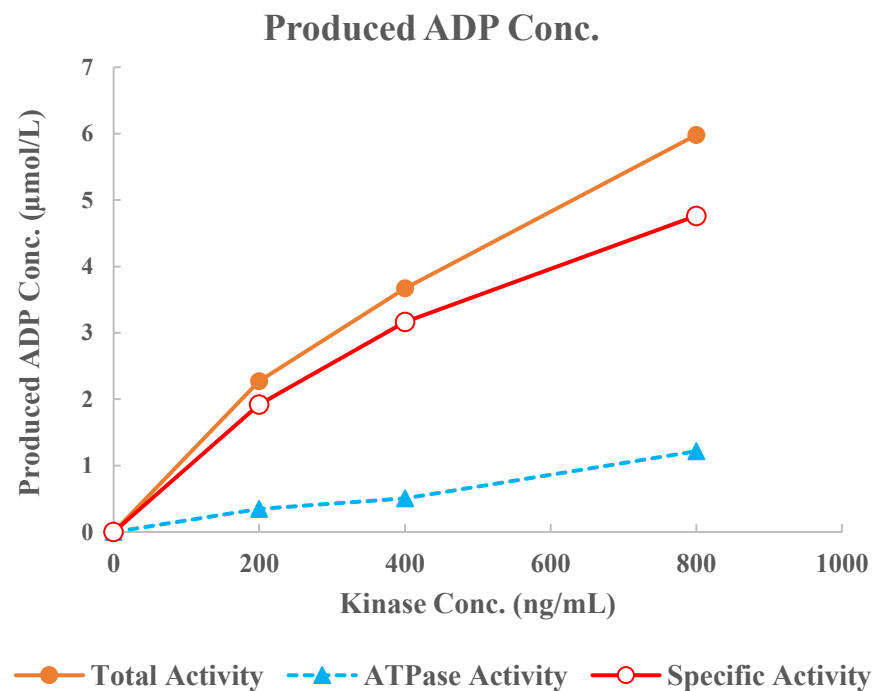
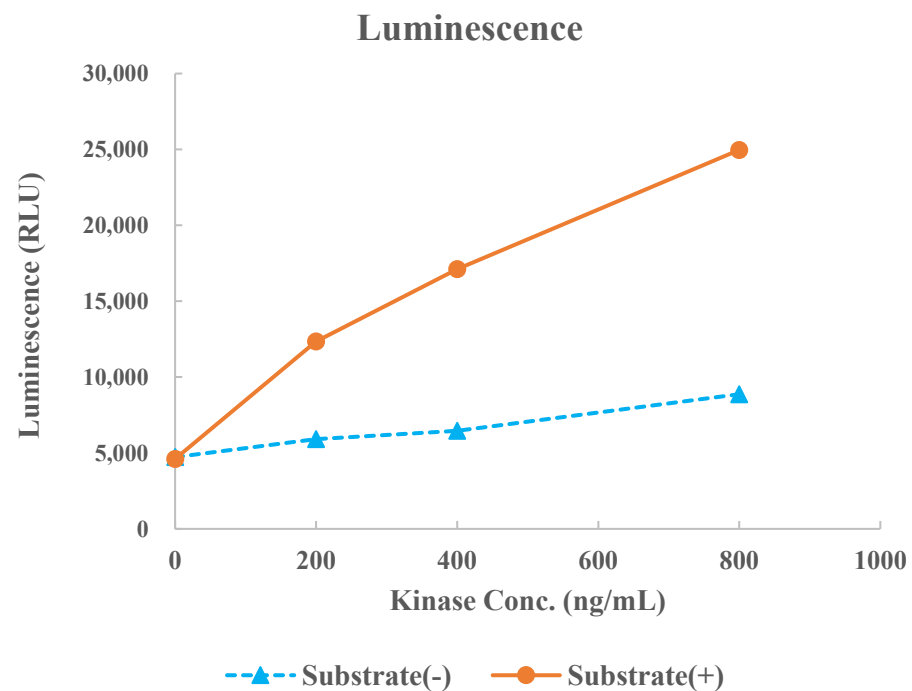


AMPK α 1/ β 1/ γ 1(PRKAA1/B1/G1)



Kinase	AMPK α 1/ β 1/ γ 1(PRKAA1/B1/G1), 02-113
Substrate	Biotin-SAMS peptide, BTSU26
Substrate Conc.	10 μ mol/L
ATP Conc.	150 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour