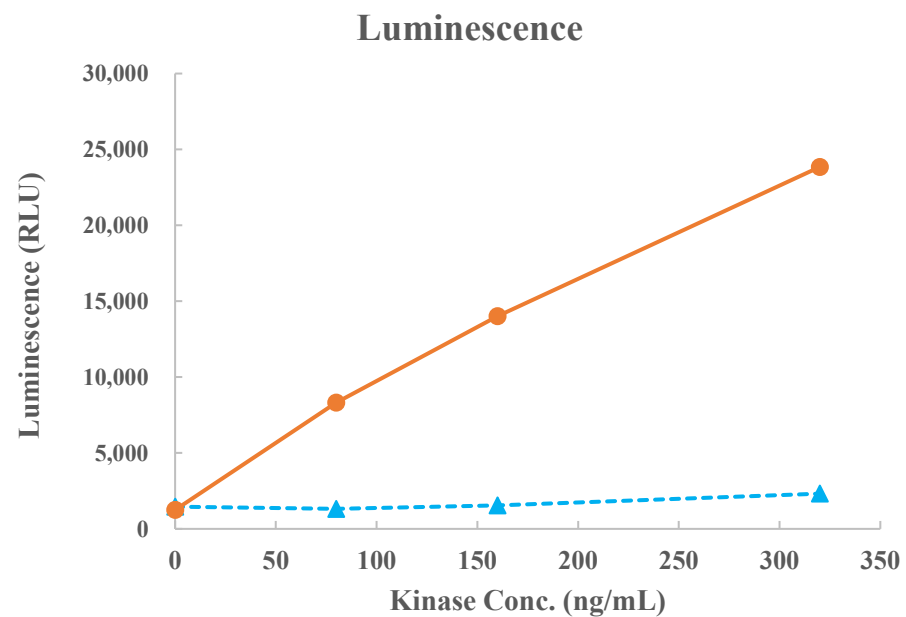
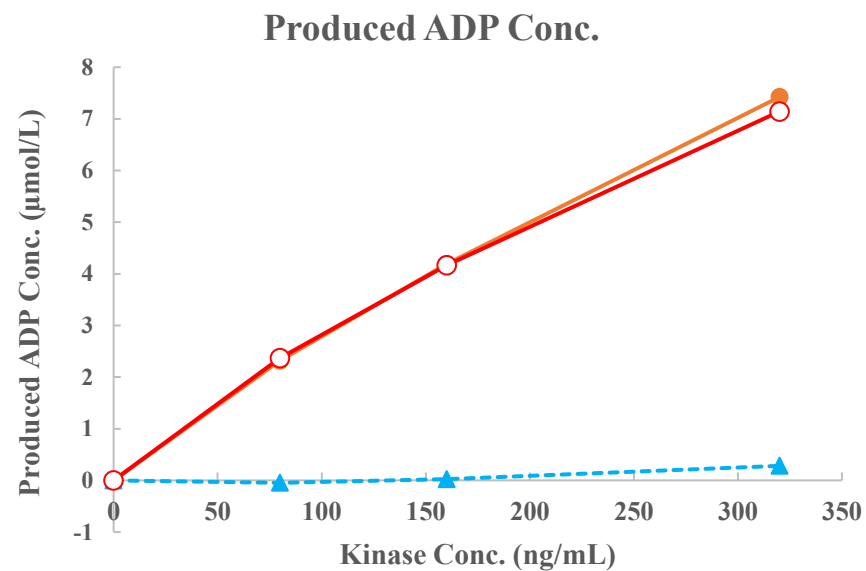


HER4(ERBB4)



--▲-- Substrate(-) —●— Substrate(+)



—●— Total Activity --▲-- ATPase Activity —○— Specific Activity

Kinase	HER4(ERBB4), 08-118
Substrate	Biotin-Srcptide, BTSU01
Substrate Conc.	10 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour