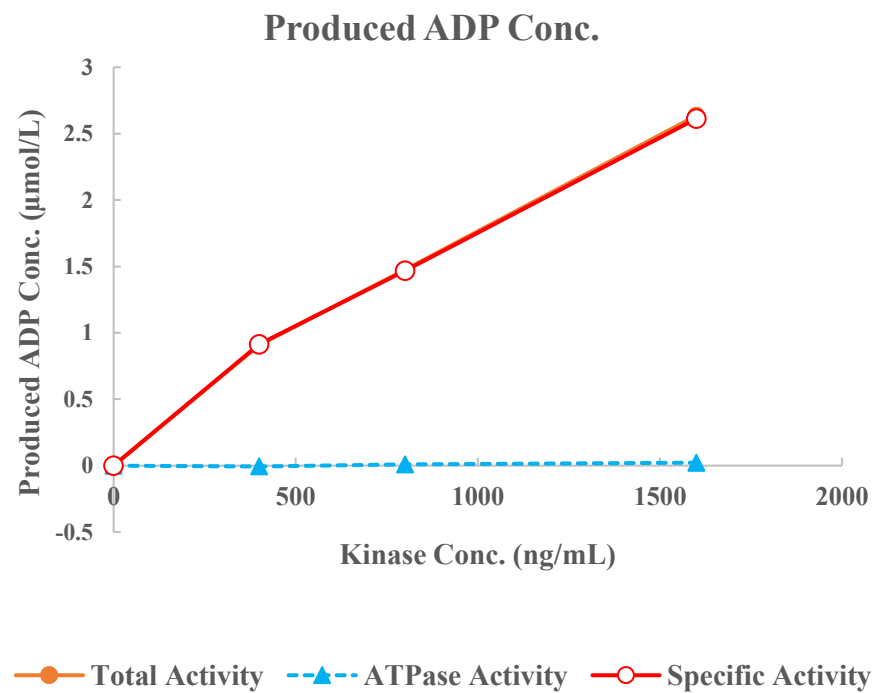
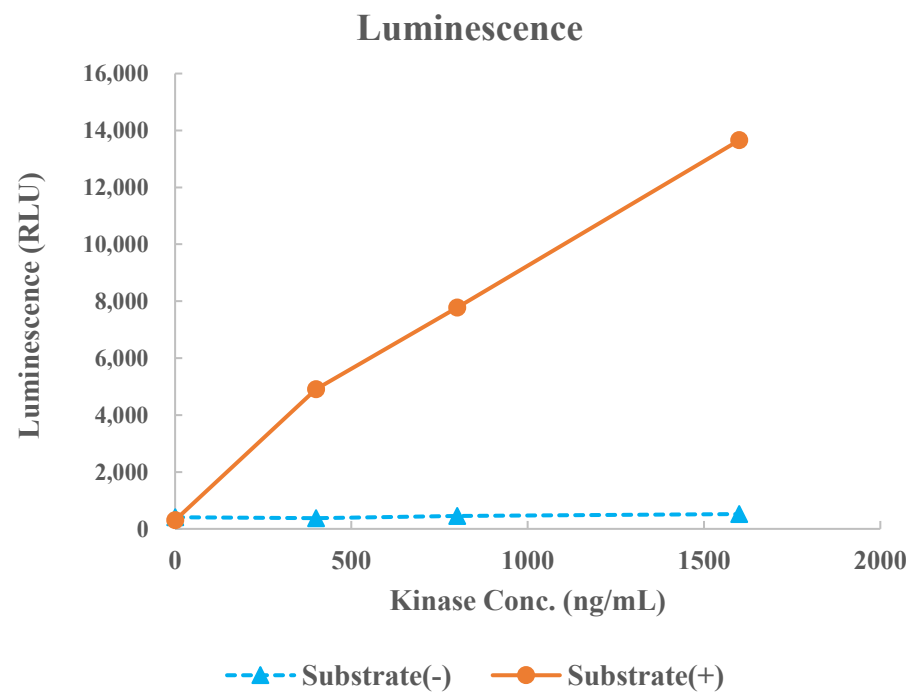


CSK



Kinase	CSK, 08-111
Substrate	Biotin-Srcptide, BTSU01
Substrate Conc.	50 μmol/L
ATP Conc.	5 μmol/L
Metal	Mg, Mn
Metal Conc.	5, 1 mmol/L
Additives	-
Reaction Time	1 hour