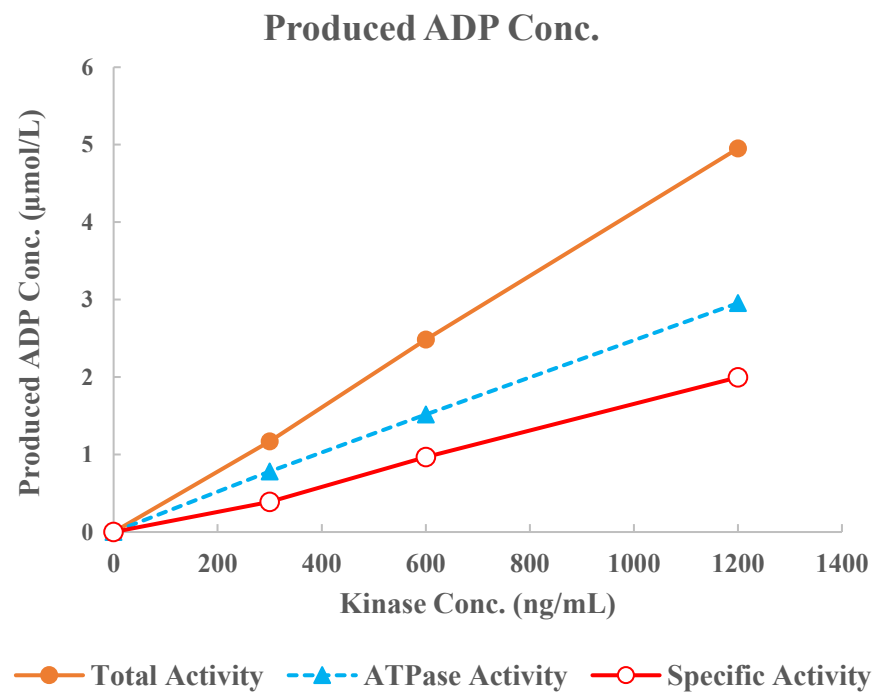
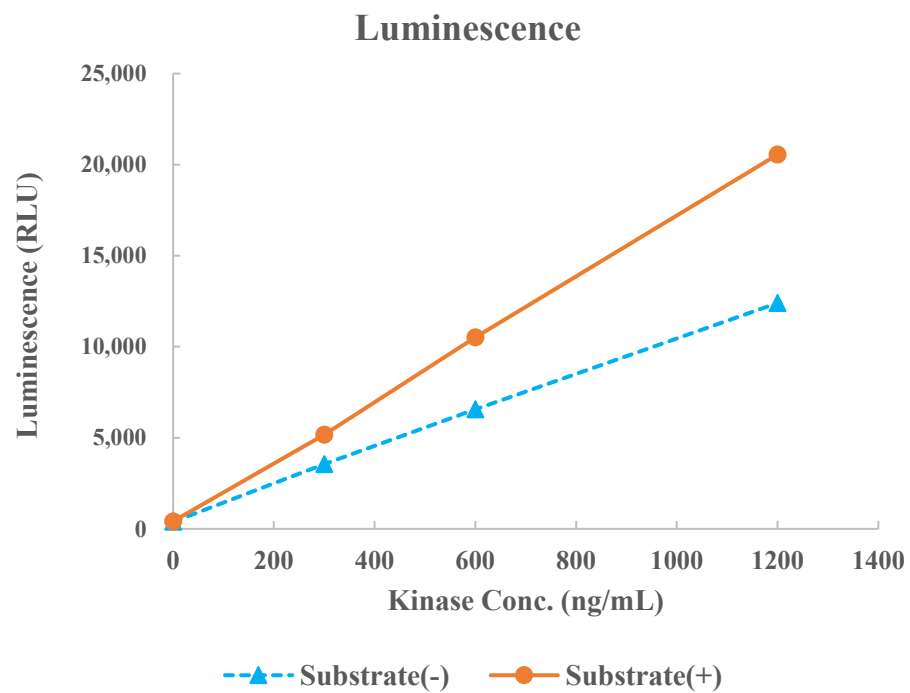


IKK ϵ (IKBKE)



Kinase	IKK ϵ (IKBKE), 05-114
Substrate	Biotin-I κ B α peptide, BTSU42
Substrate Conc.	50 μ mol/L
ATP Conc.	10 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour