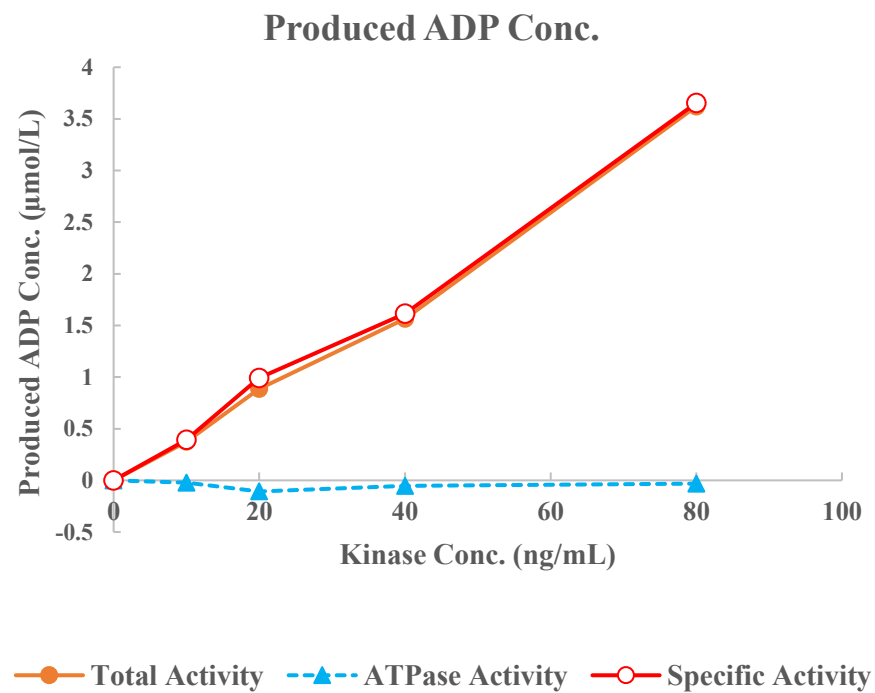
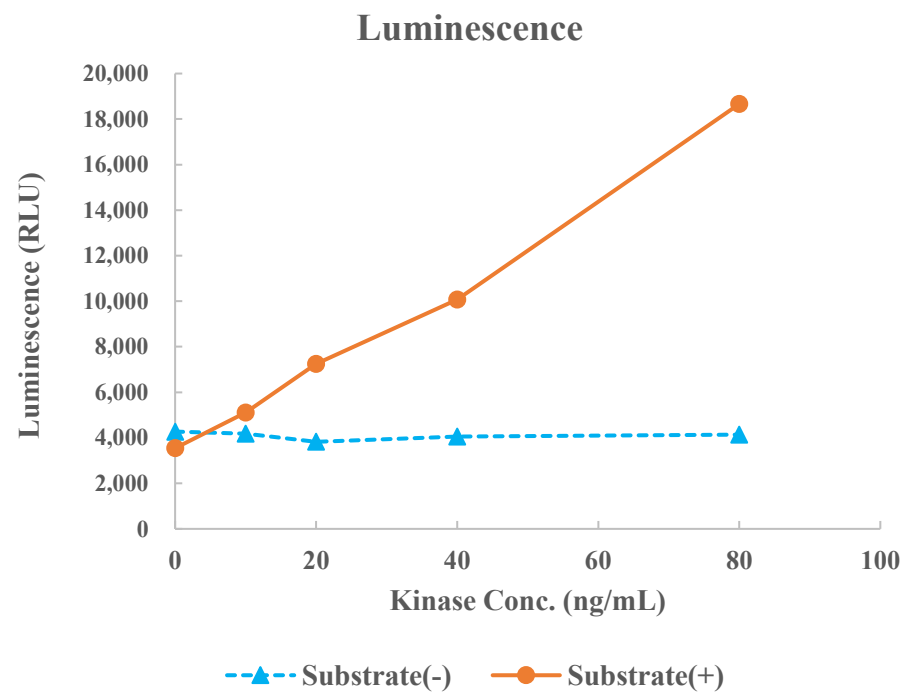


Haspin(GSG2)



Kinase	Haspin(GSG2), 05-111
Substrate	Biotin-Histone H3 peptide, BTSU23
Substrate Conc.	10 μmol/L
ATP Conc.	150 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour