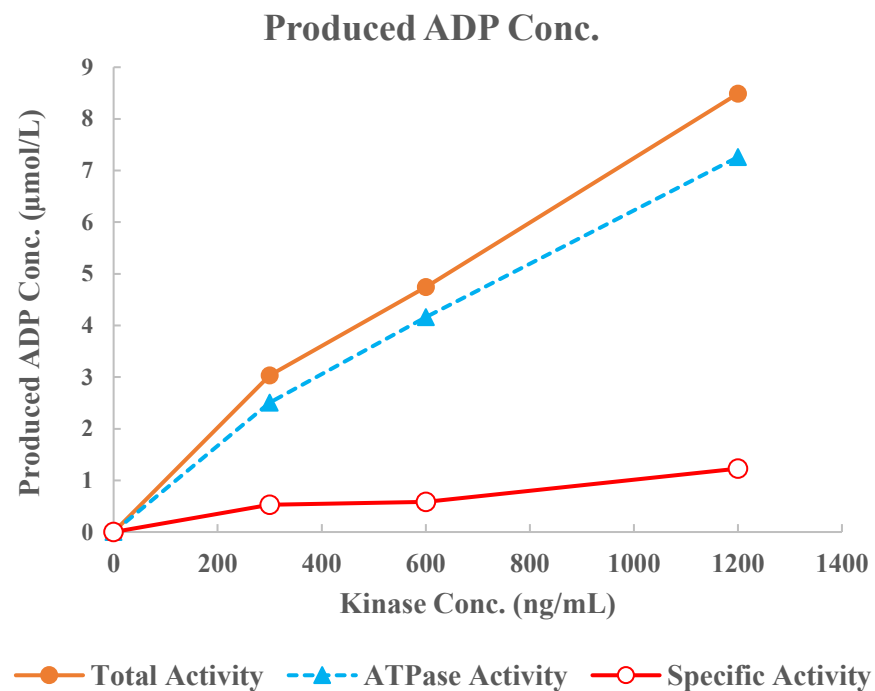
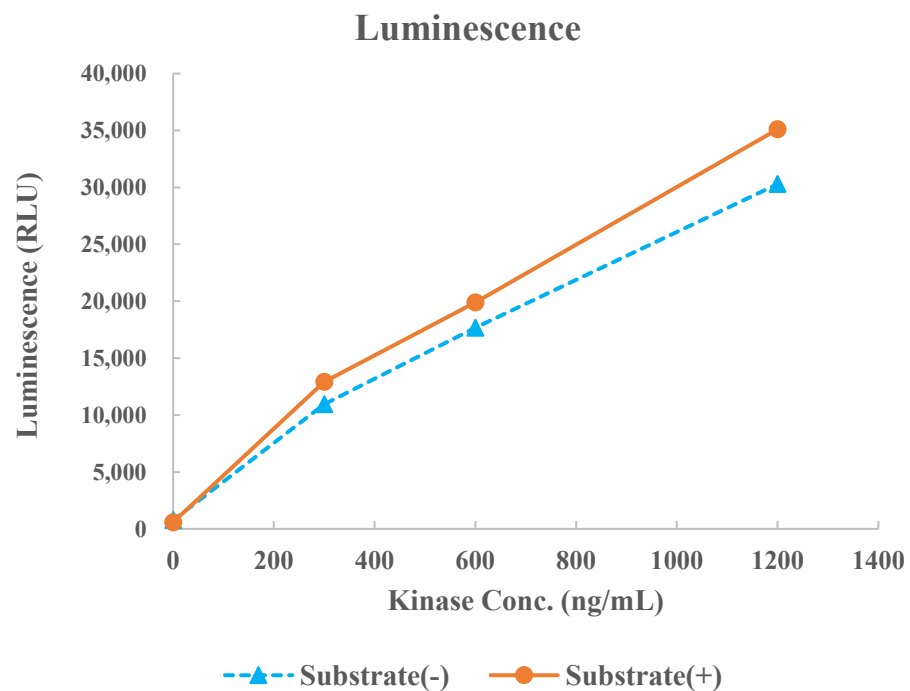


IKK β (IKBKB)



Kinase	IKK β (IKBKB), 05-084
Substrate	Biotin-Modified I κ B α -derived peptide, BTSU34
Substrate Conc.	50 μ mol/L
ATP Conc.	25 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour