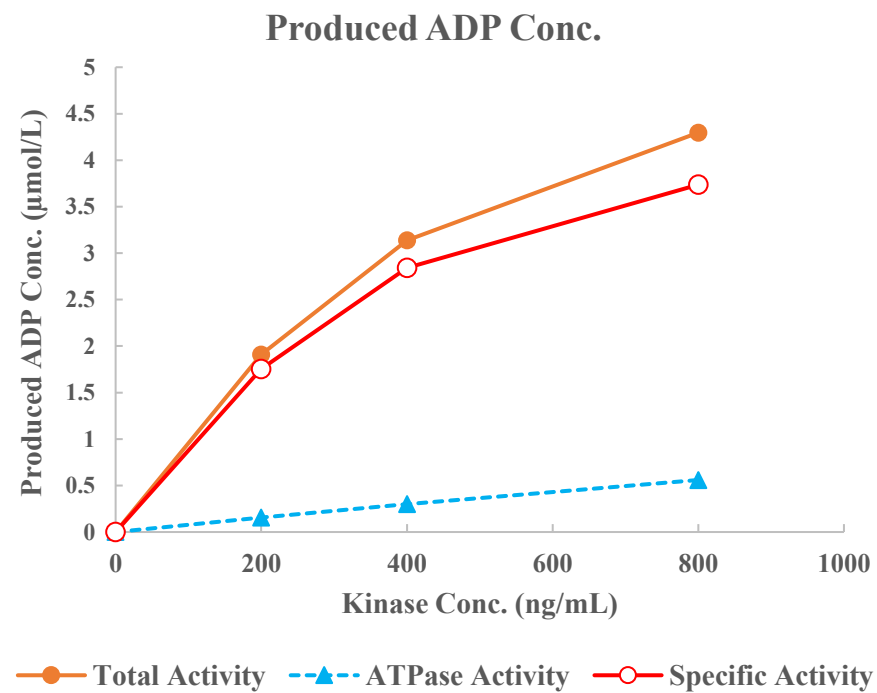
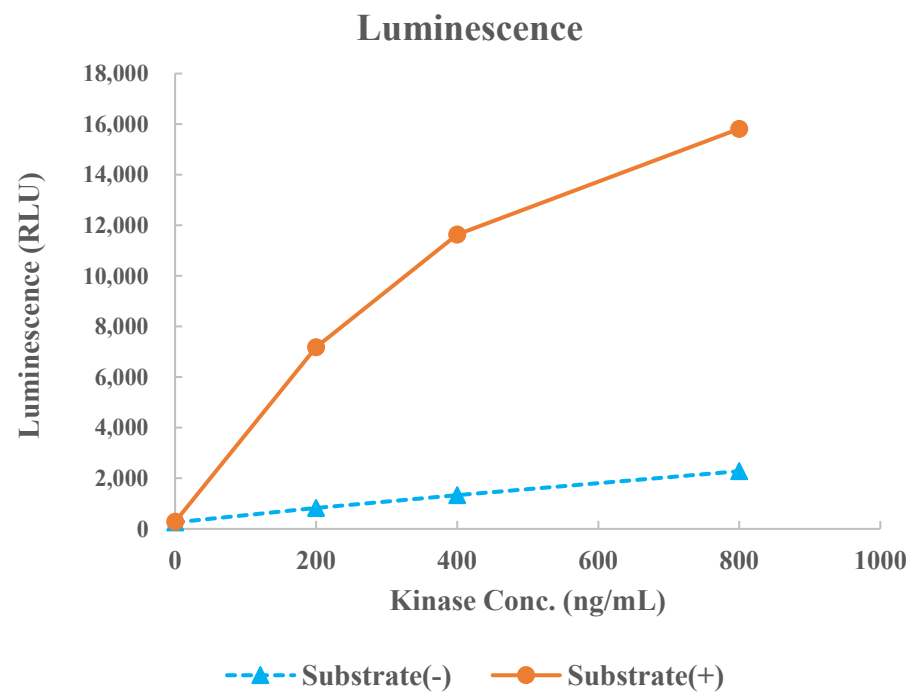


ROCK1



Kinase	ROCK1, 01-109
Substrate	Biotin-LIMKtide, BTSU24
Substrate Conc.	10 μmol/L
ATP Conc.	5 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour