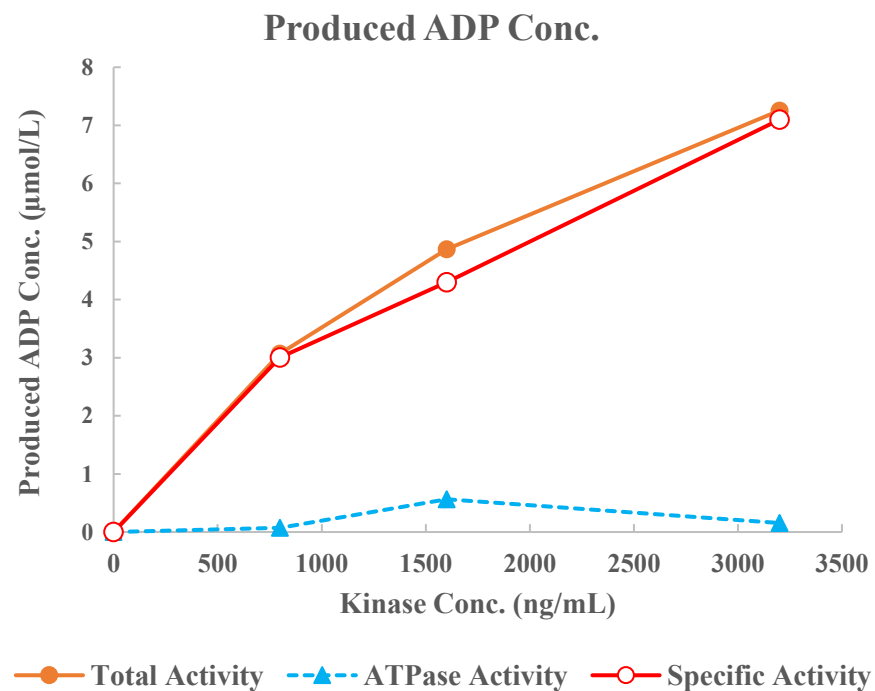
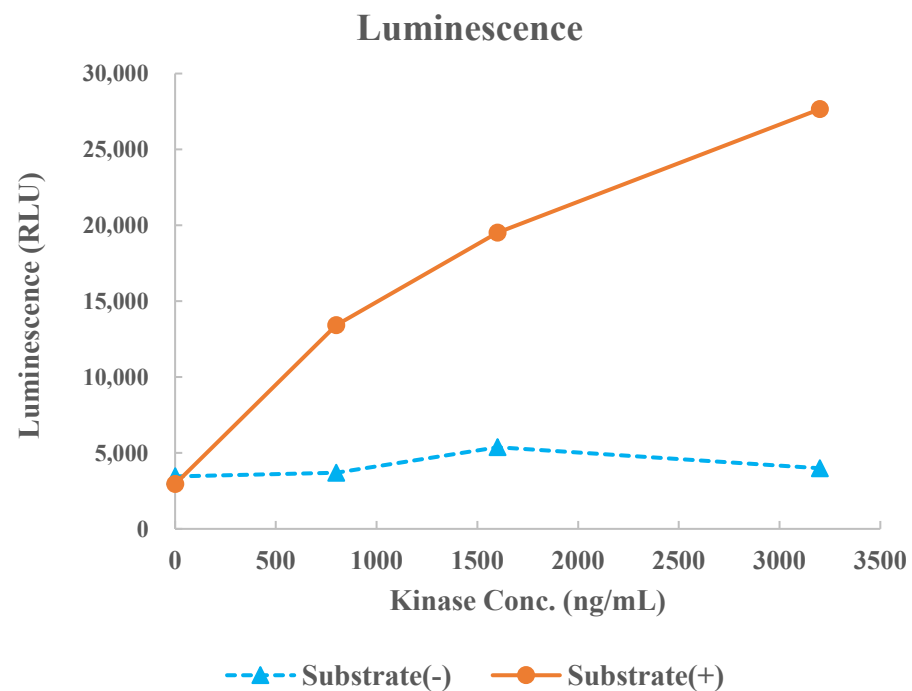


AMPK α 2/ β 1/ γ 1(PRKAA2/B1/G1)



Kinase	AMPK α 2/ β 1/ γ 1(PRKAA2/B1/G1), 02-114
Substrate	Biotin-SAMS peptide, BTSU26
Substrate Conc.	10 μ mol/L
ATP Conc.	100 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour