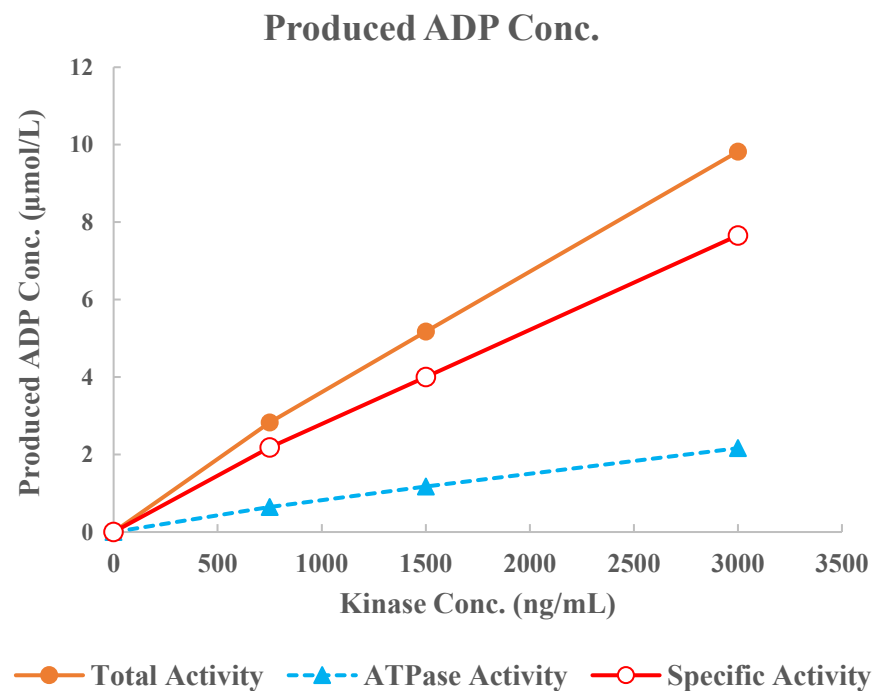
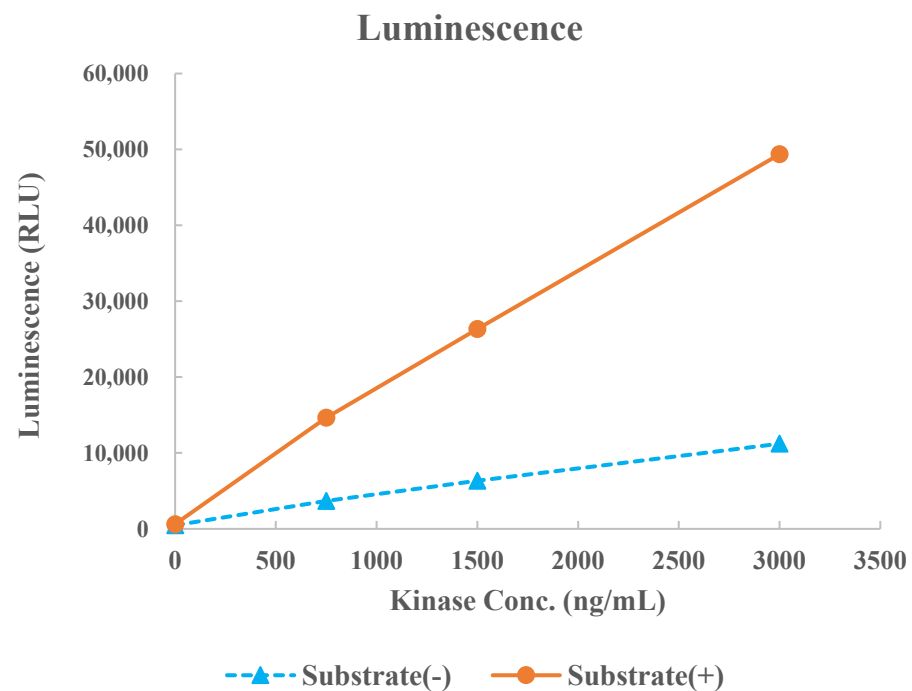


CK2 α 2/ β (CSNK2A2/B)



Kinase	CK2 α 2/ β (CSNK2A2/B), 05-185
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