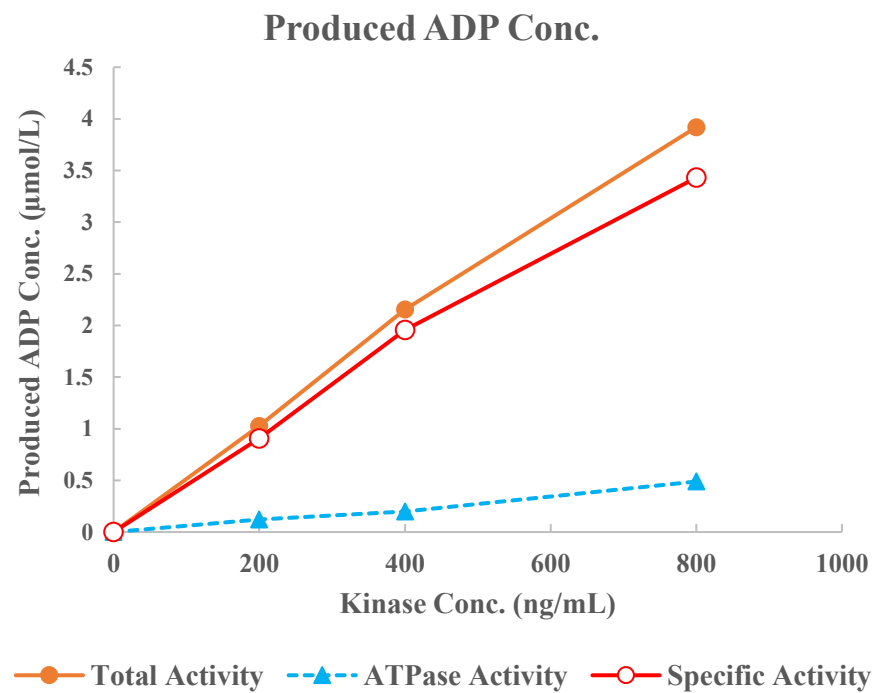
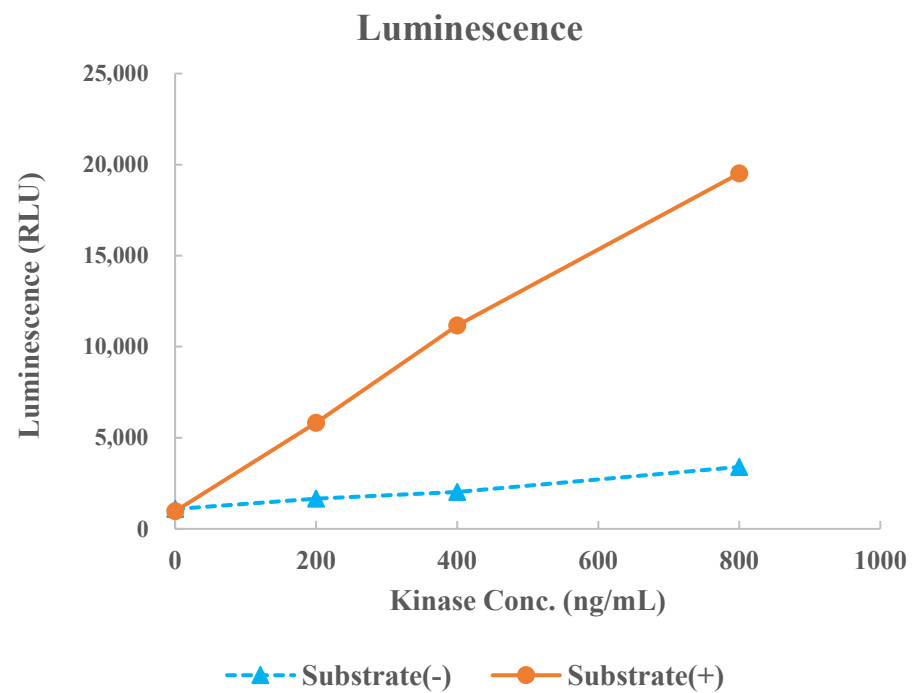


PDGFR α (PDGFRA)



Kinase	PDGFR α (PDGFRA), 08-157
Substrate	Biotin-CSKtide, BTSU02
Substrate Conc.	10 μ mol/L
ATP Conc.	25 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour