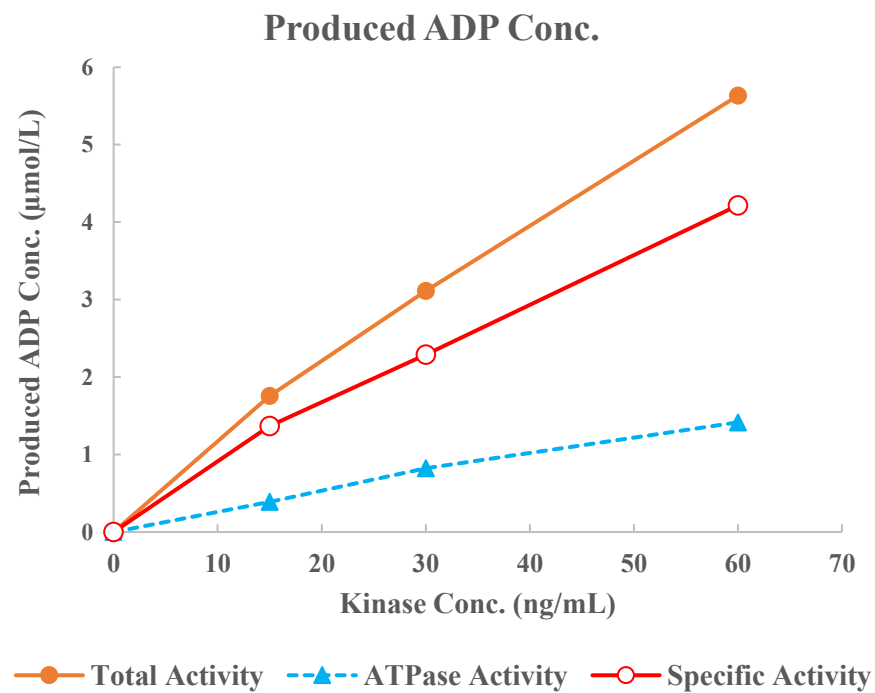
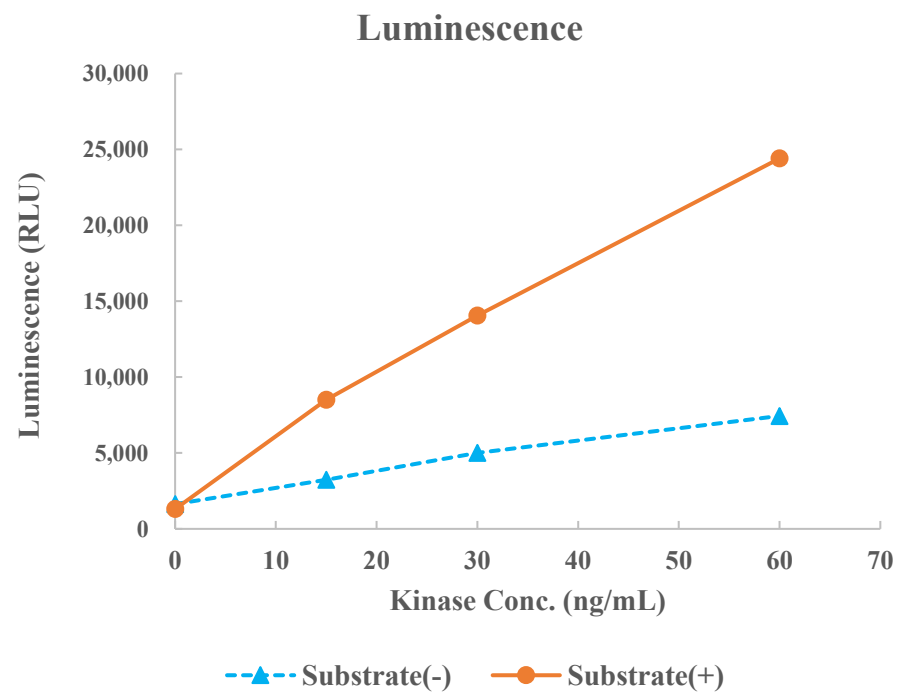


INSR



Kinase	INSR, 08-142
Substrate	Biotin-IRS1, BTSU07
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour