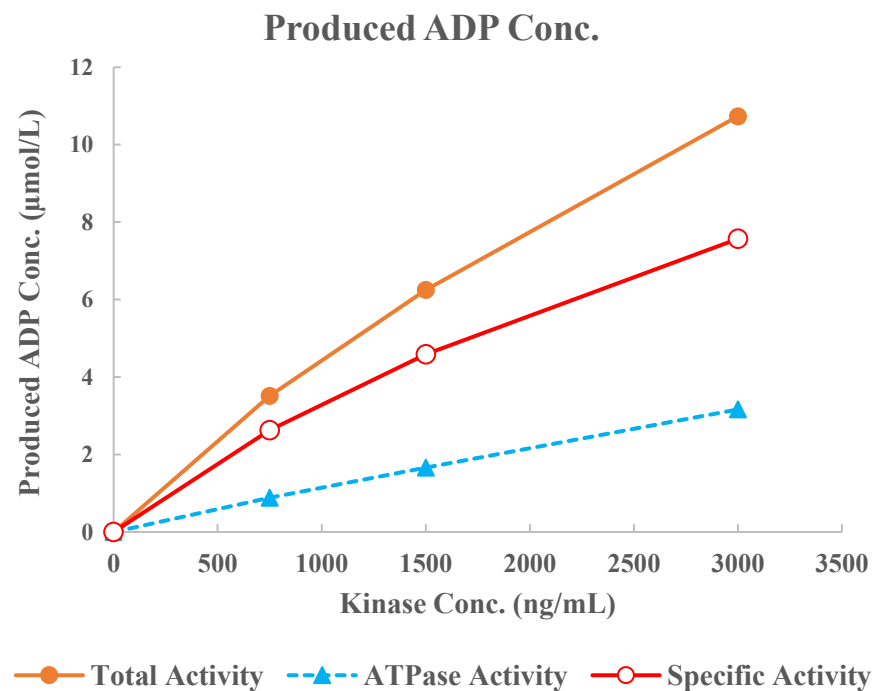
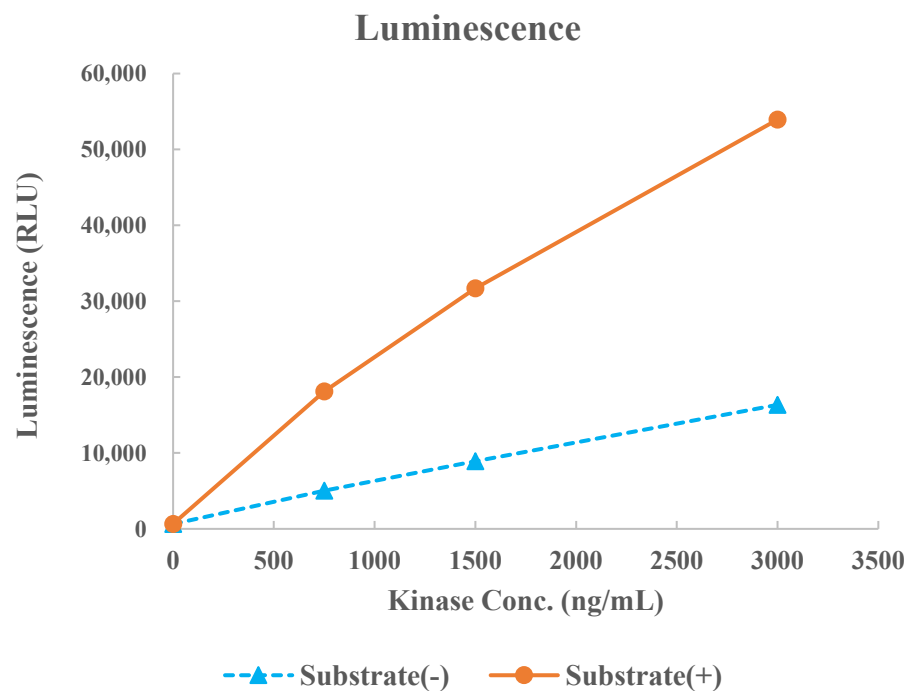


CK2 α 1/ β (CSNK2A1/B)



Kinase	CK2 α 1/ β (CSNK2A1/B), 05-184
Substrate	Biotin-CK2tide, BTSU25
Substrate Conc.	50 μ mol/L
ATP Conc.	10 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour