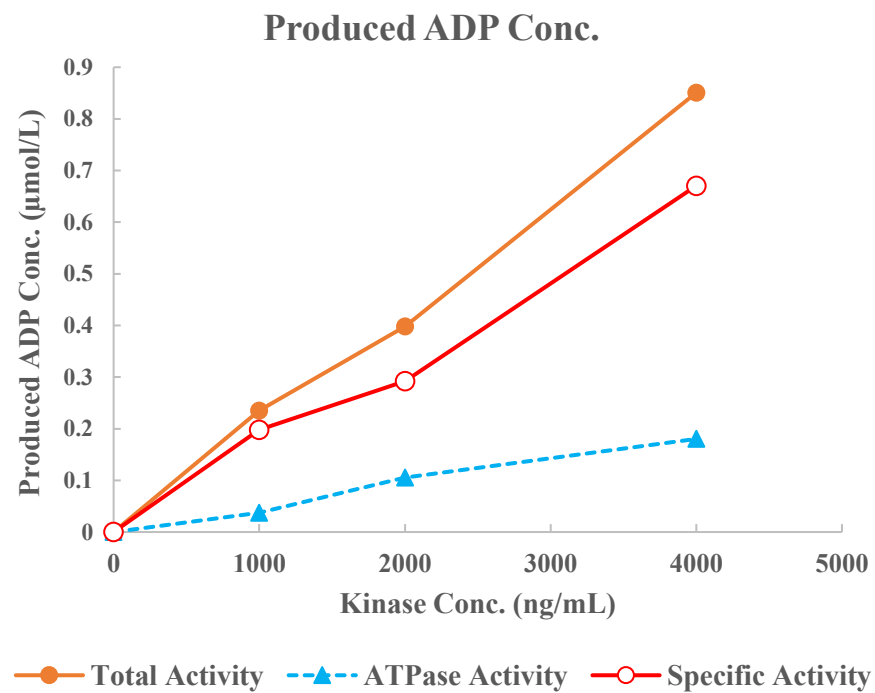
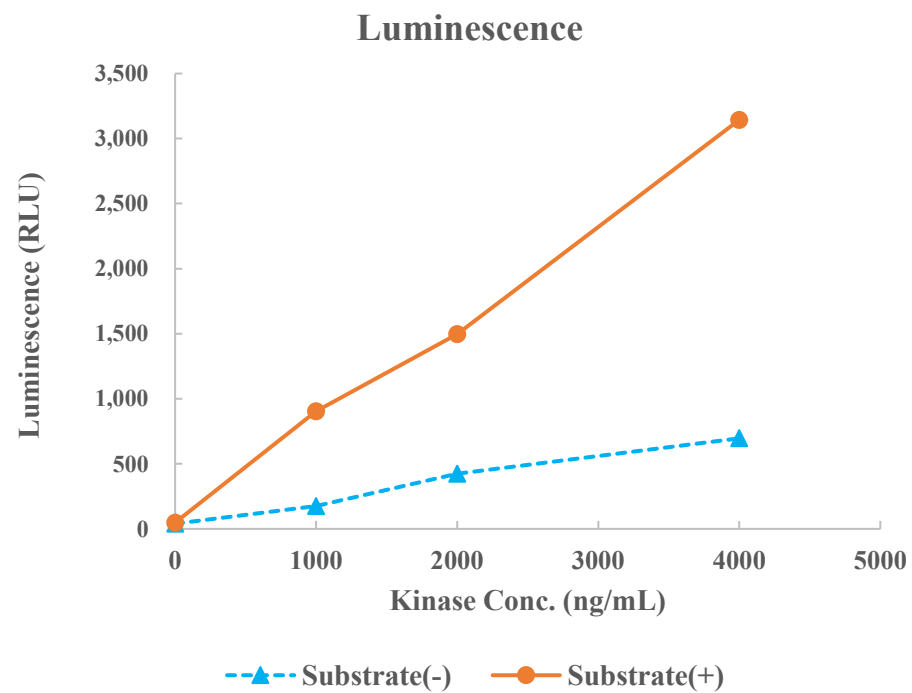


HER2(ERBB2)



Kinase	HER2(ERBB2), 08-016
Substrate	Biotin-Srcptide, BTSU01
Substrate Conc.	10 μmol/L
ATP Conc.	5 μmol/L
Metal	Mn
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour