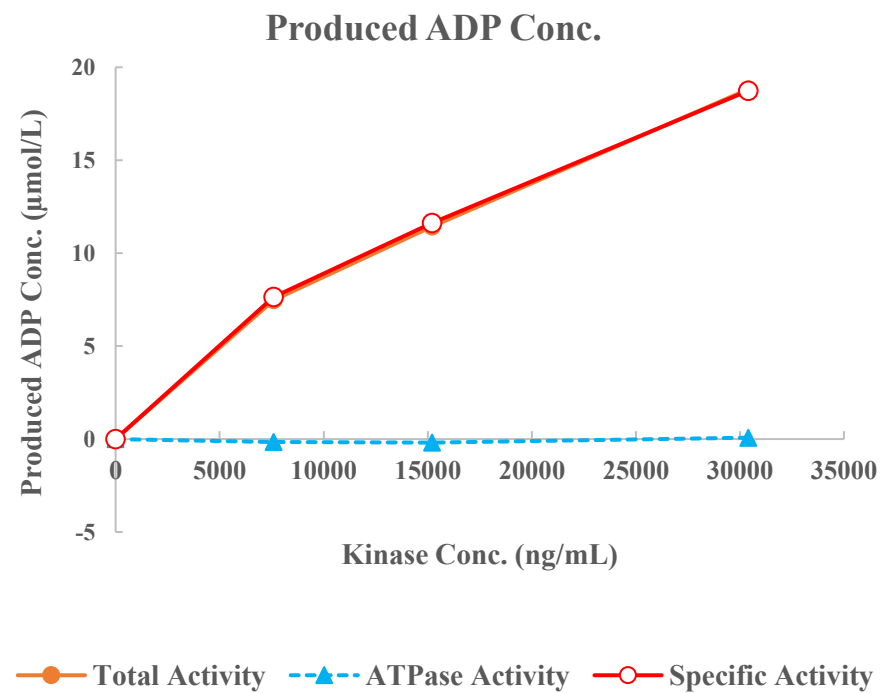
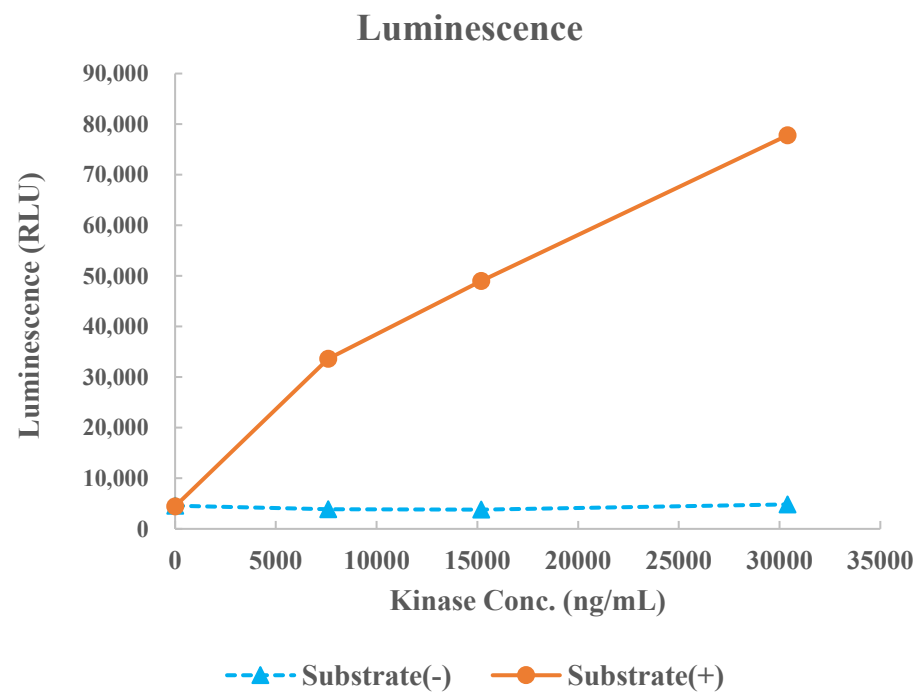


DDR1



Kinase	DDR1, 08-113
Substrate	Biotin-IRS1, BTSU07
Substrate Conc.	50 μmol/L
ATP Conc.	100 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour