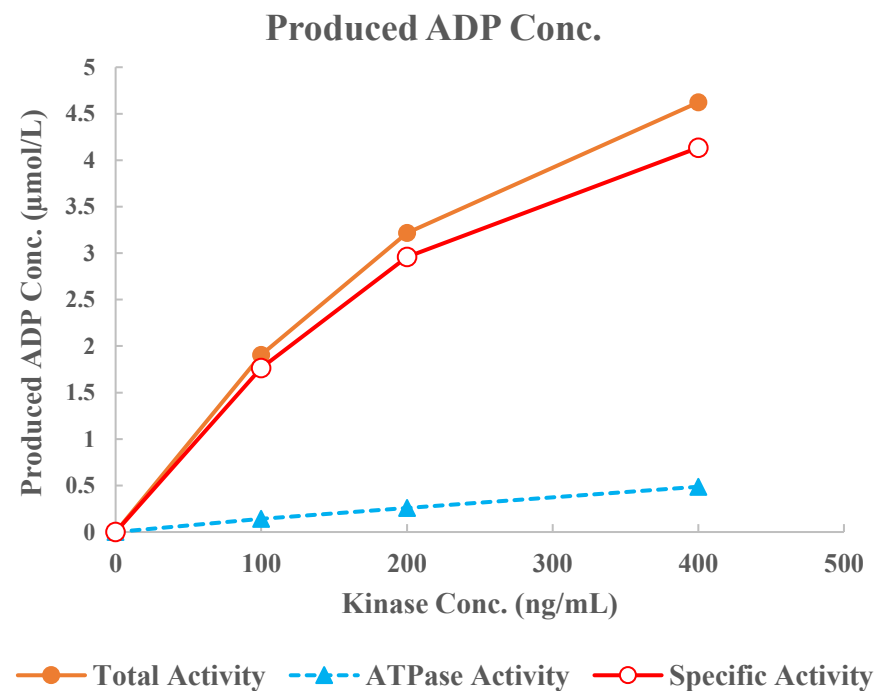
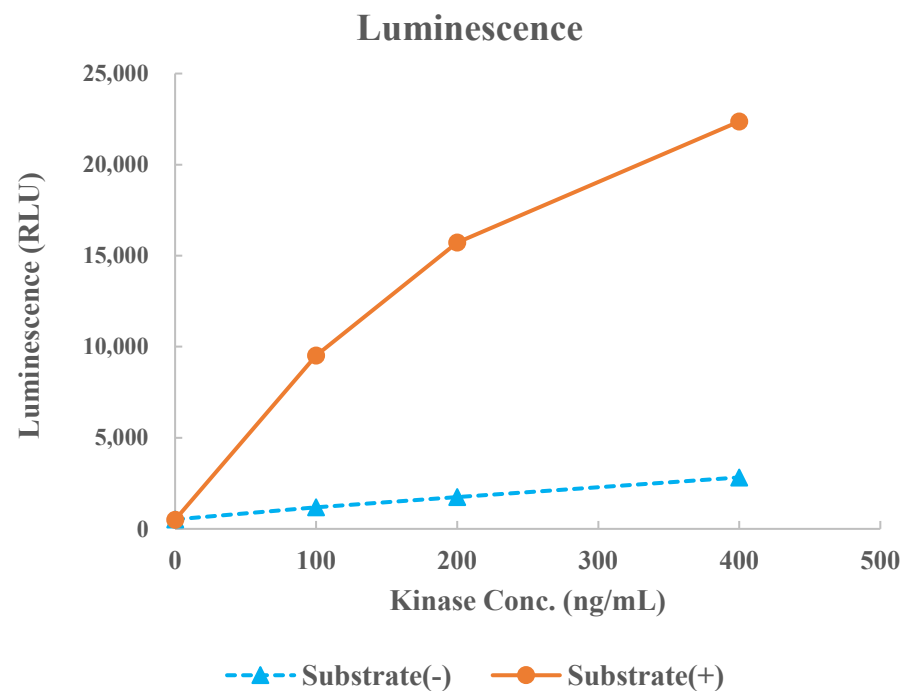


RET [S891A]



Kinase	RET [S891A], 08-523
Substrate	Biotin-CSKtide, BTSU02
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour